



Activated Bone Marrow-Derived Macrophages Eradicate Alzheimer's-Related A β ₄₂ Oligomers and Protect Synapses

Songlin Li^{1,2,3}, Eric Y. Hayden⁴, Veronica J. Garcia^{5,6}, Dieu-Trang Fuchs³, Julia Sheyn³, David A. Daley³, Altan Rentsendorj³, Tania Torbati⁷, Keith L. Black³, Ueli Rutishauser^{3,6,8}, David B. Teplow⁴, Yosef Koronyo³ and Maya Koronyo-Hamaoui^{3,6*}

¹ Institute of Neuroscience and Chemistry, Wenzhou University, Wenzhou, China, ² Institute of Life Sciences, Wenzhou University, Wenzhou, China, ³ Department of Neurosurgery, Cedars-Sinai Medical Center, Maxine-Dunitz Neurosurgical Institute, Los Angeles, CA, United States, ⁴ Department of Neurology, David Geffen School of Medicine at UCLA, Mary S. Easton Center for Alzheimer's Disease Research at UCLA, Brain Research Institute, Molecular Biology Institute, University of California, Los Angeles, Los Angeles, CA, United States, ⁵ Board of Governors Regenerative Medicine Institute, Cedars-Sinai Medical Center, Los Angeles, CA, United States, ⁶ Department of Biomedical Sciences, Cedars-Sinai Medical Center, Los Angeles, CA, United States, ⁷ College of Osteopathic Medicine of the Pacific, Western University of Health Sciences, Pomona, CA, United States, ⁸ Department of Neurology, Cedars-Sinai Medical Center, Los Angeles, CA, United States

OPEN ACCESS

Edited by:

Tanuja Chitnis,
Brigham and Women's Hospital and
Harvard Medical School,
United States

Reviewed by:

Enrico Castroflorio,
Medical Research Council Harwell
(MRC), United Kingdom
Robert Adam Harris,
Karolinska Institutet (KI), Sweden

*Correspondence:

Maya Koronyo-Hamaoui
maya.koronyo@cshs.org

Specialty section:

This article was submitted to
Multiple Sclerosis and
Neuroimmunology,
a section of the journal
Frontiers in Immunology

Received: 15 October 2019

Accepted: 09 January 2020

Published: 31 January 2020

Citation:

Li S, Hayden EY, Garcia VJ, Fuchs D-T, Sheyn J, Daley DA, Rentsendorj A, Torbati T, Black KL, Rutishauser U, Teplow DB, Koronyo Y and Koronyo-Hamaoui M (2020) Activated Bone Marrow-Derived Macrophages Eradicate Alzheimer's-Related A β ₄₂ Oligomers and Protect Synapses. *Front. Immunol.* 11:49. doi: 10.3389/fimmu.2020.00049

Impaired synaptic integrity and function due to accumulation of amyloid β -protein (A β ₄₂) oligomers is thought to be a major contributor to cognitive decline in Alzheimer's disease (AD). However, the exact role of A β ₄₂ oligomers in synaptotoxicity and the ability of peripheral innate immune cells to rescue synapses remain poorly understood due to the metastable nature of oligomers. Here, we utilized photo-induced cross-linking to stabilize pure oligomers and study their effects vs. fibrils on synapses and protection by A β -phagocytic macrophages. We found that cortical neurons were more susceptible to A β ₄₂ oligomers than fibrils, triggering additional neuritic arborization retraction, functional alterations (hyperactivity and spike waveform), and loss of VGluT1- and PSD95-excitatory synapses. Co-culturing neurons with bone marrow-derived macrophages protected synapses against A β ₄₂ fibrils; moreover, immune activation with glatiramer acetate (GA) conferred further protection against oligomers. Mechanisms involved increased A β ₄₂ removal by macrophages, amplified by GA stimulation: fibrils were largely cleared through intracellular CD36/EEA1⁺-early endosomal proteolysis, while oligomers were primarily removed via extracellular/MMP-9 enzymatic degradation. *In vivo* studies in GA-immunized or CD115⁺-monocyte-grafted APP_{SWE}/PS1 Δ E9-transgenic mice followed by pre- and postsynaptic analyses of entorhinal cortex and hippocampal substructures corroborated our *in vitro* findings of macrophage-mediated synaptic preservation. Together, our data demonstrate that activated macrophages effectively clear A β ₄₂ oligomers and rescue VGluT1/PSD95 synapses, providing rationale for harnessing macrophages to treat AD.

Keywords: Alzheimer's disease, neurodegeneration, immunomodulation therapy, amyloid-beta, regeneration, synaptogenesis

INTRODUCTION

Alzheimer's disease (AD) is a progressive, incurable and fatal neurodegenerative disorder and the most frequent cause of senile dementia (1, 2). AD is characterized by severe synaptic and neuronal loss resulting in gradual behavioral and cognitive decline (3, 4). Early pathological changes in AD involve accumulation of amyloid β -protein ($A\beta$) in the brain and retina (5–11), with the amyloidogenic 42-residue long alloform ($A\beta_{42}$) viewed as the most pathognomonic (5, 12). These $A\beta_{42}$ peptides have been shown to rapidly aggregate into insoluble fibrils that form the plaques in AD brains, but also to assemble into non-fibrillar soluble oligomers, believed to be highly neurotoxic (8, 13–20). While there is a weak correlation between cerebral plaque density and severity of AD, $A\beta_{42}$ oligomers appear to closely associate with synaptopathy and cognitive impairment (21, 22). Specifically, growing evidence demonstrates the detrimental effects of $A\beta_{42}$ oligomers on synaptic plasticity, spine morphology and density (13, 23), axonal transport (24–26), and cognition (8, 15, 17, 22, 27–29). These findings incentivize strategies that eradicate $A\beta_{42}$ oligomers in order to preserve synapses and cognitive function (30–34).

