

NEURONAL HYPERACTIVITY CAUSES Na^+/H^+ EXCHANGER-INDUCED EXTRACELLULAR ACIDIFICATION AT ACTIVE SYNAPSES

Martina Chiacchiaretta^{1,2*}, Shahrzad Latifi^{1##}, Mattia Bramini¹, Manuela Fadda², Anna Fassio^{1,2}, Fabio Benfenati^{1,2*} and Fabrizia Cesca^{1*}

*equal contribution

¹Center for Synaptic Neuroscience and Technology, Fondazione Istituto Italiano di Tecnologia, L.go Rosanna Benzi 10, 16132 Genova, Italy; ²Department of Experimental Medicine, University of Genova, Viale Benedetto XV 3, 16132, Genova, Italy

#Present address: Department of Neurology, David Geffen School of Medicine at UCLA, 710 Westwood Plaza, Los Angeles, CA 90095

Key Words: acidic pH, neural hyperexcitability, active synapses, experimental epilepsy, NHE, pH-sensitive GFP.

Corresponding author:

Fabrizia Cesca

Center for Synaptic Neuroscience

Fondazione Istituto Italiano di Tecnologia

L.go Rosanna Benzi, 10

16132 Genova - Italy

phone: +39-010-555-8382

fax: +39-010-555-8491

email: fabrizia.cesca@iit.it

SUMMARY STATEMENT

Using a novel membrane-targeted extracellular pH-indicator exquisitely sensitive to acidic shifts, we found that neuronal hyperactivity causes acidic shifts localized to active synapses and generated by Na⁺/H⁺ exchanger activity.

ABSTRACT

Extracellular pH impacts on neuronal activity, which is in turn an important determinant of extracellular H^+ concentration. The aim of this study is to describe the spatio-temporal dynamics of extracellular pH at synaptic sites during neuronal hyperexcitability. To address this issue we created ex.E²GFP, a membrane-targeted extracellular ratiometric pH indicator exquisitely sensitive to acidic shifts. By monitoring ex.E²GFP fluorescence in real time in primary cortical neurons we were able to quantify pH fluctuations during network hyperexcitability induced by convulsant drugs or high frequency electrical stimulation. Sustained hyperactivity caused a pH decrease that was reversible upon silencing of neuronal activity and localized to active synapses. This acidic shift was not attributable to the outflow of synaptic vesicle protons into the cleft nor to the activity of membrane-exposed H^+ -vATPase, but rather to the activity of the Na^+/H^+ -exchanger. Our data demonstrate that extracellular synaptic pH shifts take place during epileptic-like activity of neural cultures, underlying the strict links existing between synaptic activity and synaptic pH. This evidence may contribute to the understanding of the physio-pathological mechanisms associated with hyperexcitability in the epileptic brain.

INTRODUCTION

In biological systems, the H^+ concentration gradient between the intracellular and the extracellular environments plays a crucial role in several physiological processes (Chesler, 2003). In the central nervous system (CNS), the control of pH homeostasis is essential for the regulation of neurotransmission and neuronal excitability. While pH affects neural activity, neural activity causes pH alterations that vary in location and time (Chen and Chesler, 1992; Dulla et al., 2005; Makani and Chesler, 2007). Local pH changes occur in the brain during normal function (Chesler and Kaila, 1992; Magnotta et al., 2012), and neuronal firing causes both intracellular and extracellular pH shifts (Chesler, 2003; Chesler and Kaila, 1992; Kim and Trussell, 2009; Trapp et al., 1996). Rapid changes in intracellular pH have been observed at *Drosophila* nerve terminals during high-frequency *in vivo* activity (Caldwell et al., 2013; Rossano et al., 2013), while in the mammalian CNS intense and synchronous activity of neural networks is accompanied by an initial alkalization followed by a slower wave of acidification of the extracellular environment (DeVries, 2001; Du et al., 2014). A variety of ion channels, such as voltage-gated Ca^{2+} channels (Iijima et al., 1986), AMPA- and NMDA-type glutamate receptors (Lei et al., 2001; Tang et al., 1990; Traynelis and Cull-Candy, 1990) and $GABA_A$ receptors (Dietrich and Morad, 2010; Kaila, 1994) display sharp pH-sensitive conductances, indicating that pH *per se* can strongly influence neuronal excitability, neurotransmitter release and postsynaptic responses (Chen and Chesler, 1992; Dietrich and Morad, 2010; Makani and Chesler, 2007).

pH shifts are also associated with a number of pathological processes. Interstitial acidosis is present in several neuropathologies such as epilepsy (Gottfried and Chesler, 1996; Inamura et al., 1989; Krishtal et al., 1987; Somjen, 1984) and brain ischemia (Makani and Chesler, 2010; McCormick and Contreras, 2001; Nedergaard et al., 1991; Siesjo et al., 1996; Xiong and Stringer, 2000). The correlation between pH shifts and neuronal hyperexcitability is well established. However, because of the lack of accurate systems to measure the extracellular H^+ concentration, addressing the detailed events underlying pH-induced modulation of neural excitability is still a major challenge.

pH-sensitive microelectrodes generated some of the earliest measurements of both extracellular and intracellular pH, although their size and effects on tissue integrity strongly limited their widespread application. The use of fluorescent pH indicators opened new avenues to the study of pH fluctuations. The use of the extracellular ratiometric pH sensor SNARF allowed to measure the local acidification in PC12 cells stimulated by exposure to

high K^+ (Shuba et al., 2008) and the stimulus-evoked non-quantal release of protons from vestibular hair cells at the cell-calyx afferent synapse (Highstein et al., 2014). More recently, genetically encoded fluorescent pH reporters targeted to specific cellular compartments have been employed. One of such indicators is a variant of the green fluorescent protein (GFP) obtained via specific mutagenesis, called "ratiometric GFP", which is characterized by multiple excitation and emission maxima at distinct pH values (Bizzarri et al., 2006; Kuner and Augustine, 2000; Llopis et al., 1998). The absorption spectrum of GFP is widely used to measure the H^+ concentration in living organisms and optimized versions of GFP for pH measurement have been generated (Kneen et al., 1998). Among various forms of GFP, ecliptic and super ecliptic pHluorins display no fluorescence when protonated, which allows sensitive detection of biological processes associated with pH increases (De Angelis et al., 1998; Miesenbock et al., 1998). GFP-based pH indicators have been extensively employed to study alterations of intracellular pH related to network activity (Raimondo et al., 2012) and as markers of synaptic release (Miesenbock et al., 1998). However only a few studies employed them to monitor extracellular H^+ fluxes associated with neural excitability or under pathological conditions. Notably, very recently the superecliptic pHluorin has been targeted to synaptic compartments to detect release of protons associated with synaptic activity, suggesting the potential role of H^+ as extracellular messenger (Du et al., 2014; Wang et al., 2014). However, the use of superecliptic pHluorin, whose fluorescence is rapidly quenched in the acidic environment, makes it more sensitive to detect increases, rather than decreases, of pH.

The aim of the present study is to describe the spatio-temporal dynamics of extracellular pH fluctuations at synaptic sites during neuronal hyperexcitability. We employed E^2GFP to create a membrane-targeted extracellular ratiometric pH indicator. In contrast to other fluorescent pH sensors that are quenched at acidic pH, E^2GFP represents a very effective and exquisitely sensitive ratiometric pH indicator for acidic shifts. By monitoring variations of E^2GFP fluorescence in real time in primary hippocampal neurons, we were able to quantify pH fluctuations during network hyperactivity induced by convulsant drugs or high frequency stimulation. Sustained neural hyperactivity caused a pH decrease that was specifically localized at active synapses and was not attributable to the outflow of intravesicular protons into the cleft during intense exocytosis, but rather to the activation of the Na^+/H^+ exchanger (NHE). Our data may contribute to the understanding of the pathophysiological mechanisms associated with hyperexcitability in the epileptic brain.

RESULTS

Engineering and membrane targeting of a novel fluorescent sensor for extracellular pH

To engineer a genetically encoded indicator specifically sensitive to acidic pH shifts and targeted to the extracellular side of the plasma membrane, we inserted the E²GFP sequence (Bizzarri et al., 2006) into the mammalian expression vector pDisplayTM, which allows the exposure of proteins to the cell surface. The resulting construct was named “ex.E²GFP” (**Fig. 1A**). The construct was subsequently cloned into a lentiviral vector under the control of the ubiquitous phosphoglycerate kinase (PGK) promoter to allow effective transduction of primary neurons. The subcellular localization of the probe was studied in the HEK293 cell line and primary cortical neurons. Membrane localization was observed in both cell types (**Fig. 1B**). To quantify the extent of membrane targeting of the probe in primary neurons, we labeled membranes with the lipophilic dye FM4-64fx and performed 3D reconstruction by confocal microscopy. This analysis revealed that the vast majority of the probe was indeed localized to the plasma membrane, while the small pool of intracellular ex.E²GFP likely corresponded to the fraction of the protein that was being synthesized and translocated to the cell surface (**Fig. 1C,D**). To further demonstrate the plasma membrane expression of the probe, we performed subcellular fractionation of 17 DIV neurons infected with ex.E²GFP, lysed and analyzed by western blotting using anti-GFP antibodies (**Fig. 1E**). ex.E²GFP immunoreactivity was predominantly associated with the plasma membrane-containing fraction P2, compared to the cytosolic/microsomal fraction S2.

Characterization of the ex.E²GFP sensor

E²GFP is ideally suited for the ratiometric operation of the biosensor as its emission can increase, decrease or be independent of pH based on the excitation wavelength, and its pK_a in the cytosolic form is about 6.8 (Bizzarri et al., 2006), thus very effective to detect pH changes just below the physiological range. We asked whether the optical properties of E²GFP were affected by the exposure to the extracellular environment. Primary cortical neurons were infected with ex.E²GFP and fluorescence emission was evaluated at two extracellular pH values (5.8 and 7.4) using 405 and 488 nm as excitation wavelengths (Bizzarri et al., 2006; Bizzarri et al., 2009) (**Fig. 2A**). This analysis revealed a much higher fluorescence intensity at acidic pH compared to physiological pH when ex.E²GFP was excited at 405 nm, whereas an opposite behavior (albeit within a smaller dynamic range) was observed when 488 nm

excitation (**Fig. 2B**). Thus, the extracellular targeting of E²GFP does not alter its optical properties or its pH sensitivity.

Next, a calibration curve was made by analyzing ex.E²GFP fluorescence obtained by switching the excitation wavelength between 405 and 488 nm and by systematically changing the extracellular pH from 5.7 to 7.8. To render the signal independent of the expression level of the protein, ex.E²GFP fluorescence ratio ($\lambda_{\text{ex}405}/\lambda_{\text{ex}488}$) was calculated and found to be strongly dependent on extracellular pH (**Fig. 2C**). Data were fitted according to a logistic sigmoidal function, with a pK_a value of 6.9 ± 0.1 that was very similar to that for soluble E²GFP. This calibration curve was used to convert fluorescence ratios into absolute pH values throughout.

Neuronal hyperactivity causes localized extracellular acidic shifts in primary cultures

Several studies have addressed the question of how neuronal activity elicits changes in intracellular and extracellular H⁺ concentration, sometimes yielding contrasting results (Caldwell et al., 2013; Chen and Chesler, 2015; Cichy et al., 2015; de Curtis et al., 1998; DeVries, 2001; Lu et al., 2012; Palmer et al., 2003; Xiong et al., 2000; Xiong and Stringer, 2000; Zhang et al., 2010; Zhao et al., 2011). We studied the effects of sustained heightened neural activity on the extracellular H⁺ concentration by using *in vitro* models of epileptic-like network states characterized by synchronized neuronal discharges. Changes in fluorescence were monitored in primary cortical neurons infected with lentiviruses encoding ex.E²GFP by live imaging (0.4 Hz frame rate) and pH values were subsequently extrapolated from ratio values using the calibration curve shown in Fig. 2C. A scheme of the time course of the experiments is shown in **Fig. 3A**: pH was first measured under baseline conditions (2 min), then network hyperactivity was induced by blocking GABA_A receptors with bicuculline (30 μ M) and cultures were imaged for 10 min. Thereafter, the selective blocker of voltage-gated Na⁺ channels tetrodotoxin (TTX; 300 nM) was added to suppress firing and live imaging was continued for further 10 min. During the time frame of the experiments, ex.E²GFP fluorescence showed marked activity-dependent fluctuations, with no significant photobleaching (**Fig. S1**).

We found that *in vitro* seizure-like activity was associated with marked acidic pH transients in neurons. Interestingly, such fluorescence variations were observed both at the cell body level and along neurites of infected cells, where they mostly localized at synapse-like bouton structures (**Fig. 3B, Suppl. Movie 1**). We thus monitored extracellular pH

variations around these boutons and neuronal cell bodies. At both locations, acidic pH transients typically began at the onset of the hyperactivity triggered by the convulsant. After an initial delay, extracellular pH slowly but progressively decreased, reached a minimum acidic plateau 4-5 min after the beginning of the ‘seizure-like’ period and recovered towards baseline levels 4 min after termination of the seizure with TTX (**Fig. 3C,D**).

We subsequently asked how the time scale of increased neuronal firing compared to that of the acidic shift after addition of bicuculline. We cultured cortical neurons on 12-well multielectrode arrays (MEA) plates. Neurons were infected at day *in vitro* (DIV) 10 and analyzed at DIV15. The spiking rate was measured for 2 min under basal conditions, then bicuculline was added and the recording was continued for 10 min (**Fig. 3E**). In parallel, fluorescence was monitored in the same culture preparation. The bicuculline-induced increase in spiking rate was already observed 10 s after drug addition, while the pH decrease took about 4 min to reach the plateau. Thus, the increase in spiking activity happens first, followed by the pH decrease that appears to be a much slower process, likely due to the progressive extracellular accumulation of H⁺ in the extracellular/synaptic space.

High-frequency electrical stimulation also triggers acidosis localized at synaptic sites

To determine the variations of extracellular pH in response to a physiological electrical stimulation, cortical neurons expressing ex.E²GFP were challenged with a stimulation of 1200 action potentials (APs) at 10 Hz (**Fig. 4A, Suppl. Movie 2**). Electrical stimulation also triggered a fluorescence increase that was localized to discrete synapse-like boutons along the neurites. The acidic pH transient developed slowly during the train stimulation, and continued to reach an acidic plateau of pH 6.9 ± 0.1 about 50 sec after the end of the stimulation (**Fig. 4B,C**). pH levels remained acidic until the end of the recording, showing only a slow and modest recovery. At the cell body level the effects of electrical stimulation were much lighter and non statistically significant (**Fig. 4B,C**).

To assess whether synaptic acidification is present also under basal conditions, we analyzed the live imaging data corresponding to the ‘basal’ conditions of the various experiments. We identified single synapses and followed the behavior of ex.E²GFP fluorescence over time in the corresponding ROIs. Interestingly, we found that 27% of the analyzed synapses did not show appreciable pH shifts (> 0.1 pH units from basal), while 73% of synapses displayed spontaneous pH oscillations that reached in a small percentage of synapses values between 6.6 and 6.9 (**Fig. 4D-F**). Thus, the majority of synapses under basal conditions display spontaneous activity levels of sufficient intensity to generate synaptic pH shifts.

