

Journal Pre-proof



16pdel lipid changes in iPSC-derived neurons and function of FAM57B in lipid metabolism and synaptogenesis

Danielle L. Tomasello, Jiyeon L. Kim, Yara Khodour, Jasmine M. McCammon, Maya Mitalipova, Rudolf Jaenisch, Anthony H. Futerman, Hazel Sive

PII: S2589-0042(21)01521-2

DOI: <https://doi.org/10.1016/j.isci.2021.103551>

Reference: ISCI 103551

To appear in: *ISCIENCE*

Received Date: 1 April 2021

Revised Date: 23 September 2021

Accepted Date: 26 November 2021

Please cite this article as: Tomasello, D.L., Kim, J.L., Khodour, Y., McCammon, J.M., Mitalipova, M., Jaenisch, R., Futerman, A.H., Sive, H., 16pdel lipid changes in iPSC-derived neurons and function of FAM57B in lipid metabolism and synaptogenesis, *ISCIENCE* (2022), doi: <https://doi.org/10.1016/j.isci.2021.103551>.

This is a PDF file of an article that has undergone enhancements after acceptance, such as the addition of a cover page and metadata, and formatting for readability, but it is not yet the definitive version of record. This version will undergo additional copyediting, typesetting and review before it is published in its final form, but we are providing this version to give early visibility of the article. Please note that, during the production process, errors may be discovered which could affect the content, and all legal disclaimers that apply to the journal pertain.

© 2021

16pdel lipid changes in iPSC-derived neurons and function of FAM57B in lipid metabolism and synaptogenesis

Danielle L. Tomasello¹, Jiyeon L. Kim², Yara Khodour², Jasmine M. McCammon¹, Maya Mitalipova¹, Rudolf Jaenisch^{1,3}, Anthony H. Futerman² and Hazel Sive^{1,3*#}

¹Whitehead Institute for Biomedical Research, Cambridge, MA 02142, USA.

²Department of Biomolecular Sciences, Weizmann Institute of Science, Rehovot, 76100, Israel.

³Department of Biology, Massachusetts Institute of Technology, Cambridge, MA 02139, USA.

*Present address: Northeastern University College of Science, Boston, MA 02115, USA.

#To whom lead contact should be addressed at: 115 Richards Hall, 360 Huntington Ave, Boston MA 02115, USA. Tel: (617) 373-5085; Email: h.sive@northeastern.edu

Summary

The complex 16p11.2 Deletion Syndrome (16pdel) is accompanied by neurological disorders, including epilepsy, autism spectrum disorder and intellectual disability. We demonstrated that 16pdel iPSC differentiated neurons from affected people show augmented local field potential activity and altered ceramide-related lipid species relative to unaffected. *FAM57B*, a poorly characterized gene in the 16p11.2 interval, has emerged as a candidate tied to symptomatology. We found that *FAM57B* modulates ceramide synthase (CerS) activity, but is not a CerS per se. In *FAM57B* mutant human neuronal cells and zebrafish brain, composition and levels of sphingolipids and glycerolipids associated with cellular membranes are disrupted. Consistently, we observed aberrant plasma membrane architecture and synaptic protein mislocalization, which were accompanied by depressed brain and behavioral activity. Together, these results suggest that haploinsufficiency of *FAM57B* contributes to changes in neuronal activity and function in 16pdel Syndrome, through a crucial role for the gene in lipid metabolism.

Introduction

16p11.2 Deletion (16pdel) Syndrome, a severe and prevalent neurodevelopmental disorder, is a copy number variant with deletion of ~600 kb from chromosome 16, encompassing 25 core protein-coding genes. This haploinsufficiency syndrome is estimated to affect ~1 in 2500 worldwide and is tightly associated with autism spectrum disorder (ASD), language and intellectual disability, seizures, attention-deficit/hyperactivity disorder, macrocephaly, hypotonia and obesity (D'Angelo et al., 2016, Hanson et al., 2015, Zufferey et al., 2012, Egolf et al., 2019, Maillard et al., 2015). Strong indications of synaptic defects are associated with 16pdel symptoms, particularly epilepsy (Fetit et al., 2020, Kleinendorst et al., 2020) and ASD (Fetit et al., 2020, Kim et al., 2020, Maillard et al., 2015, Ouellette et al., 2020, Sebat et al., 2007, Pinto et al., 2010, Zufferey et al., 2012), as well as links to metabolic defects (Hoytema van Konijnenburg et al., 2020).

37 Previously, analysis in the zebrafish model suggested that *FAM57B* is a pivotal hub gene in the
38 16p11.2 interval, that encodes a protein proposed to be a ceramide synthase (Yamashita-Sugahara et al.,
39 2013). Using a pairwise partial loss of function screen for zebrafish embryonic brain morphology, we found that
40 *fam57b* interacted with numerous other 16p11.2 interval genes, suggesting haploinsufficiency of *FAM57B* is
41 critical in 16pdel Syndrome etiology (McCammon et al., 2017). *FAM57B* (family with sequence similarity 57,
42 member B) is a Tram-Lag-CLN8 (TLC) family member, containing a domain of roughly 200 amino acids found
43 in several other proteins, including ceramide synthases (CerS, the Lag of TLC) (Pewzner-Jung et al., 2006).
44 Ceramides are sphingolipids (SLs) which are key membrane components and also act as signaling molecules
45 to modulate proliferation, apoptosis, inflammation, cell cycle arrest and ER stress (Grosch et al., 2012). In
46 humans, mutations in some of the 6 known CerS are associated with autism, epilepsy and intellectual disability
47 (Vanni et al., 2014, Egawa et al., 2015, Ranta et al., 1999). In this study, to further assess the predicted
48 connection with 16pdel Syndrome, we examined *FAM57B* function through a multidisciplinary approach,
49 across human cells and the zebrafish system.

50

51 Results

52 Augmented network activity in 16pdel neuron cultures

53 Based on previous data, we hypothesized that 16pdel neurons would show an altered lipid profile due
54 to contributions of *FAM57B* and possibly other 16pdel genes with predicted roles in metabolism (McCammon
55 et al., 2017). To test this, we prepared neurons from 16pdel carrier induced-pluripotent stem cells (iPSC), part
56 of the Simons VIP Consortium, and unaffected control iPSC in culture (Simons Vip, 2012) (**Supp Table 1**).
57 Neural progenitor cells were differentiated into cortical neurons, since the cortex has consistently shown
58 anatomical differences in 16pdel affected individuals (Martin-Brevet et al., 2018, Maillard et al., 2015, Hinkley
59 et al., 2019, Lin et al., 2015, Qureshi et al., 2014, Blackmon et al., 2018). After one month in culture,
60 immunocytochemistry (ICC) indicated mature neurons by presence of vesicular glutamate 1 and 2 receptors
61 (VGlut1/2) (**Supp Figs. 1a,b**), the synaptic markers PSD95 (**Supp Fig. 1b**) and Synaptotagmin-1 (Sy1).
62 Cultures of control and 16pdel (proband) differentiated neurons showed similar percentages of mature neurons
63 by these criteria. Quantitative analysis was performed to determine equivalent maturation between control and
64 16pdel differentiated neurons. Production of synaptic proteins was measured by colocalization analysis,
65 indicating neuronal somas that were Synaptotagmin-1 positive additionally colocalized with PSD-95 (**Supp**
66 **Table 2**).

67 To further characterize these neurons, we probed network electrical activity by multi-electrode array
68 (MEA). Spontaneous activity of differentiated neurons was measured over thirty minutes; first in culture media,
69 followed by physiological solution and last in high potassium chloride solution. Comparing grouped genotypes,
70 we recorded an increased frequency of Local Field Potentials (LFPs) in 16pdel proband neuron cultures
71 relative to controls, indicating 16pdel neurons display heightened spontaneous and evoked activity compared
72 to unaffected control (**Fig. 1a**). Examining individual patient cell lines, we observed relatively similar MEA

73 activity in controls (black), and increased electrical activity in 16pdel neuron cultures (grey) (**Supp Fig. 2a**).
74 Interestingly, female 16pdel neuron cultures showed statistically increased LFP frequency compared to male
75 16pdel neuron cultures when measured in media and High KCl solution (**Fig. 1b**). Sex differences were also
76 observed in LFP firing and bursting properties, with increased burst frequency of female 16pdel neurons
77 compared to male in media (**Fig. 1c**). While behavioral deficits have been widely observed among male mouse
78 16pdel models, a recent report found stress-related sex differences in a female 16p11.2 deletion mouse model.
79 Thus, Giovanniello *et al.* discovered an increase in activity in central amygdala neurons projecting to the
80 globus pallidus in female, but not male, 16pdel model mice (Giovanniello, 2021). Our data consistently suggest
81 sex-specific differences between activity of 16pdel and control neurons. These findings expand previous
82 observations that demonstrated larger cell size and deficits in synaptic density in 16pdel neurons compared to
83 control (Deshpande *et al.*, 2017).

84

85 **Significant lipid class and individual species changes indicates complexity of 16pdel Syndrome.**

86 Using differentiated 16pdel and control neurons, we compared their lipid cohorts using untargeted
87 lipidomics (**Table S1 - iPSC Neuron Lipidomics**). Consistent with predictions, many significant changes were
88 observed in total lipid classes between 16pdel and control neurons (**Fig. 2a**). Levels of SLs (ceramide (Cer)
89 and sphingomyelin (SM)) and glycerolipids (GL) (lysophosphatidylethanolamine (LPE),
90 phosphatidylethanolamine (PE), monoacylglycerol (MG)) were significantly decreased, while GL (triacylglycerol
91 (TG)) levels significantly increased. Analyzing lipid composition, we found similar levels of unsaturated and
92 saturated species between 16pdel and control neurons, but differences in saturation of acyl carnitine (AcCa)
93 (unsaturated control 33.78% vs 16pdel 55.16%, and saturated control 66.22% and 16pdel 44.84%) and TG
94 (unsaturated control 71.15% vs 16pdel 80.28%, and saturated control 28.85% and 16pdel 19.72%) (**Fig. 2b**).
95 Polyunsaturated fatty acids (PUFAs) are important in the brain, where they are essential for signaling and
96 membrane structure (Bazinet and Laye, 2014). Chain length analysis indicated large differences in
97 lysophosphatidylglycerol (LPG) (long chain control 85.76% vs 16pdel 75.55%) and phosphatidylglycerol (PG)
98 (long chain control 71.78% vs 16pdel 51.88%). While having a similar ratio of long and very long chain PE
99 species (**Fig. 2b**), analysis of individual lipid species demonstrated significantly decreased levels of several
100 PE(18:22) species in 16pdel neurons relative to control (**Fig. 2c**). Additionally, decreased Cer(18) species were
101 observed in 16pdel (**Fig. 2d**). Comparing MG and TG, MG(18:0) decreased while many TG(18:1,18:2,18:3)
102 increased (**Fig. 2e,f**). Together, this analysis identifies differences in metabolism of ceramides and GLs in
103 16pdel neurons that are critical for function of the ER, mitochondria and plasma membrane (Flis and Daum,
104 2013). The shift in saturation and tail length of GLs between 16pdel and control neurons suggests a
105 dysfunctional neuronal membrane.

106

107 **FAM57B functions as a ceramide synthase modulator**

108 We considered that the extensive lipid differences between 16pdel and control neurons may partly
109 result from FAM57B activity. The function of this protein is not clear, although a single report suggests that
110 FAM57B has ceramide synthase activity (Yamashita-Sugahara et al., 2013). However, sequence analysis
111 indicates that while FAM57B is part of the TLC protein family, including ceramide synthases (CerS), FAM57B
112 has little sequence similarity to CerS, even in the TLC domain (**Supp Fig. 3**). To assess whether FAM57B is a
113 *bone fide* CerS, it was expressed in *CerS2^{-/-}* (KO) HEK293T cells, which lack endogenous CerS2 activity
114 (Tidhar et al., 2012) (**Fig. 3a**). No CerS2 activity was detected in *CerS2* KO cells upon transfection of FAM57B
115 alone. However, co-transfection of FAM57B with CerS2 resulted in a significant increase in CerS2 activity
116 compared to transfection of CerS2 alone (**Fig. 3a**), suggesting that FAM57B might modulate CerS2 activity.
117 There are six CerS isoforms in mammals, where each uses a restricted subset of acyl CoAs of defined chain
118 length for ceramide synthesis (Levy and Futerman, 2010). To assess whether FAM57B can modulate other
119 members of the mammalian CerS family; we expressed CerS5 and CerS6 with or without co-transfection of
120 FAM57B in wildtype HEK293T cells. Upon co-transfection of CerS2 with FAM57B in wildtype HEK293T cells,
121 levels of CerS2 activity and expression were significantly increased compared to CerS2 alone (**Fig. 3b**). While
122 co-transfection of FAM57B with CerS5 did not alter expression nor activity of this CerS (**Fig. 3c**), an opposite
123 trend was seen upon co-transfection of FAM57B with CerS6, whose activity decreased upon co-transfection
124 with FAM57B (**Fig. 3c**). These results suggest that FAM57B affect protein levels and activity of certain CerS
125 isoforms, and may do so by an indirect mechanism, dependent on interaction of the two proteins. This
126 hypothesis was confirmed by immunoprecipitation, in which Flag-tagged FAM57B was able to interact with all
127 three HA-tagged CerS isoforms (**Fig. 3d**). These data newly implicate FAM57B as a modulator of CerS, but
128 refute a previous report that this protein functions as a CerS (Yamashita-Sugahara et al., 2013).

129 130 **FAM57B modulates lipid cohorts and synaptic proteins in human cells**

131 The intriguing functional differences between 16pdel and control neurons raises the question of
132 whether *FAM57B* haploinsufficiency contributes to these differences. To address this, we used the human
133 neuroblastoma line SH-SY5Y to create knockout (*FAM57B* KO) and *FAM57B* heterozygote (*FAM57B* HET)
134 lines, using CRISPR-Cas9 editing. SH-SY5Y cells have proven useful for studying neuronal properties and
135 function (Kovalevich J, 2013). After confirmation of CRISPR induced genome editing by next generation
136 sequencing, FAM57B protein depletion was confirmed by western analysis (**Supp Fig. 4**). For our studies, SH-
137 SY5Y cells were differentiated into neurons after incubation in media containing retinoic acid. Overall, total lipid
138 classes showed significant differences between *FAM57B* KO and WT (wildtype), specifically, increased ChE
139 and MG (**Fig. 4a, Table S2 - SH-SY5Y Lipidomics**). Comparing *FAM57B* HET to WT, we observed increased
140 LPC (**Fig. 4b**). Additionally, relative to *FAM57B* HET, we found HexCer and PG significantly decreased while
141 ChE increased in *FAM57B* KO cells (**Fig. 4c**).

142 Notably, lipid class differences observed between *FAM57B* KO, *FAM57B* HET and WT were similarly
143 altered in 16pdel patient neurons compared to controls (**Fig. 2a**). This, in both *FAM57B* KO and *FAM57B* HET

144 relative to WT, we observed increased abundance of PE(18:0,18:1,22:4,22:5) (**Fig. 4d**), Cer(d18:1) (**Fig. 4e**),
145 MG(18:0) (**Fig. 4f**), and with decreased abundance of TG(16:0,16:1,18:0,18:1,18:2,22:6) (**Fig. 4g**). These
146 differences are similar to those seen in 16pdel relative to control neurons (**Fig. 2c-f**). In *FAM57B KO* relative to
147 WT, there was increased abundance of MG and decreased abundance TG (**Fig. 4f,g**), as for 16pdel neurons
148 compared to control (**Fig. 2a,e,f**). The alterations in lipid cohorts between *FAM57B KO* and *FAM57B HET*
149 human neurons is similar to lipid changes in 16pdel neurons compared to control, and consistent with a role for
150 *FAM57B* in dosage-sensitive lipid regulation. The similarities in lipid cohort alterations between *FAM57B KO*
151 compared to *FAM57B HET* in SH-SY5Y human neurons and 16pdel neurons, is consistent with a role for
152 *FAM57B* in dosage-sensitive, lipid regulation, including tight regulation of HexCer (**Fig. 4c**).

153 To understand the consequence of *FAM57B loss of function* on neuronal maturation and function, we
154 probed synaptic composition. Synaptosomes, comprising the pre- and postsynaptic membranes and
155 postsynaptic density, were isolated from unfixed cells and processed by MS/MS (**Table S3 - SH-SY5Y**
156 **Synaptosome MSMS**). Synaptosomes from *FAM57B KO* cells showed significantly decreased abundance of
157 over 100 proteins relative to WT (**Fig. 5a**). In contrast, *FAM57B HET* showed no statistically significant
158 changes in synaptosome protein composition relative to WT (**Fig. 5b**), with the exception of Dopamine Beta-
159 Hydroxylase (DBH). Among the top 20 significantly decreased proteins in the *FAM57B KO* synaptosomes were
160 those associated with protein trafficking, localization and stabilization (**Fig. 5c,d**). Additionally, we observed an
161 overall decrease in levels of hallmark synaptic proteins in *FAM57B KO* compared to *FAM57B HET* or WT (**Fig.**
162 **5e**). These decreases included α -internexin INA, small GTPase vesicle recycling RAB11B, SNARE protein
163 syntaxin STXBP1 and scaffolding protein YWHAZ. INA is a neurofilament subunit protein important for
164 neuronal cytoskeletal assembly and synaptogenesis localized to the post-synaptic terminal (Yuan and Nixon,
165 2016).

166 Separately, synaptosomes were isolated from differentiated SH-SY5Y cells and processed for lipidomic
167 analysis (**Table S4 - SH-SY5Y Synaptosome Lipidomics**). Comparing lipids localized to the pre- and post-
168 synaptic terminals between *FAM57B KO* to WT neurons, significant changes in lipid abundances were
169 observed, notably in hexosylceramide (HexCer) and monoacylglycerol (MG) (**Fig. 5f**). Relative to WT, Hex2Cer
170 and MG abundances increased, while Hex1Cer and phosphatidic acid (PA) abundances decreased. No
171 statistically significant changes in lipid group abundances were observed between *FAM57B HET* relative to WT
172 or *FAM57B KO*. These results suggest altered lipid composition, due to loss of *FAM57B* function at developing
173 synapses impacts localization and/or trafficking of synaptic proteins in *FAM57B* mutant neurons. Together, the
174 data indicate that in human neurons mutant or heterozygous for *FAM57B*, there are significant changes in lipid
175 composition and regional synaptic protein abundance. The data are consistent with the suggestion that a
176 deficit in *FAM57B* function partly contributes to 16pdel neuronal anomalies relative to control. The smaller
177 changes observed in *FAM57B HET* relative to *FAM57B KO* suggests that other 16pdel genes contribute to
178 phenotypes in the haploinsufficient syndrome.

179

FAM57B is essential for Sphingolipid (SL) and Glycerolipid (GL) homeostasis in the developing brain.

16pdel alters brain structure and function, including neuroanatomical abnormalities and increased risk of psychiatric and other brain disorders (Niarchou et al., 2019, Owen et al., 2018). To understand how *FAM57B* contributes to brain development, we analyzed zebrafish, *Danio rerio*, a powerful system for analysis of neural development and neurodevelopmental disorders (Kalueff et al., 2014, Stewart et al., 2014, Xi et al., 2011, McCammon and Sive, 2015). The zebrafish genome includes two copies of the *fam57b* gene, *fam57ba* and *fam57bb*. We used CRISPR to build double (null) mutants, *fam57ba*^{-/-};*fam57bb*^{-/-} (*fam57b mut*), and heterozygotes *fam57ba*^{+/-};*fam57bb*^{+/-} (*fam57b het*), to assess dosage effects of FAM57B.

To determine whether *fam57b* regulates lipid metabolites, we performed untargeted lipid profiling on *fam57b mut* and *fam57b het* zebrafish brain tissue at 7 days post-fertilization (7 dpf), an optimal timepoint for molecular and behavioral studies of a developing yet complex brain (**Table S5 - Zebrafish Larvae Brain Lipidomics**) (Tomasello and Sive, 2020). Striking differences in SL and GL lipid abundances were present in *fam57b mut* and *fam57b het* compared to wildtype (AB) zebrafish brain (**Fig. 6a,b**). By lipid class, there was a significant increase in Cer, LPE, MG and SM along with phosphatidylinositol (PI) and cardiolipin (CL) and decreased PS in *fam57b mut* compared to AB (**Fig. 6a**). A similar trend to *fam57b mut*, with increased hexosylceramide (HexCer) and decreased PS lipid classes, were defined in *fam57b het* brains compared to AB control. An overlap in lipid differences were observed between *fam57b mut* brains and FAM57B KO human neurons, with increased abundance of MG class, and Cer(d18:1) and MG(18:0) species (**Figs. 6a, 4a, 6d,e, 4d,e**). Many PE species similarly increased, PE(18:0,18:1,20:4), comparing *fam57b mut* and *FAM57B KO* to controls. An important finding across all systems compared, including 16pdel syndrome patient neurons, heterozygous and mutant *FAM57B* cells and larvae brains, is a change in ether-linked PE (**Figs. 2c, 4d, 6c,g**). Ether GLs differ in phase-transition temperature from gel to liquid crystalline and from lamellar to hexagonal phases, and are proposed to regulate properties of neuronal membranes (Paltauf, 1994, Lohner, 1996). In addition to PE, ceramides were altered in *fam57b mut* and *fam57b het* brain tissue compared to AB (**Fig. 4d,h**). Lipidomics resolved predominantly Cer(d18:1) species in zebrafish brains, which agrees with previously published ceramide composition at 7 dpf (Zhang et al., 2019). These findings suggest a key role for Fam57b in SL and GL regulation during brain development.

These data demonstrate that *fam57b* is crucial for regulation of SL and GL classes in the larval zebrafish brain, and that there is a gene dosage-dependent effect. The data in zebrafish brain are quite similar to changes seen in human neurons after FAM57B knockout (**Fig. 4**). These changes affect comparable lipid groups to those altered in 16pdel neurons relative to control, although they are not the same. For example, LPE significantly decreased in the 16pdel neurons (**Fig. 2a**), while the class remained unchanged or significantly increased in the *fam57b het* and null SH-SY5Y cells (**Fig. 4a-c**) and zebrafish brain (**Fig. 6a,b**), suggesting additional genes regulate 16pdel lipid metabolites or that these result from other differences between the tissue being compared.