Multiple mechanisms mediate cerebral $A\beta$ clearance, including phagocytosis and enzymatic degradation by innate immune cells such as microglia and bone marrow (BM)-derived monocytes (Mo) and macrophages ($M\Phi$) (35, 36). However, under a chronic neuroinflammatory condition typical for AD, microglia exhibit a compromised phenotype, with diminished ability to clear $A\beta$ (37, 38), further contributing to damaging inflammatory processes, and excessively pruning synapses (39–46). Nevertheless, unlike brain-resident microglia, BM-derived Mo/ $M\Phi$ appear to have an increased capacity to remove $A\beta$ fibrils, as shown by studies in both cell cultures and animal models (47–64). Indeed, various pre-clinical studies from our group and others have supported an emerging paradigm shift for the therapeutic potential of immunomodulation via enhanced cerebral recruitment of BM-derived $A\beta$ -clearing macrophages (9, 35, 50, 52, 58, 62–70). In particular, we demonstrated that successful delivery of therapeutic Mo/ $M\Phi$ to the brains of APP_{SWE}/PS1_{DE9} double-transgenic (ADtg) mice can be achieved via immunization with altered myelin-derived antigens (e.g., glatiramer acetate – GA; also known as Copaxone[®]) or peripheral-blood enrichment with BM-derived CD115⁺-monocytes (Mo^{BM}) (50, 52, 58, 63, 70). In the immunized mice, cerebral $M\Phi$ were found to be highly-phagocytic, pro-healing, and anti-inflammatory, profoundly alleviating cerebral $A\beta_{40}$ and $A\beta_{42}$ burden, reducing microgliosis and astrocytosis, and ultimately improving cognitive functions. Further, GA induced macrophage-mediated recognition and phagocytosis of fibrillar $A\beta_{42}$ (63, 70). However, the potential of clearing the non-fibrillar, soluble $A\beta_{42}$ oligomeric forms by macrophages requires further investigation.

Experimental and clinical studies have indicated that the cognitive decline seen in AD patients and animal models is related to early pathological processes occurring in entorhinal cortex layers 2 and 3 (71–75). These layers play a crucial role in connecting brain cortical regions to the hippocampus.

Our immunomodulation approach of recruiting Mo/ $M\Phi$ to brain sites of $A\beta$ deposits substantially reduced AD-related pathology within these regions (63, 70). Importantly, selective ablation of peripheral monocytes, either by diphtheria toxin-targeted depletion or Ccr2-mediated inhibition of cerebral monocyte infiltration, exacerbated cortical and hippocampal $A\beta$ pathology in ADtg mice (52, 53). In addition, our preliminary data suggested a presynaptic preservation concomitant with $A\beta$ -plaque reduction in hippocampi of these murine models (63). Yet, the underinvestigate role of peripheral Mo/ $M\Phi$ in preservation of synaptic and neurite integrity, especially postsynaptic terminals within AD-relevant cortical regions, calls for broader exploration due to promising clinical implications of immunomodulation intervention strategies for AD.

Here, we applied photo-induced cross-linking of unmodified proteins (PICUP) to assess the impact of intrinsically metastable and polydisperse $A\beta_{42}$ oligomers on neurons and to test the ability of macrophages to recognize, bind to, and remove them (76, 77). This technique allowed us to produce pure, stabilized assemblies of low-n $A\beta_{42}$ oligomers (XL-o $A\beta_{42}$). The precise detrimental impact of these oligomeric conformers, as compared with preformed fibrils (f $A\beta_{42}$), was determined on synaptic density, neuritic arborization length, and neuronal function in primary (P1) cortical neurons. Furthermore, synaptic and neurite protection against $A\beta_{42}$ oligomers was explored in co-cultures of cortical neurons accompanied with BM-derived $M\Phi$ or GA-activated $M\Phi$. Next, the innate immune mechanisms of eliminating pathogenic $A\beta_{42}$ assemblies and subsequently attaining synaptic rescue were studied. Finally, in ADtg mice, we determined synaptic loss as well as preservation of pre- and postsynaptic terminals following macrophage-mediated immunomodulation interventions in predefined hippocampal and cortical subregions.

MATERIALS AND METHODS

Mice

Alzheimer's disease double-transgenic (ADtg) B6.Cg-Tg (APP_{swe}, PSEN1 $_{\Delta E9}$) 85Dbo/J mice and their age-matched wildtype (WT) C57BL/6J littermates were purchased from Jackson Laboratories (MMRRC stock 34832-JAX) and then bred and maintained at Cedars-Sinai Medical Center. All animals in this study have a C57BL/6 congenic background. Two cohorts of 10-month-old males, WT and ADtg (n = 6 per experimental group), were used. The first cohort underwent weekly s.c. GA immunization or monthly i.v. CD115⁺-Mo^{BM} injections for a duration of 2 months. Control groups were comprised of either naïve WT mice or monthly i.v. PBS injected ADtg mice. At the completion of the experiment, mice underwent behavioral testing at 12 months of age, then were perfused, under deep anesthesia, with ice-cold 0.9% saline solution containing 0.5 mM EDTA, and brains were collected for analyses. Harvested brains were cut in half. One hemisphere was post-fixed overnight in phosphate-buffered saline (PBS) containing 2.5% paraformaldehyde. Tissue was cryoprotected and stored in PBS containing 30% sucrose and 0.1% sodium azide. Brains were cut coronally in 30 μ m serial sections for histology. The second hemisphere was further

snap-frozen for biochemical and molecular assays as described previously (63). The second cohort underwent perfusion as described above for tissue collection.

Culture and Co-culture of Primary Cortical Neurons and Macrophages

Cortical neuronal cultures were prepared from postnatal day 1 C57BL/6 mice (78, 79). Neonates were decapitated, and their brains were removed. The cerebral cortex was dissected, stripped of meninges, dissociated with a combination of calcium and magnesium free Hank's balanced salt solution (HBSS; Life Technologies) containing 0.2% w/v papain suspension (Worthington Biochemical), and digested for 12 min at 37°C. The triturated cells were passed through a 70 µm strainer and counted. The cells were plated in laminin and poly-D-lysine-coated coverslips (BD Biosciences) at a density of 8×10^4 cells per ml (in 24-well plates) in NbActiv4 (BrainBits), then supplemented with 100 units/ml penicillin and 100 µg/ml streptomycin. The purity of primary neurons was about 92–94% (78). Bone marrow-derived macrophages were obtained as described previously (63). Briefly, bone marrow cells were isolated from the femurs and tibiae of 8- to 16-week-old C57BL/6 mice, differentiated into macrophages by incubation in complete RPMI-1640 (Life Technologies), and supplemented with 10% fetal bovine serum (Life Technologies) and 20 ng/ml macrophage colony stimulating factor (M-CSF; PeproTech) for 5 days. Primary cultures of macrophages were then suspended in NbActiv4 and co-cultured with primary cortical neurons (ratio of 1:1) per well in 24-well tissue-culture plates for 2 days. Primary neurons and/or macrophages were incubated overnight with 100nM of scrambled (s), fibrillar (f), or oligomeric (o) Aβ₄₂ (see below "Preparation of Aβ_{1–42} oligomers, fibrils and Scrambled Aβ₄₂"). Additional groups included those with macrophages treated with glatiramer acetate (GA; 30 µg/ml, TEVA Neuroscience) for 24 h. Experiments were carried out at day 9 of seeding cortical neurons starting with synapse formation and maturation (78). Cultures were washed in PBS and then fixed in 4% formaldehyde for 20 min at room temperature for further experimental procedures.