Hyperactivity-induced pH shifts occur at synaptic sites

The punctate pattern of fluorescence observed in the neuropil upon induction of network hyperactivity was reminiscent of synaptic contacts. To investigate the subcellular distribution of ex.E²GFP and unambiguously identify the sites of pH change, we quantified the relative abundance of the probe in cell bodies and synaptic boutons labeled by the lipophilic dye FM4-64 upon bicuculline stimulation. Interestingly, the vast majority of ex.E²GFP codistributed with FM4-64-labeled puncta along neurites (**Fig. 5A,B**). However it was not clear whether this was a specific activity-dependent synaptic acidification, or the mere consequence of a non-specific accumulation of the sensor to synaptic boutons. To demonstrate whether this was the case, primary cortical neurons were co-infected with lentiviruses expressing mCherry-synapsin I (syn I), used as marker of presynaptic terminals (Cesca et al., 2010) and ex.E²GFP. Comparable ex.E²GFP intensity levels were observed under basal conditions at synaptic (i.e. synapsin-positive) and extra-synaptic (i.e. synapsin-negative) areas, thus confirming the ubiquitous targeting of the probe and the synaptic localization of the pH shifts (**Fig. 5C-E**). Because the expression of ex.E²GFP is driven by the constitutive PGK promoter, the probe was expressed in both excitatory and inhibitory neurons. In fact, when primary cultures at various stages of maturation (between DIV 10 and 18) were double stained for GFP and markers of excitatory neurons (VGLUT1), inhibitory neurons (VGAT) and astrocytes (GFAP), ex.E²GFP was expressed in all three cell populations, although neuronal expression was predominant due to the paucity of glial cells under our culture conditions (**Fig. S2**).

Next, we investigated the specificity of the pH shifts for synaptic sites upon hyperactivity. A digital mask corresponding to presynaptic sites was built using the mCherry fluorescence, which was constant under resting and stimulated conditions (**Fig. S3A**), and the percent overlap of the ex.E²GFP fluorescence was calculated under basal conditions and upon bicuculline-induced hyperactivity. Stimulation caused an increase in ex.E²GFP colocalization with mCherry-syn I-positive puncta up to about 25-30% (**Fig. 5F,G**), demonstrating a massive recruitment of synaptic sites in the acidic response to the convulsant. On the other hand, when the mask was built using the ex.E²GFP fluorescence, which substantially increases in intensity and area upon stimulation (**Fig. S3A**), the ex.E²GFP overlap on syn I-labeled synaptic sites did not change (\approx 60%) (**Fig. S3B**), confirming the synaptic localization of ex.E²GFP before and after stimulation and the synaptic localization of the pH shifts.

Not all synapses labeled by synaptic markers such as syn I are functionally active (silent synapses). To relate ex.E²GFP-positive puncta to active synapses we co-infected hippocampal neurons with lentiviruses expressing ex.E²GFP and synaptobrevin-pHluorin mOrange2 (Syb2O), a marker of SV exo-endocytosis at active synaptic sites with a non-overlapping emission spectrum (Ramirez et al., 2012). Upon electrical stimulation with 1200 AP@10 Hz, ex.E²GFP fluorescence and Syb2O fluorescence overlapped by approximately 60%, confirming the localization of the probe to active synapses (**Fig. S3C,D**).

We then asked whether the observed activity-dependent synaptic pH shifts were present in the whole synapse population, or whether they were preferentially localized to excitatory or inhibitory synapses. Primary cortical neurons were infected with ex.E²GFP and live-labeled with VGAT-oyster650, a specific marker of active inhibitory synapses (**Fig. 5H**). A digital mask corresponding to the active GABAergic boutons was built using the loaded VGAT-oyster650 fluorescence, which remained constant under resting and stimulated conditions (**Fig. S3E**), and the percent overlap of the ex.E²GFP fluorescence was calculated under basal conditions and upon bicuculline-induced hyperactivity. Stimulation caused an increase in ex.E²GFP colocalization up to 35% of the active GABAergic synapses (**Fig. 5I**). On the other hand, when the mask was built using the ex.E²GFP fluorescence, which increases in intensity and area upon stimulation, the ex.E²GFP overlap on VGAT-labeled synaptic sites was roughly constant ($\approx 30\%$)(**Fig. S3E-G**).

Hyperactivity-induced pH shifts at synapses are not contributed by synaptic vesicle exocytosis

Neuronal activity triggers the mobilization of synaptic vesicles (SVs), which are characterized by an acidic luminal pH due to the activity of the active proton transporter vATPase (Sudhof, 2013). ex.E²GFP, initially exposed on the plasma membrane at synapses, could get internalized in the acidic environment of endocytosed SVs. If this were the case, then the increase in ex.E²GFP fluorescence would reflect the re-acidification of the endocytosed SVs, rather than a drop in the extracellular pH. To sort out this issue, we treated primary neurons with bafilomycin, a specific inhibitor of vATPase, which prevents re-acidification of SVs after exocytosis (Forgac, 1989). Pretreatment of neurons with bafilomycin before the addition of bicuculline did not affect either the bicuculline-induced acidification, or the temporal profile of the pH decrease (**Fig. 6A,B**). These data demonstrate that acidification reflects a variation of the extracellular environment, and that the H⁺ release

of protons from exocytosed SVs and/or the incorporation of the vATPase on the plasma membrane during SV exocytosis are not primarily involved.

Hyperactivity-induced pH shifts at synapses are due to the activity of the Na⁺/H⁺ exchanger

Since the pH shifts at synaptic sites are not attributable to release of the content of SV protons in the synaptic cleft, we addressed the possibility that the secretion of H⁺ during hyperactivity occurs via the activation of the NHE exchanger. We pretreated primary neurons with the NHE inhibitor amiloride before the addition of bicuculline (**Fig. 7A,B**). The treatment with amiloride alone induced a slight, but significant, extracellular alkalinization, confirming the presence of a constitutive activity of NHE under basal conditions. Interestingly, when amiloride was associated with bicuculline stimulation, the extent of hyperactivity-induced extracellular acidification was significantly decreased, demonstrating the involvement of NHEs in mediating the activity-induced synaptic pH shifts (**Fig. 7C,D**). As amiloride also inhibits Na⁺/Ca²⁺ exchangers, we performed bicuculline treatment in the presence of 1 μM cariporide, a specific inhibitor of NHE isoform 1, one of the most expressed NHE neuronal isoforms (Harguindey et al., 2013; Luo et al., 2005). We found that synaptic acidification was strongly inhibited under this experimental condition, indicating a major role of NHE1 in the observed hyperactivity-induced synaptic acidification (**Fig. 7E-H**).

We also tested whether the blockade of bicuculline-induced extracellular acidification by amiloride was the result of an inhibition of network firing activity by the drug. To this aim, primary neurons were cultured on multi-well microelectrode arrays and the network activity analyzed in terms of firing rate, burst frequency and burst duration in the presence of the various treatments by following the same experimental protocol described in Fig. 7A (**Fig. 8A**). Notably, amiloride alone, bicuculline alone and amiloride+bicuculline all increased both the firing and the bursting frequencies (**Fig. 8B-E**), although acidification occurred only with bicuculline alone. The excitatory effect of amiloride on network activity confirms previous observations reporting a reduction of the sensitivity of GABA_A receptors toward GABA by amiloride (Fisher, 2002). Altogether, our results demonstrate the key role of NHE in coupling firing activity with a shift in extracellular pH.

DISCUSSION

Extracellular pH has important effects on neuronal activity, and in turn neuronal activity is an important determinant of the extracellular H^+ concentration. Protons released by the SVs during sustained activity act as a pseudo-transmitter, modifying several actors governing excitatory and inhibitory synaptic transmission (Cho and von Gersdorff, 2014; Du et al., 2014; Wang et al., 2014). Extracellular acidification occurs in the brain with elevated neural activity and neuronal injury (Wemmie et al., 2013). In turn, several studies demonstrated pH sensitivity of synaptic components like voltage-gated Ca^{2+} channels (VGCCs) (Iijima et al., 1986), ligand-gated AMPA and NMDA receptors (Traynelis and Cull-Candy, 1990) and GABA receptors (GABARs) (Kaila, 1994). In addition, acid sensing ion channels (ASICs) are activated by extracellular low pH and are associated with neuronal dysfunction or brain injury (Friese et al., 2007; Yin et al., 2013). However, most of the studies were done on brain slices applying high-speed pH microelectrodes (Fedirko et al., 2006) or dyes (Shuba et al., 2008), which lack accurate spatial resolution to localize pH changes to restricted neuronal domains.

In this paper we analyzed the dynamics of extracellular pH changes accompanying epileptic-like activity in primary neuronal networks. We engineered a genetically encoded ratiometric GFP-based pH sensor that provided high spatial and temporal resolution to quantify variations in extracellular pH. Superecliptic GFP (pHluorin) has been employed as an extracellular pH sensor when targeted to the synaptic cleft (Du et al., 2014; Wang et al., 2014). However, pHluorin is an excellent reporter of alkalinization but quenches very rapidly upon acidic pH shifts. We therefore focused our attention on E^2GFP (Bizzarri et al., 2006), a ratiometric GFP derivative that successfully allowed the live monitoring of the changes of intracellular pH towards acidosis (Bizzarri et al., 2006; Raimondo et al., 2012). We fused E^2GFP to the N-terminus of the PDGF receptor transmembrane domain that exposes it to the extracellular side of the plasma membrane (ex. E^2GFP). We demonstrated that the exposure to the extracellular environment does not affect E^2GFP optical properties and pH sensitivity, and that the targeted pH sensor shows high photostability with negligible photobleaching. Such chimera had several advantages, namely: (i) it is ubiquitously targeted to the plasma membrane, thus being potentially able to map acidification in any region of the neuronal network; (ii) it is a ratiometric sensor with opposite pH sensitivity depending on the excitation wavelength, thus allowing to report pH changes independently of the expression

level; and (iii) when excited at 405 nm, its fluorescence increases as the pH drops, making it more precise and sensitive in evaluating acidification than the previously used pHluorins.

We characterized hyperactivity-induced acidification by chemical stimulation with bicuculline and by high frequency electrical stimulation. Using chemical stimulation, we observed a relatively slow build-up of extracellular acidification that occurred at both cell bodies and synaptic sites and slowly returned to neutral pH upon silencing of electrical activity with TTX, with no apparent pre- or post-alkalinization phases. On the other hand, the more physiological electrical stimulation showed a slower and milder synaptic acidification specifically localized to synaptic boutons. Interestingly, our probe was sensitive enough to detect also smaller acidic shifts associated with spontaneous activity. While ex.E²GFP, exquisitely sensitive to the acidic shifts, may not be ideal to monitor alkaline shifts, the novelty of our results lays in the following points: (i) the increased network firing causes acidification at cell bodies and synapses, in both excitatory and inhibitory neurons; (ii) acidification does not result from the endocytotic capture of the membrane-exposed probe by SVs; (iii) acidification is mostly contributed by the NHE.

The acidic shifts in synaptic pH occurring during intense neuronal activity imply quantal/ non-quantal release of protons from neurons. In principle, this can occur through (i) exocytosis of SVs; (ii) exposure of vATPase on the plasma membrane; (iii) increased production of CO₂ and metabolic acids such as lactate that is externalized by monocarboxylate transporters; (iv) activation of NHE (mostly the NHE1/SLC9A1) that transports Na⁺ into the cells and extrudes H⁺; and (v) activation of Na⁺-dependent anion (Cl⁻/HCO₃⁻) exchangers (NDAE: SLC4A8 and SLC4A10) that transport Na⁺ and HCO₃⁻ into the cell and extrude protons and Cl⁻ (Sinning and Hubner, 2013). Moreover, a glial source of protons cannot be excluded (Grichtchenko and Chesler, 1994), although this is unlikely the case in our cultures, which are virtually devoid of glial cells.

The demonstration that reduction of extracellular pH occurs in synaptic regions is consistent with recent studies (Cho and von Gersdorff, 2014; Du et al., 2014; Highstein et al., 2014; Wang et al., 2014). The synaptic distribution of acidic shifts emphasizes that an alteration of synaptic transmission is the *primum movens* for the initiation and maintenance of ictal activity (Lazarevic et al., 2013). Release of SV protons in the synaptic cleft upon intense exocytosis has been recently proposed as a mechanism of protonergic transmission (Du et al., 2014; Wang et al., 2014). However, this was not the case under our experimental conditions, as the long delay between the onset of the hyperactivity state and the acidic shift makes it unlikely that the un-buffering of SV protons upon SV fusion (DeVries, 2001)

significantly contributes to the acidic shift. Moreover, blockade of SV acidification does not affect the kinetics or the extent of acidification, ruling out proton release by SVs or sustained exposure of SV vATPase on the plasma membrane. This discrepancy is likely due to the different experimental models used in the various studies, i.e. *in vitro* cultures of mouse cortical neurons (present study), zebrafish retinal neurons (Wang et al., 2014), and mouse amygdala neurons in acute brain slices (Du et al., 2014). In addition, we specifically study the pH shifts caused by epileptic-like activity, which likely rely on different mechanisms with respect to physiological network activity.

Amongst the other potential mechanisms of extracellular acidification, we focused on the NHEs, as several members of this family are known to contribute to the onset and maintenance of epilepsy and other developmental brain disorders (recently reviewed in (Zhao et al., 2016)). Our data indicate that acidification during ictal activity is mostly due to NHE, as the addition of the NHE blocker amiloride strongly reduced the bicuculline-induced synaptic acidification. NHE1 is one of the main NHE isoforms expressed in the CNS, and its mutations are known to cause epilepsy in mice (Bell et al., 1999; Cox et al., 1997). NHE1 is activated in response to intracellular acidification, and in neurons, the consequent extrusion of H⁺ into the confined space of the synaptic cleft could generate the acidic pH shifts that we observed in our cultures. The other main neuronal NHE isoform is NHE5 (Diering et al., 2011; Diering and Numata, 2014). However, treatment with the specific NHE1 inhibitor cariporide (Harguindey et al., 2013; Luo et al., 2005) virtually abolished the bicuculline induced pH shifts, thus suggesting that NHE1 is the main isoform involved in the observed synaptic acidification.

In summary, we describe a novel sensor for extracellular pH and use it to study the wave of extracellular acidification during hyperactivity of neuronal network, which is independent of protons released by SVs, and mostly contributed by the activity of NHE exchangers. These results underline the strict links between synaptic activity and synaptic pH. The extracellular sensor engineered in this study could be employed for better understanding of ionic dynamics under physiological and pathological conditions with high spatio-temporal precision.