215

216 **Changes in plasma membrane and associated proteins in *fam57b mut* and *fam57b het* brains**

217 Ceramide, hexosylceramide and GL species are integral to membrane composition, are differentially
218 distributed across inner and outer leaflets of the plasma membrane and contribute to lipid rafts (Kraft, 2016).
219 Given the changes in these lipids observed in *fam57b mut* brains, we predicted that plasma membrane
220 structure would also be altered. To assess lipid raft organization, fluorophore-conjugated Cholera Toxin subunit
221 B (CT-B), which binds ganglioside GM1 found in lipid rafts (Fishman et al., 1978), was injected into the
222 hindbrain ventricle of *fam57b mut* and AB zebrafish embryos at 24 hours post fertilization (hpf), when ventricles
223 are accessible for injection (Worstell et al., 2016) (**Fig. 7a**). Embryos fixed after 1-hour incubation
224 demonstrated a significant increase in punctate GM1 labeling in neural progenitor cells of *fam57b mut* brains
225 compared to AB (**Fig. 7b**). To assess changes in glycerophospholipid species in plasma membranes, we
226 stained with duramycin, a label for membrane PE (Marconescu and Thorpe, 2008). Mutant progenitors showed
227 statistically increased punctate PE staining, indicating altered PE localization that could impact membrane
228 architecture (**Fig. 7c**). The duramycin puncta may indicate exosomes or extracellular vesicles containing PE
229 (Beer et al., 2018). At 24 hpf, we did not observe changes in cell proliferation or cell death between *fam57b*
230 *mut* and AB (**Supp. Fig 5**). The data suggest that there is alteration in the plasma membrane of *fam57b mut*
231 brains relative to AB.

232 The changes in membranes of neural progenitor cells suggested that membrane protein localization
233 may also be altered. We therefore examined localization of membrane by biotinylation analysis. Freshly
234 dissected larval brains from *fam57b mut* or AB at 7 dpf were incubated with membrane impermeable biotin.
235 Surface proteins were affinity-purified and quantified by MS/MS (**Fig. 7d, Table S6 - Zebrafish Larvae Brain**
236 **Biotinylation MSMS**). MS/MS analysis indicated that membrane-associated protein cohorts were similar
237 between *fam57b mut* and AB brains (**Fig. 7e**), however, a small group of proteins showed altered abundance.
238 In *fam57b mut* brains, the protein whose levels most significantly decreased (2-fold) relative to AB was
239 Synaptotagmin-1a (Syt1) (**Fig. 7e**). Syt1a, homologous to human SYT1, is a vesicle membrane protein that
240 acts as a calcium sensor and regulates synaptic and endocrine vesicle exocytosis (Fernandez et al., 2001,
241 Sorensen et al., 2003, Xu et al., 2007, Gustavsson and Han, 2009, Schonn et al., 2008). Syt1 protein domains
242 interact with the lipid bilayer, including GL PS that are altered in *fam57b mut* and *fam57b het* (**Fig. 5a,b**).
243 Mammalian Syt1 can modify PS, and is able to alter curvature strain on the membrane (Lai et al., 2011).

244 To investigate the decreased membrane abundance of Syt1 in the biotinylation assay, we performed
245 immunostaining on 7 dpf larvae brains (**Fig. 7f**). Whole larval brains were cleared and tertiary structure was
246 protected using SHIELD protocols. Slice imaging of dorsal brain view showed that Syt1 protein was largely
247 confined to projections of neurons throughout the AB brain, both GABAergic (GAD65/67) and non-GABAergic,
248 while *fam57b mut* brains showed ectopic expression throughout the brain (**Fig. 7f**). By western blot, we found
249 no change in total brain Syt1 between *fam57b mut* and AB, suggesting that immunostaining demonstrates
250 Syt1a mislocalization (**Fig. 7g, Supp Fig. 6**). Imaging also revealed anatomical changes in the larval *fam57b*
251 *mut* brain, including tectum and corpus cerebelli (**Fig. 7f**).

252 Together, these results indicate that in the brain, relative to wildtype, *fam57b mut* animals show
253 changes in lipids, membrane structure and membrane protein association, including the synaptic regulator
254 Syt1 and others functioning at synapses. The data indicate that *fam57b* is required for membrane structure and
255 neuronal architecture.

256

257 **Pre- and post-synaptic proteins depleted after loss of *FAM57B***

258 To examine the implications of Syt1 mis-localization on synaptic composition, we isolated
259 synaptosomes from freshly dissected and unfixed brains of *fam57b mut* or AB larvae (**Table S7 - Zebrafish**
260 **Larvae Brain Synaptosome MSMS**). Proteomic profiling indicated a group of proteins whose representation
261 significantly increased, and another larger group whose representation significantly decreased in *fam57b mut*
262 compared to AB (**Fig. 7h**). Interestingly, we observed a decrease of the Synaptotagmin family member Syt2a-
263 like protein, similar to human SYT2, with analogous function to SYT1. Gene ontology (GO-Slim and Panther
264 Protein Class ontology) defined synaptic protein groups found only in the decreased synaptosome protein
265 group (**Fig. 7i, 7j**). Annotations in the decreased group included synapse and synapse part components,
266 cytoskeletal and membrane traffic proteins, biological adhesion, development and signaling, and numerous
267 implicated pathways including synaptic vesicle trafficking. These results indicate that synaptic protein levels
268 were significantly altered in synaptosomes from *fam57b mut* larval brain synapses relative to AB.

269 We separately examined levels of synaptotagmin family members in brain synaptosome profiles and
270 found decreased Syt1 and Syt2a protein levels in *fam57b mut* compared to AB synaptosomes (**Fig. 7k**). This
271 interesting association between FAM57B regulation and Synaptotagmin expression (**Fig. 7e,f,h**) led us to
272 analyze further human SH-SY5Y isolated synaptosomes. We found a significant decrease in elongated SYT1
273 (ESYT1) in *FAM57B KO* compared to *FAM57B HET* and WT (**Fig. 7l**), a calcium activated synaptic protein
274 found to bind GLs (Yu et al., 2016). We then characterized hallmark proteins that function at the synapse from
275 brain synaptosomes, including synaptic vesicle fusion and tethering proteins. Bayés *et al.* previously examined
276 complexity of the adult zebrafish synapse proteome relative to adult mouse synapse proteome (Bayes et al.,
277 2017). As expected, not all synaptic proteins were detected at this immature stage of zebrafish development
278 compared to the adult brain. Enrichment of synaptic vesicle proteins including Syntaxins, Slc neurotransmitter
279 transporters, SNAPs, Stx/Vps, Synaptotagmins and membrane budding proteins including Dynamins and Rabs
280 verify synaptosome isolation and give new data regarding neuronal maturation in the zebrafish larval brain
281 (**Fig. 7j**). Comparison of synaptic protein profiles between genotypes demonstrated decreased vesicle fusion
282 and transport protein Nsfa, ligand-gated ion channel Si:ch211-251b21.1, and SNARE complex proteins Snphb,
283 Stx1b, Stxbp1a and Vamp3 (**Fig. 7j**). Together, these data show that synaptic proteins essential for vesicle
284 docking, exo- and endocytosis, including synaptotagmin family members, are diminished in synaptosomes
285 isolated from *fam57b mut* brains relative to AB, suggesting *Fam57b* is essential for synapse integrity.

286

287 **Depressed spontaneous electrical activity and response to stimuli in *fam57b* mutants**

288 To understand how changes in *fam57b* gene dosage impact neuronal activity, we tested brain activity
289 by electrophysiological analysis. We previously described a noninvasive electrophysiology technique that can
290 be used in live larvae to measure spontaneous activity in the brain and spinal cord (Tomasello and Sive, 2020).
291 Using a multielectrode array (MEA), we measured local field potential (LFP) parameters and relative
292 coordinated (network) activity in the brain of 7 dpf larvae (**Figs. 8a-c**). Larva were individually immersed in
293 precooled 1.5% low-melt agarose in E3 solution and mounted in a 64-electrode containing well. We measured
294 spontaneous brain activity over a 10-minute period, comparing *fam57b mut* to AB controls. Only electrodes in
295 contact with the larval head were analyzed, ~6 to 8 electrodes, whose signal was pooled. *fam57b mut* larvae
296 had slightly smaller heads than AB at 7 dpf (**Supp Fig. 7, Table S8 - Zebrafish Larvae Head and Body**
297 **Measurements**), however these changes do not impact electrophysiological studies. Overall, *fam57b mut*
298 spontaneous brain activity was severely diminished relative to ABs. This included significant decrease in
299 number of LFPs, mean LFP rate, and inter-LFP-interval coefficient of variation measurements, indicating
300 decreased spontaneous brain activity with reduced kinetics (**Fig. 8a**). However, the decreased coefficient of
301 variation in the *fam57b mut* suggests LFP interval distributions are detected at a more regular rate. Measuring
302 electrographic bursts, at least 5 LFPs per 100 ms, we were unable to detect any burst activity under these
303 settings in the *fam57b mut*, while bursts were detected in ABs. To increase sensitivity for detection of burst
304 activity, we lowered the detection parameters to at least 3 LFPs per 200 ms (right column), and found
305 decreased electrographic burst duration, number of LFPs per burst, burst frequency and percentage in *fam57b*
306 *mut* relative to AB (**Fig. 8a**). In addition, we examined relative network activity, as defined by at least 3 LFPs
307 detected simultaneously between a minimum of two electrodes. Relative network activity was also significantly
308 decreased in *fam57b mut* compared to AB. Synchrony index of bursts did not change between the two
309 genotypes, indicating coordination of network activity did not differ. While LFP waveforms could not be
310 quantified due to small distance variations when mounting individual larva, we observed smaller relative
311 waveforms in *fam57b mut* compared to AB (**Fig.8c**), consistent with overall decreased brain activity in *fam57b*
312 *mut* larvae. A representative raster plot of LFP activity in the head region over the 10-minute recording period
313 illustrates the relative decrease in LFP propagation, burst and network detection measured (**Fig. 8d**). A
314 representative image of a mounted larva immersed in agarose on a 12-well 64 electrode plate is shown in **Fig.**
315 **8e**. These data demonstrate severely diminished spontaneous brain activity in *fam57b mut* relative to AB
316 wildtype larvae, and highlight a role for Fam57b in regulating brain function.

317 After identifying significantly diminished spontaneous brain activity in *fam57b mut* larvae, we examined
318 correlations with behavioral activity. We first tested light-responsive sensorimotor startle behavior (**Table S9 -**
319 **Zebrafish Larvae Light Startle Response Data**). Startle response, as indicated by distance traveled, was
320 measured over a 70-min time-frame with light extinguished every 10 minutes for 5 seconds (Tomasello and
321 Sive, 2020). The startle response window was in total 30 seconds, including the stimulus. We found a
322 considerable decrease in response to each light stimulus in *fam57b mut* compared to AB (**Fig. 8f**). However,
323 movement measured previous to startle (first 10 min) and relative habituation after startle cue did not overall

324 differ between the genotypes, indicating there is no alteration in movement outside of the light stimulus, and no
325 visual deficit in *fam57b mut* larvae.

326 To examine brain specific activity, we investigated seizure susceptibility (**Table S10 - Zebrafish Larvae**
327 **Seizure Assay Data**). Seizures are prevalent in individuals affected with 16pdel syndrome, and may result
328 from processes involving several neurotransmitter systems, including glutamatergic, cholinergic and
329 GABAergic (Mefford et al., 2011). To measure seizure propensity, larvae were immersed in pentylenetetrazol
330 (PTZ), a GABA_A antagonist, well characterized for use in zebrafish (Baraban et al., 2005). After a 10-minute
331 baseline movement recording, two different concentrations of PTZ, or E3 media only control, were applied to
332 the individual well of each larva and recorded over 10 minutes. There was no significant change in normalized
333 movement, compared to baseline recording, in the absence of stimulus after addition of E3 (0 mM PTZ) control
334 between the *fam57b mut* and AB (**Fig. 8g**). Increasing the PTZ dose increased normalized distance traveled
335 for both genotypes, but the increase was not significant for AB between 0 and 0.5 mM PTZ as observed in
336 *fam57b mut*. At 5 mM, we observed significantly less distance traveled in the *fam57b mut* compared to AB.
337 However, the relative fold change between 0 and 5 mM was much higher in the *fam57b mut* (roughly 32-fold)
338 compared to AB (roughly 11-fold). To understand the contribution of Fam57ba and Fam57bb to the seizure
339 phenotype, we incrossed *fam57ba^{-/-};fam57bb^{+/-}* animals and measured seizure propensity in resulting
340 genotypes (**Supp Fig. 8**). We found no statistical differences in baseline movement between AB control
341 compared to *fam57ba^{-/-};fam57bb^{+/-}*, *fam57ba^{-/-};fam57bb^{-/+}* nor *fam57ba^{+/-};fam57bb^{+/+}*. The data are consistent
342 with increased movement after addition of 5 mM PTZ in *fam57ba^{-/-};fam57bb^{+/+}* 7 dpf larvae compared to AB
343 control (McCammon et al., 2017). The enhanced PTZ-responsiveness of *fam57b mut* relative to WT suggests
344 that Fam57ba and Fam57bb function in GABA-mediated signaling with synergistic effects after the loss of
345 Fam57bb in combination with Fam57ba. These analyses indicate neuronal specific changes after loss of
346 *fam57b mut*, however, FAM57B is also expressed in muscle. We assessed neuromuscular junction
347 contribution by immunostaining, but did not observe differences between *fam57b mut* and AB (**Supp. Fig. 9**).
348 Together, altered brain activity, response to multiple stimuli including GABA_A antagonist, and gross anatomical
349 differences including the corpus cerebelli (**Fig. 7f**), suggest changes to GABAergic network activity in the
350 developing *fam57b mut* brain.

351 In sum, there are significant behavioral changes in zebrafish larvae after *fam57b* loss of function. *fam*
352 *mut* larvae move similarly to AB over time without a stimulus. However, with a stimulus – either dark or PTZ
353 application, there is altered behavioral responsiveness relative to AB. These findings are consistent with
354 changes in brain activity in *fam57b mut* relative to AB controls.

355

356 Discussion

357 This study has uncovered alterations of lipid metabolism in iPSC differentiated cortical neurons derived
358 from people affected with 16pdel Syndrome, relative to unaffected. These 16pdel neurons display increased
359 excitability relative to controls, and show a sex-linked difference. Among the set of twenty-five 16pdel genes,

360 *FAM57B* is a key candidate gene linked to symptomatology, and implicated in lipid metabolism. Consistently,
361 we find changes in 16pdel neuronal lipids, as well as those in *FAM57B* mutants in a human neuronal cell line
362 and zebrafish model. In *FAM57B* mutants, lipid alterations occur concomitantly with alterations in membrane
363 architecture, synapse lipid composition synapse-associated proteins, and in zebrafish, altered brain activity
364 and behavior. Specifically, we identified altered abundance of the monoacylglycerol (MG) group in all
365 experimental conditions, including 16p11.2 deletion neurons and SH-SY5Y *FAM57B* KO intact neurons and
366 isolated synaptosomes (**Figs. 2a, 4a, 5f, 6a**). We do not know how these outcomes are linked, but it is
367 plausible that *FAM57B* acts through sphingolipid (SL) and glycerolipid (GL) regulation as the starting point for a
368 cascade of effects after loss of function. Dysregulated lipid metabolism has a multifaceted effect on neurons,
369 for example, increased lipid energy consumption escalates oxidative stress, promoting inflammation,
370 mitochondrial and metabolic dysfunction and excitotoxicity (Tracey et al., 2018). The saturation and length of
371 individual lipids affects their intracellular localization, impacting the cytoskeleton and lipid raft composition, so
372 disrupting signaling processes that regulate neurotransmitter synthesis and release, cytoskeletal integrity,
373 myelination and intracellular transport (Pike, 2003, Tracey et al., 2018). Abnormal cholesterol metabolism has
374 been observed in patients with Asperger syndrome and other ASDs, suggesting a correlation between lipid raft
375 formation and ASD (Dziobek et al., 2007, Tierney et al., 2006).

376 In contrast to a previous report (Yamashita-Sugahara et al., 2013), we find that *FAM57B* is not a
377 ceramide synthase (CerS) but rather is a modulator of CerS activity. Supporting a functional interaction
378 between *FAM57B* with CerS2 and CerS6, lipidomic profiling uncovered altered SLs and GLs integral to the
379 lipid membrane in *FAM57B* mutants of both zebrafish brain and a human neuronal cell line, also indicating
380 consistent activity of this gene across species. Comparing these mutants with 16pdel iPSC differentiated
381 patient neurons, we identified a consistent change in ether-linked phosphatidylethanolamine (PE) species,
382 supporting a role for *FAM57B* in the altered 16pdel lipidome relative to unaffected. However, some lipid groups
383 altered in mutant human neuronal cells, zebrafish brain and 16pdel neurons relative to controls are do not
384 overlap, suggesting additional genes regulate 16pdel lipid metabolites or that these result from other
385 differences between tissues being compared. Beyond *FAM57B*, multiple genes in the 16p11.2 interval encode
386 enzymes with predicted roles in metabolic processing or interconversion including *ALDOA*, *CDIPT*, *GDPD3*,
387 *BOLA2*, *SULT1A3*, *SULT1A4* and *YPEL3* (Giannuzzi et al., 2019, Arbogast et al., 2016). Together with
388 *FAM57B* activity, this set of genes may function to modulate lipid metabolism.

389 Early in brain development, *fam57b* mutants displayed altered plasma membrane architecture, while in
390 more mature neurons, synaptic proteins were present at significantly diminished levels in synaptosomes
391 prepared from mutants relative to controls. One important affected protein in zebrafish brain was Syt1a, a
392 member of the Synaptotagmin family and calcium sensor SNARE binding complex protein that contributes to
393 synchronous synaptic vesicle release (Li et al., 2017). Baker-Gordon Syndrome, a SYT1-associated
394 neurodevelopmental disorder, maps to an autosomal dominant heterozygous mutation of *SYT1*, that is
395 associated with reduced neurotransmitter release (Baker et al., 2018). This association supports our previous

396 findings of a genetic interaction between *fam57ba* and *doc2a*, another 16p11.2 interval gene, encoding a
397 calcium sensor SNARE binding complex protein for spontaneous vesicle release, where double heterozygotes
398 showed hyperactivity and increased seizure propensity (McCammon et al., 2017). In further connections,
399 ESYT1 is a related Synaptotagmin family member whose synaptic levels were diminished in human neuronal
400 cell lines, that may play a role in cellular transport of PC, PE, PI, and translocates to sites of contact between
401 the presynaptic endoplasmic reticulum and the cell membrane in response to increased cytosolic calcium
402 levels (Yu et al., 2016). Neurotransmission is decreased in *Esy1 D.melanogaster* mutants, with a proposed role
403 in synapse extension, highlighting the essential homeostasis of lipids at the synapse (Kikuma et al., 2017).

404 Diminished spontaneous brain activity and altered behavioral response after stimulation seen in *fam57b*
405 zebrafish mutants is consistent with alteration of brain synaptic composition. These studies suggest Fam57b is
406 essential early in development of the brain, and loss of *fam57b* leads to linked events starting with changes in
407 plasma membrane architecture followed by disturbance in protein organization at the membrane and
408 detriments to basic neuronal function that impacts brain activity and behavior. Together, we propose a model
409 whereby Fam57b functions to maintain normal plasma membrane physiology, necessary for proper formation
410 and function of neurons (**Fig. 8h**).

411 In a recent study that analyzed the largest ASD-associated exome sequence to date (Satterstrom et al.,
412 2019), 102 high risk genes were identified as tightly associated with ASD. In the analysis, a rare G:A mutation
413 was discovered in *FAM57B*, located in the 5'UTR of one *FAM57B* transcript isoform, and residing in the
414 promoter/enhancer region of the five other *FAM57B* isoforms. This synonymous mutation is predicted to create
415 binding sites for several transcription factors and may impact enhancer activity in neurons, affecting gene
416 expression (prediction with information from dbSNP, JASPAR, GTE_x). The association of a single gene *FAM57B*
417 mutation with ASD outside of the complex CNV 16pdel gene cohort, encourages further evaluation of *FAM57B*
418 in brain development and function.

419 Correlating the multitude of symptoms to specific genes associated with a multigenic copy number
420 variant region is extremely challenging. Variation in clinical phenotypes of 16p11.2 deletion syndrome patients
421 further indicates the need to understand the biology of this CNV (Fetit et al., 2020). Previously, we identified
422 gene interactions among the 16p11.2 interval (McCammon et al., 2017) that do not converge on functional
423 networks predicted by ASD gene-set enrichment analysis by Pinto *et al.* (Pinto et al., 2010), suggesting indirect
424 mechanisms of genetic interaction. We investigated the interaction between *doc2a*, encoding for a synaptic
425 vesicle-associated calcium-binding protein, and *fam57ba*. Haploinsufficiency of this genetic interaction
426 identified both a body and central nervous system phenotype, including seizure activity. No evidence for
427 physical interactions between proteins encoded by 16p11.2 genes has been found (Lin et al., 2015). Along with
428 *DOC2A*, several genes within the interval are candidate contributors to neurodevelopment and
429 neuropsychiatric phenotypes, including *KCTD13*, *SEZ6L2*, *KIF22*, *MVP*, *TAOK2* and *QPR1*. In a functional
430 screen, we defined genetic interaction between *fam57ba*, *kctd13*, *sez6l2* and *kif22* (McCammon et al., 2017).
431 *KCTD13* encodes a ubiquitin ligase adaptor with cognitive defects identified in mice heterozygous for the gene

432 (Chen et al., 2009, Golzio et al., 2012, Martin Lorenzo et al., 2021). KIF22 is a kinesin-like protein necessary
433 for embryonic chromosome segregation and axonal branching patterns (Park et al., 2016, Antonio et al., 2000,
434 Ohsugi et al., 2008). SEZ6L2 has been connected to seizure activity and modulates neurite outgrowth (Boonen
435 et al., 2016). MVP can function as a regulator of the homeostatic component of experience-dependent
436 behavior (Ip et al., 2018). *TAOK2* encodes a serine/threonine kinase that can play a role in dendrite formation
437 (de Anda et al., 2012). *QPRT* encodes quinolinate phosphoribosyltransferase that catabolizes quinolinic acid,
438 whose elevation has been linked to epilepsy and has also showed regulation by ASD candidate genes
439 (Chiocchetti et al., 2016, Haslinger et al., 2018). The strong genetic interaction between *FAM57B* and most of
440 the above-described genes highlights the connection between lipid regulation and brain development or
441 maturation. Although *FAM57B* haploinsufficiency alone cannot account for the multitude of disrupted
442 biochemical and cellular properties in 16pdel affected neurons, disrupted lipid metabolism is tightly correlated
443 to 16pdel Syndrome. The insight into lipid alterations and a potential role for *FAM57B* in mediating these
444 changes, gives a new view of mechanisms underlying 16pdel Syndrome, and holds promise for new
445 therapeutic directions.