Microelectrode Array Recording

Cortical neuronal cultures (as described above) were plated in a M768-GLx 12-, 24-, or 48-well plate (Axion Biosystems). Preformed Aβ₄₂ (100 nM) were added to the culture on day 9. Simultaneous recordings from 64 extracellular electrodes per well were made using the Maestro (Axion BioSystems) microelectrode array (MEA) system at a constant temperature of 37°C. Data were sampled at 12.5 kHz, digitized and analyzed using Axion Integrated Studio software (Axion BioSystems) where spike events were detected using an adaptive spike detection threshold of 5.5 SD for each electrode with 1 s binning. Detected waveforms were then further sorted using Offline Sorter v4 software (Plexon) and exported through NeuroExplorer (Nex Technologies) into MatLab for waveform analysis using custom code. All data is presented as means ± s.e.m. and include recordings from 5 independent experiments.

Amyloid-β₄₂ Determination by Sandwich ELISA

At day 7, macrophages were incubated for 30 min or 24 h with fibrillar (f) or oligomeric (o) Aβ₄₂ and thereafter culture medium was collected. Total protein levels in the supernatant were determined using the Pierce BCA Protein Assay Kit (#23227; Thermo Scientific). The Aβ₄₂ levels in the supernatant were assessed using a sandwich ELISA for an anti-human Aβ₄₂ end-specific kit (KHB3442, Invitrogen; which does not recognize mouse Aβ nor human Aβ₄₀/Aβ₄₃). The kit was used according to the manufacturer's instructions. This assay is based on a combination of two antibodies specific for the N- and the COOH-termini of Aβ₄₂ sequences. The bound rabbit anti-COOH-terminus was detected with a horseradish peroxidase-labeled anti-rabbit antibody and was read at 450 nm using a microplate reader (Spectra Max 384 plus, Molecular Devices). Concentrations detected at 24 h incubation were normalized to basal levels at 30 min incubation.

Immunocytochemistry and Quantification of Synapses and Neurites

Primary neurons/macrophages in coverslips and brain coronal cryosections at bregma –2.50, –2.65, and –2.80 mm (80) per animal were washed in PBS and then treated with a permeabilization/blocking solution containing 20% normal horse serum (Invitrogen) and 0.05% Triton X-100 (Sigma-Aldrich). Cells or tissue sections were stained overnight at 4°C with combinations of presynaptic marker VGluT1 (Chemicon, 1:6,000), postsynaptic marker PSD95 (Abcam, 1:600; for synapses), or neuron-specific class III beta-tubulin/Tuj1 (Abcam) or Tubb3 (BioLegend; for neurites; **Supplementary Table 1**) in 2% blocking solution in PBS. Secondary antibodies of Cy3 donkey anti-guinea pig or anti-mouse or Cy5 donkey anti-rabbit (**Supplementary Table 1**) were incubated for 1 h at room temperature. The samples were washed in PBS and mounted using ProLong® Gold with DAPI (Life Technologies). All groups were immunostained using identical procedures. Negative controls were processed using the same protocol with the omission of the primary or secondary antibodies to assess nonspecific labeling. There was no specific staining with these negative controls. Microphotographs were shot using a Carl Zeiss Axio Imager Z1 fluorescence microscope equipped with ApoTome (AxioVision 4.6.3, Carl Zeiss). The parameters for scanning were consistent across all groups.

Synaptic and neuritic quantification of primary cultures *in vitro* was carried out from 16 images, each coverslipped at a 40× objective lens. At least 2 coverslips, 32 images, and 150 neurons for each condition were analyzed. For synaptic analysis *in vivo* and to cover the hippocampal area, 3 of the same rectangular fields (90 × 70 µm) under 100× oil objective lens were precisely selected in the lateral and medial blade molecular layer (ML) of the dentate gyrus (DG), the stratum lacunosum-moleculare (SLM), the stratum radium (SR) and the stratum oriens (SO) of cornu ammonis 1 (CA1) in each condition, respectively. In addition, 2 of the same fields were carefully chosen in layers 2 and 3 of the entorhinal

cortex (ENT). Fifteen optical sections per field, 15 fields per hippocampal area, 4 fields per entorhinal cortex per section, and 855 total images per brain were analyzed. Single optical section images at 0.25 μm intervals and 3.75 μm Zeiss ApoTome high-resolution scans were performed. Synaptic puncta number and synaptic immunoreactive (IR) area were quantified using Puncta Analyzer (81, 82) and ImageJ (NIH) macro and batch process. Total neurite length was measured using the NeuriteTracer program (83). Briefly, the cultures were immunostained with Tuj1 for neurite and NeuN for the neuronal nucleus. For each condition, at least 150 primary neurons, 32 images in random fields from 2 coverslips in 2 independent experiments were analyzed. The NeuriteTracer was utilized to detect the neurites strongly stained for Tuj1. Following optimization of parameters to separate neurites from the neuronal cell body and tracing the neurite through skeletonization, positively labeled neurites and respective lengths were quantified (Figure 3B). The observer was blind to the treatment conditions. Average puncta number, synaptic area, and percentage of the area per image or per neuron were calculated for each condition.

Primary Cultures of Bone Marrow-Derived Macrophages and the Phagocytosis Assay

To test amyloid- β phagocytosis by macrophages, in repeated experiments monocytes were isolated from the bone marrow of wild-type mice ($n = 18$ mice) and differentiated into macrophages by 7-day cultivation in complete RPMI-1640 medium (#21870; Life Technologies) with 10% serum and 20 ng/ml MCSF (#315-02; PeproTech). Macrophage primary cultures were then plated at 1.2×10^5 cells per well (3–4 wells for each condition) in 24-well tissue-culture plates on glass coverslips overnight. Next, macrophages were either treated with 30 $\mu\text{g}/\text{ml}$ GA (Copaxone[®]; TEVA Neuroscience) for the duration of 1, 3, or 24 h, or not treated (control group). Before addition of *f*/*o*/sA β_{42} or vehicle, the cells were chilled in a 4°C ice bath for 5 min; immediately after addition of the preformed A β_{42} (100 nM), the plates were centrifuged at 515g in 25°C followed by incubation at 37°C for 30 or 60 min. The cells were then rinsed with A β -free medium to remove non-incorporated A β and later washed twice with PBS. Methanol (99.8%) at –20°C for 20 min or 4% paraformaldehyde at room temperature for 12 min were used for fixation of the cells followed by repeated washes with PBS. For immunostaining, the cells were first stained using the mouse anti-human amyloid- β mAb clone 6E10 (1:100; SIG-39320; Covance), rat anti-CD36 mAb clone MF3 (1:200; ab80080; Abcam), rat anti-CD204 scavenger receptor type I/II (SCARA1) mAb (1:100; MCA1322; AbD Serotec), rabbit anti-EEA1 pAb (1:100; Millipore #07-1820) and goat anti-MMP-9 pAb (1:100; AF909; R&D systems; **Supplementary Table 1**). Secondary polyclonal antibodies included donkey anti-mouse, anti-rat, anti-rabbit, and anti-goat conjugated with Cy2, Cy3, or Cy5 (1:200; Jackson ImmunoResearch Laboratories). The cells were mounted using ProLong[®] Gold with DAPI (Molecular Probes, Life Technologies; **Supplementary Table 1**). Several fields (minimum $n = 5$ randomly selected per group) were obtained from each well using a Carl Zeiss Axio Imager Z1 ApoTome-equipped microscope (an average of 120 cells in