MATERIALS AND METHODS

Reagents. All biochemical reagents and drugs were purchased from Sigma-Aldrich (St. Louis, MO), unless otherwise specified. Tissue culture reagents and media were from Gibco/Thermo-Fischer Scientific (Waltham, MA, USA) or Sigma-Aldrich. The following primary antibodies were used: anti-GFP (Thermo Fisher Scientific, #A-11122, dilution 1:500; Synaptic System, #132004, dilution 1:500 for triple staining with anti-homer and anti-bassoon antibodies), anti-glial fibrillary acidic protein (GFAP; Sigma-Aldrich, #G3893, dilution 1:500), anti-vesicular glutamate transporter-1 (VGLUT1, Synaptic Systems, Goettingen, Germany; #1350304, dilution 1:500), anti-mCherry (Clontech, #632543, dilution 1:500), anti-Homer (Synaptic Systems, #16011, dilution 1:500), anti-Bassoon (Synaptic System; #141002, dilution 1:500), anti-PSD95 (Synaptic System, #124011, dilution 1:1000) and anti-actin (Sigma-Aldrich, #A4700, dilution 1:1000). Fluorophore-conjugated secondary antibodies were Alexa Fluor® 488 anti-rabbit (#A1108; dilution 1:1000), Alexa Fluor® 488 anti-guinea pig Alexa Fluor® 568 anti-mouse (#A-11004; dilution 1:1000), Alexa Fluor® 546 anti-rabbit (#A-11035 dilution 1:1000), Alexa Fluor® 647 anti-mouse (#A-21236; dilution 1:1000) and Alexa Fluor® 647 anti-guinea pig (#A-2145; dilution 1:1000; all Thermo Fisher Scientific).

Plasmids and transfection. The ex.E²GFP sequence was amplified from the pcDNA3.1 + E²GFP plasmid (kindly provided by dr. Bizzarri, CNR Pisa, Italy) (Bizzarri et al., 2006) and inserted into the pDisplay™ vector (Thermo Fisher Scientific) between the BglII and PstI sites, to obtain pDisplay-E²GFP (ex.E²GFP). The ex.E²GFP sequence was subsequently cloned into the pLenti-hPGK viral vector (kind gift of drs. Amendola and Naldini, Tiget, Milano, Italy) between the AgeI and XhoI sites. The sequence of all constructs was verified by direct sequencing. Vectors were transiently transfected into cultured cells using Lipofectamine 2000 (Thermo-Fischer Scientific).

Cell culture procedures. All experiments were carried out in accordance with the guidelines established by the European Community Council (Directive 2010/63/EU of 22 September 2010) and approved by the Italian Ministry of Health. Human embryonic kidney cells (HEK293, purchased from ATCC, Manassas, VA, USA) and primary cortical cultures from wild-type C57Bl/6 mice (Charles River, Calco, Italy) were prepared and maintained following standard procedures (Baldelli et al., 2007; Cesca et al., 2015).

Viral transduction procedures. The production of VSV-pseudotyped third-generation lentiviruses was performed as previously described (De Palma and Naldini, 2002). Primary neurons were infected with ex.E²GFP alone or co-infected with mCherry-syn I (Verstegen et al., 2014) at 10DIV at 10 multiplicity of infection and the experiments were performed 7-8 days post-infection (17-18DIV). For MEA experiments, neurons were infected at 10DIV and recordings performed at 17-18DIV. For pHluorin experiments, neurons were co-infected with Synaptobrevin-pHluorin m-Orange2 (Syb2O) (Ramirez et al., 2012) and ex.E²GFP lentiviruses at 13DIV, and optical recordings performed at 17-18DIV. Subcellular fractionation of transduced 17DIV neurons was performed as previously described (Huttner et al., 1983).

Immunofluorescence and confocal microscopy. Immunohistochemistry on fixed cells was performed following standard procedures (Cesca et al., 2015). Coverslips were imaged by confocal microscopy (SP8, Leica Microsystems GmbH, Wetzlar, Germany) using a 63× (1.4 NA) magnification lens. Confocal images were analyzed with the Leica LAS AF software (Leica Application Suite Advance Fluorescence, version 3.3, Leica Microsystems). Z-stack images for 3D reconstruction were acquired every 300 nm steps and offline analysis for colocalization quantification and intensity profiles determination was performed using the JACoP plugin of the ImageJ software (Bolte and Cordelieres, 2006).

Time-lapse live cell imaging. For live cell imaging experiments, an SP8 confocal laser-scanning microscope was used. The sequential excitation of ex.E²GFP at 405 nm and 488 nm was achieved with a multiline argon laser. Emitted fluorescence was collected between 510 and 560 nm using a single photomultiplier tube (PMT) at a constant voltage. To verify fluctuations in laser intensity and bleaching of the probe, fluorescence was recorded in the absence of stimulation for the whole duration of the typical experiments. Neurons and cells were grown and maintained in 25 mm cover glasses and placed in a chamber with Tyrode solution (140 mM NaCl, 4 mM KCl, 1 mM MgCl₂, 2 mM CaCl₂, 10 mM glucose and 10 mM HEPES at pH 7.4). To build the calibration curve, neurons were incubated in solutions at controlled pH. For the acidic buffer (pH 5.8-6.2), HEPES was replaced with MES. Solutions were applied with a valve control perfusion system (Warner Instruments, Hamden, CT, USA) after 2 min of baseline acquisition in physiological solution; 30 μM of bicuculline were

added, followed by 300 μM of TTX. For amiloride experiments, after 2 min of baseline acquisition, 20 μM amiloride were added, followed by 30 μM bicuculline. For cariporide (Tocris, Avonmouth, Bristol, UK) experiments, after 2 min of baseline acquisition, 1 μM cariporide was added, followed by 30 μM bicuculline. The acquisition rate was 0.4 Hz for each excitation wavelength. For field stimulation, time-lapse images were acquired every 2.58 s for 360 sec. Action potentials were evoked by passing 1 ms current pulses, yielding fields of 10 V/cm, through platinum-iridium electrodes using an AM 2100 stimulator (AM Systems, Carlsborg, WA, USA). After 30 s of baseline acquisition, neurons were stimulated with a train of 1200 APs at 10 Hz.

Image analysis. Confocal images were analyzed with the Leica LAS X software. First, regions of interest (ROIs) were chosen for each image. For cell bodies (CBs), the area of each ROI was 5 μm^2 , while for nerve terminals (NTs) was 2 μm^2 , located on the center of mass of each NT. Fluorescence intensity ratios (R) were calculated according to the equation below:

$$R = \frac{MV(\lambda_x 405 - \lambda_e 510 - 560) - MV(\text{background})}{MV(\lambda_x 488 - \lambda_e 510 - 560) - MV(\text{background})}$$

where MV (mean value) is the averaging pixel intensity for each ROI and $\lambda_x - \lambda_e$ is the excitation (collection) wavelength. pH values were calculated by fitting each average fluorescence intensity trace with the sigmoidal dose-response equation below:

$$y = -0,307 + \frac{(4,83 - 0,372)}{1 + 10^{6,9-x}}$$

Images were acquired with constant gain and exposure times across all experiments.

MEA recordings. Dissociated cortical neurons were plated onto 12-well planar Multi Electrode Arrays (768-GL1-30Au200 from Axion BioSystems, Atlanta, GA, USA), consisting of 768 electrodes in 12 wells (64 electrodes/well). Extracellular spontaneous activity was recorded throughout the experiment and analyzed as previously described (Cesca et al., 2015; Chiappalone et al., 2009). Data were analyzed using the Neural metric tool (Axion Biosystems, Atlanta, GA, USA).

Quantitative live cell imaging for mCherry-syn I, VGAT-Oyster650 and pHluorin experiments. Co-infected mCherry-syn I/ex.E²GFP-expressing neurons were kept in a stimulation chamber (Warner Instruments, Hamden, CT, USA) and imaged in Tyrode solution. Bicuculline (30 μ M) was added to the medium and images were acquired after 10 min incubation using an SP8 fluorescence confocal laser-scanning microscope as described above. Five fields were analyzed using the JaCoP plug-in of ImageJ (Bolte and Cordelieres, 2006).

For pHluorin experiments, Syb2O/ex.E²GFP-expressing neurons were kept in a stimulation chamber and imaged in physiological solution. The chamber was positioned on the stage of an IX-81 motorized inverted epifluorescence microscope (Olympus, Tokio, Japan). An MT20 Hg-Xe lamp (Olympus, Tokio, Giappone) was used as a light source with 480 \pm 20 nm excitation, 495 nm dichroic and 525 \pm 50 nm emission filters to detect ex.E²GFP signal. 560 \pm 40 nm excitation, 585 nm dichroic and 630 \pm 75 nm emission filters were instead used to detect the Syb2O signal. An electrical stimulation of 400 APs at 20 Hz was delivered to visualize active synapses. Images were acquired with a Hamamatsu Orca-ER CCD camera (Hamamatsu Photonics, Hamamatsu City, Japan) using an UplanSapo 60X1.35 NA oil immersion objective (Olympus, Tokio, Japan) and the software Excellence RT (Olympus, Tokio, Japan). Fields expressing both signals were analyzed to evaluate the co-localization under basal and stimulated conditions; 11 fields were analyzed using the JaCoP plugin of ImageJ (Bolte and Cordelieres, 2006).

Labeling with VGAT-Oyster650 (Synaptic Systems, Goettingen, Germany; #131103; dilution 1:200) was performed by incubation of ex.E²GFP-infected neurons for 20 min at 37 $^{\circ}$ C, 5% CO₂, to load the dye into inhibitory presynaptic terminals by spontaneous activity. For basal conditions, images were acquired in the Tyrode solution. For the hyperexcited state, 30 μ M bicuculline was added to the medium and images were acquired after 10 min of incubation. Experiments were performed at an SP8 fluorescence confocal laser-scanning microscope, as described above. For all the experiments, colocalization was analyzed by generating Mander's colocalization coefficients (Manders et al., 1992).

Statistical Analysis. Data are expressed as means \pm sem for number of cells (n) and mouse preparations as detailed in the figure legends. Normal distribution of data was assessed using D'Agostino-Pearson's normality test (for n>8 values/experimental group) or the

Kolmogorov-Smirnov test ($n < 8$ values/experimental group). To compare two normally distributed sample groups, the paired or unpaired Student's *t*-test was used. To compare more than two normally distributed sample groups, one- or two-way ANOVA, followed by the Bonferroni's multiple comparison test was used. Alpha levels for all tests were 0.05% (95% confidence intervals). Statistical analysis was carried out using Origin Pro 9.1 (OriginLab Corp., Northampton, MA) and GraphPad Prism 5 (GraphPad Software, Inc., La Jolla, CA) softwares.

Acknowledgements. We thank dr. Kavalali (Dept. Neuroscience, The University of Texas Southwestern Medical Center, Dallas, TX) for the SynaptobrevinOrange2 constructs, dr. Naldini (Tiget, Milano, Italy) for lentiviral constructs and protocols, and drs. Nanni and Mehili (Istituto Italiano di Tecnologia, Genova, Italy) for assistance in the preparation of primary cultures. We are also grateful to Ms Donato for helping in the initial phases of the project and to drs. Moretti and Pecoraro for helping with the MEA experiments.

Conflict of interest. No competing interests declared.

Author Contributions. MC and SL performed the cell and molecular biology studies and the live imaging experiments and analyzed the data. MB performed image analysis and quantification. MF performed the SybO2 experiments. FC and AF supervised the live imaging and cell biology experiments, analyzed data and contributed to paper writing. FB and FC designed and supervised research, wrote the paper and supported the study.

Funding. This study was supported by research grants from the European Commission (FP7 Integrating Project "Desire" n. 602531 and ITN "ECMED" n. 642881 to FB). The support of Telethon-Italy (Grant GGP13033 to FB), CARIPLO Foundation (Grant 2013 0879 to FB) and the Italian Ministry of Health Ricerca Finalizzata 2013 (to FB) is also acknowledged.

REFERENCES

- Baldelli, P., Fassio, A., Valtorta, F. and Benfenati, F.** (2007). Lack of synapsin I reduces the readily releasable pool of synaptic vesicles at central inhibitory synapses. *J Neurosci* **27**, 13520-31.
- Bell, S. M., Schreiner, C. M., Schultheis, P. J., Miller, M. L., Evans, R. L., Vorhees, C. V., Shull, G. E. and Scott, W. J.** (1999). Targeted disruption of the murine Nhe1 locus induces ataxia, growth retardation, and seizures. *Am J Physiol* **276**, C788-95.
- Bizzarri, R., Arcangeli, C., Arosio, D., Ricci, F., Faraci, P., Cardarelli, F. and Beltram, F.** (2006). Development of a novel GFP-based ratiometric excitation and emission pH indicator for intracellular studies. *Biophys J* **90**, 3300-14.
- Bizzarri, R., Serresi, M., Luin, S. and Beltram, F.** (2009). Green fluorescent protein based pH indicators for in vivo use: a review. *Anal Bioanal Chem* **393**, 1107-22.
- Bolte, S. and Cordelieres, F. P.** (2006). A guided tour into subcellular colocalization analysis in light microscopy. *J Microsc* **224**, 213-32.
- Caldwell, L., Harries, P., Sydlik, S. and Schwiening, C. J.** (2013). Presynaptic pH and vesicle fusion in Drosophila larvae neurones. *Synapse* **67**, 729-40.
- Cesca, F., Baldelli, P., Valtorta, F. and Benfenati, F.** (2010). The synapsins: key actors of synapse function and plasticity. *Prog Neurobiol* **91**, 313-48.
- Cesca, F., Satapathy, A., Ferrea, E., Nieu, T., Benfenati, F. and Scholz-Starke, J.** (2015). Functional Interaction between the Scaffold Protein Kidins220/ARMS and Neuronal Voltage-Gated Na⁺ Channels. *J Biol Chem* **290**, 18045-55.
- Chen, H. Y. and Chesler, M.** (2015). Autocrine boost of NMDAR current in hippocampal CA1 pyramidal neurons by a PMCA-dependent, perisynaptic, extracellular pH shift. *J Neurosci* **35**, 873-7.
- Chen, J. C. and Chesler, M.** (1992). pH transients evoked by excitatory synaptic transmission are increased by inhibition of extracellular carbonic anhydrase. *Proc Natl Acad Sci U S A* **89**, 7786-90.
- Chesler, M.** (2003). Regulation and modulation of pH in the brain. *Physiol Rev* **83**, 1183-221.
- Chesler, M. and Kaila, K.** (1992). Modulation of pH by neuronal activity. *Trends Neurosci* **15**, 396-402.
- Chiappalone, M., Casagrande, S., Tedesco, M., Valtorta, F., Baldelli, P., Martinoia, S. and Benfenati, F.** (2009). Opposite changes in glutamatergic and GABAergic transmission underlie the diffuse hyperexcitability of synapsin I-deficient cortical networks. *Cereb Cortex* **19**, 1422-39.
- Cho, S. and von Gersdorff, H.** (2014). Proton-mediated block of Ca²⁺ channels during multivesicular release regulates short-term plasticity at an auditory hair cell synapse. *J Neurosci* **34**, 15877-87.
- Cichy, A., Ackels, T., Tsitoura, C., Kahan, A., Gronloh, N., Sochtig, M., Engelhardt, C. H., Ben-Shaul, Y., Muller, F., Spehr, J. et al.** (2015). Extracellular pH Regulates Excitability of Vomeronasal Sensory Neurons. *J Neurosci* **35**, 4025-39.