446

447 **Limitations of the Study**

448 In the large 16p11.2 copy number variant interval, haploinsufficient symptoms do not resolve to a single causal
449 gene. *FAM57B* interacts genetically with many other 16p11.2 interval genes and is a strong candidate for
450 contribution to symptomatology. We studied the poorly defined function of *FAM57B* to understand its role in
451 lipid regulation and function in the brain. *FAM57B* mutant neuronal lines do not recapitulate all phenotypes
452 associated with 16p11.2 deletion syndrome neurons. For instance, the *FAM57B* heterozygote SH-SY5Y
453 differentiated neurons and zebrafish brains do not show exactly the same lipid profiles as neurons derived from
454 16p11.2 deletion iPSC, although there is overlap. Since 16pdel syndrome results from changes in gene
455 dosage through haploinsufficiency, rescue of *FAM57B* levels must be precise to match endogenous levels to
456 avoid spurious gene dosage effects. However, in the 16p11.2 chromosomal deletion, the *FAM57B* promoter
457 and enhancer regions are also deleted, it is not an easy task to rescue *FAM57B* expression to endogenous
458 levels. The nature of the copy number variant phenotype indicates that exogenous expression to rescue
459 *FAM57B* could result in outcomes modulated by dosage, and that the entire gene with its regulatory regions
460 would have to be used for rescue. The *FAM57B* regulatory regions have not been mapped, and so are not
461 available for rescue constructs.

462

463 **Acknowledgements**

464 We thank Olivier Paugois for excellent fish care and lab management. Special thank you to the Whitehead
465 Institute core facilities as follows: Wendy Salmon (Keck Imaging Facility) for help with confocal imaging, Eric
466 Spooner (MS Core) for proteomic work, Dr. Caroline Lewis (Metabolomics Core) for lipidomics analysis, Dr.
467 George Bell (Bioinformatics and Research Computing) for analysis and help in compiling MS/MS data and Dr.

468 Heather Keys (Functional Genomics Platform) for aid in CRISPR-induced *FAM57B* mutations. We thank Dr.
469 Shifra Ben-Dor, Bioinformatics Unit, Life Sciences Core Facilities, Weizmann Institute of Science, for sequence
470 alignment in Supplemental Fig 3. We thank Dr. Mike Gallagher for help analyzing the published FAM57B
471 mutation from Satterstrom *et al.* 2019 (Satterstrom et al., 2019), Dr. Monika Saxena for constructive and
472 helpful edits to the manuscript, Master of Science student Adrianna Vandeuren for help maintaining the fish
473 lines, and Undergraduate Research Opportunity student Allysa Allen for help in expanding iPSCs. Finally, we
474 would like to thank colleagues at the Simons Center for the Social Brain at MIT for collegial interactions. This
475 work was supported by a grant from the Simons Foundation to the Simons Center for the Social Brain at MIT,
476 National Institute of Mental Health (NIMH) of the National Institutes of Health (1R01MH119173-01A1), a
477 Balkin-Weinberg-Markell Fellowship, the Israel Science Foundation Grant #3172/19 and additional support
478 from Jim and Pat Poitras. RJ was supported by NIMH grant 2R01MH104610-20.

479

480 **Author Contributions**

481 DLT and HS designed the study, interpreted results and wrote the manuscript. AHF helped conceive,
482 supervise, interpret and write the CerS study (**Fig. 3**). JK and YK performed the experiments in **Fig. 3**. JMM
483 created the *fam57b* mutant zebrafish line. MM and RJ instructed on stem cell culture work and provided
484 unaffected control iPSC (originally from Coriell Institute Biobank). DLT performed all other studies.

485

486 **Declaration of Interests**

487 The authors declared no potential conflicts of interest with respect to the research, authorship, and/or
488 publication of this article.

489

490 **STAR Methods**

491

492 **Resource Availability**

493 *Lead Contact*

494 Further information and requests for resources and reagents should be directed to and will be fulfilled by the
495 lead contact, Hazel Sive (h.sive@northeastern.edu).

496 *Materials Availability*

497 pcDNA3.1 FAM57B and CerS constructs, SH-SY5Y *FAM57B* HET and *FAM57B* KO, and Zebrafish lines
498 *fam57ba*^{+/-};*fam57bb*^{+/-} and *fam57ba*^{-/-};*fam57bb*^{-/-} created in this manuscript are available upon request from
499 lead contact.

500 *Data Availability*

- 501 • The authors declare that all data supporting the findings of this study are available within the article and
502 its supplementary information.
- 503 • Raw data, including proteomics and lipidomics results, are available in the Supplemental Excel
504 spreadsheet. Additional data can be requested from lead contact.
- 505 • Any additional information required to reanalyze the data reported in this paper is available from the
506 lead contact upon request.

507 **Experimental model and subject details**

508

509 **Animal Model**

510 Adult zebrafish of the wildtype AB strain were maintained at 28°C on 12h/12h light/dark cycle. Embryos were
511 obtained from natural spawning and staged as previously described by Kimmel *et al.* (Kimmel et al., 1995).
512 Due to the polygenic nature of sex determination and timing of gonadal development in zebrafish, we are
513 unable to determine the sexes of the embryos and larvae for our assays. However, because our assays
514 utilized large numbers of embryos and larvae, both sexes should be adequately represented. Embryos were
515 obtained from separate crosses of *fam57b mut* mutant fish. *fam57b het* fish were generated by crossing
516 *fam57b mut* to AB fish. The MIT Committee on Animal Care approved animal experimentation under protocol
517 0417-036-20. The Whitehead Institute Biosafety Committee approved of all materials under protocol HS001.
518 All experiments conform to the relevant regulatory standards.

519

520 *fam57ba*^{-/-} mutants were injected with *fam57bb* targeted sgRNA at 1 - 4 cell stage, previously described in
521 McCammon *et al.* CRISPR/Cas9 induced mutation resulted in 17 bp deletion and early stop codon.
522 Experiments were performed after 4 generations of crosses with AB controls.

523 *fam57bb* 5' to 3' TAGGTGATGTCCTGGCAGGAAG

524 *fam57bb* 3' to 5' AACCTTCCTGCCAGGACATCA

525 For genotyping, PCR amplified region of in/del. PCR was digested with Earl restriction enzyme, with
526 homozygous mutation detected by loss of Earl restriction site. *fam57b mut* line was outcrossed with AB
527 periodically to avoid chromosomal abnormalities.

528

529 **Generation and characterization of iPSC lines**

530 Unaffected control male and female iPSC lines, 599 and 657, were a generous gift from Rudolf Jaenisch,
531 originally obtained as fibroblasts from Coriell Institute Biobank. iPSC of 16p11.2 deletion carriers were obtained
532 from Simons Variation in Individuals Project (Supplemental Table 1) (Simons Vip, 2012). Cell line corresponds
533 to subjects with abbreviated ID from RUCDR. Acquisition of lines were in consideration of potential sex
534 differences in 16pdel syndrome. All iPSCs were tested for negative mycoplasma and normal karyotype.

535 Cytogenetic analysis was performed on twenty G-banded metaphase cells at Cell Line Genetics. All
536 experiments involving cells from human donors were performed in compliance with established IRB protocols
537 at the Whitehead Institute. The Whitehead Institute for Biomedical Research and MIT Biosafety Committees
538 approved safety considerations around the experiments performed.

539

540 *iPSCs* - Cells were cultured on plates coated with Matrigel (Corning #CB-40234A) in mTeSR+ media
541 (STEMCELL Technologies #85850) with pen/strep. Y27632 (STEMCELL Tech #72302) was added to cells
542 prior to passaging (single colonies), then passaged with ReLeSR or Accutase for single colonies (STEMCELL
543 Technologies #05872 and #07922). Cells were maintained at 37°C with 5% O₂.

544

545 *Generation of Cortical Neurons* - iPSCs were differentiated into neural progenitor cells (NPCs) by FGF
546 exchange. FGF was slowly removed by exchanging mTeSR+ with -FGF media (DMEM/F12/HEPES (Thermo
547 Fisher Scientific #12400024), Neurobasal (Thermo Fisher Scientific #21103049), N2 (Gibco #17502048),
548 Gem21 (GeminiBio #400-160) MEM non-essential amino acids (NEAA) (Thermo Fisher Scientific #11140050),
549 GlutaMAX (Gibco #35050061), pen/strep, D(+) Glucose and NaCl) every day over 2 - 3 weeks. When rosettes
550 were present, media was exchanged with +FGF media (DMEM/F12/HEPES, Neurobasal, N2, Gem21 -
551 Vitamin A (GeminiBio #400161), MEM NEAA, GlutaMAX, pen/strep, Beta-Mercaptoethanol (Sigma-Aldrich
552 #M3148) and 4 ng/mL FGF (Peprotech #100-18B)) plus 2.5 µM/mL dorsomorphin (Tocris #3093). Cells were
553 incubated with Y27632 before passaging with Accutase, expanded and passaged at least 3 times until
554 homogeneous NPC culture. NPCs were passaged on poly-D-lysine (Thermo Fisher Scientific #A3890401) and
555 laminin (Sigma-Aldrich #L2020) coated plates for cortical neuron differentiation. NPCs media was exchanged
556 with Neuronal Differentiation media (Neurobasal, GlutaMAX, NEAA, D(+) Glucose, Gem21, Culture One
557 (Gibco #A3320201), 5 µg/mL BDNF and GDNF (Peprotech #450-02 & #450-10), pen/strep) for 1 month,
558 changing media every 2 to 3 days. Cells were maintained at 37°C under normoxic conditions.

559

560 **Generation and characterization of SH-SY5Y Neuroblastoma cell line**

561 SH-SY5Y cells, originally from ATCC, were a kind gift from David Bartel, Whitehead Institute for Biomedical
562 Research. Cells were maintained in EMEM (ATCC # 30-2003), F12 (ATCC # 30-2006) media supplemented
563 with fetal bovine serum (FBS Sigma-Aldrich #12306C) and pen/strep in a 37°C incubator with 5% CO₂.
564 Differentiation of cells to neuronal model were induced with media containing Neurobasal, Gem21, GlutaMAX,
565 All-trans-retinoic acid (Sigma-Aldrich # R2625) and pen/strep, for 4 days in dark to prevent retinoic acid
566 degradation from light exposure (Kovalevich J, 2013).

567

568 CRISPR sgRNA designs were identified from Target Guide Sequence Cloning Protocol, Zhang lab, with
569 sequence overlapping the TLC domain of FAM57B (Cong et al., 2013). 10 targeted guides to FAM57B
570 sequence were individually transformed in pLC OPTI-Stuffer plasmid, a kind gift from David Sabatini,

571 Whitehead Institute, and lentivirus was grown in HEK293T cells. Generation of CRISPR/Cas9 induced
572 mutations via lentiviral transduction was performed according to Wang *et al.* protocols (Wang et al., 2014,
573 Wiles et al., 2015). After puromycin selection, cells were gently triturated and diluted to approximate 1 cell per
574 well in 96 well plate. Wildtype cells were simultaneously single cell diluted and sorted to serve as additional
575 control for experiments. Incorporation of mutation was determined by Next Generation Sequencing.

576

577 FAM57B Homozygote deletion (KO)

578 sgFAM57B1 5' to 3' - GGTGCTCCACCATGCCGCCA

579 Mutation resulted in frameshift with 111 and 121 bp deletion on either strand, resulting in early stop codon.

580

581 FAM57B Heterozygote deletion (HET)

582 sgFAM57B2 5' to 3' - GGGCACAGCAAATTGCGTGT

583 Mutation resulted in frameshift with 20 bp deletion on one strand, resulting in early stop codon.

584

585 Adeno-Associated Virus Integration Site 1 (AAVS1) targeted control

586 sgAAVS1 5' to 3' - CACCGGGGCCACTAGGGACAGGAT

587 Mutation resulted in frameshift and 51 bp and 1 bp deletion on either stand, resulting in early stop codon. The

588 AAVS1 served as a control for all SH-SY5Y experiments. WT was compared to AAVS1 to determine

589 confidence of statistical significance when compared to *FAM57B HET* and *KO*.

590 **Method Details**

591 **HEK293 Cell Culture and Co-Immunoprecipitation**

592 HEK293T cells were cultured in Dulbecco's modified Eagle's medium (Corning # MT15017CV) supplemented
593 with 10% fetal bovine serum, 100 IU/ml penicillin, 100 µg/ml streptomycin, and 110 µg/ml sodium pyruvate.

594 Transfections were performed with the polyethylenimine reagent (Sigma-Aldrich # 08719) using 8 µg of
595 plasmid per 10 cm culture dish for 36–48 h; medium was exchanged after 6 hours. pcDNA3.1 was used as a
596 control. DYKDDDDK (Flag)-tagged human FAM57B plasmid (sequence sent to Genscript and available upon
597 request). Hemagglutinin (HA)-tagged human CerS plasmids were generated as described (Laviad et al., 2012).

598 Co-immunoprecipitation was performed using cells transfected with a variety of plasmids in pcDNA3.1-C-DYK
599 or pcDNA3.1-C-HA (sequences sent to Genscript and available upon request). HA-tagged CerSs were used to
600 confirm non-specific binding to Flag affinity resins. Cells were washed twice with cold PBS and lysed in lysis
601 buffer (20 mM Tris (pH 7.4), 150 mM NaCl, 1 mM EDTA, 1% NP-40, 5% glycerol and protease inhibitor
602 (Sigma-Aldrich #200-664-3)). Lysates were incubated on ice for 10-15 min. Protein was determined using the
603 BCA reagent. FLAG-tagged human FAM57B, using an anti-FLAG affinity resin (Genscript #L00432). Lysates

604 were incubated with 40 μ l of beads overnight at 4°C with rotation. The resin was washed three times in 1 ml of
605 lysis buffer at 4°C with rotation. Proteins were eluted using 4X SDS sample buffer (BioRad #161-0747). Eluted
606 proteins were analyzed by Western blotting for detection of HA-tagged interacting proteins.

607

608 **Ceramide Synthase Assays**

609 Cell homogenates were prepared in 20 mM HEPES-KOH, pH 7.2, 25 mM KCl, 250 mM sucrose, and 2 mM
610 MgCl₂ containing a protease inhibitor mixture. Protein was determined using the BCA reagent (Thermo Fisher
611 Scientific). Samples were incubated with 15 μ M NBD-sphinganine (Avanti Polar Lipids # 810206P), 20 μ M
612 defatted BSA (Sigma-Aldrich #10775835001), and 50 μ M 16- or 24-fatty acyl-CoA (Avanti Polar Lipids 870743
613 & 870725) in a 20 μ l reaction volume. CerS (40 μ g protein, 25 min reaction time) was assayed using C24.1-
614 CoA and Cer5/6 (5 μ g protein, 5 min reaction time) assayed using C16-CoA. Reactions were terminated by
615 chloroform/methanol (1:2, v/v) and lipids extracted. Lipids were dried under N₂, resuspended in
616 chloroform/methanol (9:1, v/v), and separated by thin layer chromatography using chloroform/methanol, 2M
617 NH₄OH (40:10:1, v/v/v) as the developing solvent. NBD-labeled lipids were visualized using an Amersham
618 Typhoon5 imager and quantified by ImageQuantTL (GE Healthcare, Chalfont St Giles, UK). All solvents were
619 of analytical grade and were purchased from Bio-Lab (Jerusalem, Israel).

620

621 **Sample Collection for Lipidomics**

622 *iPSC Differentiated Neurons* - 2 x 10⁶ NPCs were plated in 6 well plate, at least 3 wells per genotype.
623 Differentiation to cortical neurons was performed as stated above. Cells were washed with phosphate buffered
624 saline solution (PBS). Cells were scraped in LC grade methanol and homogenized in eppendorf tube
625 containing water and LC grade chloroform with pestle mixer, followed by vortexing for 10 minutes at 4°C. Lipids
626 were separated by centrifuging top speed at 4°C. This was repeated three times with all samples run together
627 in positive ion mode for lipidomic analysis. Raw data are provided in the Supplemental Excel spreadsheet.

628

629 *Differentiated SH-SY5Y Cells* - 1 x 10⁶ cells were plated per well in 6 well plate, 3 wells per genotype. SH cells
630 were differentiated over 4 days in media containing retinoic acid. Cells were washed with PBS. Cells were
631 scraped in LC grade methanol and homogenized in eppendorf tube containing water and LC grade chloroform
632 with pestle mixer, followed by vortexing for 10 minutes at 4°C. Lipids were separated by centrifuging top speed
633 at 4°C. This was repeated twice with all samples run together. Raw data are provided in the Supplemental
634 Excel spreadsheet.

635

636 *Zebrafish* - At 7 dpf, larvae were deeply anesthetized with tricaine. Brains with surrounding epidermal layer
637 were dissected and flash frozen on dry ice. Collections were pooled at 20 brains per sample. Brains were
638 homogenized in an eppendorf tube with a pestle mixer in LC grade methanol, LC grade chloroform and water,
639 followed by vortexing for 10 minutes at 4°C. Lipids were separated by centrifuging top speed at 4°C. Brains

640 were collected and stored in -80°C over many dissections to acquire adequate tissue for analysis. This was
641 repeated twice per genotype, with the *fam57b mut* and AB cohort run at different times (different normalization)
642 while the *fam57b het* and AB cohort were run at the same time. Raw data are provided in the Supplemental
643 Excel spreadsheet.

644

645 **Untargeted Lipidomics**

646 Lipids were separated on an Ascentis Express C18 2.1 x 150 mm 2.7 μm column (Sigma-Aldrich) connected to
647 a Vanquish Horizon UPLC system and an ID-X tribrid mass spectrometer (Thermo Fisher Scientific) equipped
648 with a heated electrospray ionization (HESI) probe. External mass calibration was performed using the
649 standard calibration mixture every seven days. Dried lipid extracts were reconstituted in 50 μL 65:30:5
650 acetonitrile: isopropanol: water (v/v/v). Typically, 2 μL of sample were injected onto the column, with separate
651 injections for positive and negative ionization modes. Mobile phase A in the chromatographic method consisted
652 of 60:40 water: acetonitrile with 10 mM ammonium formate and 0.1% formic acid, and mobile phase B
653 consisted of 90:10 isopropanol: acetonitrile, with 10 mM ammonium formate and 0.1% formic acid. The
654 chromatographic gradient was adapted from Hu *et al.* 2008 (Hu et al., 2008) and Bird *et al.* 2011 (Bird et al.,
655 2011). Briefly, the elution was performed with a gradient of 40 min; during 0–1.5 min isocratic elution with 32%
656 B; from 1.5 to 4 min increase to 45% B, from 4 to 5 min increase to 52% B, from 5 to 8 min to 58% B, from 8 to
657 11 min to 66% B, from 11 to 14 min to 70% B, from 14 to 18 min to 75% B, from 18 to 21 min to 97% B, during
658 21 to 35 min 97% B is maintained; from 35–35.1 min solvent B was decreased to 32% and then maintained for
659 another 4.9 min for column re-equilibration. The flow rate was set to 0.260 mL/min. The column oven and
660 autosampler were held at 55°C and 15°C , respectively. The mass spectrometer parameters were as follows:
661 The spray voltage was set to 3.25 kV in positive mode and 3.0 kV in negative mode, and the heated capillary
662 and the HESI were held at 300°C and 375°C , respectively. The S-lens RF level was set to 45, and the sheath
663 and auxiliary gas were set to 40 and 10 units, respectively. These conditions were held constant for both
664 positive and negative ionization mode acquisitions. The mass spectrometer was operated in full-scan-
665 ddMS/MS mode with an orbitrap resolution of 120,000 (MS1) and 30,000 (MS/MS). Internal calibration using
666 Easy IC was enabled. Quadrupole isolation was enabled, the AGC target was 1×10^5 , the maximum injection
667 time was 50 msec, and the scan range was $m/z = 200\text{--}2000$. For data-dependent MS/MS, the cycle time as 1.5
668 sec, the isolation window was 1, and an intensity threshold of 1×10^3 was used. HCD fragmentation was
669 achieved using a step-wise collision energy of 15, 25, and 35 units, and detected in the orbitrap with an AGC
670 target of 5×10^4 and a maximum injection time of 54 msec. Isotopic exclusion was on, a dynamic exclusion
671 window of 2.5 sec was used, and an exclusion list was generated using a solvent bank.

672

673 High-throughput annotation and relative quantification of lipids was performed using LipidSearch v4.2.21
674 (Thermo Fisher Scientific/ Mitsui Knowledge Industries) using the HCD database (Taguchi and Ishikawa, 2010,
675 Yamada et al., 2013). LipidSearch matches MS/MS data in the experimental data with spectral data in the

676 HCD database. Precursor ion tolerance was set to 5 ppm, product ion tolerance was set to 10 ppm.
677 LipidSearch nomenclature uses underscores to separate the fatty acyl chains to indicate the lack of *sn*
678 positional information. In cases where there is insufficient MS/MS data to identify all acyl chains, only the sum
679 of the chains is displayed. Following the peak search, positive and negative mode data were aligned together
680 where possible and raw peak areas for all annotated lipids were exported to Microsoft Excel and filtered
681 according to the following predetermined quality control criteria: Rej (“Reject” parameter calculated by
682 LipidSearch) equal to 0; PQ (“Peak Quality” parameter calculated by LipidSearch software) greater than 0.75;
683 CV (standard deviation/ mean peak area across triplicate injections of a represented (pooled) biological
684 sample) below 0.4; R (linear correlation across a three-point dilution series of the representative (pooled)
685 biological sample) greater than 0.9. Typically, ~70% of annotated lipids passed all four quality control criteria.
686 Redundant lipid ions (those with identical retention times and multiple adducts) were removed such that only
687 one lipid ion per species/ per unique retention time is reported in merged alignments. For data where positive
688 and negative mode data were aligned separately some redundancies may still exist. Raw peak areas of the
689 filtered lipids were normalized to total lipid signal (positive or negative ionization mode) in each sample to
690 control for sample loading. Data presented are shown as Log₂FC compared to wildtype/control samples.
691 Statistics were performed in Prism, with each run analyzed separately.

692

693 **Zebrafish brain staining and imaging**

694 For 24 hours post-fertilization staining, embryos were deeply anesthetized in tricaine after being
695 dechorionated. Embryos were placed into wells in 1% Agarose dishes. 1 ng Cholera Toxin subunit B (CT-B)
696 (Recombinant Alexa Fluor 488 conjugate, Invitrogen #C34775) was injected into the hindbrain ventricle
697 (Gutzman and Sive, 2010). Embryos were washed with E3 and incubated for 1 hour to allow CT-B binding.
698 Embryos were then fixed in fresh 4% PFA in phosphate buffered solution (PBS) overnight at 4°C. Embryos
699 were washed in PBS + Tween-20 (PBT) and incubated with 555-Phalloidin (Invitrogen #A34055) for 1 hour.
700 Alternatively, embryos were incubated with Duramycin-Cy3 conjugate (Molecular Targeting Technologies #D-
701 1006) PE stain for 45 min. Embryos were washed in PBT and mounted in DAPI Antifade (Thermo Fisher
702 Scientific #P36931) overnight. Imaging was performed on an inverted Zeiss LSM700 Laser Scanning Confocal
703 and processed on Fiji (ImageJ). CT-B and Duramycin images were processed on ImageJ to measure relative
704 puncta from staining. Particles were measured after drawing a size circle in each hemisphere comparing AB to
705 *fam57b mut* embryos. Threshold was set to intermodes to assume for bimodal histogram, particle size set
706 between 0 – 2 μm^2 .