each field). Images were obtained using the same exposure time on each occasion. The fluorescent signal and its total area were determined and quantified by the conversion of individual images to greyscale and standardizing to baseline using histogram-based thresholds with NIH ImageJ software. The “mean area per cell” was a result of a numerical average of the individual cell’s immunoreactive area per field. The “area/cell” measures the total fluorescent signal (area) divided by the total number of cells (DAPI count) of the same field (image). For all experiments, the investigators were blinded to the treatment condition. Colocalization of EEA1-6E10 (puncta analysis) was performed as synaptic puncta number analysis described above (81, 82).

Preparation of A β_{1-42} Oligomers, Fibrils, and Scrambled A β_{42}

Filtration was used to prepare LMW A β_{42} as described (84). Microcon YM-30 filters (EMD Millipore) were washed in 200 μl of distilled deionized water. A β was dissolved in 10% (v/v) 60 mM NaOH and 90% (v/v) 10 mM phosphate buffer, pH 7.4, at a concentration of 1 mg/ml. The solution was sonicated for 1 min and then placed into the washed filter. Following centrifugation at $14,000 \times g$ for 20 min, the filtrate, LMW A β_{42} , was collected. This LMW A β was stabilized using photo-induced cross-linking of unmodified proteins (PICUP) (76, 85–87) to yield a distribution of “oligomeric A β .” Briefly, 2 mM Tris(2,2’-bipyridyl)di-chlororuthenium(II) hexahydrate (Ru(Bpy); Aldrich) and 40 mM ammonium persulfate (APS; Sigma) were prepared in distilled deionized water. An 18 μl aliquot of 80 μM LMW A β_{42} was placed in a PCR tube, followed by 1 μl of Ru(Bpy) and 1 μl of APS. The sample was irradiated (150 W incandescent lamp) for 1 s and the reaction was quenched immediately with 1 M dithiothreitol. Cross-linking reagents were removed by dialysis using 3.5 kDa MWCO Slide-A-Lyzer cassettes (Pierce) against 10 mM sodium phosphate pH 7.4. More than 5 changes of buffer were completed, and protein purity and concentration were confirmed with optical absorption spectroscopy and SDS-PAGE.

For preparation of fibrillar A β_{42} , LMW A β_{42} was prepared as described above, and the peptide preparation was incubated at 100 μM with orbital shaking at 37°C until fibril formation occurred, approximately 2 weeks. The presence of fibrils was confirmed by electron microscopy. We used a previously designed “scrambled” A β_{42} peptide that: (1) had an amino acid composition identical to that of wild type A β_{42} , and (2) did not display amphipathicity (as does the wild type peptide). We accomplished this by using word scrambler software (for example: <https://www.wordunscrambler.net/word-scrambler.aspx>) to randomly permute the A β amino acid sequence. We then used Kyte-Doolittle analysis (88) to determine the hydrophathy profiles of the permuted sequences. This sequence met the 2 criteria for scrambled A β_{42} : YHAGVDKEVVFDEGAGAEHGLAQK - IVRGFGVSDVSMIHINLF.

Both WT and scrambled A β_{42} were synthesized using 9-fluorenylmethoxycarbonyl (Fmoc) chemistry, purified by reverse phase-high performance liquid chromatography, and

characterized by mass spectrometry and amino acid analysis, as described previously (89). Quantitative amino acid analysis and mass spectrometry yielded the expected compositions and molecular weights for each peptide. Purified peptides were stored as lyophilizates at -20°C . During fibril assembly studies, scrambled $\text{A}\beta_{42}$ has been shown to maintain a statistical coil secondary structure, bind thioflavin T poorly, and exhibit oligomerization characteristics distinct from those of wild type $\text{A}\beta_{42}$ (90). Scrambled $\text{A}\beta_{42}$ was dissolved in 10% (v/v) 60 mM NaOH and 90% (v/v) 10 mM phosphate buffer, pH 7.4, at a concentration of 1 mg/ml.

Transmission Electron Microscopy

Ten microliters of $\text{A}\beta$ were spotted onto 400 mesh carbon-coated Formvar grids (Electron Microscopy Sciences) and incubated for 2 min. Each grid was then negatively stained with 1% (v/v) filtered (0.2 μm) uranyl acetate (Ted Pella), which was immediately wicked off. Electron microscopy analysis was performed using JEOL 1200 EX at 80 KV (91).

All other experimental protocols, including genotyping, GA immunization, isolation and adoptive transfer of bone-marrow CD115^{+} monocytes, quantification and stereological counting, and Barnes maze behavioral tests were previously described (63).

Statistics

GraphPad Prism 5.0b (GraphPad Software) was used to analyze the data. A comparison of three or more groups was performed using two-way or one-way ANOVA followed by the Tukey's or Bonferonni's (92) *post hoc* multiple comparison test for paired groups. Two-group comparisons were analyzed using two-tailed paired *t*-tests (StatPlus). Correlation analyses were performed using the Pearson's coefficient (*r*) tests (Prism). Results are expressed as mean \pm standard error of the mean (s.e.m.). A *p* value less than 0.05 was considered significant.

RESULTS

Excitatory Synapses Are More Vulnerable to $\text{A}\beta_{42}$ Oligomers Than Fibrils