Cox, G. A., Lutz, C. M., Yang, C. L., Biemesderfer, D., Bronson, R. T., Fu, A., Aronson, P. S., Noebels, J. L. and Frankel, W. N. (1997). Sodium/hydrogen exchanger gene defect in slow-wave epilepsy mutant mice. *Cell* **91**, 139-48.

De Angelis, D. A., Miesenbock, G., Zemelman, B. V. and Rothman, J. E. (1998). PRIM: proximity imaging of green fluorescent protein-tagged polypeptides. *Proc Natl Acad Sci U S A* **95**, 12312-6.

de Curtis, M., Manfredi, A. and Biella, G. (1998). Activity-dependent pH shifts and periodic recurrence of spontaneous interictal spikes in a model of focal epileptogenesis. *J Neurosci* **18**, 7543-51.

De Palma, M. and Naldini, L. (2002). Transduction of a gene expression cassette using advanced generation lentiviral vectors. *Methods Enzymol* **346**, 514-29.

DeVries, S. H. (2001). Exocytosed protons feedback to suppress the Ca²⁺ current in mammalian cone photoreceptors. *Neuron* **32**, 1107-17.

Diering, G. H., Mills, F., Bamji, S. X. and Numata, M. (2011). Regulation of dendritic spine growth through activity-dependent recruitment of the brain-enriched Na⁽⁺⁾/H⁽⁺⁾ exchanger NHE5. *Mol Biol Cell* **22**, 2246-57.

Diering, G. H. and Numata, M. (2014). Endosomal pH in neuronal signaling and synaptic transmission: role of Na⁽⁺⁾/H⁽⁺⁾ exchanger NHE5. *Front Physiol* **4**, 412.

Dietrich, C. J. and Morad, M. (2010). Synaptic acidification enhances GABAA signaling. *J Neurosci* **30**, 16044-52.

Du, J. Y., Reznikov, L. R., Price, M. P., Zha, X. M., Lu, Y., Moninger, T. O., Wemmie, J. A. and Welsh, M. J. (2014). Protons are a neurotransmitter that regulates synaptic plasticity in the lateral amygdala. *Proceedings of the National Academy of Sciences of the United States of America* **111**, 8961-8966.

Dulla, C. G., Dobelis, P., Pearson, T., Frenguelli, B. G., Staley, K. J. and Masino, S. A. (2005). Adenosine and ATP link PCO₂ to cortical excitability via pH. *Neuron* **48**, 1011-23.

Fedirko, N., Svichar, N. and Chesler, M. (2006). Fabrication and use of high-speed, concentric h⁺- and Ca²⁺-selective microelectrodes suitable for in vitro extracellular recording. *J Neurophysiol* **96**, 919-24.

Fisher, J. L. (2002). Amiloride inhibition of gamma-aminobutyric acid(A) receptors depends upon the alpha subunit subtype. *Mol Pharmacol* **61**, 1322-8.

Forgac, M. (1989). Structure and function of vacuolar class of ATP-driven proton pumps. *Physiol Rev* **69**, 765-96.

Friese, M. A., Craner, M. J., Etzensperger, R., Vergo, S., Wemmie, J. A., Welsh, M. J., Vincent, A. and Fugger, L. (2007). Acid-sensing ion channel-1 contributes to axonal degeneration in autoimmune inflammation of the central nervous system. *Nat Med* **13**, 1483-9.

Gottfried, J. A. and Chesler, M. (1996). Temporal resolution of activity-dependent pH shifts in rat hippocampal slices. *J Neurophysiol* **76**, 2804-7.

Grichtchenko, I. and Chesler, M. (1994). Depolarization-induced acid secretion in gliotic hippocampal slices. *Neuroscience* **62**, 1057-70.

Harguindey, S., Arranz, J. L., Polo Orozco, J. D., Rauch, C., Fais, S., Cardone, R. A. and Reshkin, S. J. (2013). Cariporide and other new and powerful NHE1 inhibitors as potentially selective anticancer drugs--an integral molecular/biochemical/metabolic/clinical approach after one hundred years of cancer research. *J Transl Med* **11**, 282.

Highstein, S. M., Holstein, G. R., Mann, M. A. and Rabbitt, R. D. (2014). Evidence that protons act as neurotransmitters at vestibular hair cell-calyx afferent synapses. *Proc Natl Acad Sci U S A* **111**, 5421-6.

Huttner, W. B., Schiebler, W., Greengard, P. and De Camilli, P. (1983). Synapsin I (protein I), a nerve terminal-specific phosphoprotein. III. Its association with synaptic vesicles studied in a highly purified synaptic vesicle preparation. *J Cell Biol* **96**, 1374-88.

Iijima, T., Ciani, S. and Hagiwara, S. (1986). Effects of the external pH on Ca channels: experimental studies and theoretical considerations using a two-site, two-ion model. *Proc Natl Acad Sci U S A* **83**, 654-8.

Inamura, K., Smith, M. L., Hansen, A. J. and Siesjo, B. K. (1989). Seizure-induced damage to substantia nigra and globus pallidus is accompanied by pronounced intra- and extracellular acidosis. *J Cereb Blood Flow Metab* **9**, 821-9.

Kaila, K. (1994). Ionic basis of GABAA receptor channel function in the nervous system. *Prog Neurobiol* **42**, 489-537.

Kim, Y. and Trussell, L. O. (2009). Negative shift in the glycine reversal potential mediated by a Ca²⁺- and pH-dependent mechanism in interneurons. *J Neurosci* **29**, 11495-510.

Kneen, M., Farinas, J., Li, Y. and Verkman, A. S. (1998). Green fluorescent protein as a noninvasive intracellular pH indicator. *Biophys J* **74**, 1591-9.

Krishtal, O. A., Osipchuk, Y. V., Shelest, T. N. and Smirnov, S. V. (1987). Rapid extracellular pH transients related to synaptic transmission in rat hippocampal slices. *Brain Res* **436**, 352-6.

Kuner, T. and Augustine, G. J. (2000). A genetically encoded ratiometric indicator for chloride: capturing chloride transients in cultured hippocampal neurons. *Neuron* **27**, 447-59.

Lazarevic, V., Pothula, S., Andres-Alonso, M. and Fejtova, A. (2013). Molecular mechanisms driving homeostatic plasticity of neurotransmitter release. *Front Cell Neurosci* **7**, 244.

Lei, S., Orser, B. A., Thatcher, G. R., Reynolds, J. N. and MacDonald, J. F. (2001). Positive allosteric modulators of AMPA receptors reduce proton-induced receptor desensitization in rat hippocampal neurons. *J Neurophysiol* **85**, 2030-8.

Llopis, J., McCaffery, J. M., Miyawaki, A., Farquhar, M. G. and Tsien, R. Y. (1998). Measurement of cytosolic, mitochondrial, and Golgi pH in single living cells with green fluorescent proteins. *Proc Natl Acad Sci U S A* **95**, 6803-8.

Lu, Y., Yi, L., Liu, D., Li, J., Sun, L. and Zhang, Z. (2012). Alkalosis leads to the over-activity of cortical principal neurons. *Neurosci Lett* **525**, 117-22.

Luo, J., Chen, H., Kintner, D. B., Shull, G. E. and Sun, D. (2005). Decreased neuronal death in Na⁺/H⁺ exchanger isoform 1-null mice after in vitro and in vivo ischemia. *J Neurosci* **25**, 11256-68.

Magnotta, V. A., Heo, H. Y., Dlouhy, B. J., Dahdaleh, N. S., Follmer, R. L., Thedens, D. R., Welsh, M. J. and Wemmie, J. A. (2012). Detecting activity-evoked pH changes in human brain. *Proc Natl Acad Sci U S A* **109**, 8270-3.

Makani, S. and Chesler, M. (2007). Endogenous alkaline transients boost postsynaptic NMDA receptor responses in hippocampal CA1 pyramidal neurons. *J Neurosci* **27**, 7438-46.

Makani, S. and Chesler, M. (2010). Rapid rise of extracellular pH evoked by neural activity is generated by the plasma membrane calcium ATPase. *J Neurophysiol* **103**, 667-76.

Manders, E. M., Stap, J., Brakenhoff, G. J., van Driel, R. and Aten, J. A. (1992). Dynamics of three-dimensional replication patterns during the S-phase, analysed by double labelling of DNA and confocal microscopy. *J Cell Sci* **103 (Pt 3)**, 857-62.

McCormick, D. A. and Contreras, D. (2001). On the cellular and network bases of epileptic seizures. *Annu Rev Physiol* **63**, 815-46.

Miesenbock, G., De Angelis, D. A. and Rothman, J. E. (1998). Visualizing secretion and synaptic transmission with pH-sensitive green fluorescent proteins. *Nature* **394**, 192-5.

Nedergaard, M., Goldman, S. A., Desai, S. and Pulsinelli, W. A. (1991). Acid-induced death in neurons and glia. *J Neurosci* **11**, 2489-97.

Palmer, M. J., Hull, C., Vigh, J. and von Gersdorff, H. (2003). Synaptic cleft acidification and modulation of short-term depression by exocytosed protons in retinal bipolar cells. *J Neurosci* **23**, 11332-41.

Raimondo, J. V., Irkle, A., Wefelmeyer, W., Newey, S. E. and Akerman, C. J. (2012). Genetically encoded proton sensors reveal activity-dependent pH changes in neurons. *Front Mol Neurosci* **5**, 68.

Ramirez, D. M., Khvotchev, M., Trauterman, B. and Kavalali, E. T. (2012). Vt1a identifies a vesicle pool that preferentially recycles at rest and maintains spontaneous neurotransmission. *Neuron* **73**, 121-34.

Rossano, A. J., Chouhan, A. K. and Macleod, G. T. (2013). Genetically encoded pH-indicators reveal activity-dependent cytosolic acidification of *Drosophila* motor nerve termini in vivo. *J Physiol* **591**, 1691-706.

Shuba, Y. M., Dietrich, C. J., Oermann, E., Cleemann, L. and Morad, M. (2008). Local extracellular acidification caused by Ca²⁺-dependent exocytosis in PC12 cells. *Cell Calcium* **44**, 220-9.

Siesjo, B. K., Katsura, K. and Kristian, T. (1996). Acidosis-related damage. *Adv Neurol* **71**, 209-33; discussion 234-6.

Sinning, A. and Hubner, C. A. (2013). Minireview: pH and synaptic transmission. *FEBS Lett* **587**, 1923-8.

Somjen, G. G. (1984). Acidification of interstitial fluid in hippocampal formation caused by seizures and by spreading depression. *Brain Res* **311**, 186-8.

Sudhof, T. C. (2013). Neurotransmitter release: the last millisecond in the life of a synaptic vesicle. *Neuron* **80**, 675-90.

Tang, C. M., Dichter, M. and Morad, M. (1990). Modulation of the N-methyl-D-aspartate channel by extracellular H⁺. *Proc Natl Acad Sci U S A* **87**, 6445-9.

Trapp, S., Luckermann, M., Kaila, K. and Ballanyi, K. (1996). Acidosis of hippocampal neurones mediated by a plasmalemmal Ca²⁺/H⁺ pump. *Neuroreport* **7**, 2000-4.

Traynelis, S. F. and Cull-Candy, S. G. (1990). Proton inhibition of N-methyl-D-aspartate receptors in cerebellar neurons. *Nature* **345**, 347-50.

Verstegen, A. M., Tagliatti, E., Lignani, G., Marte, A., Stoloro, T., Atias, M., Corradi, A., Valtorta, F., Gitler, D., Onofri, F. et al. (2014). Phosphorylation of synapsin I by cyclin-dependent kinase-5 sets the ratio between the resting and recycling pools of synaptic vesicles at hippocampal synapses. *J Neurosci* **34**, 7266-80.

Wang, T. M., Holzhausen, L. C. and Kramer, R. H. (2014). Imaging an optogenetic pH sensor reveals that protons mediate lateral inhibition in the retina. *Nat Neurosci* **17**, 262-8.

Wemmie, J. A., Taugher, R. J. and Kreple, C. J. (2013). Acid-sensing ion channels in pain and disease. *Nat Rev Neurosci* **14**, 461-71.

Xiong, Z. Q., Saggau, P. and Stringer, J. L. (2000). Activity-dependent intracellular acidification correlates with the duration of seizure activity. *J Neurosci* **20**, 1290-6.

Xiong, Z. Q. and Stringer, J. L. (2000). Extracellular pH responses in CA1 and the dentate gyrus during electrical stimulation, seizure discharges, and spreading depression. *J Neurophysiol* **83**, 3519-24.

Yin, T., Lindley, T. E., Albert, G. W., Ahmed, R., Schmeiser, P. B., Grady, M. S., Howard, M. A. and Welsh, M. J. (2013). Loss of Acid sensing ion channel-1a and bicarbonate administration attenuate the severity of traumatic brain injury. *PLoS One* **8**, e72379.

Zhang, Z., Nguyen, K. T., Barrett, E. F. and David, G. (2010). Vesicular ATPase inserted into the plasma membrane of motor terminals by exocytosis alkalizes cytosolic pH and facilitates endocytosis. *Neuron* **68**, 1097-108.

Zhao, H., Cai, Y., Yang, Z., He, D. and Shen, B. (2011). Acidosis leads to neurological disorders through overexciting cortical pyramidal neurons. *Biochem Biophys Res Commun* **415**, 224-8.

Zhao, H., Carney, K. E., Falgoust, L., Pan, J. W., Sun, D. and Zhang, Z. (2016). Emerging roles of Na⁽⁺⁾/H⁽⁺⁾ exchangers in epilepsy and developmental brain disorders. *Prog Neurobiol* **138-140**, 19-35.