707

708 For 7 dpf larvae, the following protocol was adapted from the mouse protocol provided by LifeCanvas
709 Technologies (SHIELD kit, LifeCanvas Technologies). At 7 dpf, zebrafish were collected into Eppendorf tubes,
710 25 zebrafish per tube and anesthetized on ice briefly. Embryo buffer E3 was removed and replaced with 1 mL
711 SHIELD Perfusion Solution with diluted 4% paraformaldehyde (PFA) (Electron Microscopy Sciences # 50-980-

712 495), shaking overnight at 4°C. Whole zebrafish brains were dissected the next day and placed into tubes with
713 fresh SHIELD Perfusion Solution, shaking overnight at 4°C. Tissue was placed into 1 mL SHIELD OFF
714 solution, shaking overnight at 4°C. Tissue was transferred into SHIELD ON Buffer, shaking overnight at 37°C
715 in MaxQ 4450 (ThermoFisher Scientific). Tissue was then cleared with 1 mL passive clearing protocol using
716 SDS Clearing Solution, shaking for 5 days at 45°C. Clearing solution was washed off with 1 mL PBS + 1%
717 Triton-X (PBT) with 0.02% Sodium Azide 3 times over 24 hours shaking at 37°C. Tissue was blocked in 1 mL
718 PBT + 1% BSA for 2 hours shaking at room temperature, then incubated in primary antibody, shaking
719 overnight at 4°C. Antibodies: 1:100 Synaptotagmin-1 (Lifespan Bioscience # LS-B12889), GAD65 + GAD67
720 (Abcam #ab11070), Beta-Actin (Proteintech 60008-1) and 1:500 DAPI (Life Technologies # D1306), in 0.5 mL
721 PBT + 1% BSA. Primary antibody was washed off 3 times in PBT and incubated in secondary antibody
722 shaking overnight at 4 °C (1:500 488- 555- 680-conjugated antibodies (Jackson ImmunoResearch, 488 anti-
723 goat #805-545-180, 488 anti-mouse #715-545-151, 594 anti-mouse #715-585-150, 594 anti-rabbit #711-585-
724 152, anti-mouse 680 #715-625-150, anti-rabbit 680 #711-625-152) in PBT + 1% BSA). Secondary was washed
725 off 3 times in PBT, then 1 mL EasyIndex was added to tissue, shaking overnight at room temperature. Whole
726 brains were mounted in fresh EasyIndex on slides, placing coverslip with vacuum grease. Imaging was
727 performed on an inverted Zeiss LSM700 Laser Scanning Confocal and processed on Fiji (ImageJ).

728

729 **PH3 Staining**

730 At 24 hpf, *fam57b mut* and AB embryos were dechorionated and fixed overnight at 4°C in paraformaldehyde.
731 Embryos were washed with phosphate buffered saline with Tween-20 (PBT) and yolk sac was removed.
732 Embryos were incubated with 10% H₂O₂ for 1.5 hrs, then washed in PBT. Embryos were blocked in PBT with
733 bovine serum albumin at room temp for 4 hrs, then incubated with α-PH3 antibody (1:1000, Upstate
734 Biotechnology #06-570) overnight at room temp. Embryos were washed with PBT and incubated with
735 secondary antibody (1:500 goat α-rabbit IgG HRP, Invitrogen #31460) in PBT overnight at room temp.
736 Embryos were washed in PBT and flat mounted on glass slide with propidium iodide in glycerol. Imaging was
737 performed on a confocal microscope.

738

739 **TUNEL Staining**

740 Embryos were collected, fixed and processed as PH3 staining. Embryos were then dehydrated then
741 rehydrated interchanging ethanol and PBT. Proteinase K (Invitrogen # EO0491) was incubated in PBT on
742 neutator, then rinsed in PBT. TdT labeling was followed per manufacturer's instructions, ApopTag kit
743 (Chemicon # S7101). α-DIG (1:100, Thermo Fisher Scientific #700772) was used to detect the DIG labeled
744 ends. Embryos were washed in PBT and flat mounted on glass slide with propidium iodide in glycerol. Imaging
745 was performed on a confocal microscope.

746

747 **Immunocytochemistry**

748 Patient derived neurons were washed with PBS and fixed in fresh 4% paraformaldehyde in PBS overnight
749 rocking at 4°C. Cells were washed with PBT and blocked in PBT + BSA for 1 hour at room temperature.
750 Primary antibody was added to PBT overnight rocking at 4°C. Antibodies: 1:100 Syt-1, Vesicular Glutamate 1
751 and 2 (VGlut1/2, Synaptic Systems #135503) or Postsynaptic Density 95 (PSD95, Abcam #ab18258),
752 Acetylated-Tubulin (Ac-Tubulin, Abcam #ab179513). Primary antibody was washed off 3 times in PBT and
753 incubated in secondary antibody shaking overnight at 4°C (1:500 488- 555- 680- (Jackson ImmunoResearch,
754 see above) in PBT + BSA). Secondary was washed off 3 times in PBT, rocking for 2 hours at room
755 temperature. Cells were washed 3 times in PBT and mounted on slides with DAPI (Prolong Gold Antifade with
756 DAPI (Life Technologies #P36935). Imaging was performed on an inverted Zeiss LSM700 Laser Scanning
757 Confocal and processed on Fiji (ImageJ).

758

759 SH-SY5Y cells were plated on coverslips and differentiated over 4 days with retinoic acid medium. The same
760 imaging protocol was performed as above. Antibody: 1:200 Beta-Actin.

761

762 **Western blot**

763 HEK293T studies - Proteins were separated by SDS-PAGE and transferred to nitrocellulose membranes. HA-
764 tagged constructs were identified using antibodies against HA or Flag peptides (1:5,000, Abcam #ab18181,
765 #ab1162), and goat anti-rabbit or mouse horseradish peroxidase (1:10,000, #323-001-021, #223-005-024)
766 were used as secondary antibodies (Jackson). Equal loading was confirmed using a mouse anti-GAPDH
767 (Abcam #ab8245). Detection was performed using the ECL detection system.

768

769 Larvae brain tissue (25 larvae brains pooled per sample) or differentiated SH-SY5Y cells (1 x 10⁶ cells per
770 sample) were washed with PBS then lysed in RIPA buffer (Thermo Fisher Scientific #89900) with protease
771 inhibitor cocktail with a pestle homogenizer. Tissue/cells were rotate at 4°C for 30 min, then spun full speed 10
772 min. The supernatant was removed containing proteins, with denature in laemmli buffer for 1 hr at RT. Protein
773 was separated on 10-40% gel and transferred PVDF by wet transfer. Membranes were blocked in 5% dry milk
774 in TBS + Tween 20. Primary antibody was incubated overnight. Same antibodies were used for
775 immunofluorescence and western analysis. Antibodies: Syt-1, Beta-Actin, GAPDH), FAM57B (Proteintech
776 20760-1-AP). Secondary antibodies 1:2000 IRDye (Li-Cor 800CW Rabbit #92632211, 680RD Mouse
777 #92668070) were incubated for 1 hour at room temperature. Blots were imaged and quantified on a Li-Cor
778 Odyssey.

779

780 **Biotinylation and MS/MS**

781 At 7 dpf, larvae were deeply anesthetized with tricaine. Larvae were dissected in PBS with protease inhibitor
782 cocktail on ice, pooling 20 brains per genotype per sample. Assay was performed according to protocol
783 utilizing Pierce Cell Surface Protein Isolation Kit (Thermo Fisher Scientific #89881) with the following

784 modifications. 1 vial of biotin was resuspended in 2 mL PBS and fresh brains were incubated with 1 mL
785 biotin/PBS solution rotating for 45 minutes at 4°C. Elution of biotin-bound proteins in water + DTT 1 hour at
786 room temperature. Eluates were reduced, alkylated and digested with trypsin at 37°C overnight. This solution
787 was subjected to solid phase extraction to concentrate the peptides and remove unwanted reagents. Solution
788 was injected onto a Waters NanoAcquity HPLC equipped with a self-packed Aeris 3.6 µm C18 analytical
789 column 0.075 mm by 20 cm, (Phenomenex). Peptides were eluted using standard reverse-phase gradients.
790 The effluent from the column was analyzed using a Thermo Orbitrap Elite mass spectrometer (nanospray
791 configuration) operated in a data dependent manner for 54 minutes. The resulting fragmentation spectra were
792 correlated against the known database using Mascot (Matrix Science). Scaffold Q+S (Proteome Software) was
793 used to provide consensus reports for the identified proteins. PEAKS Studio 8.5 was used for data analysis as
794 a supplement to Mascot. Raw data are provided in the Supplemental Excel spreadsheet.

795

796 **Synaptosome Isolation**

797 7 dpf zebrafish larvae were anesthetized in Tricaine with larvae buffer E3. Whole brains were dissected from
798 the larvae and placed into a 1.5 mL Eppendorf tube on ice, pooling 20 brains per genotype. The tissue was
799 centrifuged at top speed and excess liquid was removed. 200 µl of SYN-Per Reagent (Thermo Fisher
800 Scientific #87793) plus protease cocktail inhibitor was added to the eppendorf on ice. With a pestle, the brains
801 were homogenized with 10 strokes and the tube was gently turned 3 times to dissociate cells. The tissue was
802 centrifuged at 1200 x G for 10 min at 4°C. The supernatant was collected and added to a new Eppendorf tube.
803 The sample was centrifuged at 15,000 x G for 20 min at 4°C. The supernatant was removed and the pellet was
804 gently resuspended in 100 µl of SYN-Per Reagent plus protease cocktail inhibitor. The samples were flash
805 frozen and processed for tandem mass spectrometry (MS/MS) after reducing alkylating and digesting with
806 trypsin as indicated above.

807

808 SH-SY5Y cells were plated at 1×10^6 per well in 6 well plates, with 1 well per independent sample for all
809 genotypes. After 4 days differentiation, cells were washed with PBS plus protease inhibitor cocktail. Cells were
810 scraped in the same solution and centrifuges in 1.5 mL Eppendorf tube at top speed for 30 seconds at 4°C.
811 The wash was removed and the same protocol was used as for larval brain synaptosome isolation. After
812 samples were flash frozen, they were then processed for MS/MS or LC/MS analysis. Raw data are provided in
813 the Supplemental Excel spreadsheet.

814

815 **Electrophysiology**

816 iPSC Differentiated Neurons - 1×10^4 NPCs were plated and matured over 1 month in a PDL and Laminin
817 coated 48-well CytoView plate (Axion Biosystems # M768-tMEA-48B). Recordings of spontaneous activity
818 were taken over 10-minute periods on the Maestro system (Axion Biosystems). AxIS software compiled the
819 data collected from recordings. Data were collected for LFPs (firing frequency in Hz), electrographic burst

820 events (minimum 5 LFPs/100 ms) and relative network activity (minimum 3 LFPs detected simultaneously
821 between a minimum of two electrodes). LFP detection was filtered at $6 \times$ standard deviation to remove potential
822 artifacts. The external physiological solution contained (in mM) 128 NaCl, 5 KCl, 2 CaCl₂, 1 MgCl₂, 25 HEPES
823 and 30 glucose, pH 7.3, Osmolarity 315 - 325. The High KCl solution contained (in mM) 63 NaCl, 70 KCl, 2
824 CaCl₂, 1 MgCl₂, 25 HEPES and 30 glucose, pH 7.3, Osmolarity 315 - 325.

825

826 Live larvae MEA recordings were performed as detailed in Tomasello and Sive 2020 (Tomasello and Sive,
827 2020). For these recordings, larva was immersed in low-melt agarose in 12-well 64 electrode Cytoview plates
828 (Axion Biosystems (discontinued, recommend 6-well plate)). LFP activity was recorded for 10 minutes, noting
829 the electrodes in contact with larva head region. Larva were immediately sacrificed after recordings. Data was
830 processed with AxIS and Axion Neural Metric Tool (Axion Biosystems).

831

832 **Larval Behavior**

833 At 7 dpf, dishes containing larvae were moved to the bench to allow acclimation to RT. For experimentation,
834 only larvae with an inflated swim bladder and no other morphology phenotypes, such as a crooked tail, were
835 selected. With a cut 200 μ l tip, larvae were individually pipetted into 96-well plates with 200 μ l E3 media and
836 moved to the Noldus Daniovision for 10 min habituation period. The larvae were exposed to a testing period of
837 70 minutes, with light (at 10%) extinguished for 5 seconds at 10-minute intervals. Point tracking collected
838 distance and velocity traveled. Distance moved was calculated using the Ethovision XT 11 software from
839 Noldus. Raw data are provided in the Supplemental Excel spreadsheet.

840

841 The same method is performed as above up to habituation. Baseline activity was then recorded for 10 minutes,
842 followed by exchange of 100 μ l E3 from each well with 100 μ l of varying concentrations of PTZ to test a range
843 of doses. Plates were immediately placed back on the Daniovision system for another 10 min recording. Point
844 tracking collected distance and velocity traveled. Distance moved was calculated using the Ethovision XT 11
845 software from Noldus, normalizing to habituation time. Raw data are provided in the Supplemental Excel
846 spreadsheet.

847

848 **Neuromuscular Junction Staining**

849 Larvae were fixed in 4% PFA. Alpha-Bungarotoxin AlexaFluor 488 conjugate (Invitrogen B13422) was used at
850 1:500, znp1 (anti-SYT2, Abcam ab154035) monoclonal antibody was used at 1:200, with secondary antibody
851 594 anti-mouse (Jackson #715-585-150) incubated at 1:500.

852

853 **Larval Head Measurements**

854 Larvae were deeply anesthetized in tricaine and immersed in methylcellulose for brightfield imaging on Leica
855 microscope. Larvae were oriented for dorsal measurements of dorsal head length, hindbrain head length, inter-

856 eye width and forebrain head length, and oriented for lateral measurements of head height, lateral head length,
 857 eye width, eye height and lateral length. Schematic of measurements can be found in McCammon *et al.* 2017
 858 (McCammon *et al.*, 2017). Raw data and conversion measurements are provided in the Supplemental Excel
 859 spreadsheet.

860

861 **Quantification and statistical analysis**

862 With exception to proteomic analysis in Figures 5C, 5D, 7E, 7H, all statistical analysis was performed in Prism.
 863 Statistical test is denoted in the figure legend corresponding to the appropriate figure. Error bars represent
 864 standard error of the mean. Asterisks are defined in each relevant figure legend. On the lipidomic analysis of
 865 individual species, to correct p-values we had utilized Sidak's multiple comparisons post hoc test rather than
 866 FDR. Note, we tried FDR for multiple hypothesis correction, and that method produced similar results. GO
 867 SLIM analysis was performed with PANTHER Classification System (www.pantherdb.org) that combines
 868 genomes, gene function classifications, pathways and statistical analysis tools to enable biologists to analyze
 869 large-scale genome-wide experimental data (Mi *et al.*, 2019). For the Proteomics analysis, the peptide
 870 intensities of biological samples were analyzed with MSstats, an R package for statistical analysis, using the
 871 options FDR=0.05 or 0.01, 'removeProtein_with1Feature=TRUE' and 'fewMeasurements="remove"'. The
 872 labeled genes indicate lowest p-values. For post hoc colocalization analysis of iPSC differentiated neurons
 873 (Table 2), individual cell somas were outlined by freehand in FIJI, followed by Coloc 2 colocalization analysis
 874 between Synaptotagmin-1 (Channel 1) and PSD-95 (Channel 2). Neurons were isolated between 3 images per
 875 genotype of representative images from Supplemental Fig. 1. Similar statistics are indicated between control
 876 16pdel neurons.

877

878 **Additional Resources**

REAGENT or RESOURCE	SOURCE	IDENTIFIER
Antibodies		
Synaptotagmin-1	Lifespan Bioscience	LS-B12889
GAD65 + 67	Abcam	ab11070
Beta-Actin	Proteintech	60008-1
DAPI	Life Technologies	D1306
555-Phalloidin	Invitrogen	A34055
488 anti-goat	Jackson	805-545-180,
488 anti-mouse	Jackson	715-545-151
594 anti-mouse	Jackson	715-585-150
594 anti-rabbit	Jackson	711-585-152
anti-mouse 680	Jackson	715-625-150
anti-rabbit 680	Jackson	711-625-152
VGlut1/2	Synaptic Systems	135503
PSD95	Abcam	ab18258
Ac-Tubulin	Abcam	ab179513
Prolong Gold Antifade with DAPI	LifeTech	P36935
HA	Abcam	ab18181

FLAG	Abcam	ab1162
GAPDH	Abcam	ab8245
Li-Cor 800CW Rabbit	Li-Cor	92632211
Li-Cor 680RD Mouse	Li-Cor	92668070
Chemicals, peptides, and recombinant proteins		
Matrigel	Corning	CB-40234A
mTeSR+	STEMCELL Tech	85850
Y27632	STEMCELL Tech	72302
ReLeSR	STEMCELL Tech	05872
Accutase	STEMCELL Tech	07922
DMEM/F12/HEPES	Thermo	12400024
Neurobasal	Thermo	21103049
N2	Gibco	17502048
Gem21	GeminiBio	400-160
MEM NEAA	Thermo	11140050
GlutaMAX	Gibco	35050061
Gem21 - Vit A	GeminiBio	400161
Beta-Mercaptoethanol	Sigma	M3148
FGF	Peprtech	100-18B
BDNF	Peprtech	450-02
GDNF	Peprtech	450-10
Dorsomorphin	Tocis	3093
Poly-D-Lysine	Thermo	A3890401
Laminin	Sigma	L2020
EMEM	ATCC	30-2003
F12	ATCC	30-2006
FBS	Sigma	12306C
All-trans-RA	Sigma	R2625
DMEM	Corning	MT15017CV
Polyethylenimine	Sigma	08719
Protease Inhibitor Cocktail	Sigma	200-664-3
NBD-spinganine	Avanti Polar	810206P
BSA -FA	Sigma	10775835001
16.1 Coenzyme A	Avanti Polar	870743
24.1 Coenzyme 1	Avanti Polar	870725
CT-B	Invitrogen	C34775
Duramycin-Cy3	Molecular Targeting	D-1006
SHIELD	LifeCanvas Tech	https://lifecanvastech.com/products/shield
PFA	EMS	50-970-495
RIPA	Thermo	89900
SYN-Per	Thermo	87793
Critical commercial assays		
Pierce Cell Surface Protein Isolation Kit	Thermo	89881
Deposited data		
Lipidomics	This Paper	Supplemental Excel
Biotinylation Proteomics	This Paper	Supplemental Excel
Synaptosome Proteomics	This Paper	Supplemental Excel
Experimental models: Cell lines		
SH-SY5Y	ATCC	CRL-2266
HEK293T	ATCC	CRL-3216

IPSC 16pdel 1453	Simons VIP	SV0001453
IPSC 16pdel 1455	Simons VIP	SV0001455
IPSC 16pdel 1459	Simons VIP	SV0001459
IPSC 16pdel 1473	Simons VIP	SV0001473
IPSC 16pdel 1481	Simons VIP	SV0001481
IPSC 16pdel 1495	Simons VIP	SV0001495
IPSC 16pdel 3104	Simons VIP	SS0013104
IPSC CTR 599	Corriell Institute Biobank	AG07599
IPSC CTR 675	Corriell Institute Biobank	AG07657
Experimental models: Organisms/strains		
<i>fam57ba</i> ^{-/-} ; <i>fam57bb</i> ^{-/-} Zebrafish	This Paper	
FAM57B HET SH-SY5Y	This Paper	
FAM57B KO SH-SY5Y	This Paper	
FAM57B AAVS1 SH-SY5Y	This Paper	
Oligonucleotides		
fam57bb 5' to 3' TAGGTGATGTCCTGGCAGGAAG	This Paper	
fam57bb 3' to 5' AAACCTTCCTGCCAGGACATCA	This Paper	
sgFAM57B1 5' to 3' - GGTGCTCCACCATGCCGCCA	This Paper	
sgFAM57B2 5' to 3' - GGGCACAGCAAATTGCGTGT	This Paper	
sgAAVS1 5' to 3' - CACCGGGGCCACTAGGGACAGGAT	This Paper	
Recombinant DNA		
FAM57B-Flag	This Paper	
CerS2-HA	This Paper	
CerS5-HA	This Paper	
CerS6-HA	This Paper	
Software and algorithms		
AxIS and Neural Metric Tool	Axion Biosystems	Axionbiosystems.com
FIJI ImageJ	ImageJ	imagej.net
GO SLIM analysis	PANTHER	pantherdb.org
LipidSearch	Thermo Scientific	IQLAAEGABSFAPC MBFK
Mascot	Matrix Science	Matrixscience.com
Scaffold Q+S	Proteome Software	Proteomesoftware.com
PEAKS Studio 8.5	Bioinformatics Solutions Inc.	Bioinfor.com
EthoVision XT	Noldus	Noldus.com

879

880 **Main Figure Titles and Legends**881 **Figure 1. Augmented local field potential activity in 16pdel syndrome differentiated neuronal culture.**

882 **a)** Local Field Potential (LFP) summary analyzed by log₂ fold change between control and 16pdel patient
883 differentiated neurons. MEA activity was recorded over 30 min starting in media, followed by physiological and
884 high potassium chloride (KCl) solution. Data was summarized and pooled from 3 experiments. Control n = 9
885 (media), n = 17 (Physiological Solution), n = 13 (High KCl Solution). 16pdel n = 35 (media), n = 76

886 (Physiological Solution), n = 35 (High KCl Solution). Violin plot group analysis: Control - 16pdel 2-way ANOVA.
 887 *p ≤ 0.05. Technical experimental replicates n = 3.

888 **b)** Increased sex specific activity in female 16pdel probands drives overall increased LFPs, compared to
 889 unaffected controls. Media unaffected neurons (Control) male (♂) n = 3, Control female (♀) n = 6, 16pdel
 890 neurons (Proband) ♂ n = 21, Proband ♀ n = 14, Physiological Solution Control ♂ n = 7, Control ♀ n = 10,
 891 Proband ♂ n = 38, Proband ♀ n = 38, High KCl Solution Control ♂ n = 6, Control ♀ n = 7, Proband ♂ n = 18,
 892 Proband ♀ n = 17. Violin plot analysis: male vs female T-Test. ****p ≤ 0.0001.