To assess the effect of $\text{A}\beta_{42}$ oligomers vs. fibrils on synaptic integrity and neuronal structure, postnatal day 1 (P1) primary cortical neurons (CN) were incubated for 12 h with 100 nM of either defined and stabilized oligomers (XL- $\text{oA}\beta_{42}$), preformed fibrils (f $\text{A}\beta_{42}$), or non-aggregated scrambled $\text{A}\beta$ control peptides (s $\text{A}\beta_{42}$) (90), compared to medium alone without $\text{A}\beta$ (vehicle; **Figures 1A–D**). The pure populations of $\text{A}\beta_{42}$ fibrils and stabilized low-n oligomers were verified by SDS-PAGE gel and transmission electron microscopy (**Figures 1A–C**). We observed no statistical difference in synaptic puncta number [analyzed by colocalized presynaptic VGluT1 (vesicular glutamate transporter 1) and PSD95 (postsynaptic density protein 95)] or neurite length between vehicle and s $\text{A}\beta_{42}$ groups (**Supplementary Figures 1A,B**). There was also no significant difference between CN co-cultured for 24 h with BM-derived M Φ for 24 h supplemented with either medium alone or s $\text{A}\beta_{42}$ (**Supplementary Figures 1A,B**). Interestingly, there was a small but significant increase in VGluT1/PSD95 synaptic number in

cortical neurons following 24 h co-culture with GA-treated M Φ (**Supplementary Figure 1A**). While s $\text{A}\beta_{42}$ did not induce loss of synapses or neuritic length retraction in CN, incubation with f $\text{A}\beta_{42}$, and moreover, with XL- $\text{oA}\beta_{42}$ (**Figure 1D**), decreased presynaptic VGluT1 $^{+}$ number by 63 and 81%, respectively, as compared to vehicle (**Figure 1E**; $P < 0.001$). Similarly, but to a lesser extent, postsynaptic PSD95 $^{+}$ puncta numbers were significantly reduced by 48 and 62% after exposure to f $\text{A}\beta_{42}$ or XL- $\text{oA}\beta_{42}$, respectively (**Figure 1E**; $P < 0.001$). Analysis of co-localized VGluT1 and PSD95 puncta number indicated a 50 and 70% reduction in synaptic density after 12 h exposure to f $\text{A}\beta_{42}$ or XL- $\text{oA}\beta_{42}$, respectively (**Figure 1F**; $P < 0.001$). Importantly, cortical excitatory synapses were substantially more vulnerable to low-n $\text{A}\beta_{42}$ oligomers than to their fibril counterparts (**Figure 1F**; $P < 0.001$).

Spontaneous Neuronal Hyperactivity and Altered Spike Waveforms Induced by $\text{A}\beta_{42}$ Oligomers

To examine the functional consequences of exposing neurons to $\text{A}\beta_{42}$ oligomers vs. fibrils, we recorded the spontaneous extracellular activity of P1 cortical neurons using Microelectrode arrays (MEA) after incubation with vehicle, f $\text{A}\beta_{42}$, or XL- $\text{oA}\beta_{42}$ for either 24 or 48 h (**Figures 2A–G** and **Supplementary Figures 2A–D**). We first compared the spontaneous activity levels across all neurons with a level of activity of at least 0.2 Hz. We found that after incubation with $\text{A}\beta_{42}$ oligomers for 24 h, spontaneous activity levels were significantly larger compared to incubation with vehicle or fibrils (**Figure 2A**; $P < 0.05$), whereas activity levels for incubation with fibrils was not significantly different from incubation with vehicle (**Figure 2A**; $P < 0.05$). After 48 h, neurons incubated with either f $\text{A}\beta_{42}$ or XL- $\text{oA}\beta_{42}$ had significantly elevated levels of spontaneous activity compared to vehicle (**Figure 2B**; $P < 0.001–0.05$). This shows that incubation with $\text{A}\beta_{42}$ oligomers is accompanied by hyperactivity as early as 24 h. This was also true when examining all neurons regardless of spontaneous activity levels and when only included neurons recorded at both time points (see **Supplementary Figures 2A–C**).

Did the presence of oligomers change the biophysical properties of neurons? To address this question, we next assessed whether the extracellular waveforms of the recorded action potentials differed between neurons exposed to fibrillar and oligomeric forms (**Figures 2D–G** and **Supplementary Figure 2D**). We found that two critical features of the extracellular waveform changed after incubation with oligomeric $\text{A}\beta_{42}$. First, the trough-to-peak width of waveforms (**Figure 2D**) recorded from neurons incubated with oligomeric $\text{A}\beta_{42}$ was significantly decreased relative to controls after 48 h (**Figure 2E**; $P < 0.01$); this effect was not apparent after 24 h (**Supplementary Figure 1D**). Second, the repolarization slope significantly decreased for neurons incubated with $\text{A}\beta_{42}$ oligomers, but not fibrils, relative to controls after 24 h (**Figure 2F**; $P = 0.0103$), an effect which persisted through 48 h (**Figure 2G**; $P < 0.05$). Together, this shows that incubation