Figures

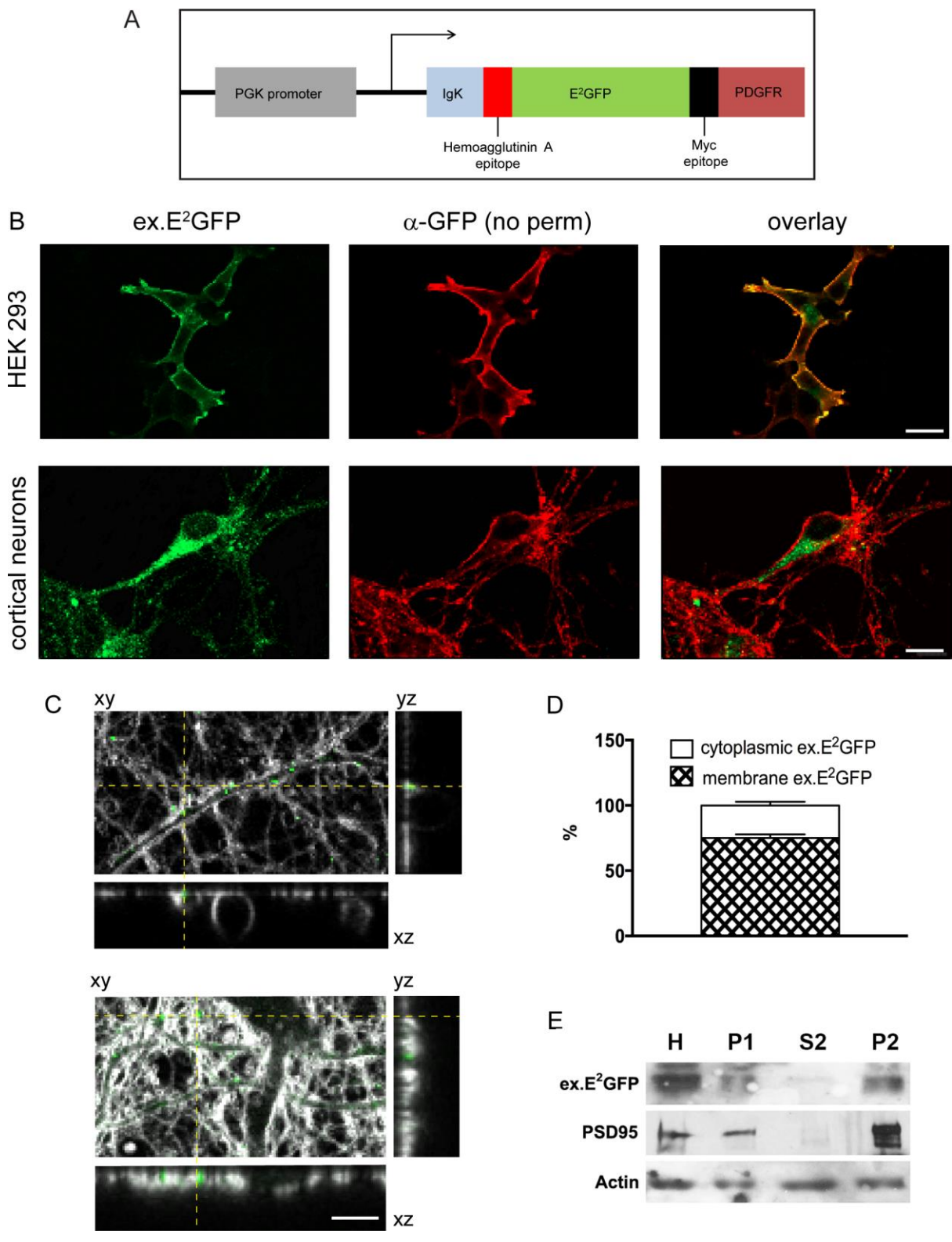


Figure 1. Engineering and cellular localization of ex.E²GFP.

A. Schematic illustration of the ex.E²GFP lentiviral construct, showing: the phosphoglycerate kinase (PGK) promoter, the murine I μ chain signal peptide, the *myc* epitope and the platelet-derived growth factor receptor (PDGFR) transmembrane domain. **B.** Plasma membrane expression of ex.E²GFP in HEK cells and primary cortical neurons. ex.E²GFP expression in green; anti-GFP in red, overlay in yellow. Scale bars, 10 μ m. **C, D.** Representative images of primary cortical neurons expressing ex.E²GFP (green) and stained with FM4-64fx (grey) for cell membrane visualization. Z-stack images were acquired to quantify the percentage of ex.E²GFP targeted to the plasma membrane (**D**). **E.** Subcellular fractionation of 17 DIV neurons infected with ex.E²GFP, lysed and analyzed by western blotting using anti-GFP antibodies. PSD95 was used as marker of membrane distribution, and actin to verify equal loading. Immunoreactive bands corresponding to ex.E²GFP are mostly present in the plasma membrane-containing section P2.

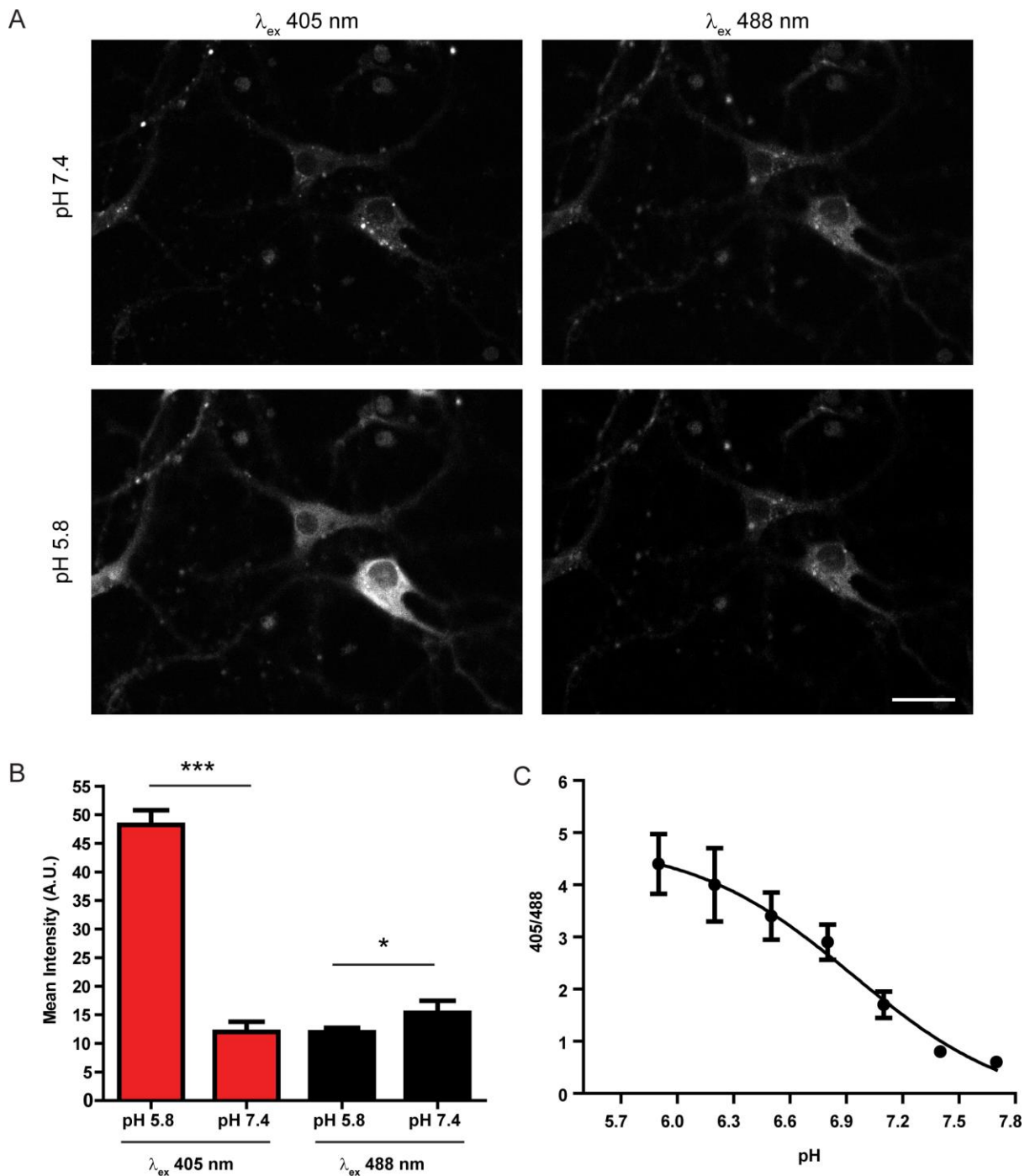


Figure 2. Characterization of ex.E²GFP.

A, B. Representative images and quantitative evaluation of ex.E²GFP fluorescence intensity in cortical neurons at various extracellular pH values. The fluorophore was excited at 405 and 488 nm (λ_{ex}) and the emitted photons were collected by a single window (510-560 nm). Scale bar, 20 μ m. The histograms show the mean \pm sem values of fluorescence intensity under the various λ_{ex} and pH conditions (* $p < 0.05$, *** $p < 0.001$, Student's t -test; $n = 15$ cells from 3 independent experiments). **C.** Calibration curve relating fluorescence intensity

ratios to extracellular pH. Data (means \pm sem of $n = 15$ cells from 3 independent experiments) were best fitted according to a sigmoidal dose-response equation: $y = -0,307 + (4,83 - 0,372) / (1 + 10^{6,9-x})$, $R^2 = 0.99$.

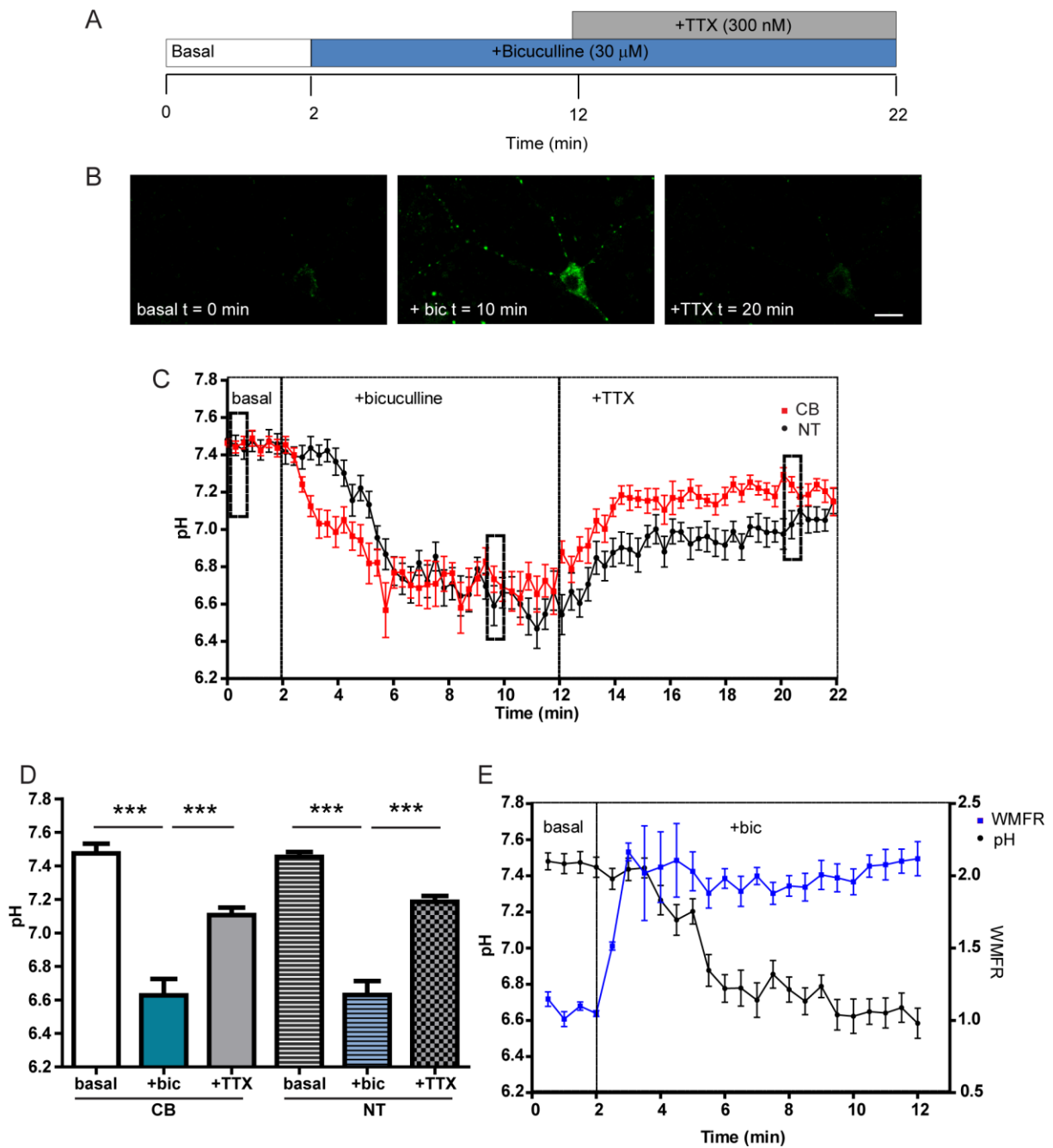


Figure 3. Neural hyperactivity causes localized acidic shifts in primary neurons.

A. Schematic representation of the time course of a typical experiment with the addition of bicuculline (30 μ M) and TTX (300 nM). **B.** Representative images showing the fluorescence intensity under basal conditions (t = 0 min), after bicuculline treatment (t = 10 min), and after TTX addition (t = 20 min). Scale bar, 20 μ m. **C.** pH fluctuations as a function of time in cell bodies (CB, red squares; n = 32) and nerve terminals (NT, black circles; n = 20). Data are means \pm sem from at least 3 independent experiments. Values are shown every 6 time-points for clarity. Rectangles indicate values binned for the statistical analysis. **D.** Quantification of

pH shifts observed under basal conditions (*basal*), after treatment with bicuculline (*+bic*) and after TTX addition (*+TTX*) in nerve terminals (*NT*) and cell bodies (*CB*). Mean \pm sem values were calculated by binning data of each independent trace as follows: “*basal*”, from 0 to 60 s; “*+bic*”, from 560 to 600 s; “*+TTX*”, from 1250 to 1318 s. *** $p < 0.001$, one-way ANOVA followed by Bonferroni’s multiple comparison test. **E.** Network firing rate (weighted mean firing rate, WMFR, Hz; blue squares) and pH changes (black circles) were recorded under basal conditions (2 min) and after administration of bicuculline (10 min). Data are calculated from 8 MEA dishes from 3 independent experiments.