893 **c)** Increased sex specific female electrogenic burst frequency analyzed by log2 fold change between 16pdel
 894 male and female Media ♂ n = 21, ♀ n = 14, Physiological Solution ♂ n = 38, ♀ n = 38, High KCl Solution ♂ =
 895 18, ♀ n = 17. Violin plot analysis: male vs female T-Test. *p ≤ 0.05. Technical experimental replicates n = 3.

896

897 **Figure 2. Significant lipid changes between control and 16pdel differentiated neurons.**

898 **a)** Total log2 fold change from normalized peak area of lipid class analysis from untargeted lipidomics. Bolded
 899 and colored indicate statistically significant changes by T-Test, p ≤ 0.05 - 0.0001.

900 AcCa acyl carnitine, AEA N-arachidonylethanolamine, Cer ceramide, ChE cholesterol ester, Co coenzyme,
 901 DG diacylglycerol, Hex1Cer hexosylceramide, LPC lysophosphatidylcholine, LPE lysophosphatidylethanolamine,
 902 LPG lysophosphatidylglycerol, MG monoacylglycerol, PA phosphatidic acid, PC phosphatidylcholine, PE
 903 phosphatidylethanolamine, PG phosphatidylglycerol, PI phosphatidylinositol, PS phosphatidylserine, SM
 904 sphingomyelin, SPH sphingosine, TG triacylglycerol.

905 **b)** Total lipid composition analysis from untargeted lipidomics between control and 16pdel neuron. Chain
 906 Length: Small 1-5, Medium 6-12, Long 13-21, Very Long 22+, and Unresolved.

907 **c – f)** Selected analysis of lipid species from untargeted lipidomics classes. Lipid Class specified for each
 908 histogram (c - phosphatidylethanolamine, d - ceramide, e - monoacylglycerol, f – triacylglycerol) normalized
 909 peak area between control (grey) and 16pdel (orange). Statistical analysis by 2-Way ANOVA, *p ≤ 0.05 **p ≤
 910 0.01, ***p ≤ 0.001, ****p ≤ 0.0001. Control n = 10, 16pdel n = 69, error bars represent SEM. TG and PE long and
 911 very long chain species not shown as no significant differences were found by ANOVA. Technical experimental
 912 replicates n = 3.

913

914 **Figure 3. FAM57B interacts with CerS but does not have CerS activity.**

915 **a)** CerS2 activity assayed using C24:1-CoA in CerS2 KO HEK293T cells. Statistical analysis by T-test *p ≤
 916 0.05, **p ≤ 0.01, error bars SEM. Technical experimental replicates n = 3.

917 **b)** (Upper) Western blot analysis of total human FAM57B-Flag and CerS2-HA after transfection in HEK293T
 918 cells. Proteins were prepared from HEK293T cells overexpressing the indicated constructs. Anti-HA and anti-
 919 Flag are indicated. (Lower) CerS2 activity assayed using C24:1-CoA in HEK293T cells. GAPDH was used as a
 920 loading control. Statistical analysis by T-test *p ≤ 0.05, **p ≤ 0.01, error bars SEM. Technical experimental
 921 replicates n = 3.

922 **c)** (Upper) Western blot analysis of total human FAM57B-Flag, CerS5-HA and CerS6-HA after transfection in
 923 HEK293T cells. Proteins were prepared from cells overexpressing the indicated constructs. Anti-HA and anti-
 924 Flag are indicated. (Lower) CerS5 and CerS6 activity was assayed using C16:0-CoA in HEK293T cells. Anti-
 925 HA and anti-Flag are indicated. GAPDH was used as a loading control. Technical experimental replicates n =
 926 4.

927 **d)** Total cell lysates were prepared from the co-transfected cells with FAM57B-Flag and CerS2, 5 or 6-HA
 928 constructs and solubilized with 1% NP-40. Total lysates (input) or proteins immuno-precipitated with anti-Flag
 929 M2 agarose (IP) were subjected to immunoblotting with anti-HA or anti-Flag antibodies. GAPDH was used as a
 930 loading control. Technical experimental replicates n = 3.

931

932 **Figure 4. Significant lipid changes in sphingolipids and glycerolipids between WT and *FAM57B* mutant**
 933 **human differentiated SH-SY5Y neuronal cells.**

934 **a - c)** Total log₂ fold change from normalized peak area of lipid class analysis from untargeted lipidomics. **a)**
 935 *FAM57B* KO – WT, **b)** *FAM57B* HET – WT, **c)** *FAM57B* KO – *FAM57B* HET. Bolded and colored indicate
 936 statistically significant changes by T-Test, p ≤ 0.05 - 0.0001. AcCa acyl carnitine, Cer ceramide, ChE
 937 cholesterol ester, CL cardiolipin, Co coenzyme, DG diacylglycerol, HexCer Hexosylceramide, LPC
 938 lysophosphatidylcholine, LPE lysophosphatidylethanolamine, MG monoacylglycerol, PC phosphatidylcholine, PE
 939 phosphatidylethanolamine, PG phosphatidylglycerol, PI phosphatidylinositol, PS phosphatidylserine, SM
 940 sphingomyelin, TG triacylglycerol. Technical experimental replicates n = 3.

941 **d – g)** Selected analysis of lipid species from untargeted lipidomics classes. Lipid Class specified for each
 942 histogram, normalized peak area between WT (black) *FAM57B* HET (orange) and *FAM57B* KO (blue).
 943 Statistical analysis by 2-Way ANOVA, *p ≤ 0.05 **p ≤ 0.01, ***p ≤ 0.001, ****p ≤ 0.0001. Color of asterisks
 944 indicate comparison between WT – HET (orange), WT – KO (blue), HET – KO (black). WT n = 3, *FAM57B*
 945 *HET* n = 3, *FAM57B* KO n = 3, error bars represent SEM. Experiment repeated twice, analysis was similar
 946 between two separate runs.

947

948 **Figure 5. *FAM57B* knockout human neurons indicate altered synaptic composition.**

949 **a-b)** Isolated synaptosome protein abundance changes between **a)** *FAM57B* KO relative to WT and **b)**
 950 *FAM57B* HET relative to WT (Log₂ Fold). Labeled and colored indicating increased (purple) or decreased
 951 (green) abundance. Only the top 20 proteins of statistical significance were labeled in **c** and analyzed in **e-g**. WT
 952 n = 8, HET n = 10, KO n = 7. Technical experimental replicates n = 3.

953 **c-e)** Gene ontology analysis of statistically significant synaptosome isolated proteins (**e**) in *FAM57B* KO
 954 relative to WT. **e)** gene ontology pie graphs of top 20 decreased protein groups of cellular components,
 955 molecular function, protein classes, biological processes and pathways. **f)** gene ontology figure legend.

956 **g)** Analysis of synaptic markers from isolated synaptosomes between all 3 genotypes. Bolded are significantly
 957 decreased protein abundance of synaptic structural and maturation proteins, and vesicle regulation machinery.

958 INA – WT - HET & WT - KO, RAB11B – HET - KO, STXBP1 – WT - HET & WT - KO, YWHAZ – WT - HET &
 959 WT - KO. 2-Way ANOVA, $p \leq 0.05 - 0.0001$.

960 **f)** Isolated synaptosome lipid abundance between *FAM57B* KO relative to WT. Bolded text and color (purple
 961 increased and green decreased) indicates statistically significant changes by T-Test, $p \leq 0.05 - 0.01$. $n = 6$ per
 962 genotype, technical experimental replicates $n = 2$. No statistically significant differences observed when
 963 comparing *FAM57B* HET to WT, nor *FAM57B* KO to *FAM57B* HET.

964

965 **Figure 6. Significant lipid changes in ceramides and glycerols between AB and *fam57b mut* brain**
 966 **tissue.**

967 **a)** Total log₂ fold change from normalized peak area of lipid class analysis from untargeted lipidomics. Bolded
 968 and colored indicate statistically significant changes by T-Test, $p \leq 0.05 - 0.0001$. AcCa acyl carnitine, Cer
 969 ceramide, ChE cholesterol ester, CL cardiolipin, Co coenzyme, DG diacylglycerol, HexCer Hexosylceramide,
 970 LPC lysophosphatidylcholine, LPE lysophosphatidylethanolamine, MG monoacylglycerol, PC
 971 phosphatidylcholine, PE phosphatidylethanolamine, PG phosphatidylglycerol, PI phosphatidylinositol, PS
 972 phosphatidylserine, SM sphingomyelin, TG triacylglycerol. Technical experimental replicates $n = 3$.

973 **b – g)** Selected analysis of lipid species from untargeted lipidomics classes. Lipid Class specified for each
 974 histogram, normalized peak area between AB (grey) and *fam57b mut* (green). Statistical analysis by 2-Way
 975 ANOVA, * $p \leq 0.05$ ** $p \leq 0.01$, *** $p \leq 0.001$, **** $p \leq 0.0001$. AB $n = 3$, *fam57b mut* $n = 3$, error bars represent
 976 SEM. Experiment repeated twice, analysis was similar between two separate runs. Individual MG species
 977 between *fam57b het* to AB n.s. Not shown for space: 2-Way ANOVA analysis of TG. Increase in TG
 978 (16:0_16:0_16:1) in *fam57b het* compared to AB ($p \leq 0.01$).

979

980 **Figure 7. Loss of *fam57b* results in altered plasma membrane architecture early in development and**
 981 **decreased localization of Synaptotagmin family at the synapse later in development.**

982 **a)** Schematic of Cholera toxin-B-488 (CT-B) injection into hindbrain ventricle of embryo and flat-mounted
 983 midbrain region for imaging at 24 hpf.

984 **b)** Representative embryo midbrain imaging and quantification of CT-B labeling of AB compared to *fam57b*
 985 *mut*. Punctate CT-B labeling (arrows), actin marker phalloidin indicates labelling of CT-B at the plasma
 986 membrane, merged with DAPI. Quantification of puncta between WT (grey) and *fam57b mut* (green) CT-B ($p \leq$
 987 0.05) T-Test. Scale bar = 5 μm . AB $n = 16$, *fam57b mut* $n = 18$. Error bars SEM, statistical analysis by T-test * p
 988 ≤ 0.05 . Technical experimental replicates $n = 4$.

989 **c)** Representative embryo midbrain imaging and quantification of duramycin-488 labeling of AB compared to
 990 *fam57b mut*. Punctate duramycin labeling (arrows), actin marker phalloidin indicates labelling of duramycin at
 991 the plasma membrane, merged with DAPI. Quantification of puncta between WT (grey) and *fam57b mut*
 992 (green) Duramycin PE staining ($p \leq 0.05$) T-Test. Scale bar = 5 μm . AB $n = 8$, *fam57b mut* $n = 8$. Error bars
 993 SEM, statistical analysis by T-test * $p \leq 0.05$. Technical experimental replicates $n = 3$.

- 994 **d)** Schematic of membrane protein labeling biotinylation assay and processing for MS/MS in 7 dpf larvae
 995 brains.
- 996 **e)** Larvae brain total plasma membrane protein abundance changes between *fam57b mut* relative to AB (Log2
 997 Fold). Statistically significant $p \leq 0.05 - 0.0001$ proteins labelled, indicating increased (purple) or decreased
 998 (green) abundance. Lowest abundance membrane protein Synaptotagmin-1a (red box). $n = 3$ per genotype.
- 999 **f)** Representative slice of 7 dpf whole larva brain mount with Sytaptotagmin-1a (green), GAD65/67 (red) and
 1000 Beta-actin (magenta). Z-stack composite image merged with DAPI. Forebrain and midbrain areas of diffused
 1001 Syt-1 localization (white arrows). Anatomical differences noted throughout brain, including (1) optic tectum and
 1002 (2) corpus cerebelli (red arrows). Scale bar = 10 μm . Technical experimental replicates $n = 2$.
- 1003 **g)** Representative 7 dpf whole brain western blot indicate no significant change in total Syt-1a protein levels
 1004 between *fam57b mut* relative to AB. Zebrafish larvae brains pooled (20 per genotype). Syt-1a protein
 1005 abundance normalized to Beta-Actin loading control, repeated twice.
- 1006 **h)** Larvae brain total isolated synaptosome protein abundance changes between *fam57b mut* relative to AB
 1007 (Log2 Fold). Statistically significant $p \leq 0.05 - 0.0001$ proteins labelled, indicating increased (purple) or
 1008 decreased (green) abundance. Low abundance Synaptotagmin-2a like protein (red box). $n = 7$ per genotype.
 1009 Technical experimental replicates $n = 2$.
- 1010 **i-j)** Gene ontology analysis of statistically significant larvae synaptosome isolated proteins (h) in *fam57b mut*
 1011 relative to AB. i) Gene ontology pie graphs of increased and decreased protein groups of cellular components,
 1012 molecular function, protein classes, biological processes and pathways. j) Gene ontology figure legend.
- 1013 **k)** Analysis of Synaptotagmin family members from larvae isolated synaptosomes. Significantly decreased
 1014 protein abundance of Syt1a and Syt2a by 2-Way ANOVA, $**p \leq 0.01$, $****p \leq 0.0001$.
- 1015 **l)** Analysis of Synaptotagmin family members from differentiated SH-SY5Y isolated synaptosomes between all
 1016 3 genotypes. Significantly decreased protein abundance of elongated ESYT1 by 2-Way ANOVA, $*p \leq 0.05$, $**p$
 1017 ≤ 0.01 . Error bars SEM.
- 1018 **m)** Analysis of synaptic markers from larvae isolated synaptosomes. Significantly decreased protein
 1019 abundance (bolded) of vesicle regulation machinery and glutamate receptor activity by 2-Way ANOVA, $**p \leq$
 1020 0.01 , $****p \leq 0.0001$.

1021

1022 **Figure 8. Decreased spontaneous brain activity and diminished behavioral response after stimuli**
 1023 **presentation in *fam57b* mutants.**

- 1024 **a-b)** Local field potential (LFP) recordings in unanesthetized live larvae at 7dpf. Brain localized LFP recordings
 1025 were pooled for each larva. **a)** Decreased average number, mean rate and inter-LFP interval (ILI) coefficient of
 1026 variation of LFP in *fam57b mut* compared to AB (orange). No electrographic burst activity was identified in
 1027 *fam57b mut* at standard 5 LFPs/100 ms. Decreased electrographic burst parameters, including duration,
 1028 number of LFPs per burst, frequency and percentage at 3 LFPs/200ms (blue). **b)** Decreased average
 1029 electrographic burst network activity and frequency, defined as a minimum of 3 electrographic bursts between

1030 2 electrodes simultaneously, in *fam57b mut* compared to AB (grey). AB n = 21, *fam57b mut* n = 24 over 6
1031 experiments. Statistical significance by unpaired T-test, *p ≤ 0.05, **p ≤ 0.01, *** p ≤ 0.001, **** p ≤ 0.0001.
1032 Technical experimental replicates n = 7.

1033 **c)** Representative LFP waveform in brain region, indicating smaller relative waveform in *fam57b mut* compared
1034 to AB.

1035 **d)** Representative LFP raster plot over experimental time frame, indicating less overall activity in *fam57b mut*
1036 compared to AB.

1037 **e)** Representative image of 7 dpf immersed in cooled agarose in contact with electrodes on 12-well CytoView
1038 MEA plate.

1039 **f)** Startle response behavioral assay. Light source was removed for 5 secs at 10 min intervals. Mean distance
1040 reported from tracked movement during 70 min assay. Decreased light startle response identified in *fam57b*
1041 *mut* compared to AB. Statistical analysis of each startle response by T-test **p ≤ 0.01, *** p ≤ 0.001, **** p ≤
1042 0.0001. Error bars SEM. No overall significant change in movement outside of the startle identified. AB n =
1043 125, *fam57b mut* n = 33 over 5 experiments.

1044 **g)** Seizure response behavioral assay. Normalized (baseline recording subtracted) mean distance from tracked
1045 movement after absence or presence of pentylenetetrazol (PTZ) 0.5 mM and 5 mM. Significantly increased
1046 seizure-induced movement observed at 5mM in AB, while increased movement observed at 0.5 and 5 mM
1047 PTZ in *fam57b mut*. Diminished overall seizure-induced movement at 5 mM in *fam57b mut* compared to AB.
1048 Relative fold change compared to absence of PTZ indicated below histogram. Statistical analysis of each
1049 condition by T-test *p ≤ 0.05, *** p ≤ 0.001, **** p ≤ 0.0001. Error bars SEM. AB n = 166 (0 mM), 92 (0.5 mM),
1050 150 (5 mM), *fam57b mut* n = 91 (0 mM), 70 (0.5 mM), 56 (5mM) over 6 experiments.

1051 **h)** Model proposing role of Fam57b activity in the brain. Loss of function in *fam57b* mutants indicate significant
1052 changes in plasma membrane lipid groups alter architecture of plasma membrane early the developing brain.
1053 Architectural changes indicated by increased lipid raft abundance and aggregation. Altered plasma membrane
1054 homeostasis results in mis-localization of synaptic proteins, including synaptotagmins, after maturation.
1055 Decreased spontaneous brain and network activity suggests diminished synaptic function and developed
1056 circuits. As evidence has suggested spontaneous network activity shapes synaptic development, this cycles
1057 back to declined neuronal maturation and circuitry. Molecular changes to synaptic function and decreased
1058 spontaneous brain activity translate to altered behavioral response after stimuli presentation.

1059

1060 **Supplementary Tables**

1061 **Table S1 - iPSC Neuron Lipidomics. Related to figure 2.**

1062 **Table S2 - SH-SY5Y Lipidomics. Related to figure 4.**

1063 **Table S3 - SH-SY5Y Synaptosome MSMS. Related to figure 5.**

1064 **Table S4 - SH-SY5Y Synaptosome Lipidomics. Related to figure 5.**

1065 **Table S5 - Zebrafish Larvae Brain Lipidomics. Related to figure 7.**

- 1066 **Table S6 - Zebrafish Larvae Brain Biotinylation MSMS. Related to figure 7.**
 1067 **Table S7 - Zebrafish Larvae Brain Synaptosome MSMS. Related to figure 7.**
 1068 **Table S8 - Zebrafish Larvae Head and Body Measurements. Related to figure 8.**
 1069 **Table S9 - Zebrafish Larvae Light Startle Response Data. Related to figure 8.**
 1070 **Table S10 - Zebrafish Larvae Seizure Assay Data. Related to figure 8.**

1071

1072 **References**

- 1073 ANTONIO, C., FERBY, I., WILHELM, H., JONES, M., KARSENTI, E., NEBRED, A. R. & VERNOS, I.
 1074 2000. Xkid, a chromokinesin required for chromosome alignment on the metaphase plate. *Cell*, 102, 425-
 1075 35.
 1076 ARBOGAST, T., OUAGAZZAL, A. M., CHEVALIER, C., KOPANITSA, M., AFINOWI, N.,
 1077 MIGLIAVACCA, E., COWLING, B. S., BIRLING, M. C., CHAMPY, M. F., REYMOND, A. &
 1078 HERAULT, Y. 2016. Reciprocal Effects on Neurocognitive and Metabolic Phenotypes in Mouse Models
 1079 of 16p11.2 Deletion and Duplication Syndromes. *PLoS Genet*, 12, e1005709.
 1080 BAKER, K., GORDON, S. L., MELLAND, H., BUMBAK, F., SCOTT, D. J., JIANG, T. J., OWEN, D.,
 1081 TURNER, B. J., BOYD, S. G., ROSSI, M., AL-RAQAD, M., ELPELEG, O., PECK, D., MANCINI, G.
 1082 M. S., WILKE, M., ZOLLINO, M., MARANGI, G., WEIGAND, H., BORGGRAEFE, I., HAACK, T.,
 1083 STARK, Z., SADEDIN, S., BROAD CENTER FOR MENDELIAN, G., TAN, T. Y., JIANG, Y., GIBBS,
 1084 R. A., ELLINGWOOD, S., AMARAL, M., KELLEY, W., KURIAN, M. A., COUSIN, M. A. &
 1085 RAYMOND, F. L. 2018. SYT1-associated neurodevelopmental disorder: a case series. *Brain*, 141, 2576-
 1086 2591.
 1087 BARABAN, S. C., TAYLOR, M. R., CASTRO, P. A. & BAIER, H. 2005. Pentylentetrazole induced changes
 1088 in zebrafish behavior, neural activity and c-fos expression. *Neuroscience*, 131, 759-68.
 1089 BAYES, A., COLLINS, M. O., REIG-VIADER, R., GOU, G., GOULDING, D., IZQUIERDO, A.,
 1090 CHOUDHARY, J. S., EMES, R. D. & GRANT, S. G. 2017. Evolution of complexity in the zebrafish
 1091 synapse proteome. *Nat Commun*, 8, 14613.
 1092 BAZINET, R. P. & LAYE, S. 2014. Polyunsaturated fatty acids and their metabolites in brain function and
 1093 disease. *Nat Rev Neurosci*, 15, 771-85.
 1094 BEER, K. B., RIVAS-CASTILLO, J., KUHN, K., FAZELI, G., KARMANN, B., NANCE, J. F., STIGLOHER,
 1095 C. & WEHMAN, A. M. 2018. Extracellular vesicle budding is inhibited by redundant regulators of TAT-
 1096 5 flippase localization and phospholipid asymmetry. *Proc Natl Acad Sci U S A*, 115, E1127-E1136.
 1097 BLACKMON, K., THESEN, T., GREEN, S., BEN-AVI, E., WANG, X., FUCHS, B., KUZNIECKY, R. &
 1098 DEVINSKY, O. 2018. Focal Cortical Anomalies and Language Impairment in 16p11.2 Deletion and
 1099 Duplication Syndrome. *Cereb Cortex*, 28, 2422-2430.
 1100 BOONEN, M., STAUDT, C., GILIS, F., OORSCHOT, V., KLUMPERMAN, J. & JADOT, M. 2016. Cathepsin
 1101 D and its newly identified transport receptor SEZ6L2 can modulate neurite outgrowth. *J Cell Sci*, 129,
 1102 557-68.
 1103 CHEN, Y., YANG, Z., MENG, M., ZHAO, Y., DONG, N., YAN, H., LIU, L., DING, M., PENG, H. B. & SHAO,
 1104 F. 2009. Cullin mediates degradation of RhoA through evolutionarily conserved BTB adaptors to control
 1105 actin cytoskeleton structure and cell movement. *Mol Cell*, 35, 841-55.
 1106 CHIOCCHETTI, A. G., HASLINGER, D., STEIN, J. L., DE LA TORRE-UBIETA, L., COCCHI, E.,
 1107 ROTHAMEL, T., LINDLAR, S., WALTES, R., FULDA, S., GESCHWIND, D. H. & FREITAG, C. M.
 1108 2016. Transcriptomic signatures of neuronal differentiation and their association with risk genes for
 1109 autism spectrum and related neuropsychiatric disorders. *Transl Psychiatry*, 6, e864.
 1110 D'ANGELO, D., LEBON, S., CHEN, Q. X., MARTIN-BREVET, S., SNYDER, L. G., HIPPOLYTE, L.,
 1111 HANSON, E., MAILLARD, A. M., FAUCETT, W. A., MACE, A., PAIN, A., BERNIER, R.,