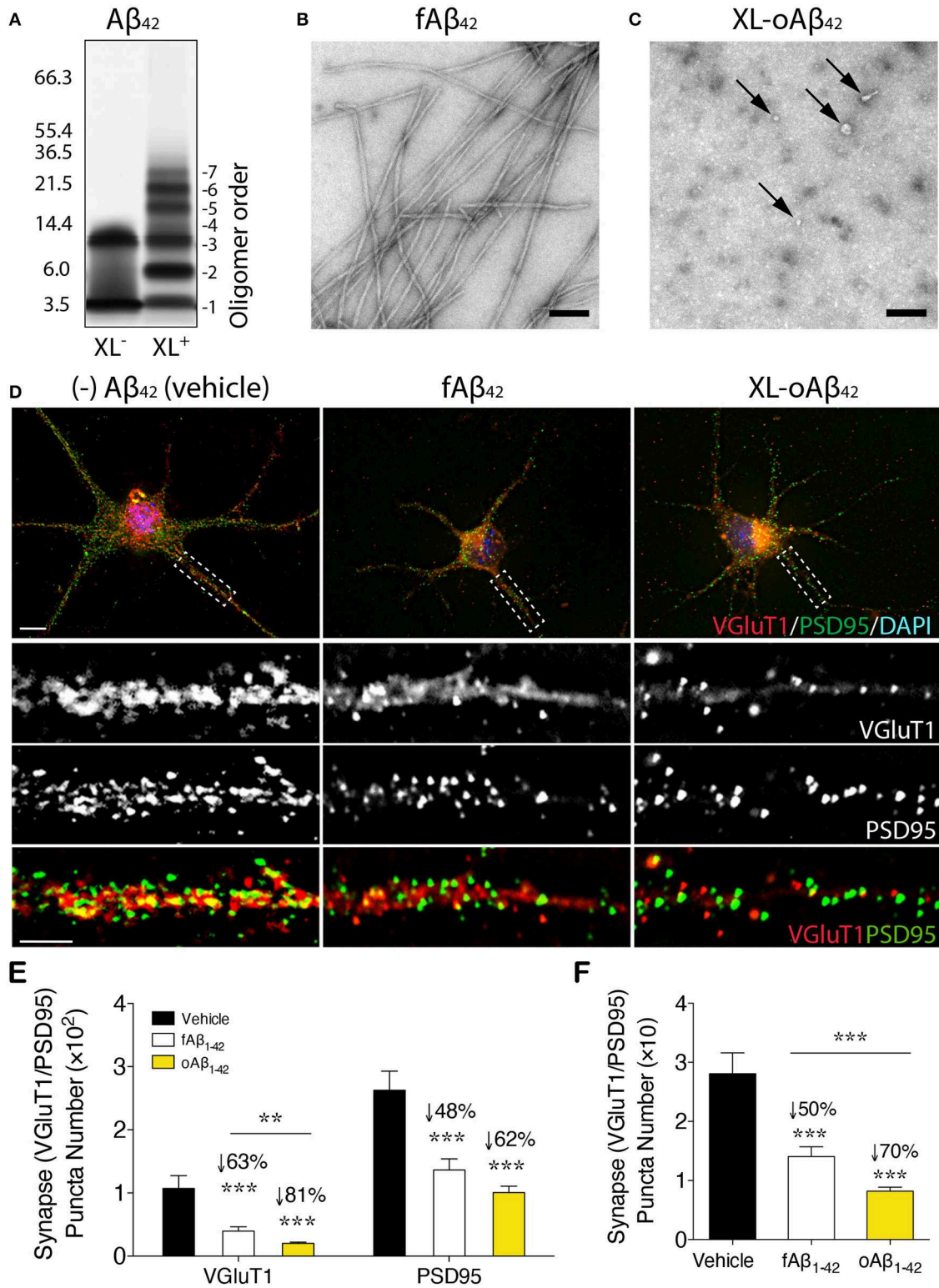
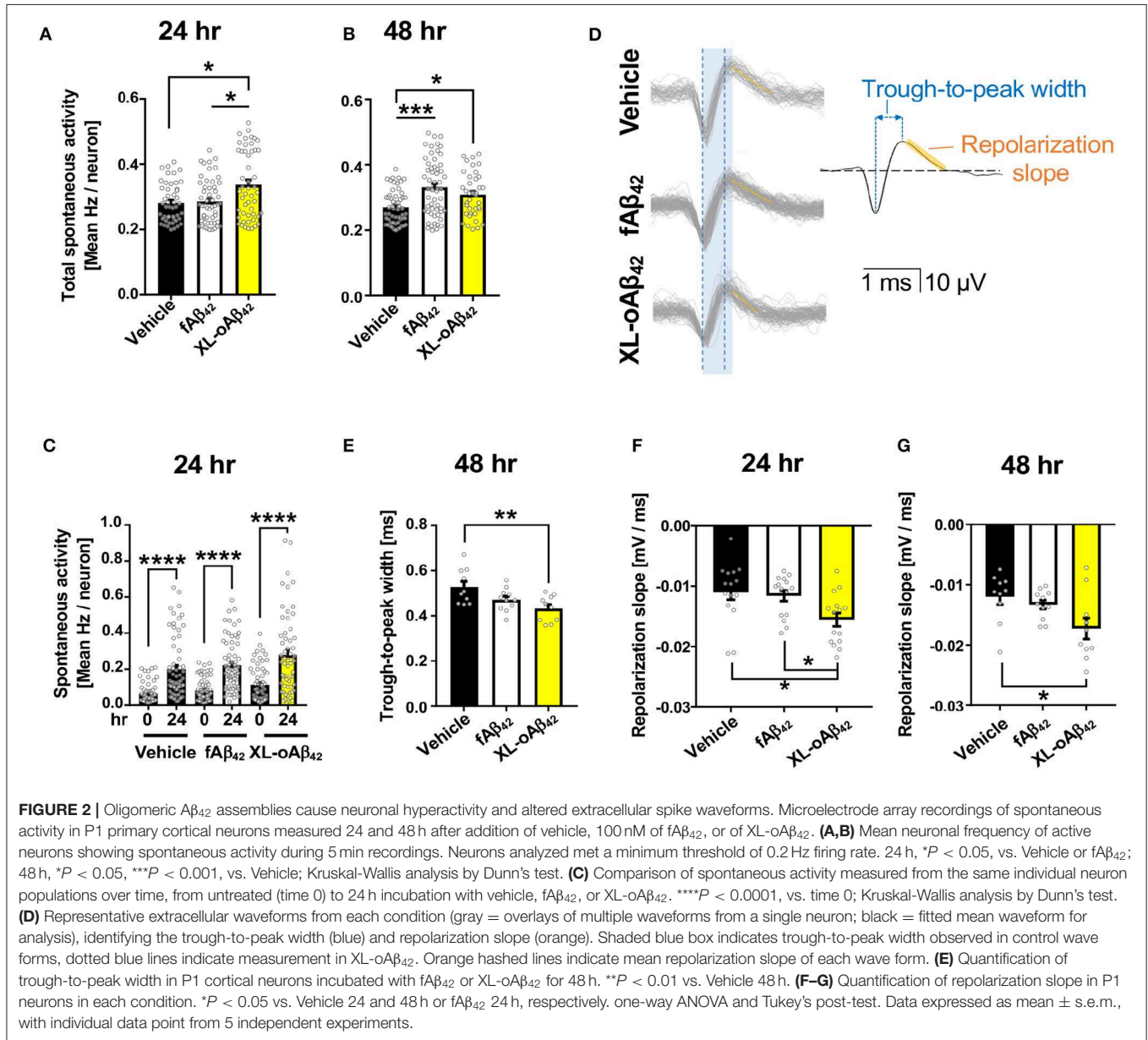


FIGURE 1 | Impact of defined and stabilized $A\beta_{42}$ oligomers vs. purified fibrils on synaptic integrity in primary cortical neurons. **(A)** SDS-PAGE image of cross-linked (XL^+) and non- XL (XL^-) $A\beta_{42}$ oligomers. **(B)** Electron microscopy (EM) of preformed fibrillar ($fA\beta_{42}$) and **(C)** $XL-oA\beta_{42}$, consisting of a defined distribution of covalently linked oligomers that were characterized by negative EM stain. Spherical (arrows) and amorphous structures were observed for $XL-oA\beta_{42}$, whereas typical amyloid (Continued)

FIGURE 1 | fibrils were seen in fA β_{42} . **(D)** Representative high-resolution images of pre-VGLuT1 and post-PSD95 synapses in primary cortical neurons, treated with no (-)A β_{42} (vehicle), fA β_{42} or XL-oA β_{42} . Areas within dashed boxes are magnified below. **(E,F)** Quantification of pre- and postsynapses in postnatal day 1 (P1) mouse primary cortical neurons. **(E)** VGLuT1- and PSD95-immunoreactive area per neuron and **(F)** colocalized VGLuT1/PSD95 synaptic puncta number in P1 cortical neurons treated with vehicle, fA β_{42} or XL-oA β_{42} . *** P < 0.001, fA β_{42} vs. vehicle or XL-oA β_{42} vs. vehicle; ** P < 0.01, fA β_{42} vs. XL-oA β_{42} , by one-way ANOVA and Tukey's post-test. Data expressed as mean \pm s.e.m.; n = 32 fields analyzed from 2 independent experiments; Scale bars = 100 nm **(B,C)** and 5 μ m **(D)**.



with A β_{42} oligomers was accompanied by spontaneous neuronal hyperactivity and altered extracellular waveforms.