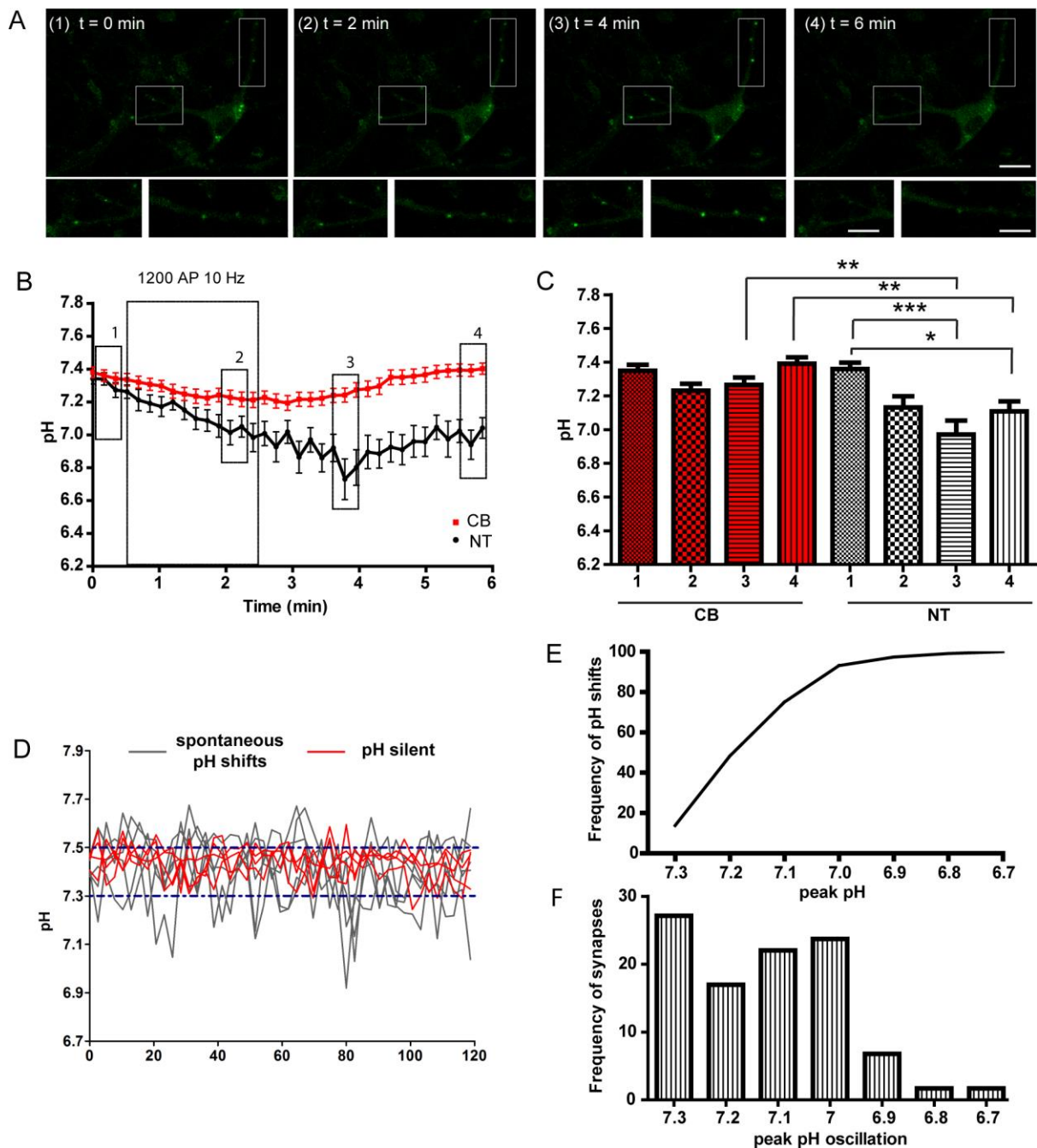


Figure 4. High frequency stimulation triggers a slowly developing acidosis.

A. Electrical field stimulation (1200 APs, 10 Hz) was applied to cortical neurons infected with ex.E²GFP. Representative fluorescence images taken at $\lambda_{\text{ex}} = 405$ nm from neurons under basal conditions ($t = 0$ min), during the stimulus ($t = 2$ min), after reaching the low pH plateau ($t = 4$ min) and during the subsequent recovery ($t = 6$ min). Scale bar, 20 μm and 10 μm for low and high (inset) magnification respectively. **B.** pH fluctuations as a function of time in nerve terminals (NT, black squares; $n = 28$) and cell bodies (CB, red circles; $n = 31$) kept under basal conditions for 30 s, stimulated for 120 s with 1200 APs at 10 Hz and left to

recover for 360 s. Data are means \pm sem from at least 3 independent experiments, with values shown every 3 time points for clarity. Rectangles indicate values binned for the statistical analysis. **C.** Quantification of the extracellular pH shifts (means \pm sem) in response to high frequency stimulation: data were calculated by binning data of each independent trace as follows: 1, 0-30 s; 2, 100-130 s; 3, 200-230 s; 4, 307-348 s. * $p < 0.05$, ** $p < 0.01$, *** $p < 0.001$, one-way ANOVA followed by Bonferroni's multiple comparison test. **D.** Representative traces of synaptic pH changes under basal conditions in 14-17 DIV cortical neurons. Two populations of synapses were identified, characterized by spontaneous acidic shifts ($\text{pH} < 7.3$; gray traces) or by substantially stable pH values ($7.3 < \text{pH} < 7.5$; red traces). **E.** Cumulative distribution of the amplitude of pH shifts. **F.** Frequency distribution of synapses as function of peak pH oscillations observed during spontaneous activity. Data are from $n = 60$ synapses from 3 independent experiments.

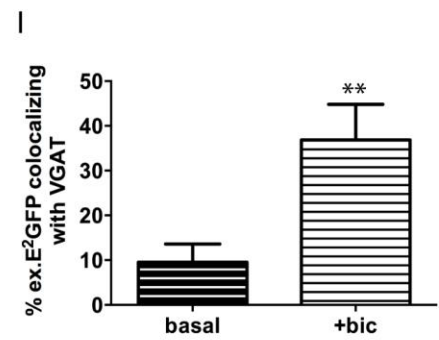
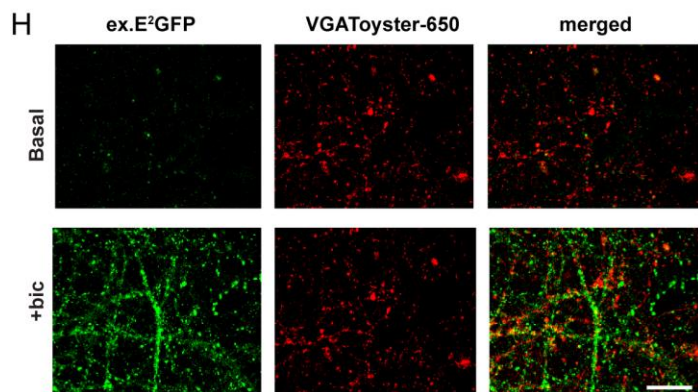
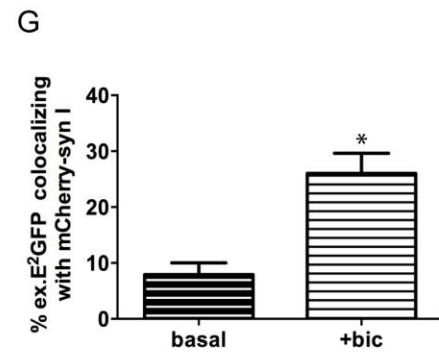
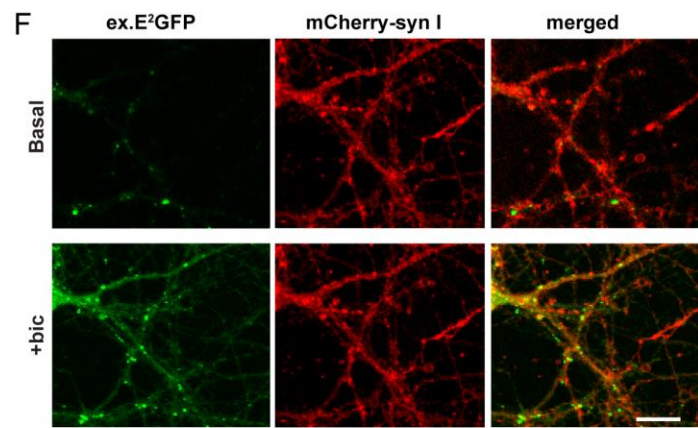
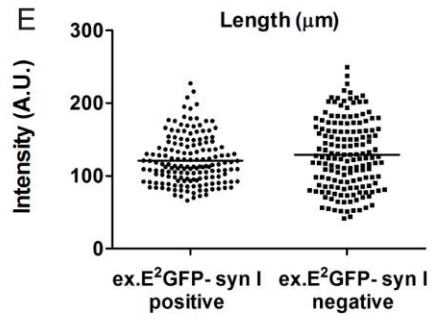
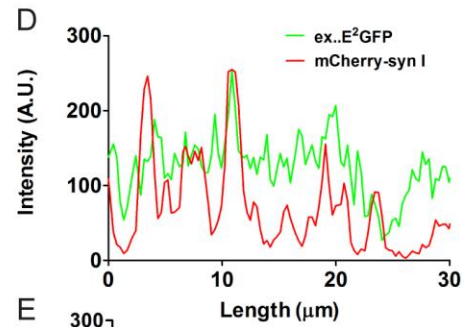
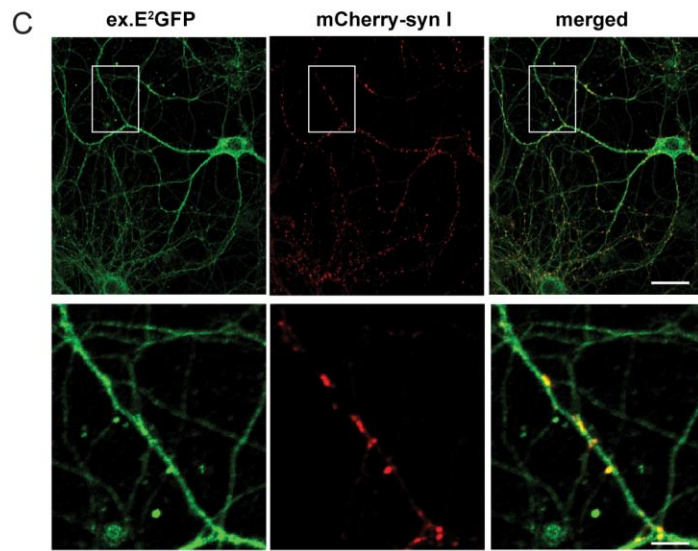
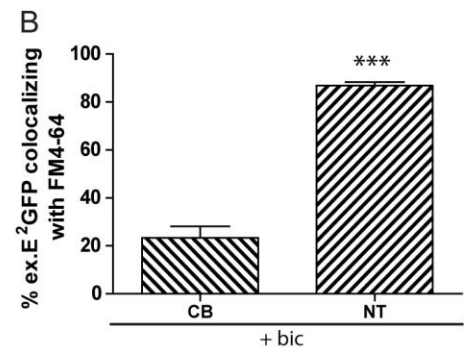
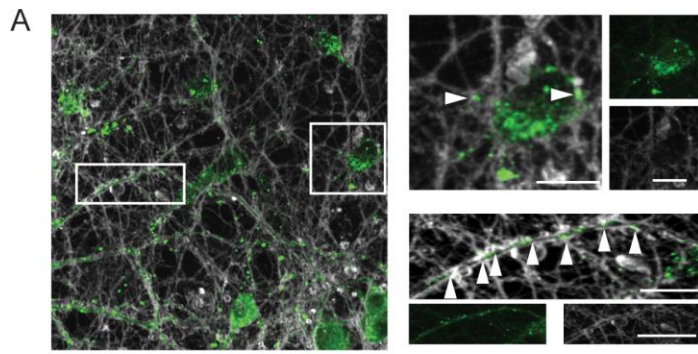


Figure 5. Acidic shifts are mainly localized to nerve terminals and hyperactivity-induced acidification occurs at active synapses.

A. Representative images taken at $\lambda_{\text{ex}} = 405$ nm show overlay of ex.E²GFP and FM4-64 at low (left panel) and high (right panel) magnification. Scale bars: 20 and 10 μm for low and high magnification respectively **B.** Quantification of ex.E²GFP positive puncta colocalization with the cell membrane. Bar graph shows cell bodies (CB) and nerve terminal (NT) colocalization of ex.E²GFP with FM4-64 fluorescence under bicuculline stimulation. Means \pm sem of $n = 5$ fields from 3 independent experiments. *** $p < 0.001$, unpaired Student's *t*-test. **C-E.** **C.** Primary cortical neurons co-infected with ex.E²GFP and mCherry-syn I expressing viruses were fixed and co-stained with anti-GFP and anti-mCherry antibodies to evaluate the synaptic *vs* extra-synaptic expression of ex.E²GFP. Scale bars, 25 and 5 μm for low (upper row) and high (lower row) magnification, respectively. **D.** Representative intensity profiles of ex.E²GFP and syn I signals along neurites. **E.** Quantification of fluorescence intensity of ex.E²GFP evaluated along neurites at synaptic (syn I-positive) and extra-synaptic (syn I-negative) regions. Data are calculated from $n = 160$ ROIs from 2 independent experiments. **F, G.** **F.** Representative images of ex.E²GFP ($\lambda_{\text{ex}} = 405$ nm) and mCherry fluorescence under basal conditions (upper panels) and upon bicuculline (+*bic*) treatment (10 min, 30 μM ; lower panels). The bicuculline-induced increase in ex.E²GFP fluorescence (green channel) is mostly localized to syn I-positive puncta (red channel). Scale bar, 10 μm . **G.** Quantification of the synaptic localization of ex.E²GFP fluorescence under basal conditions and after 10 min of bicuculline treatment. The fraction of the ex.E²GFP positive area that overlapped with mCherry-syn I fluorescence under basal and bicuculline conditions evaluated as Mander's coefficient is given as means \pm sem of $n = 5$ fields from 3 independent experiments. * $p < 0.05$, paired Student's *t*-test. **H, I.** **H.** Representative images of ex.E²GFP fluorescence and Oyster-VGAT650 live labeling in synaptic fields under basal conditions (upper panels) and upon bicuculline treatment (10 min, 30 μM ; lower panels). Bicuculline-induced increase of ex.E²GFP fluorescence (green channel) partially colocalizes with VGAT-positive puncta (red channel). Scale bar, 10 μm . **I.** Quantification of co-localization evaluated as Mander's coefficient between ex.E²GFP and Oyster-VGAT650 under basal conditions and after 10 min of bicuculline treatment ($n = 6$ fields, from 3 independent experiments) ** $p < 0.01$, paired Student's *t*-test.

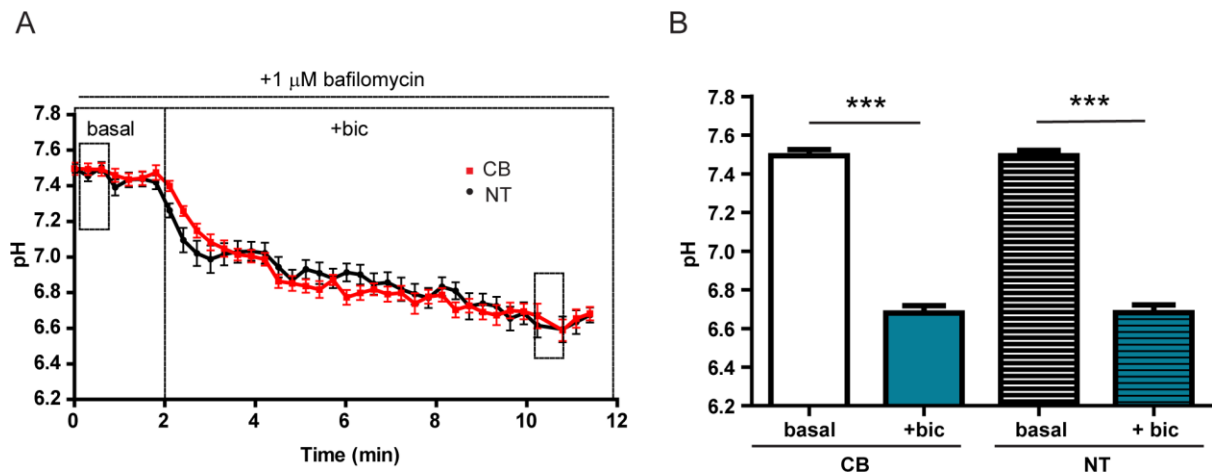


Figure 6. The hyperactivity-induced pH shifts are not due to internalization of ex.E²GFP.