- 1112 CHAWNER, S. J. R. A., DAVID, A., ANDRIEUX, J., AYLWARD, E., BAUJAT, G., CALDEIRA, I.,
 1113 CONUS, P., FERRARI, C., FORZANO, F., GERARD, M., GOIN-KOCHEL, R. P., GRANT, E.,
 1114 HUNTER, J. V., ISIDOR, B., JACQUETTE, A., JONCH, A. E., KEREN, B., LACOMBE, D., LE
 1115 CAIGNEC, C., MARTIN, C. L., MAANNIK, K., METSPALU, A., MIGNOT, C., MUKHERJEE, P.,
 1116 OWEN, M. J., PASSEGGERI, M., ROORYCK-THAMBO, C., ROSENFELD, J. A., SPENCE, S. J.,
 1117 STEINMAN, K. J., TJERNAGEL, J., VAN HAELST, M., SHEN, Y. P., DRAGANSKI, B., SHERR, E.
 1118 H., LEDBETTER, D. H., VAN DEN BREE, M. B. M., BECKMANN, J. S., SPIRO, J. E., REYMOND,
 1119 A., JACQUEMONT, S., CHUNG, W. K., CHILDREN, C. U. E., CONSORTIUM, P. E. & PROJE, S. V.
 1120 I. 2016. Defining the Effect of the 16p11.2 Duplication on Cognition, Behavior, and Medical
 1121 Comorbidities. *Jama Psychiatry*, 73, 20-30.
- 1122 DE ANDA, F. C., ROSARIO, A. L., DURAK, O., TRAN, T., GRAFF, J., MELETIS, K., REI, D., SODA, T.,
 1123 MADABHUSHI, R., GINTY, D. D., KOLODKIN, A. L. & TSAI, L. H. 2012. Autism spectrum disorder
 1124 susceptibility gene TAOK2 affects basal dendrite formation in the neocortex. *Nat Neurosci*, 15, 1022-31.
- 1125 DESHPANDE, A., YADAV, S., DAO, D. Q., WU, Z. Y., HOKANSON, K. C., CAHILL, M. K., WIITA, A. P.,
 1126 JAN, Y. N., ULLIAN, E. M. & WEISS, L. A. 2017. Cellular Phenotypes in Human iPSC-Derived Neurons
 1127 from a Genetic Model of Autism Spectrum Disorder. *Cell Rep*, 21, 2678-2687.
- 1128 DZIOBEK, I., GOLD, S. M., WOLF, O. T. & CONVIT, A. 2007. Hypercholesterolemia in Asperger syndrome:
 1129 independence from lifestyle, obsessive-compulsive behavior, and social anxiety. *Psychiatry Res*, 149,
 1130 321-4.
- 1131 EGAWA, J., WATANABE, Y., WANG, C., INOUE, E., SUGIMOTO, A., SUGIYAMA, T., IGETA, H.,
 1132 NUNOKAWA, A., SHIBUYA, M., KUSHIMA, I., ORIME, N., HAYASHI, T., OKADA, T., UNO, Y.,
 1133 OZAKI, N. & SOMEYA, T. 2015. Novel rare missense variations and risk of autism spectrum disorder:
 1134 whole-exome sequencing in two families with affected siblings and a two-stage follow-up study in a
 1135 Japanese population. *PLoS One*, 10, e0119413.
- 1136 EGOLF, L. E., VAKSMAN, Z., LOPEZ, G., ROKITA, J. L., MODI, A., BASTA, P. V., HAKONARSON, H.,
 1137 OLSHAN, A. F. & DISKIN, S. J. 2019. Germline 16p11.2 Microdeletion Predisposes to Neuroblastoma.
 1138 *Am J Hum Genet*, 105, 658-668.
- 1139 FERNANDEZ, I., ARAC, D., UBACH, J., GERBER, S. H., SHIN, O., GAO, Y., ANDERSON, R. G., SUDHOF,
 1140 T. C. & RIZO, J. 2001. Three-dimensional structure of the synaptotagmin 1 C2B-domain: synaptotagmin
 1141 1 as a phospholipid binding machine. *Neuron*, 32, 1057-69.
- 1142 FETIT, R., PRICE, D. J., LAWRIE, S. M. & JOHNSTONE, M. 2020. Understanding the clinical manifestations
 1143 of 16p11.2 deletion syndrome: a series of developmental case reports in children. *Psychiatr Genet*, 30,
 1144 136-140.
- 1145 FISHMAN, P. H., MOSS, J. & OSBORNE, J. C., JR. 1978. Interaction of cholera toxin with the oligosaccharide of
 1146 ganglioside GM1: evidence for multiple oligosaccharide binding sites. *Biochemistry*, 17, 711-6.
- 1147 FLIS, V. V. & DAUM, G. 2013. Lipid transport between the endoplasmic reticulum and mitochondria. *Cold
 1148 Spring Harb Perspect Biol*, 5.
- 1149 GIANNUZZI, G., SCHMIDT, P. J., PORCU, E., WILLEMIN, G., MUNSON, K. M., NUTTLE, X., EARL, R.,
 1150 CHRAST, J., HOEKZEMA, K., RISSO, D., MANNIK, K., DE NITTIS, P., BARATZ, E. D., P. C.,
 1151 HERAULT, Y., GAO, X., PHILPOTT, C. C., BERNIER, R. A., KUTALIK, Z., FLEMING, M. D.,
 1152 EICHLER, E. E. & REYMOND, A. 2019. The Human-Specific BOLA2 Duplication Modifies Iron
 1153 Homeostasis and Anemia Predisposition in Chromosome 16p11.2 Autism Individuals. *Am J Hum Genet*,
 1154 105, 947-958.
- 1155 GIOVANNIELLO, J. A., S.; YU, K.; LI, B. 2021. Sex-Specific Stress-Related Behavioral Phenotypes and Central
 1156 Amygdala Dysfunction in a Mouse Model of 16p11.2 Microdeletion. *Biological Psychiatry Global Open
 1157 Science*, 1, 59 - 69.
- 1158 GOLZIO, C., WILLER, J., TALKOWSKI, M. E., OH, E. C., TANIGUCHI, Y., JACQUEMONT, S.,
 1159 REYMOND, A., SUN, M., SAWA, A., GUSELLA, J. F., KAMIYA, A., BECKMANN, J. S. &
 1160 KATSANIS, N. 2012. KCTD13 is a major driver of mirrored neuroanatomical phenotypes of the 16p11.2
 1161 copy number variant. *Nature*, 485, 363-7.

- 1162 GROSCH, S., SCHIFFMANN, S. & GEISLINGER, G. 2012. Chain length-specific properties of ceramides.
1163 *Prog Lipid Res*, 51, 50-62.
- 1164 GUSTAVSSON, N. & HAN, W. 2009. Calcium-sensing beyond neurotransmitters: functions of synaptotagmins
1165 in neuroendocrine and endocrine secretion. *Biosci Rep*, 29, 245-59.
- 1166 HANSON, E., BERNIER, R., PORCHE, K., JACKSON, F. I., GOIN-KOCHEL, R. P., SNYDER, L. G., SNOW,
1167 A. V., WALLACE, A. S., CAMPE, K. L., ZHANG, Y., CHEN, Q. X., D'ANGELO, D., MORENO-DE-
1168 LUCA, A., ORR, P. T., BOOMER, K. B., EVANS, D. W., KANNE, S., BERRY, L., MILLER, F. K.,
1169 OLSON, J., SHERR, E., MARTIN, C. L., LEDBETTER, D. H., SPIRO, J. E., CHUNG, W. K. & PROJE,
1170 S. V. I. 2015. The Cognitive and Behavioral Phenotype of the 16p11.2 Deletion in a Clinically Ascertained
1171 Population. *Biological Psychiatry*, 77, 785-793.
- 1172 HASLINGER, D., WALTES, R., YOUSAF, A., LINDLAR, S., SCHNEIDER, I., LIM, C. K., TSAI, M. M.,
1173 GARVALOV, B. K., ACKER-PALMER, A., KREZDORN, N., ROTTER, B., ACKER, T.,
1174 GUILLEMIN, G. J., FULDA, S., FREITAG, C. M. & CHIOCCHETTI, A. G. 2018. Loss of the
1175 Chr16p11.2 ASD candidate gene QPRT leads to aberrant neuronal differentiation in the SH-SY5Y
1176 neuronal cell model. *Mol Autism*, 9, 56.
- 1177 HINKLEY, L. B. N., DALE, C. L., LUKS, T. L., FINDLAY, A. M., BUKSHPUN, P., POJMAN, N., THIEU,
1178 T., CHUNG, W. K., BERMAN, J., ROBERTS, T. P. L., MUKHERJEE, P., SHERR, E. H. &
1179 NAGARAJAN, S. S. 2019. Sensorimotor Cortical Oscillations during Movement Preparation in 16p11.2
1180 Deletion Carriers. *J Neurosci*, 39, 7321-7331.
- 1181 HOYTEMA VAN KONIJNENBURG, E. M. M., LUIRINK, I. K., SCHAGEN, S. E. E., ENGELEN, M.,
1182 BERENDSE, K., POLL-THE, B. T. & CHEGARY, M. 2020. Hyperinsulinism in a patient with a
1183 Zellweger Spectrum Disorder and a 16p11.2 deletion syndrome. *Mol Genet Metab Rep*, 23, 100590.
- 1184 IP, J. P. K., NAGAKURA, I., PETRAVICZ, J., LI, K., WIEMER, E. A. C. & SUR, M. 2018. Major Vault Protein,
1185 a Candidate Gene in 16p11.2 Microdeletion Syndrome, Is Required for the Homeostatic Regulation of
1186 Visual Cortical Plasticity. *J Neurosci*, 38, 3890-3900.
- 1187 KALUEFF, A. V., STEWART, A. M. & GERLAI, R. 2014. Zebrafish as an emerging model for studying
1188 complex brain disorders. *Trends Pharmacol Sci*, 35, 63-75.
- 1189 KIKUMA, K., LI, X., KIM, D., SUTTER, D. & DICKMAN, D. K. 2017. Extended Synaptotagmin Localizes to
1190 Presynaptic ER and Promotes Neurotransmission and Synaptic Growth in *Drosophila*. *Genetics*, 207, 993-
1191 1006.
- 1192 KIM, S. H., GREEN-SNYDER, L., LORD, C., BISHOP, S., STEINMAN, K. J., BERNIER, R., HANSON, E.,
1193 GOIN-KOCHEL, R. P. & CHUNG, W. K. 2020. Language characterization in 16p11.2 deletion and
1194 duplication syndromes. *Am J Med Genet B Neuropsychiatr Genet*, 183, 380-391.
- 1195 KLEINENDORST, L., VAN DEN HEUVEL, L. M., HENNEMAN, L. & VAN HAELST, M. M. 2020. Who
1196 ever heard of 16p11.2 deletion syndrome? Parents' perspectives on a susceptibility copy number variation
1197 syndrome. *Eur J Hum Genet*, 28, 1196-1204.
- 1198 KOVALEVICH J, L. D. 2013. Considerations for the Use of SH-SY5Y Neuroblastoma Cells in Neurobiology.
1199 *Methods Mol Biol*, 9-21.
- 1200 KRAFT, M. L. 2016. Sphingolipid Organization in the Plasma Membrane and the Mechanisms That Influence It.
1201 *Front Cell Dev Biol*, 4, 154.
- 1202 LAI, A. L., TAMM, L. K., ELLENA, J. F. & CAFISO, D. S. 2011. Synaptotagmin 1 modulates lipid acyl chain
1203 order in lipid bilayers by demixing phosphatidylserine. *J Biol Chem*, 286, 25291-300.
- 1204 LEVY, M. & FUTERMAN, A. H. 2010. Mammalian ceramide synthases. *IUBMB Life*, 62, 347-56.
- 1205 LI, Y. C., CHANADAY, N. L., XU, W. & KAVALALI, E. T. 2017. Synaptotagmin-1- and Synaptotagmin-7-
1206 Dependent Fusion Mechanisms Target Synaptic Vesicles to Kinetically Distinct Endocytic Pathways.
1207 *Neuron*, 93, 616-631 e3.
- 1208 LIN, G. N., COROMINAS, R., LEMMENS, I., YANG, X., TAVERNIER, J., HILL, D. E., VIDAL, M., SEBAT,
1209 J. & IAKOUCHEVA, L. M. 2015. Spatiotemporal 16p11.2 protein network implicates cortical late mid-
1210 fetal brain development and KCTD13-Cul3-RhoA pathway in psychiatric diseases. *Neuron*, 85, 742-54.

- 1211 LOHNER, K. 1996. Is the high propensity of ethanolamine plasmalogens to form non-lamellar lipid structures
 1212 manifested in the properties of biomembranes? *Chem Phys Lipids*, 81, 167-84.
- 1213 MAILLARD, A. M., RUEF, A., PIZZAGALLI, F., MIGLIAVACCA, E., HIPPOLYTE, L., ADASZEWSKI, S.,
 1214 DUKART, J., FERRARI, C., CONUS, P., MANNIK, K., ZAZHYTSKA, M., SIFFREDI, V., MAEDER,
 1215 P., KUTALIK, Z., KHERIF, F., HADJIKHANI, N., BECKMANN, J. S., REYMOND, A., DRAGANSKI,
 1216 B., JACQUEMONT, S. & P11.2 EUROPEAN, C. 2015. The 16p11.2 locus modulates brain structures
 1217 common to autism, schizophrenia and obesity. *Mol Psychiatry*, 20, 140-7.
- 1218 MARCONESCU, A. & THORPE, P. E. 2008. Coincident exposure of phosphatidylethanolamine and anionic
 1219 phospholipids on the surface of irradiated cells. *Biochim Biophys Acta*, 1778, 2217-24.
- 1220 MARTIN LORENZO, S., NALESSO, V., CHEVALIER, C., BIRLING, M. C. & HERAULT, Y. 2021. Targeting
 1221 the RHOA pathway improves learning and memory in adult Kctd13 and 16p11.2 deletion mouse models.
 1222 *Mol Autism*, 12, 1.
- 1223 MARTIN-BREVET, S., RODRIGUEZ-HERREROS, B., NIELSEN, J. A., MOREAU, C., MODENATO, C.,
 1224 MAILLARD, A. M., PAIN, A., RICHTIN, S., JONCH, A. E., QURESHI, A. Y., ZURCHER, N. R.,
 1225 CONUS, P., P11.2 EUROPEAN, C., SIMONS VARIATION IN INDIVIDUALS PROJECT, C.,
 1226 CHUNG, W. K., SHERR, E. H., SPIRO, J. E., KHERIF, F., BECKMANN, J. S., HADJIKHANI, N.,
 1227 REYMOND, A., BUCKNER, R. L., DRAGANSKI, B. & JACQUEMONT, S. 2018. Quantifying the
 1228 Effects of 16p11.2 Copy Number Variants on Brain Structure: A Multisite Genetic-First Study. *Biol*
 1229 *Psychiatry*, 84, 253-264.
- 1230 MCCAMMON, J. M., BLAKER-LEE, A., CHEN, X. & SIVE, H. 2017. The 16p11.2 homologs fam57ba and
 1231 doc2a generate certain brain and body phenotypes. *Hum Mol Genet*, 26, 3699-3712.
- 1232 MCCAMMON, J. M. & SIVE, H. 2015. Challenges in understanding psychiatric disorders and developing
 1233 therapeutics: a role for zebrafish. *Dis Model Mech*, 8, 647-56.
- 1234 MEFFORD, H. C., YENDLE, S. C., HSU, C., COOK, J., GERAGHTY, E., MCMAHON, J. M., EEG-
 1235 OLOFSSON, O., SADLEIR, L. G., GILL, D., BEN-ZEEV, B., LERMAN-SAGIE, T., MACKAY, M.,
 1236 FREEMAN, J. L., ANDERMANN, E., PELAKANOS, J. T., ANDREWS, I., WALLACE, G., EICHLER,
 1237 E. E., BERKOVIC, S. F. & SCHEFFER, I. E. 2011. Rare copy number variants are an important cause of
 1238 epileptic encephalopathies. *Ann Neurol*, 70, 974-85.
- 1239 NIARCHOU, M., CHAWNER, S., DOHERTY, J. L., MAILLARD, A. M., JACQUEMONT, S., CHUNG, W.
 1240 K., GREEN-SNYDER, L., BERNIER, R. A., GOIN-KOCHEL, R. P., HANSON, E., LINDEN, D. E. J.,
 1241 LINDEN, S. C., RAYMOND, F. L., SKUSE, D., HALL, J., OWEN, M. J. & VAN DEN BREE, M. B.
 1242 M. 2019. Correction: Psychiatric disorders in children with 16p11.2 deletion and duplication. *Transl*
 1243 *Psychiatry*, 9, 107.
- 1244 OHSUGI, M., ADACHI, K., HORAI, R., KAKUTA, S., SUDO, K., KOTAKI, H., TOKAI-NISHIZUMI, N.,
 1245 SAGARA, H., IWAKURA, Y. & YAMAMOTO, T. 2008. Kid-mediated chromosome compaction
 1246 ensures proper nuclear envelope formation. *Cell*, 132, 771-82.
- 1247 OUELLETTE, J., TOUSSAY, X., COMIN, C. H., COSTA, L. D. F., HO, M., LACALLE-AURIOLES, M.,
 1248 FREITAS-ANDRADE, M., LIU, Q. Y., LECLERC, S., PAN, Y., LIU, Z., THIBODEAU, J. F., YIN, M.,
 1249 CARRIER, M., MORSE, C. J., DYKEN, P. V., BERGIN, C. J., BAILLET, S., KENNEDY, C. R.,
 1250 TREMBLAY, M. E., BENOIT, Y. D., STANFORD, W. L., BURGER, D., STEWART, D. J. &
 1251 LACOSTE, B. 2020. Vascular contributions to 16p11.2 deletion autism syndrome modeled in mice. *Nat*
 1252 *Neurosci*, 23, 1090-1101.
- 1253 OWEN, J. P., BUKSHUPUN, P., POJMAN, N., THIEU, T., CHEN, Q., LEE, J., D'ANGELO, D., GLENN, O. A.,
 1254 HUNTER, J. V., BERMAN, J. I., ROBERTS, T. P., BUCKNER, R., NAGARAJAN, S. S.,
 1255 MUKHERJEE, P. & SHERR, E. H. 2018. Brain MR Imaging Findings and Associated Outcomes in
 1256 Carriers of the Reciprocal Copy Number Variation at 16p11.2. *Radiology*, 286, 217-226.
- 1257 PALTAUF, F. 1994. Ether lipids in biomembranes. *Chem Phys Lipids*, 74, 101-39.
- 1258 PARK, S. M., LITTLETON, J. T., PARK, H. R. & LEE, J. H. 2016. Drosophila Homolog of Human KIF22 at
 1259 the Autism-Linked 16p11.2 Loci Influences Synaptic Connectivity at Larval Neuromuscular Junctions.
 1260 *Exp Neurobiol*, 25, 33-9.

- 1261 PEWZNER-JUNG, Y., BEN-DOR, S. & FUTERMAN, A. H. 2006. When do Lasses (longevity assurance genes)
 1262 become CerS (ceramide synthases)? Insights into the regulation of ceramide synthesis. *J Biol Chem*, 281,
 1263 25001-5.
- 1264 PIKE, L. J. 2003. Lipid rafts: bringing order to chaos. *J Lipid Res*, 44, 655-67.
- 1265 PINTO, D., PAGNAMENTA, A. T., KLEI, L., ANNEY, R., MERICO, D., REGAN, R., CONROY, J.,
 1266 MAGALHAES, T. R., CORREIA, C., ABRAHAMS, B. S., ALMEIDA, J., BACCHELLI, E., BADER,
 1267 G. D., BAILEY, A. J., BAIRD, G., BATTAGLIA, A., BERNEY, T., BOLSHAKOVA, N., BOLTE, S.,
 1268 BOLTON, P. F., BOURGERON, T., BRENNAN, S., BRIAN, J., BRYSON, S. E., CARSON, A. R.,
 1269 CASALLO, G., CASEY, J., CHUNG, B. H., COCHRANE, L., CORSELLO, C., CRAWFORD, E. L.,
 1270 CROSSETT, A., CYTRYNBAUM, C., DAWSON, G., DE JONGE, M., DELORME, R., DRMIC, I.,
 1271 DUKETIS, E., DUQUE, F., ESTES, A., FARRAR, P., FERNANDEZ, B. A., FOLSTEIN, S. E.,
 1272 FOMBONNE, E., FREITAG, C. M., GILBERT, J., GILLBERG, C., GLESSNER, J. T., GOLDBERG, J.,
 1273 GREEN, A., GREEN, J., GUTER, S. J., HAKONARSON, H., HERON, E. A., HILL, M., HOLT, R.,
 1274 HOWE, J. L., HUGHES, G., HUS, V., IGLIOZZI, R., KIM, C., KLAUCK, S. M., KOLEVZON, A.,
 1275 KORVATSKA, O., KUSTANOVICH, V., LAJONCHERE, C. M., LAMB, J. A., LASKAWIEC, M.,
 1276 LEBOYER, M., LE COUTEUR, A., LEVENTHAL, B. L., LIONEL, A. C., LIU, X. Q., LORD, C.,
 1277 LOTSPEICH, L., LUND, S. C., MAESTRINI, E., MAHONEY, W., MANTOULAN, C., MARSHALL,
 1278 C. R., MCCONACHIE, H., MCDUGLE, C. J., MCGRATH, J., MCMAHON, W. M., MERIKANGAS,
 1279 A., MIGITA, O., MINSHEW, N. J., MIRZA, G. K., MUNSON, J., NELSON, S. F., NOAKES, C., NOOR,
 1280 A., NYGREN, G., OLIVEIRA, G., PAPANIKOLAOU, K., PARR, J. R., PARRINI, B., PATON, T.,
 1281 PICKLES, A., PILORGE, M., et al. 2010. Functional impact of global rare copy number variation in
 1282 autism spectrum disorders. *Nature*, 466, 368-72.
- 1283 QURESHI, A. Y., MUELLER, S., SNYDER, A. Z., MUKHERJEE, P., BERMAN, J. I., ROBERTS, T. P.,
 1284 NAGARAJAN, S. S., SPIRO, J. E., CHUNG, W. K., SHERR, E. H., BUCKNER, R. L. & SIMONS, V.
 1285 I. P. C. 2014. Opposing brain differences in 16p11.2 deletion and duplication carriers. *J Neurosci*, 34,
 1286 11199-211.
- 1287 RANTA, S., ZHANG, Y., ROSS, B., LONKA, L., TAKKUNEN, E., MESSER, A., SHARP, J., WHEELER, R.,
 1288 KUSUMI, K., MOLE, S., LIU, W., SOARES, M. B., BONALDO, M. F., HIRVASNIEMI, A., DE LA
 1289 CHAPELLE, A., GILLIAM, T. C. & LEHESJOKI, A. E. 1999. The neuronal ceroid lipofuscinoses in
 1290 human EPMP and mnd mutant mice are associated with mutations in CLN8. *Nat Genet*, 23, 233-6.
- 1291 SATTERSTROM, F. K., WALTERS, R. K., SINGH, T., WIGDOR, E. M., LESCAI, F., DEMONTIS, D.,
 1292 KOSMICKI, J. A., GROVE, J., STEVENS, C., BYBJERG-GRAUHOLM, J., BAEKVAD-HANSEN, M.,
 1293 PALMER, D. S., MALLER, J. B., I, P.-B. C., NORDENTOFT, M., MORS, O., ROBINSON, E. B.,
 1294 HOUGAARD, D. M., WERGE, T. M., BO MORTENSEN, P., NEALE, B. M., BORGLUM, A. D. &
 1295 DALY, M. J. 2019. Autism spectrum disorder and attention deficit hyperactivity disorder have a similar
 1296 burden of rare protein-truncating variants. *Nat Neurosci*, 22, 1961-1965.
- 1297 SCHONN, J. S., MAXIMOV, A., LAO, Y., SUDHOF, T. C. & SORENSEN, J. B. 2008. Synaptotagmin-1 and -
 1298 7 are functionally overlapping Ca²⁺ sensors for exocytosis in adrenal chromaffin cells. *Proc Natl Acad
 1299 Sci U S A*, 105, 3998-4003.
- 1300 SEBAT, J., LAKSHMI, B., MALHOTRA, D., TROGE, J., LESE-MARTIN, C., WALSH, T., YAMROM, B.,
 1301 YOON, S., KRASNITZ, A., KENDALL, J., LEOTTA, A., PAI, D., ZHANG, R., LEE, Y. H., HICKS, J.,
 1302 SPENCE, S. J., LEE, A. T., PUURA, K., LEHTIMAKI, T., LEDBETTER, D., GREGERSEN, P. K.,
 1303 BREGMAN, J., SUTCLIFFE, J. S., JOBANPUTRA, V., CHUNG, W., WARBURTON, D., KING, M.
 1304 C., SKUSE, D., GESCHWIND, D. H., GILLIAM, T. C., YE, K. & WIGLER, M. 2007. Strong association
 1305 of de novo copy number mutations with autism. *Science*, 316, 445-9.
- 1306 SIMONS VIP, C. 2012. Simons Variation in Individuals Project (Simons VIP): a genetics-first approach to
 1307 studying autism spectrum and related neurodevelopmental disorders. *Neuron*, 73, 1063-7.
- 1308 SORENSEN, J. B., FERNANDEZ-CHACON, R., SUDHOF, T. C. & NEHER, E. 2003. Examining
 1309 synaptotagmin 1 function in dense core vesicle exocytosis under direct control of Ca²⁺. *J Gen Physiol*,
 1310 122, 265-76.