Activated Macrophages Effectively Prevent Synaptic Loss and Neuritic Arborization Retraction Caused by A β_{42} Oligomers

To determine whether naïve BM-derived (M Φ ^{BM}) or GA-activated macrophages (GA-M Φ) can confer neuroprotection, P1

CN were co-cultured with M Φ ^{BM} or GA-M Φ in the presence of fA β_{42} or XL-oA β_{42} (**Figure 3**; experimental scheme and neurite tracing quantification method are shown in **Figures 3A,B**). Substantial synaptic loss and neuritic retraction were observed, especially after exposure to A β_{42} oligomers (**Figures 3C,D,E,G** vs. **Figure 3E**; P < 0.01; 32 and 65% VGLuT1/PSD95-synaptic loss in response to fibrils and oligomers, respectively). Importantly, reductions in pre- and postsynaptic puncta number (and area)

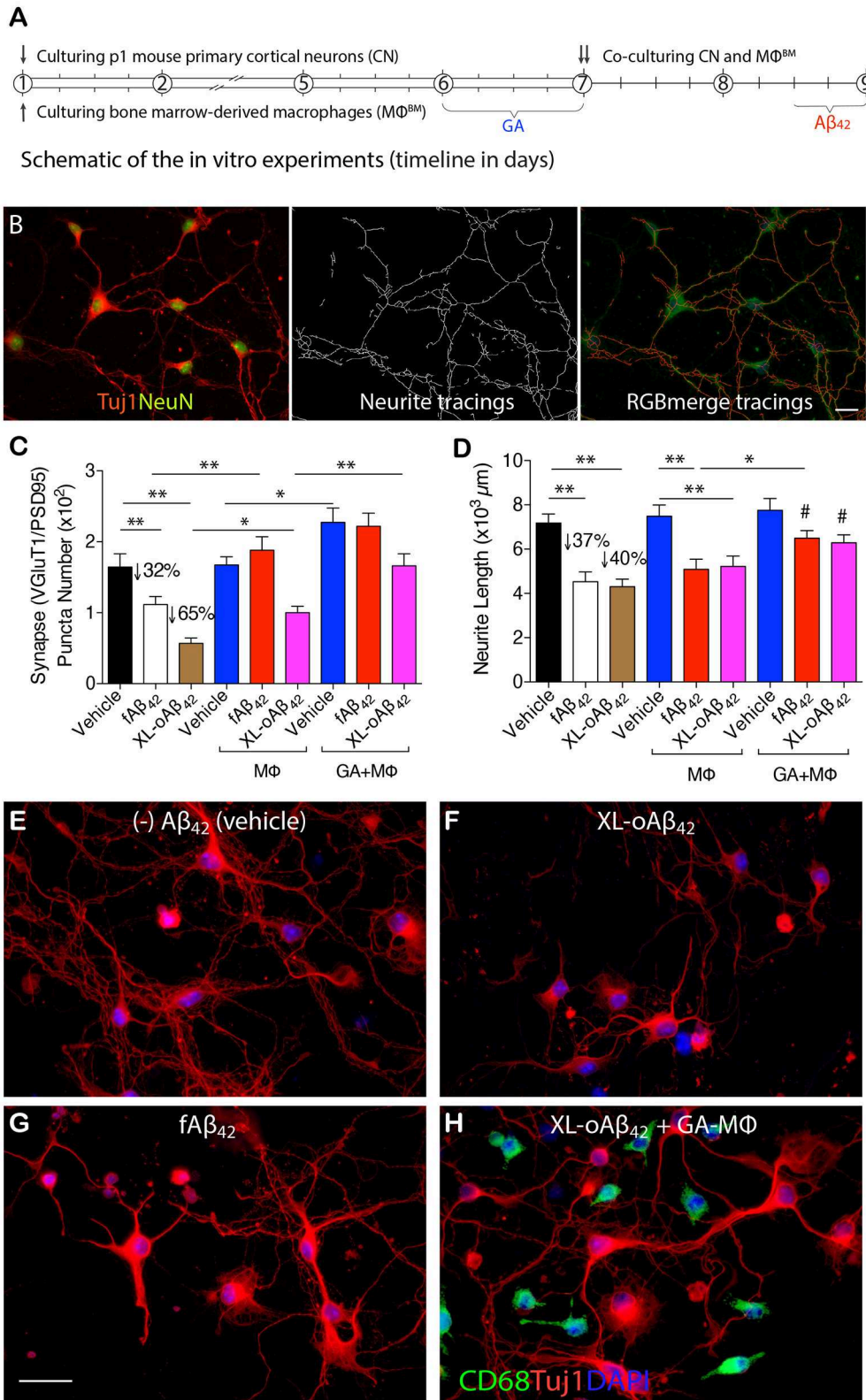


FIGURE 3 | Activated MΦ effectively protect against oligomeric Aβ₄₂-induced synaptic and neuritic arborization loss in primary cortical neurons. **(A)** Schematic of the *in vitro* experiments (timeline in days). P1 cortical neurons (treated with 100 nM XL-oAβ₄₂, fAβ₄₂, or vehicle for 12 h, respectively), bone marrow-derived MΦ (MΦ^{BM}),
 (Continued)

FIGURE 3 | and GA-activated M Φ ^{BM} (GA-M Φ) were cultured for 9 d. **(B)** Representative microphotographs of P1 neurons labeled with anti-Tuj1 and -NeuN serum (left), neurite tracings with NeuriteTracer (83) (middle), and RGB merge tracings (right). Scale bar represents 20 μ m. **(C)** Quantification of colocalized VGlut1/PSD95 synaptic puncta number in P1 neurons incubated with fA β ₄₂, XL-oA β ₄₂, or vehicle, and P1 neurons co-cultured with M Φ or with GA-M Φ . Note that fA β ₄₂ and XL-oA β ₄₂ both reduced the VGlut1/PSD95 synaptic density that was significantly preserved by co-culturing with M Φ . This effect was enhanced by co-culturing with GA-M Φ . **(D)** Quantification of neuritic length of P1 neurons incubated with fA β ₄₂, XL-oA β ₄₂, or vehicle, and P1 neurons co-cultured with M Φ or with GA-M Φ . Note that co-culturing with GA-M Φ significantly prevented decreases in neuritic length from fA β ₄₂ or XL-oA β ₄₂. Data expressed as mean \pm s.e.m.; $n = 48$ fields analyzed from 3 independent experiments; * $P < 0.05$, ** $P < 0.01$, comparisons as indicated by lines; # $P < 0.05$, vs. fA β ₄₂ or XL-oA β ₄₂ alone (no M Φ), by one-way ANOVA and Tukey's post-test. **(E-H)** Representative microphotographs of primary P1 neurons incubated with **(E)** vehicle, **(F)** XL-oA β ₄₂, **(G)** fA β ₄₂, and **(H)** co-cultured with GA-M Φ + XL-oA β ₄₂. Scale bar = 20 μ m.