A. Time-dependence of pH variations at synaptic puncta upon addition of bicuculline to neurons pretreated with 1 μ M bafilomycin. The dynamics of pH changes are superimposable to those obtained in the absence of bafilomycin (see Fig. 3). The graph displays average traces of ex.E²GFP fluorescence plotted as pH values for CB (n = 37) and NT (n = 39). Values (means \pm sem from at least 3 independent experiments) are shown every 6 time-points. Rectangles indicate values binned for the statistical analysis. **B.** Quantification of pH shifts in the presence of bafilomycin. Data (means \pm sem) were calculated by binning data of each independent trace as follows: “basal”, from 0 to 46 s; “+bic”, from 637 to 683 s; ***p < 0.001, one-way ANOVA followed by the Bonferroni’s multiple comparison test.

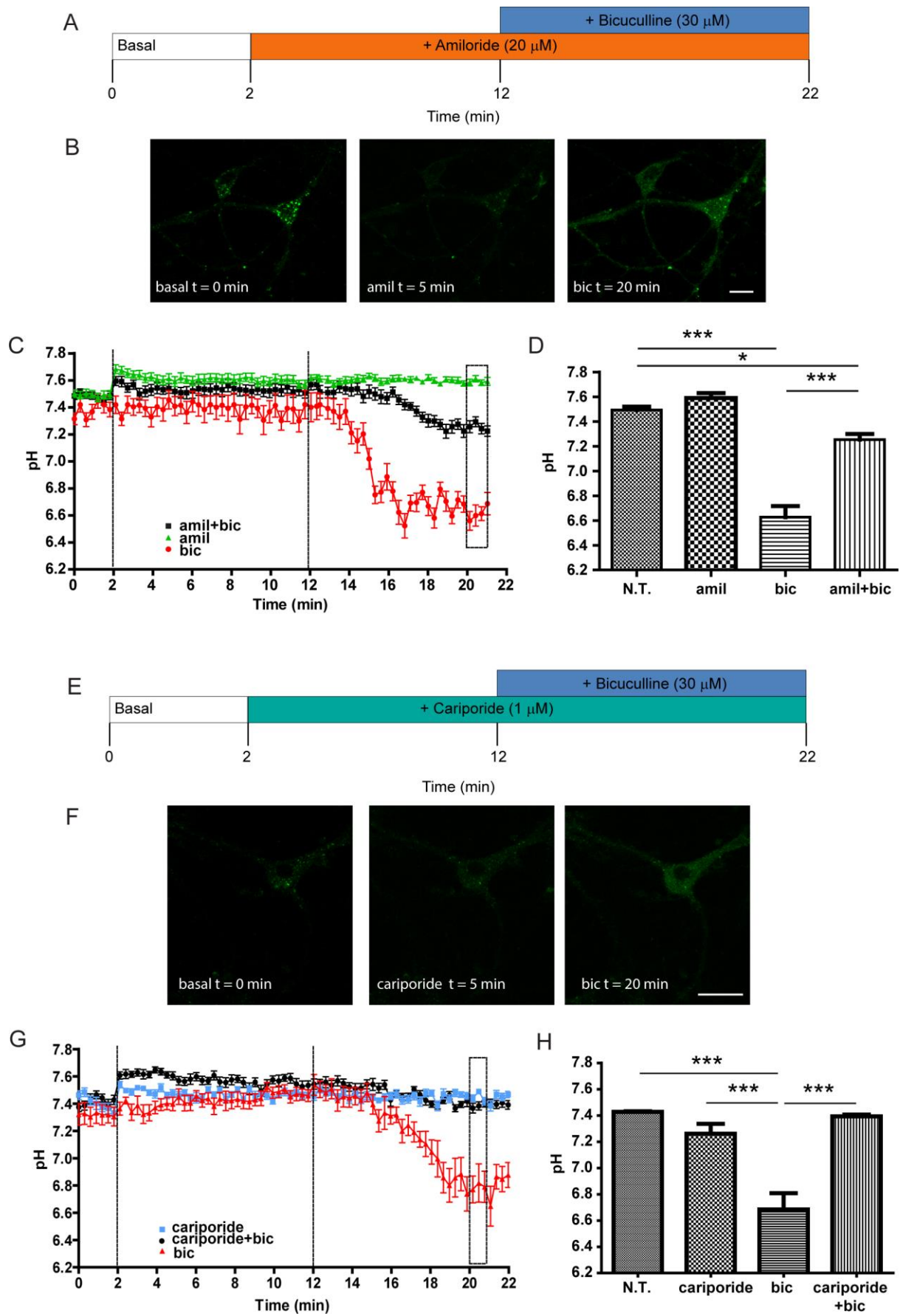


Figure 7. NHE channels modulate extracellular pH shifts.

A. Schematic representation of the time course of a typical experiment in which cells are treated with amiloride (20 μ M) for 10 min before addition of bicuculline (30 μ M). **B.** Representative images showing the fluorescence intensity under basal conditions ($t = 0$ min), after amiloride treatment ($t = 5$ min), and 10 min after bicuculline addition ($t = 20$ min). Scale bar, 20 μ m. **C.** Dynamics of pH fluctuations as a function of time in nerve terminals ($n = 20$) from neurons treated with: i) amiloride (amil, green triangles), ii) bicuculline (bic, red circles), and iii) amiloride + bicuculline (amil+bic, black squares). Non-treated (N.T.) samples were analyzed in parallel (trace not shown for clarity) and showed no appreciable pH changes over the course of the experiments. Data are means \pm sem from at least 3 independent experiments. Values are shown every 6 time-points. The rectangle indicates values binned for the statistical analysis. **D.** Quantification of pH shifts observed after drug administration at plateau values from 1266 to 1318 s and compared to untreated samples (N.T.). Data (means \pm sem) were calculated by binning data of each independent trace $***p < 0.001$, one-way ANOVA followed by Bonferroni's multiple comparison test. **E.** Schematic representation of the time course of a typical experiment in which cells are treated with cariporide (1 μ M) for 10 min before addition of bicuculline (30 μ M). **F.** Representative images showing the fluorescence intensity under basal conditions ($t = 0$ min), after cariporide treatment ($t = 5$ min), and 10 min after bicuculline addition ($t = 20$ min). Scale bar, 20 μ m. **G.** Dynamics of pH fluctuations in nerve terminals from neurons treated with: i) cariporide (blue squares, $n = 19$) bicuculline (bic, red triangles, $n = 15$), and iii) cariporide + bicuculline (black circles $n = 18$). Non-treated (N.T.) samples were analyzed in parallel (trace not shown for clarity) and showed no appreciable pH changes over the course of the experiments. Data are means \pm sem from at least 2 independent experiments. Values are shown every 6 time points. The rectangle indicates values binned for the statistical analysis. **H.** Quantification of pH shifts observed after drug administration at plateau values from 1200 to 1260 s and compared to untreated samples (N.T.). Data (means \pm sem) were calculated by binning data of each independent trace $***p < 0.001$, one-way ANOVA followed by Bonferroni's multiple comparison test.

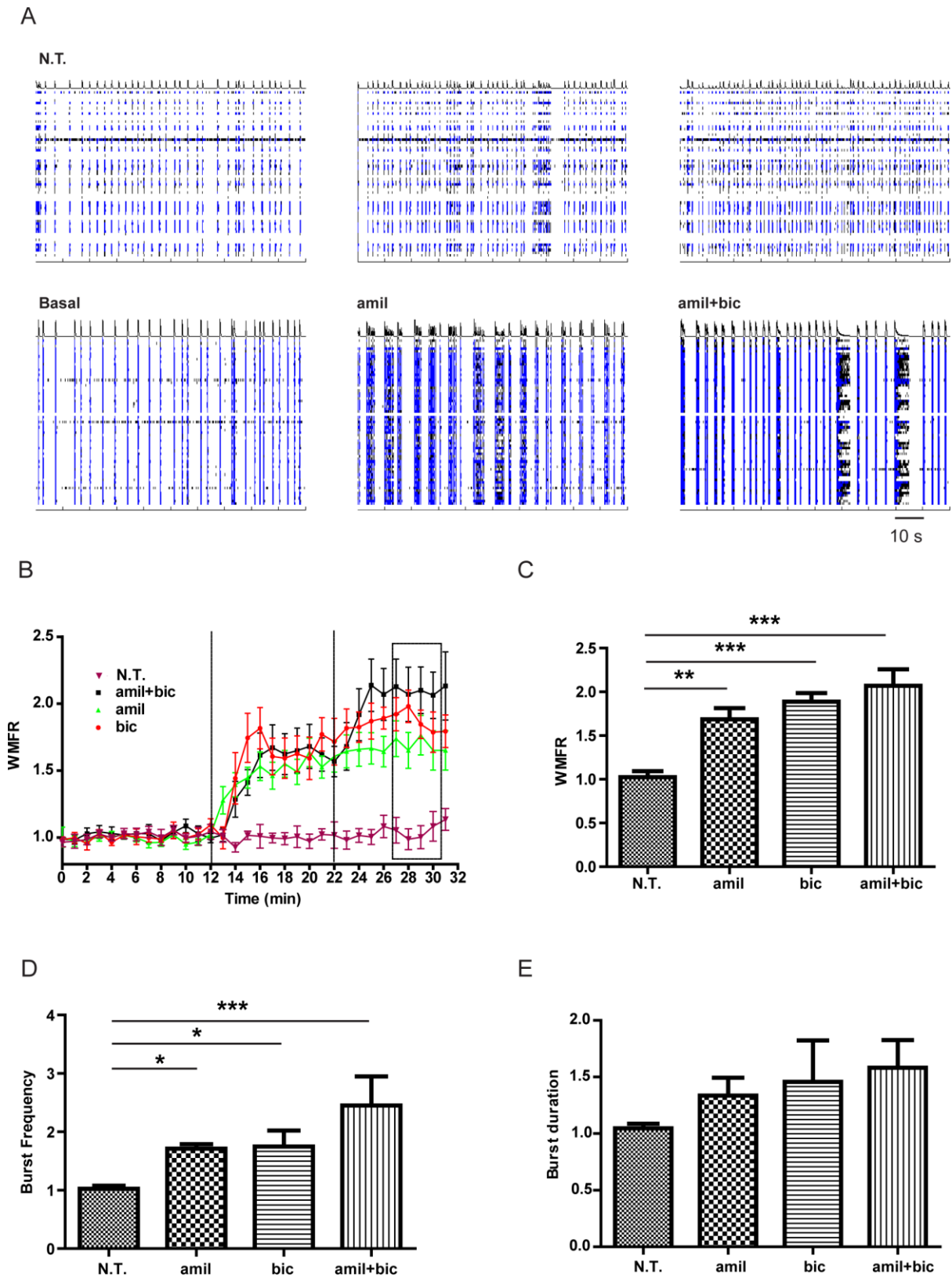


Figure 8. NHE inhibition does not affect firing rate and burst frequency.

A. Representative raster plots of network activity recorded from DIV21 cortical neurons treated with vehicle only (N.T.; upper panel) or with amiloride and amiloride+bicuculline

(lower panel). Each row along the vertical axis represents an individual microelectrode of the MEA device for 100-s time windows (horizontal axis). **B.** Dynamics of the weighted mean firing rate (WMFR) for neuronal cultures treated with: i) vehicle (N.T., pink triangles), ii) bicuculline for 20 min (bic, *red circles*), iii) amiloride for 20 min (amil, *green triangles*) and iv) amiloride for 10 min followed by bicuculline for 10 min (amil+bic *black squares*). The rectangle indicates values binned for the analysis shown in (C-E). **C.** Bar graphs for WMFR were calculated by binning data from 1560 to 1860 s for each experimental group. Data (means \pm sem) were calculated from 9 MEA dishes from 3 independent experiments. ** $p < 0.01$, *** $p < 0.001$, two-way ANOVA followed by Bonferroni's multiple comparison test. **D, E.** Bar graphs representing the changes in bursting rate and burst duration, respectively, in the same experimental groups. Data (means \pm sem) were calculated from 8 MEA dishes from 3 independent experiments. * $p < 0.05$, *** $p < 0.001$, two-way ANOVA followed by Bonferroni's multiple comparison test.

SUPPLEMENTARY MATERIALS

Fig. S1 Chiacchiaretta et al

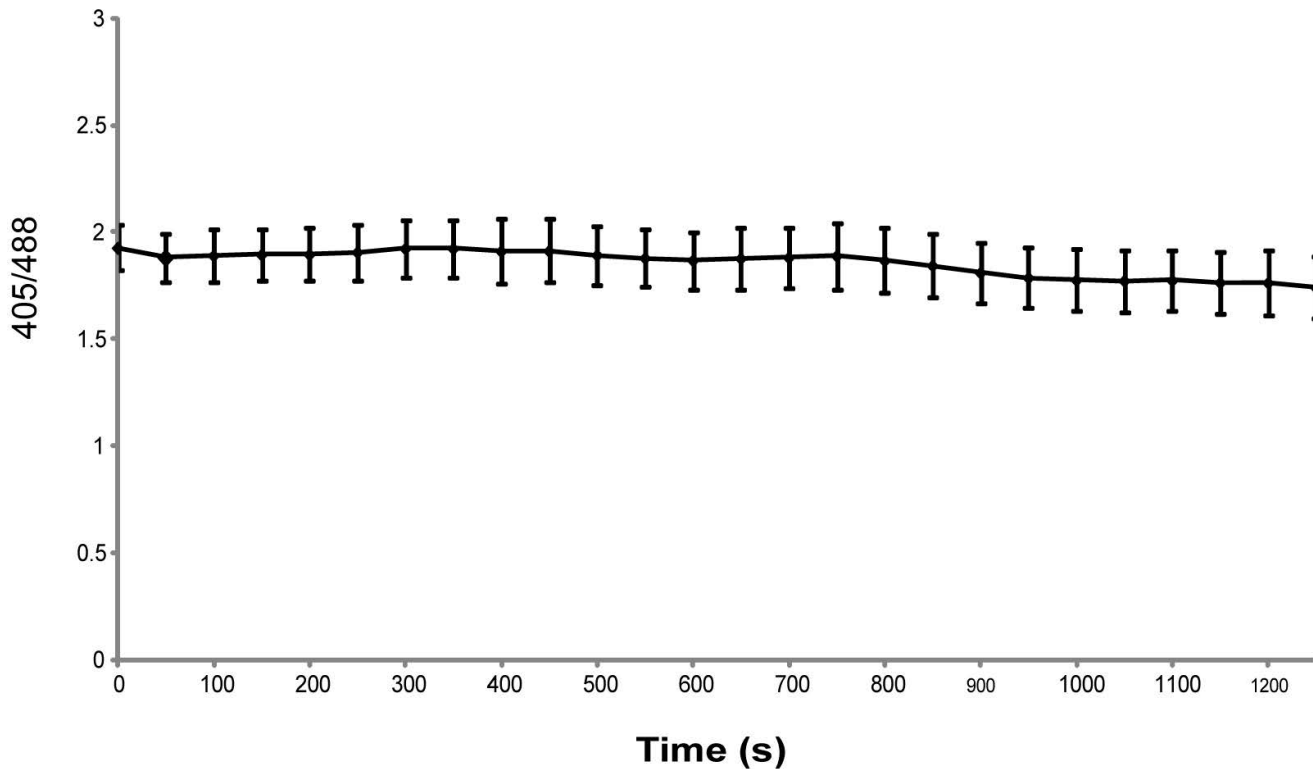


Figure S1. ex.E²GFP fluorescence is not sensitive to photobleaching.