- 1311 STEWART, A. M., NGUYEN, M., WONG, K., POUDEL, M. K. & KALUEFF, A. V. 2014. Developing
1312 zebrafish models of autism spectrum disorder (ASD). *Prog Neuropsychopharmacol Biol Psychiatry*, 50,
1313 27-36.
- 1314 TIDHAR, R., BEN-DOR, S., WANG, E., KELLY, S., MERRILL, A. H., JR. & FUTERMAN, A. H. 2012. Acyl
1315 chain specificity of ceramide synthases is determined within a region of 150 residues in the Tram-Lag-
1316 CLN8 (TLC) domain. *J Biol Chem*, 287, 3197-206.
- 1317 TIERNEY, E., BUKELIS, I., THOMPSON, R. E., AHMED, K., ANEJA, A., KRATZ, L. & KELLEY, R. I. 2006.
1318 Abnormalities of cholesterol metabolism in autism spectrum disorders. *Am J Med Genet B Neuropsychiatr*
1319 *Genet*, 141B, 666-8.
- 1320 TOMASELLO, D. L. & SIVE, H. 2020. Noninvasive Multielectrode Array for Brain and Spinal Cord Local Field
1321 Potential Recordings from Live Zebrafish Larvae. *Zebrafish*, 17, 271-277.
- 1322 TRACEY, T. J., STEYN, F. J., WOLVETANG, E. J. & NGO, S. T. 2018. Neuronal Lipid Metabolism: Multiple
1323 Pathways Driving Functional Outcomes in Health and Disease. *Front Mol Neurosci*, 11, 10.
- 1324 VANNI, N., FRUSCIONE, F., FERLAZZO, E., STRIANO, P., ROBBIANO, A., TRAVERSO, M., SANDER,
1325 T., FALACE, A., GAZZERRO, E., BRAMANTI, P., BIELAWSKI, J., FASSIO, A., MINETTI, C.,
1326 GENTON, P. & ZARA, F. 2014. Impairment of ceramide synthesis causes a novel progressive myoclonus
1327 epilepsy. *Ann Neurol*, 76, 206-12.
- 1328 WORSTELL, N. C., KRISHNAN, P., WEATHERSTON, J. D. & WU, H. J. 2016. Binding Cooperativity Matters:
1329 A GM1-Like Ganglioside-Cholera Toxin B Subunit Binding Study Using a Nanocube-Based Lipid
1330 Bilayer Array. *PLoS One*, 11, e0153265.
- 1331 XI, Y., NOBLE, S. & EKKER, M. 2011. Modeling neurodegeneration in zebrafish. *Curr Neurol Neurosci Rep*,
1332 11, 274-82.
- 1333 XU, J., MASHIMO, T. & SUDHOF, T. C. 2007. Synaptotagmin-1, -2, and -9: Ca(2+) sensors for fast release that
1334 specify distinct presynaptic properties in subsets of neurons. *Neuron*, 54, 567-81.
- 1335 YAMASHITA-SUGAHARA, Y., TOKUZAWA, Y., NAKACHI, Y., KANESAKI-YATSUKA, Y.,
1336 MATSUMOTO, M., MIZUNO, Y. & OKAZAKI, Y. 2013. Fam57b (family with sequence similarity 57,
1337 member B), a novel peroxisome proliferator-activated receptor gamma target gene that regulates
1338 adipogenesis through ceramide synthesis. *J Biol Chem*, 288, 4522-37.
- 1339 YU, H., LIU, Y., GULBRANSON, D. R., PAINE, A., RATHORE, S. S. & SHEN, J. 2016. Extended
1340 synaptotagmins are Ca²⁺-dependent lipid transfer proteins at membrane contact sites. *Proc Natl Acad Sci*
1341 *U S A*, 113, 4362-7.
- 1342 YUAN, A. & NIXON, R. A. 2016. Specialized roles of neurofilament proteins in synapses: Relevance to
1343 neuropsychiatric disorders. *Brain Res Bull*, 126, 334-346.
- 1344 ZHANG, T., TRAUGER, S. A., VIDOUDEZ, C., DOANE, K. P., PLUIMER, B. R. & PETERSON, R. T. 2019.
1345 Parallel Reaction Monitoring reveals structure-specific ceramide alterations in the zebrafish. *Sci Rep*, 9,
1346 19939.
- 1347 ZUFFEREY, F., SHERR, E. H., BECKMANN, N. D., HANSON, E., MAILLARD, A. M., HIPPOLYTE, L.,
1348 MACE, A., FERRARI, C., KUTALIK, Z., ANDRIEUX, J., AYLWARD, E., BARKER, M., BERNIER,
1349 R., BOUQUILLON, S., CONUS, P., DELOBEL, B., FAUCETT, W., GOIN-KOCHEL, R. P., GRANT,
1350 E., HAREWOOD, L., HUNTER, J. V., LEBON, S., LEDBETTER, D. H., MARTIN, C. L., MANNIK,
1351 K., MARTINET, D., MUKHERJEE, P., RAMOCKI, M. B., SPENCE, S. J., STEINMAN, K. J.,
1352 TJERNAGEL, J., SPIRO, J. E., REYMOND, A., BECKMANN, J. S., CHUNG, W. K., JACQUEMONT,
1353 S., CONSORTIUM, S. V. & CONSORTIUM, P.-E. 2012. A 600 kb deletion syndrome at 16p11.2 leads
1354 to energy imbalance and neuropsychiatric disorders. *Journal of Medical Genetics*, 49, 660-668.
1355
1356

REAGENT or RESOURCE	SOURCE	IDENTIFIER
Antibodies		
Synaptotagmin-1	Lifespan Bioscience	LS-B12889
GAD65 + 67	Abcam	ab11070
Beta-Actin	Proteintech	60008-1
DAPI	Life Technologies	D1306
555-Phalloidin	Invitrogen	A34055
488 anti-goat	Jackson	805-545-180,
488 anti-mouse	Jackson	715-545-151
594 anti-mouse	Jackson	715-585-150
594 anti-rabbit	Jackson	711-585-152
anti-mouse 680	Jackson	715-625-150
anti-rabbit 680	Jackson	711-625-152
VGlut1/2	Synaptic Systems	135503
PSD95	Abcam	ab18258
Ac-Tubulin	Abcam	ab179513
Prolong Gold Antifade with DAPI	LifeTech	P36935
HA	Abcam	ab18181
FLAG	Abcam	ab1162
GAPDH	Abcam	ab8245
Li-Cor 800CW Rabbit	Li-Cor	92632211
Li-Cor 680RD Mouse	Li-Cor	92668070
Chemicals, peptides, and recombinant proteins		
Matrigel	Corning	CB-40234A
mTeSR+	STEMCELL Tech	85850
Y27632	STEMCELL Tech	72302
ReLeSR	STEMCELL Tech	05872
Accutase	STEMCELL Tech	07922
DMEM/F12/HEPES	Thermo	12400024
Neurobasal	Thermo	21103049
N2	Gibco	17502048
Gem21	GeminiBio	400-160
MEM NEAA	Thermo	11140050
GlutaMAX	Gibco	35050061
Gem21 - Vit A	GeminiBio	400161
Beta-Mercaptoethanol	Sigma	M3148
FGF	Peprtech	100-18B
BDNF	Peprtech	450-02
GDNF	Peprtech	450-10
Dorsomorphin	Tocis	3093
Poly-D-Lysine	Thermo	A3890401
Laminin	Sigma	L2020
EMEM	ATCC	30-2003
F12	ATCC	30-2006
FBS	Sigma	12306C
All-trans-RA	Sigma	R2625
DMEM	Corning	MT15017CV
Polyethylenimine	Sigma	08719

Protease Inhibitor Cocktail	Sigma	200-664-3
NBD-spanganine	Avanti Polar	810206P
BSA -FA	Sigma	10775835001
16.1 Coenzyme A	Avanti Polar	870743
24.1 Coenzyme 1	Avanti Polar	870725
CT-B	Invitrogen	C34775
Duramycin-Cy3	Molecular Targeting	D-1006
SHIELD	LifeCanvas Tech	https://lifecanvastech.com/products/shield
PFA	EMS	50-970-495
RIPA	Thermo	89900
SYN-Per	Thermo	87793
Critical commercial assays		
Pierce Cell Surface Protein Isolation Kit	Thermo	89881
Deposited data		
Lipidomics	This Paper	Supplemental Excel
Biotinylation Proteomics	This Paper	Supplemental Excel
Synaptosome Proteomics	This Paper	Supplemental Excel
Experimental models: Cell lines		
SH-SY5Y	ATCC	CRL-2266
HEK293T	ATCC	CRL-3216
IPSC 16pdel 1453	Simons VIP	SV0001453
IPSC 16pdel 1455	Simons VIP	SV0001455
IPSC 16pdel 1459	Simons VIP	SV0001459
IPSC 16pdel 1473	Simons VIP	SV0001473
IPSC 16pdel 1481	Simons VIP	SV0001481
IPSC 16pdel 1495	Simons VIP	SV0001495
IPSC 16pdel 3104	Simons VIP	SS0013104
IPSC CTR 599	Corriell Institute Biobank	AG07599
IPSC CTR 675	Corriell Institute Biobank	AG07657
Experimental models: Organisms/strains		
<i>fam57ba^{-/-};fam57bb^{-/-}</i> Zebrafish	This Paper	
FAM57B HET SH-SY5Y	This Paper	
FAM57B KO SH-SY5Y	This Paper	
FAM57B AAVS1 SH-SY5Y	This Paper	
Oligonucleotides		
fam57bb 5' to 3' TAGGTGATGTCCTGGCAGGAAG	This Paper	
fam57bb 3' to 5' AACCTTCCTGCCAGGACATCA	This Paper	
sgFAM57B1 5' to 3' - GGTGCTCCACCATGCCGCCA	This Paper	
sgFAM57B2 5' to 3' - GGGCACAGCAAATTGCGTGT	This Paper	
sgAAVS1 5' to 3' - CACCGGGGCCACTAGGGACAGGAT	This Paper	
Recombinant DNA		
FAM57B-Flag	This Paper	
CerS2-HA	This Paper	

CerS5-HA	This Paper	
CerS6-HA	This Paper	
Software and algorithms		
AxIS and Neural Metric Tool	Axion Biosystems	Axionbiosystems.com
FIJI ImageJ	ImageJ	imagej.net
GO SLIM analysis	PANTHER	pantherdb.org
LipidSearch	Thermo Scientific	IQLAAEGABSFAPC MBFK
Mascot	Matrix Science	Matrixscience.com
Scaffold Q+S	Proteome Software	Proteomesoftware.com
PEAKS Studio 8.5	Bioinformatics Solutions Inc.	Bioinfor.com
EthoVision XT	Noldus	Noldus.com

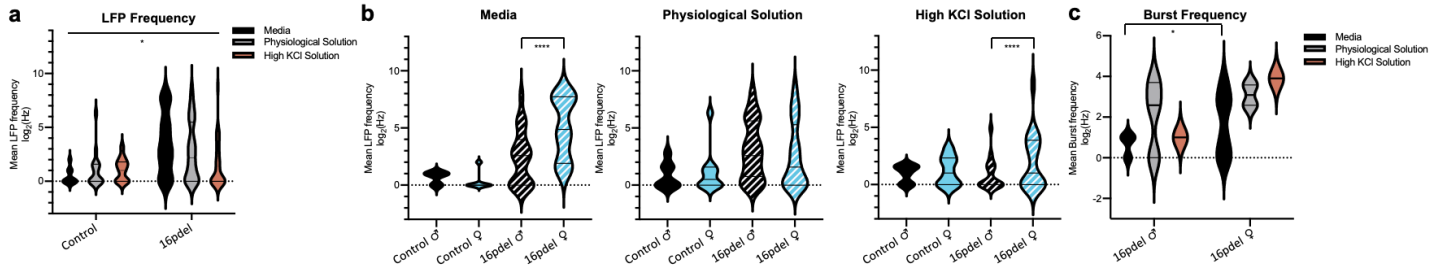


Figure 1. Augmented local field potential activity in 16pdel syndrome differentiated neuronal culture.

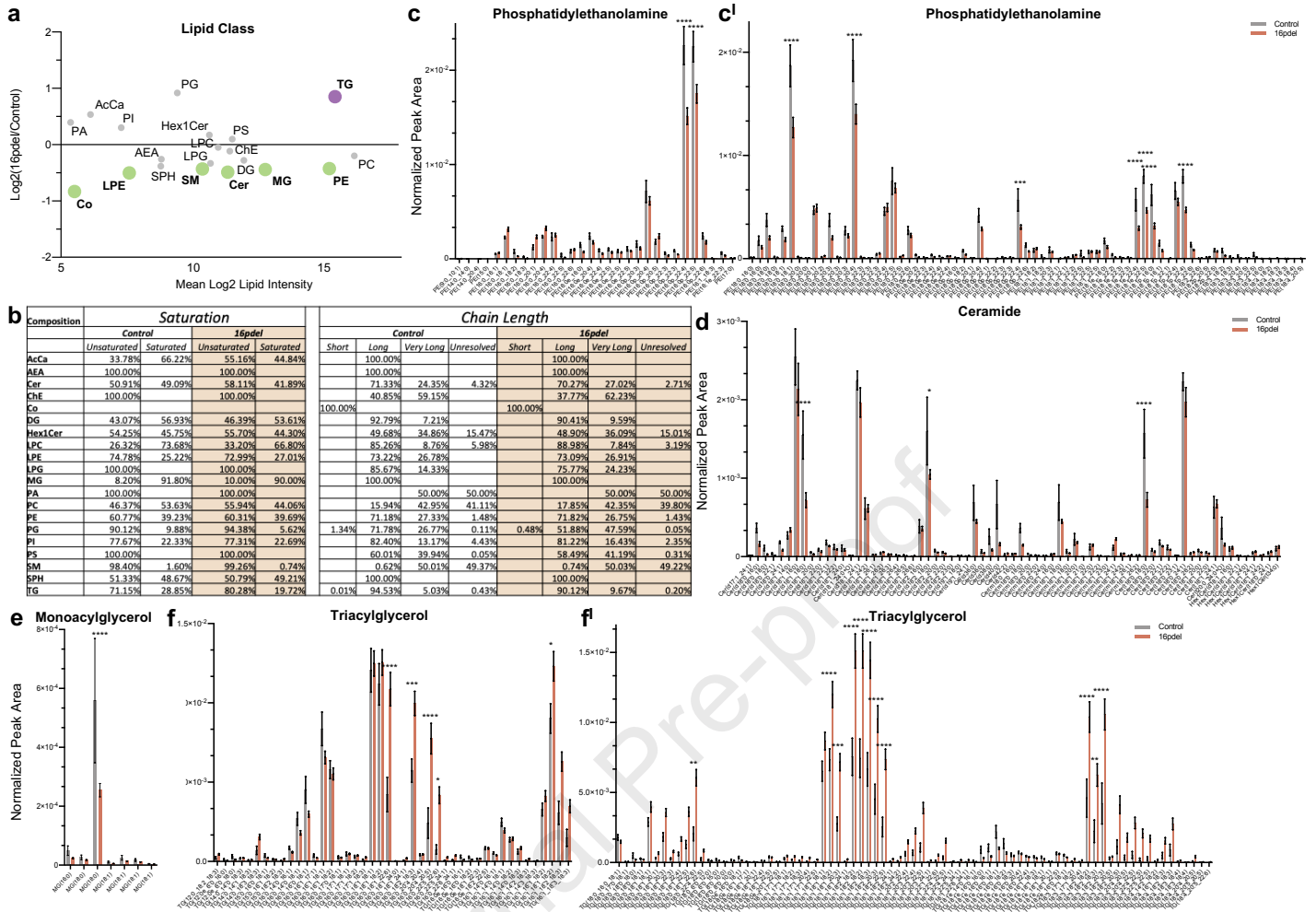


Figure 2. Significant lipid changes between control and 16pdel differentiated neurons.

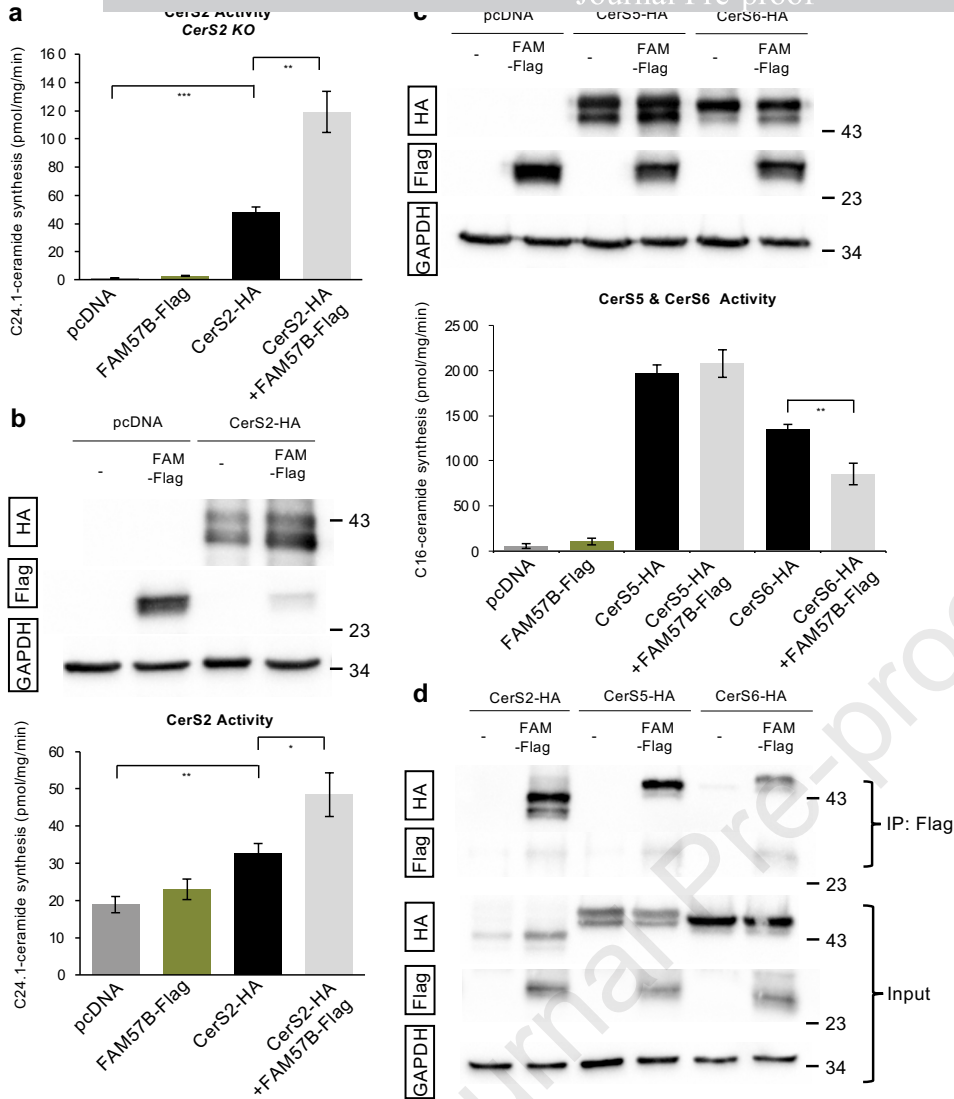


Figure 3. FAM57B interacts with CerS but does not have CerS activity.