To quantify the extent of ex.E²GFP photobleaching, changes in fluorescence intensity were followed for a period of time corresponding to our time-lapse experiments under constant pH (pH 6). Emitted photons were monitored by live imaging throughout the experiments at a frame rate of 0.1 Hz. For collecting excitation and emission, the same optical parameters were used as described for the calibration curve, and fluorescence ratios were calculated using the equation described in Materials and Methods. Values are means \pm sem. No significant photobleaching was observed during the time course of the experiment.

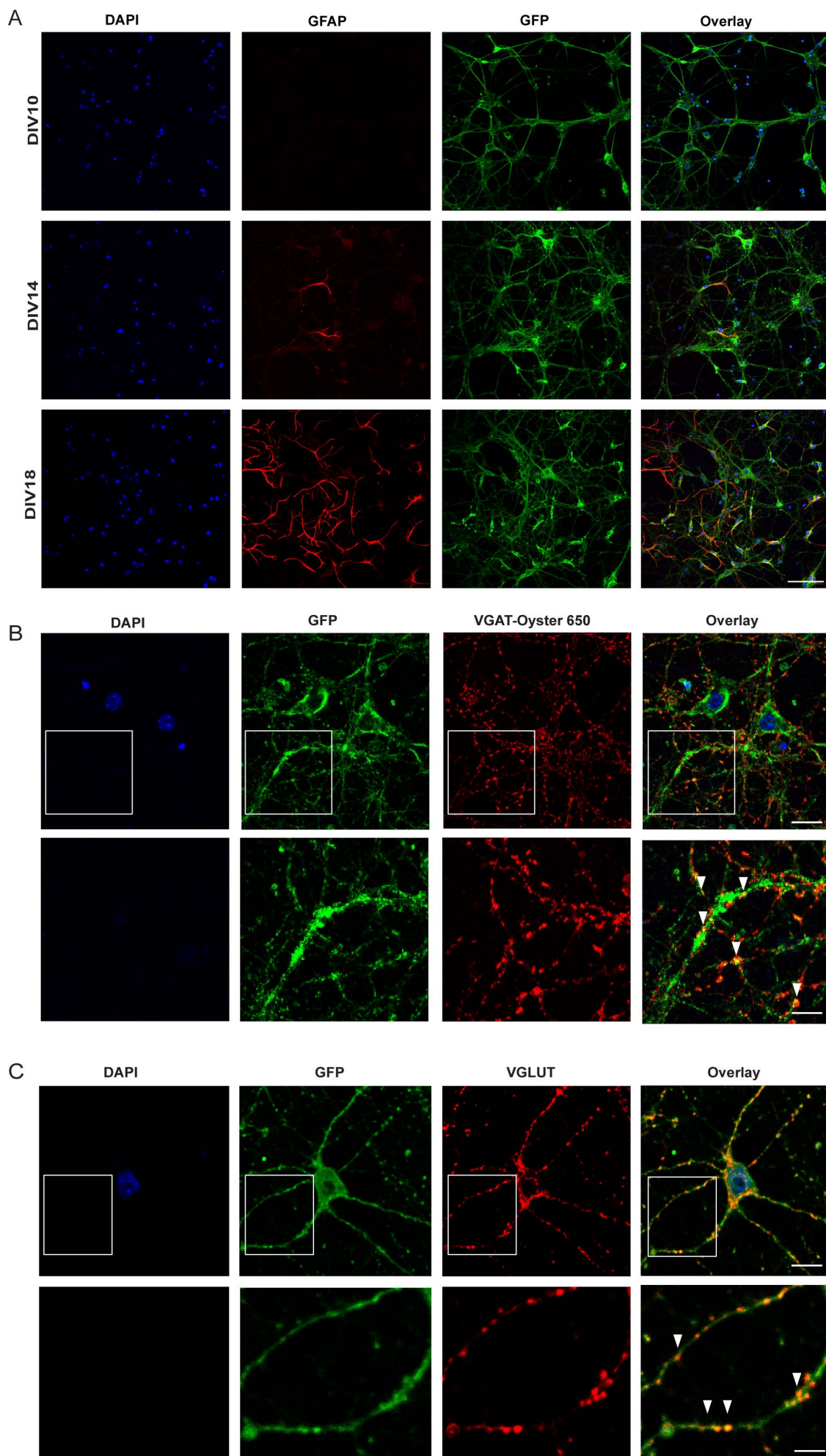


Figure S2. ex.E²GFP expression pattern in vitro. **A.** ex.E²GFP-infected neurons were fixed and stained with antibodies to GFP (green), glial fibrillary acidic protein (GFAP, red) and DAPI for nuclei. Between DIV 10-14, very few GFAP-positive cells were present. More glial cells were present at DIV 18; however also at this stage no significant expression of ex.E²GFP in GFAP-positive cells was detected. Scale bars, 20 μ m. **B.** Representative images of primary cortical neurons infected with ex.E²GFP and stained for GFP (green), VGAT-oyster 650 (red) and DAPI (blue), to monitor expression of the probe in inhibitory neurons/terminals. Scale bars, 20 μ m and 10 μ m for low (upper row) and high (lower row) magnification, respectively. **C.** Representative images of primary cortical neurons infected with ex.E²GFP and stained for GFP (green), VGLUT (red) and DAPI (blue), to monitor expression of the probe in excitatory neurons/terminals. Scale bars, 20 μ m and 5 μ m for low (upper row) and high (lower row) magnification, respectively.

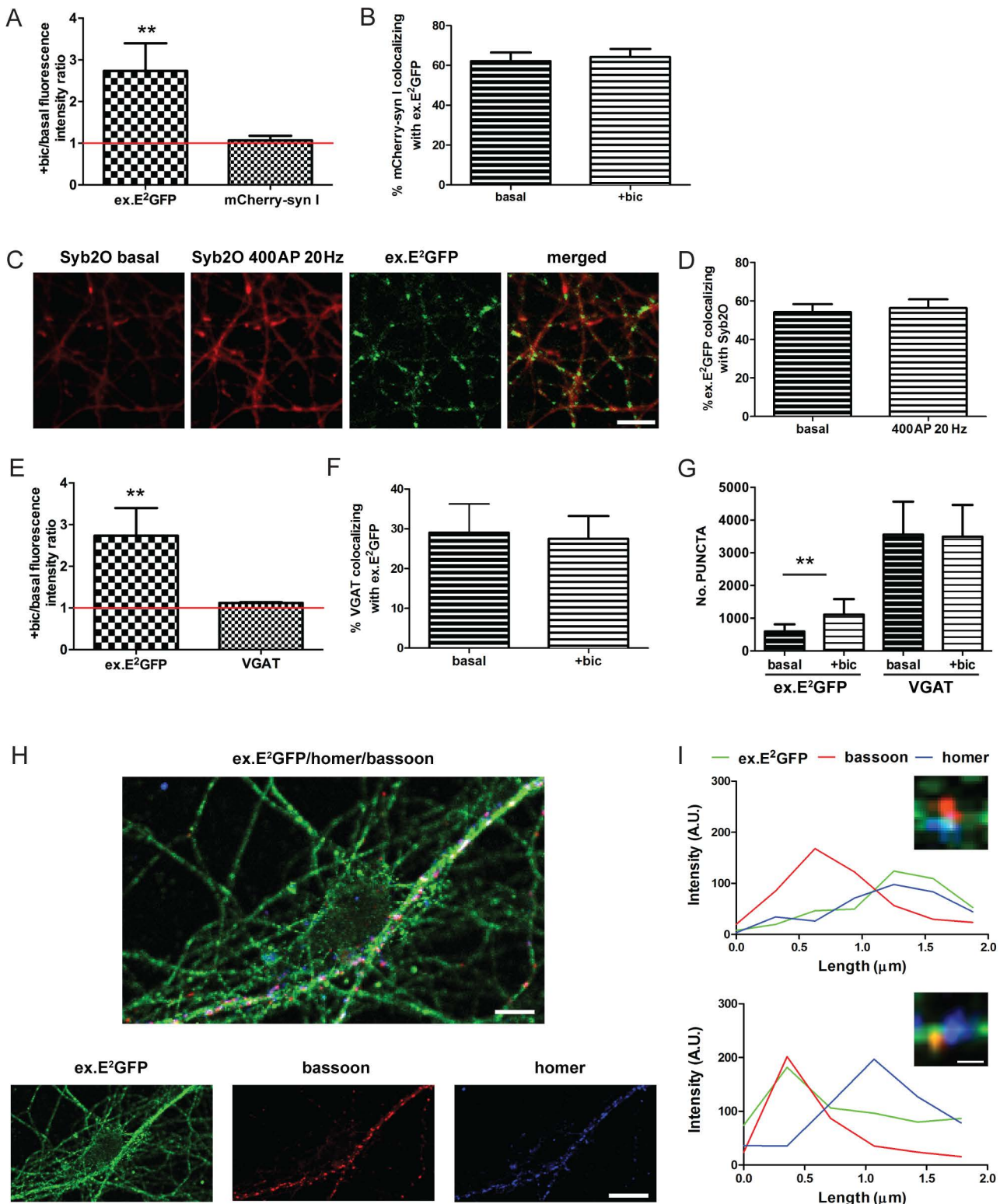
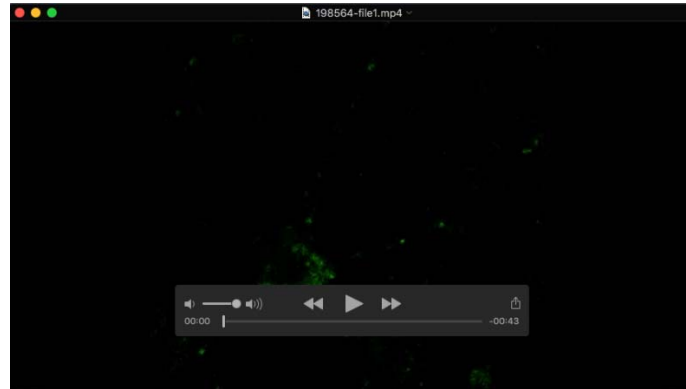


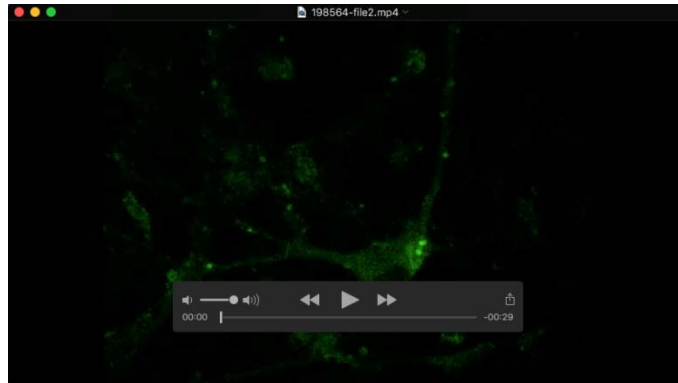
Figure S3. ex.E²GFP-positive puncta increase in number and intensity after bicuculline stimulation.

A. Primary cortical neurons were co-infected at DIV7 with ex.E²GFP and mCherry-syn I expressing viruses. Quantitative evaluation of ex.E²GFP and mCherry syn I fluorescence intensity after bicuculline treatment (+*bic*) was measured and normalized to the values of untreated samples, set to 1 (red line); n = 5 fields from 3 independent experiments. **B.** Quantification of the fraction of the mCherry-syn I area that overlapped with ex.E²GFP positive fluorescence under basal and bicuculline conditions evaluated as Mander's coefficient is given as means ± sem (n = 5 fields from 3 independent experiments). No significant differences were observed. **C,D. C.** Primary cortical neurons were co-infected at DIV13 with ex.E²GFP and synaptobrevin-pHluorin m-Orange2 (Syb2O) expressing viruses. Optical recordings were performed at DIV 17-18. Representative images of ex.E²GFP and Syb2O fluorescence under basal conditions and after electrical stimulation (1200 AP, 20 Hz). Scale bar, 5 μm. **D.** Quantification of ex.E²GFP/Syb2O colocalization evaluated as Manders's coefficient (n = 11 fields from 3 independent experiments). No significant differences were observed. **E.** ex.E²GFP infected neurons were stained with VGAT-oyster650 and quantitative evaluation of ex.E²GFP and VGAT-oyster650 fluorescence intensity after bicuculline treatment (+*bic*) was measured and normalized to the values of untreated samples, set to 1 (red line); n = 4 fields from 3 independent experiments. **F.** Quantification of the fraction of Oyster-VGAT650 puncta that overlapped with ex.E²GFP fluorescence evaluated as Mander's coefficient under basal and bicuculline conditions, respectively (means ± sem of n = 6 fields from 3 independent experiments). No significant differences were observed. **G.** Density of ex.E²GFP and VGAT-oyster650 positive puncta under basal (*basal*) conditions and after bicuculline treatment (+*bic*); n = 5 fields from 3 independent experiments. Data are means ± sem, **p < 0.01, one-way ANOVA and Bonferroni's post-hoc test. **H,I. H.** Representative images of primary neurons infected with ex.E²GFP and stained for GFP (green), bassoon (red) and homer (blue), to monitor expression of the probe at pre- or post- synaptic membranes. Scale bar, 10 μm. **I.** Intensity profiles of ex.E²GFP at single synaptic puncta in which pre- and post-synaptic signals did not completely overlap. Representative images of synaptic puncta analyzed are shown (top right). Scale bar, 1 μm.



Movie S1: pH fluctuations upon network hyperactivity

Neurons were imaged by time-lapse fluorescence microscopy at $\lambda_{\text{ex}} = 405 \text{ nm}$ (1 image/2.58 s, total time: 1320 s, movie played at 29 frames/s) under basal conditions and after administration of bicuculline (30 μM) and TTX (300 nM).



Movie S2: pH fluctuations upon high frequency stimulation

Neurons were imaged by time-lapse fluorescence microscopy, at $\lambda_{\text{ex}} = 405 \text{ nm}$ (1 image/2.58 s, total time: 350 s, movie played at 29 frames/s) under high frequency electrical stimulation (1200 AP, 10 Hz).