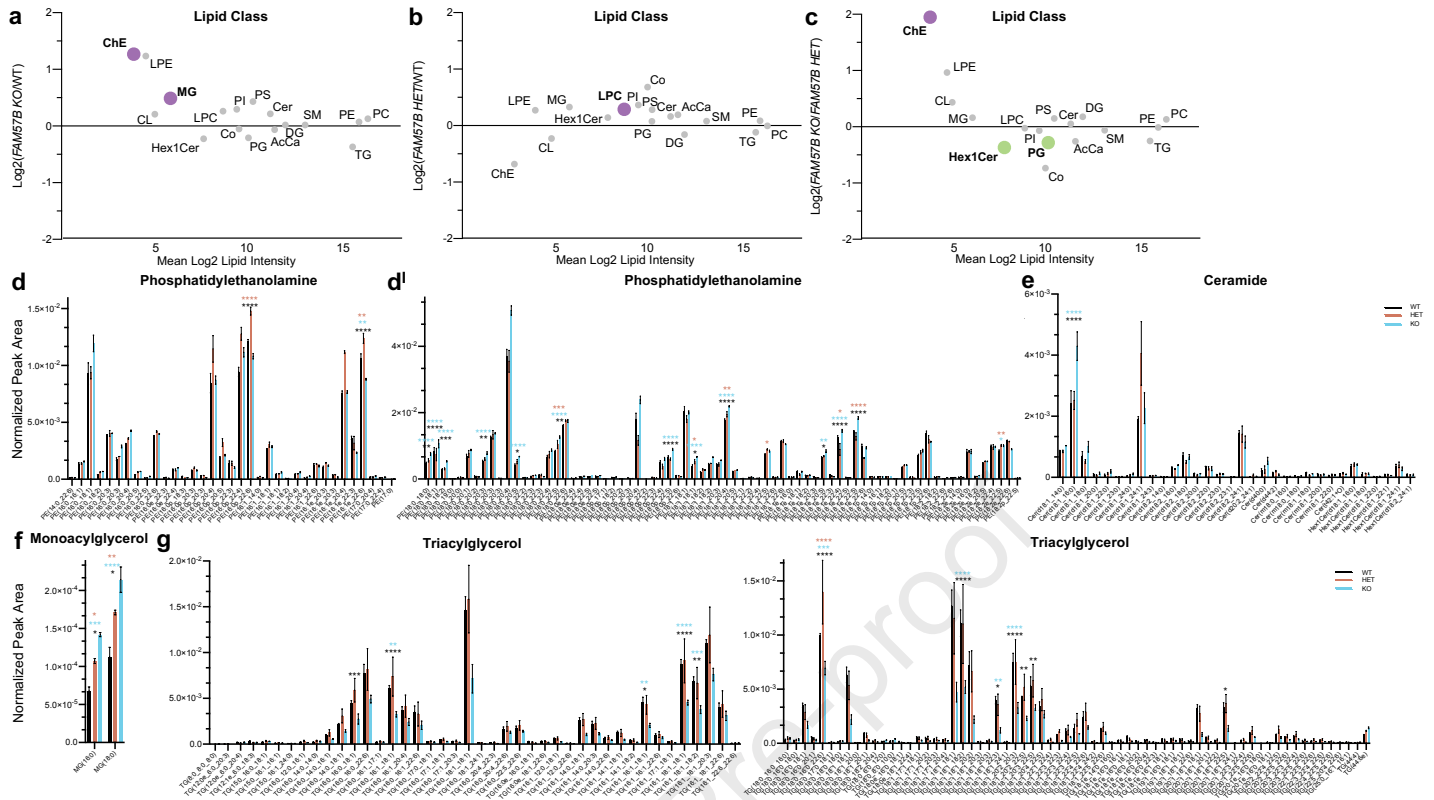


Figure 4. Significant lipid changes in sphingolipids and glycerolipids between WT and *FAM57B* mutant human differentiated SH-SY5Y neuronal cells.

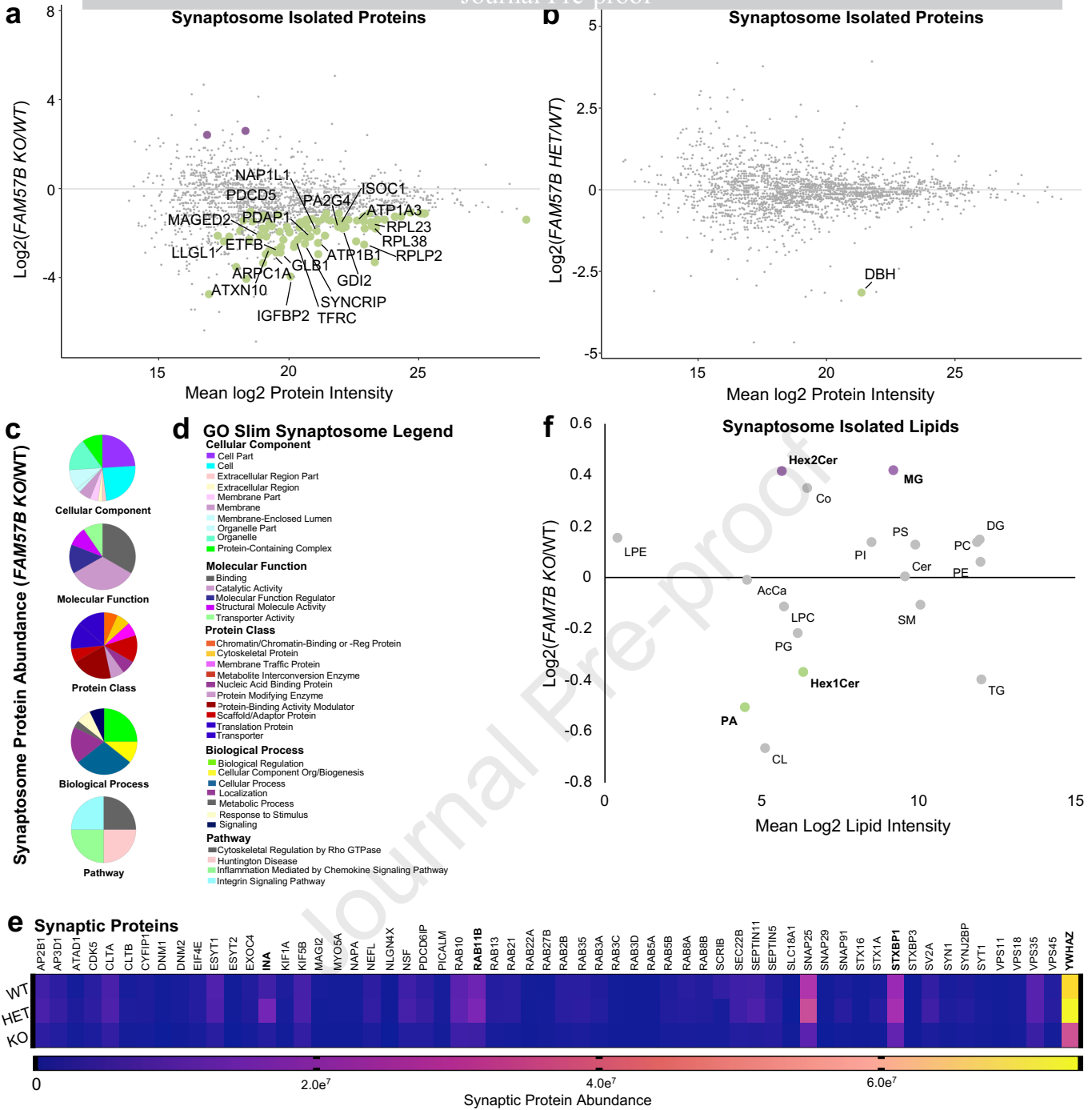


Figure 5. *FAM57B* knockout human neurons indicate altered synaptic composition.

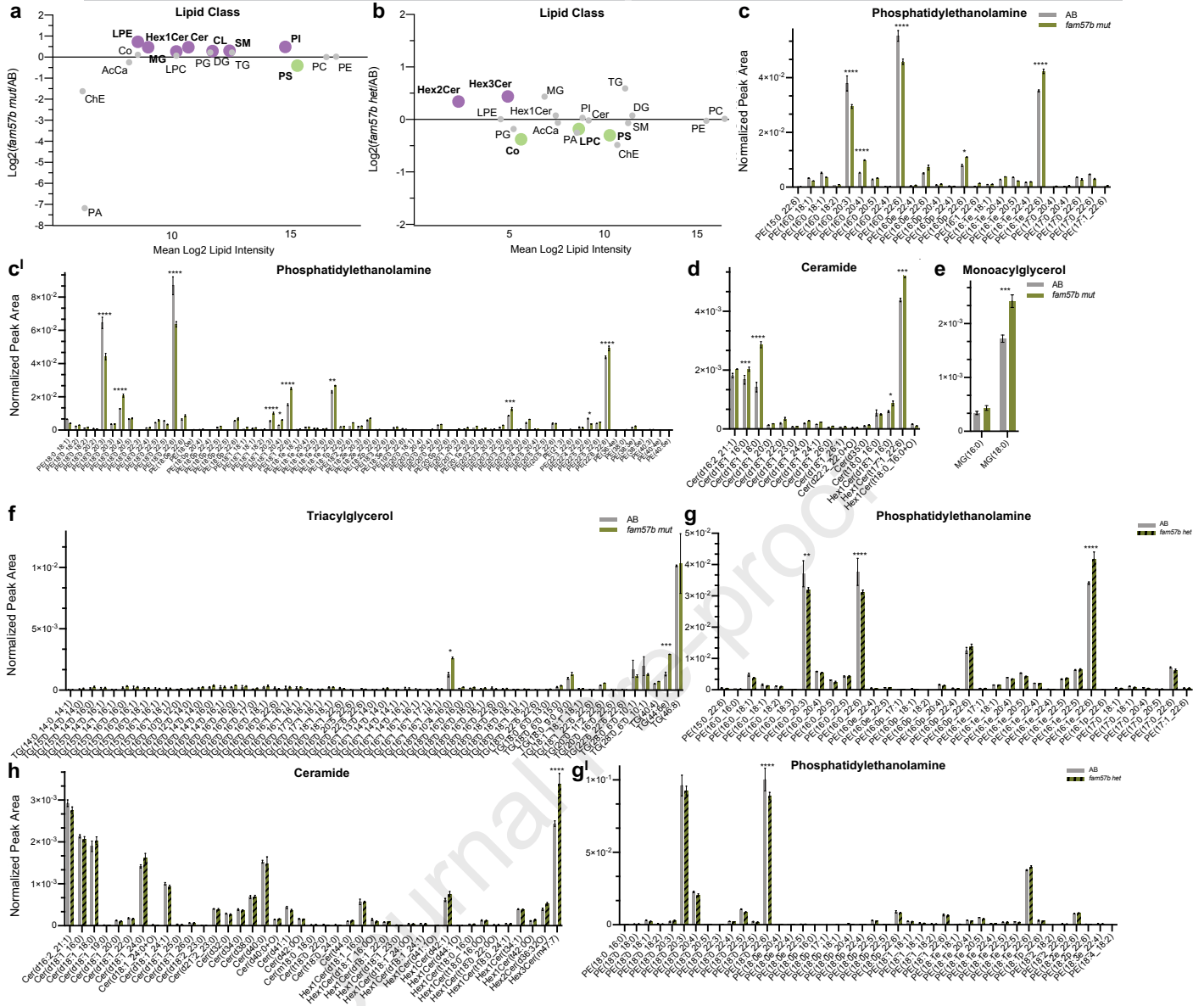


Figure 6. Significant lipid changes in ceramides and glycerols between AB and *fam57b mut* brain tissue.

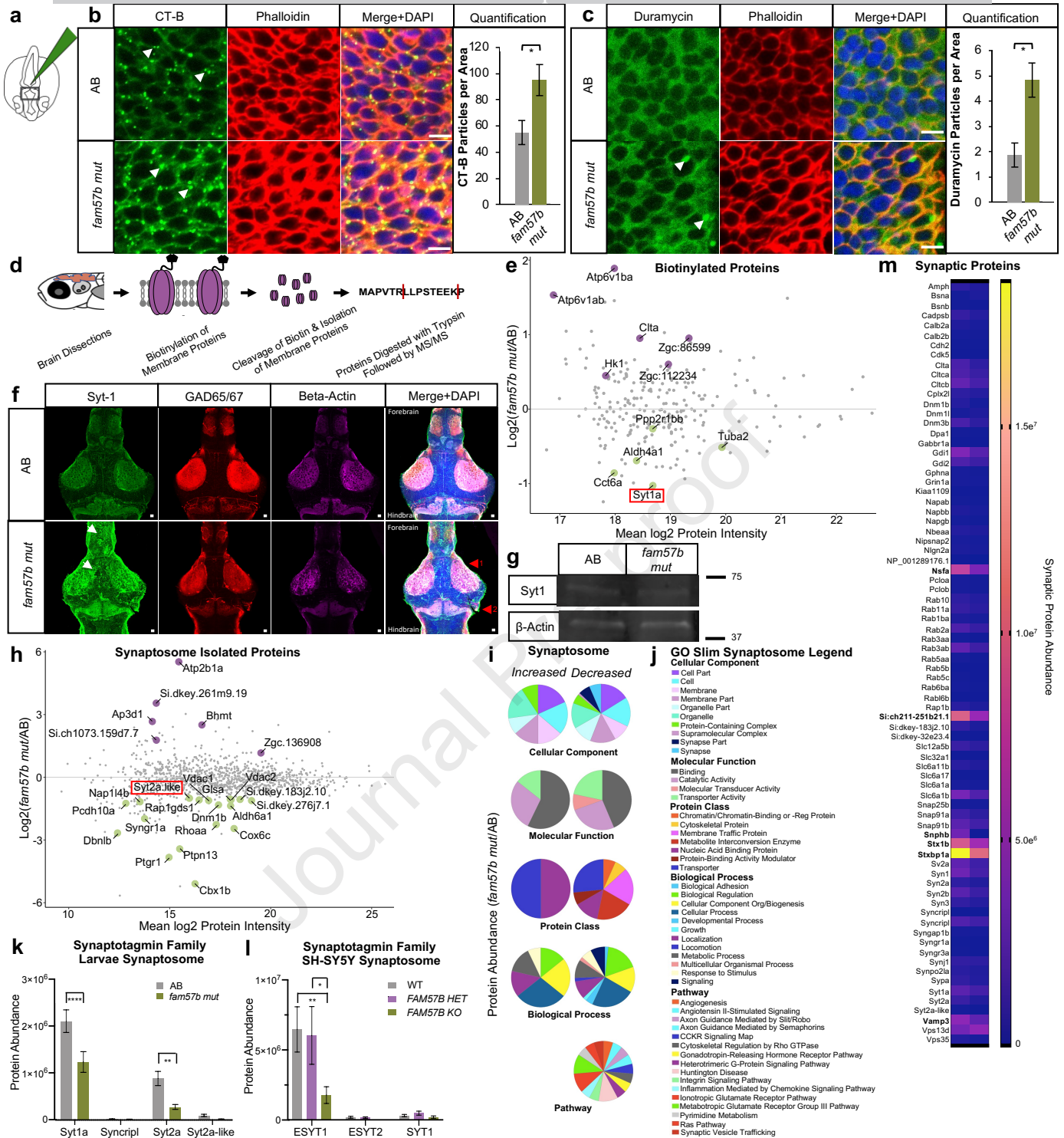


Figure 7. Loss of *fam57b* results in altered plasma membrane architecture early in development and decreased localization of Synaptotagmin family at the synapse later in development.

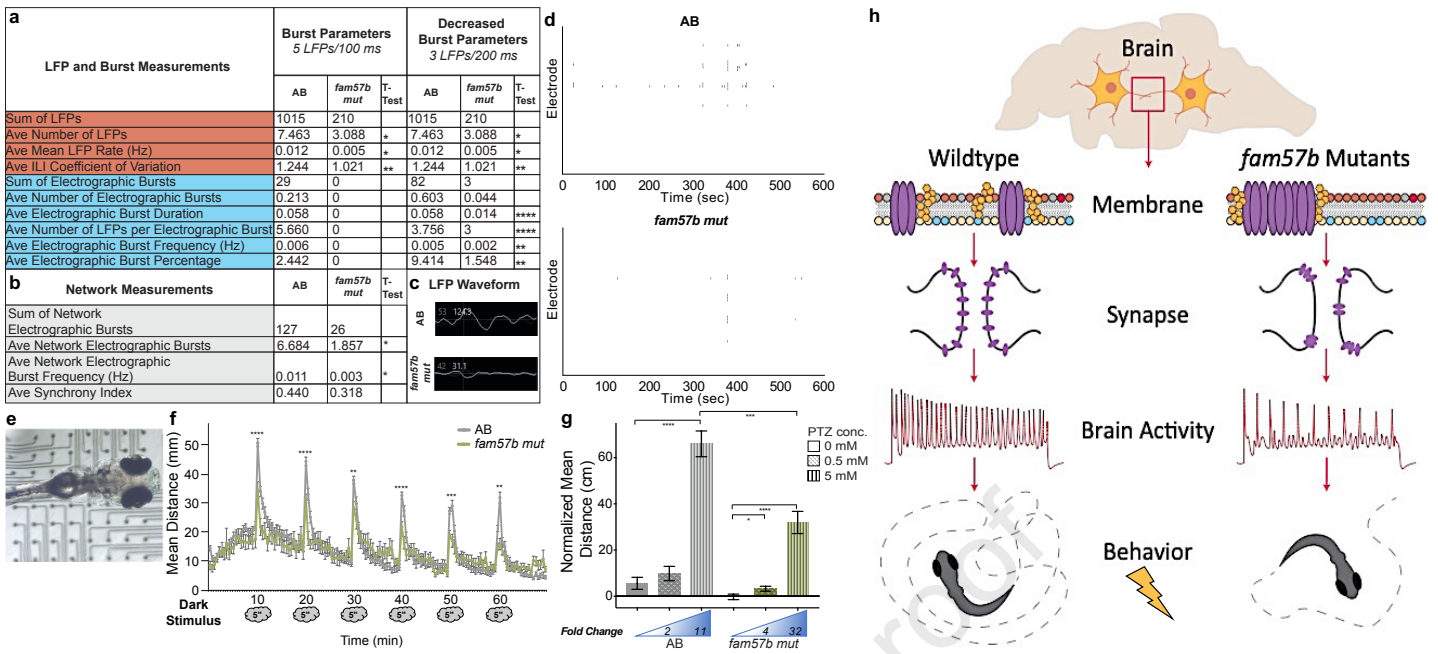
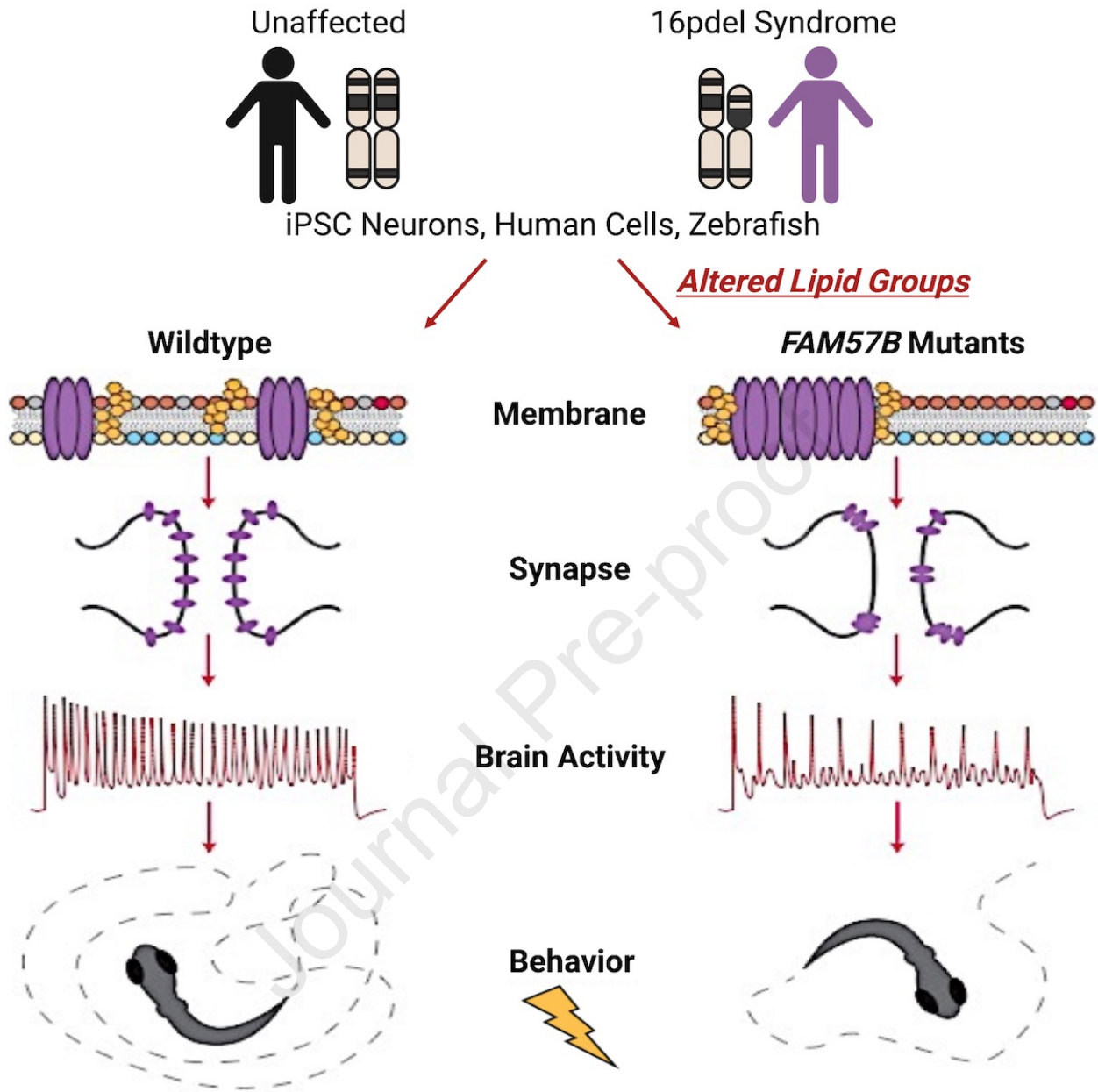


Figure 8. Decreased spontaneous brain activity and diminished behavioral response after stimuli presentation in *fam57b* mutants.



Highlights

- Augmented LFP activity and sex-specific differences in 16pdel neurons.
- 16pdel neuronal lipidome indicated altered ceramide related species.
- FAM57B is a ceramide synthase modulator essential for lipid regulation in the brain.
- FAM57B functions in synaptogenesis, synapse architecture and composition.

Journal Pre-proof