after exposure to fA β ₄₂ were fully restored when co-cultured with M Φ ^{BM} or GA-M Φ (**Figure 3C** and **Supplementary Figures 3A, B**; $P < 0.01$), suggesting that M Φ have the capacity to effectively protect synapses against A β ₄₂ fibril toxicity within this time frame. However, untreated M Φ had limited ability to protect synapses against damage induced by A β ₄₂ oligomers (**Figure 3C**; $P < 0.05$), although importantly, GA activation of M Φ promoted full protection also against these highly synaptotoxic oligomers (**Figure 3C** and **Supplementary Figures 3A,B**; $P < 0.01$). While minimal to no protection against neuritic arborization loss was observed following co-culturing cortical neurons with naïve M Φ in the presence of either fA β ₄₂ or XL-oA β ₄₂ (**Figure 3D**), significant protection to neuronal dendritic structure was obtained with GA-activated M Φ , as measured by neurite process lengths (**Figure 3D**; $P < 0.05$). Interestingly, even in the absence of A β , co-cultures of cortical neurons with GA-M Φ vs. naïve M Φ had a significant increase in synaptic density (**Figure 3C**; $P < 0.05$). Thus, in primary cortical neurons, GA-activated M Φ effectively prevented loss of synapses and defects in arborization of neurite processes induced by both A β ₄₂ fibrils and oligomers.

MMP-9/Extracellular Degradation Is Predominant for Macrophage Clearance of A β ₄₂ Oligomers

To investigate how macrophages, including GA-activated macrophages, facilitate clearance of A β ₄₂ oligomers vs. fibrils and potentially increase synaptic protection, we characterized M Φ phenotypes, with and without GA stimulation, in response to the same quantities of XL-oA β ₄₂ and pre-formed fA β ₄₂ (100 nM; **Figure 4**). First, to assess surface recognition and intracellular uptake of A β ₄₂ assemblies, M Φ were incubated with XL-oA β ₄₂ or fA β ₄₂ for 30 min and labeled for A β (6E10), early endosome antigen 1 (EEA1), and class B scavenger receptor CD36 or class A type 1 scavenger receptor SCARA1 (encoded by macrophage scavenger receptor 1 – *MSR1* gene). There was a notable increase in CD36 expression, intracellular o/fA β ₄₂, and o/fA β ₄₂ colocalized within EEA1 endosomes in GA-M Φ vs. naïve M Φ (**Figures 4A, B**). A quantitative analysis of scavenger receptors indicated a significant increase in CD36 expression (**Figure 4C**; $P < 0.01$; **Supplementary Figure 4A**; $P < 0.05$) in GA-M Φ compared to untreated M Φ , but not in SCARA1 for XL-oA β ₄₂ (**Supplementary Figure 4B**). Regardless of GA, exposure to A β ₄₂ oligomers substantially upregulated surface expression of CD36 in M Φ (**Figure 4C**; $P < 0.001$). GA-M Φ exhibited significantly more intracellular uptake of both fA β ₄₂ and XL-oA β ₄₂ vs. naïve M Φ (**Figures 4D,E**; $P < 0.0001$ and $P < 0.001$, respectively).

Both naïve and GA-M Φ were on average six times less effective in intracellular uptake of oligomeric vs. fibrillar A β ₄₂ (**Figure 4D** vs. **Figure 4E**). Strikingly, as compared with fibrils, A β ₄₂ oligomers were approximately 2,500 times less targeted into early endosomes (**Figure 4F** vs. **Figure 4G**), as measured by 6E10⁺-A β puncta number colocalized in EEA1⁺ vesicles after 30 min phagocytosis assay. Hence, these data suggest that while intracellular A β ₄₂ fibrils are predominantly targeted to the endosomal-lysosomal proteolysis pathway, intracellular oligomers mostly escape these proteolytic vesicles in M Φ . Nonetheless, GA-M Φ significantly increased the presence of both A β ₄₂ fibrils and oligomers in EEA1⁺ endosomes (**Figures 4F,G**; $P < 0.01$ – 0.0001).

Next, we evaluated the extent to which M Φ ^{BM} degrade A β ₄₂ oligomers as compared to fibrillar forms in the extracellular space. To this end, we first assessed the expression of matrix metalloproteinase 9 (MMP-9)—an A β -degrading enzyme—by analysis of MMP-9⁺-immunoreactive area in M Φ (**Figure 4H**). Then, we determined A β _{1–42} concentrations in supernatant of M Φ following incubation with 100nM of either fA β ₄₂ or XL-oA β ₄₂, using a highly sensitive ELISA assay (**Figure 4I**). We previously showed that after exposure to A β ₄₂ fibrils, GA-M Φ exhibited a significant increase in MMP-9 levels (63, 70). Yet, the response to oligomers has not been reported. Here, we found that MMP-9 expression in GA-M Φ was *markedly* increased after incubation with XL-oA β ₄₂ (**Figure 4H**; $P < 0.0001$). Examination of extracellular degradation as a function of A β ₄₂ conformation indicated significant decreases in concentration of A β ₄₂ in the media of GA-M Φ vs. naïve M Φ after 24 h incubation with either fA β ₄₂ or XL-oA β ₄₂ (**Figure 4I**; 36 or 62% reduction, respectively, $P < 0.01$). Likewise, the substantial reduction in fA β ₄₂ and XL-oA β ₄₂ concentrations in the extracellular media of GA-treated vs. untreated M Φ can be seen after 30 min of incubation with these A β species (**Supplementary Figures 4C,D**; $P < 0.001$). Notably, at the two incubation time points (30 min and 24 h incubation), A β ₄₂ oligomers were primarily degraded in the extracellular rather than the intracellular space of macrophages and were considerably more accessible for extracellular degradation than their fibril counterparts (**Figure 4I**; 32% vs. 12 and 62% vs. 36% for M Φ and GA-M Φ , respectively; **Supplementary Figures 4C,D**; $P < 0.001$). Representative microscopic images in **Figure 4J** panel illustrate the increase in oA β ₄₂ uptake and CD36 expression by GA-activated M Φ , along with a decrease in extracellular or synaptic-bound oA β ₄₂, which resulted in synaptic and neuritic outgrowth preservation in primary cortical neurons. Overall, these data suggest that GA induces CD36/EEA1-mediated macrophage

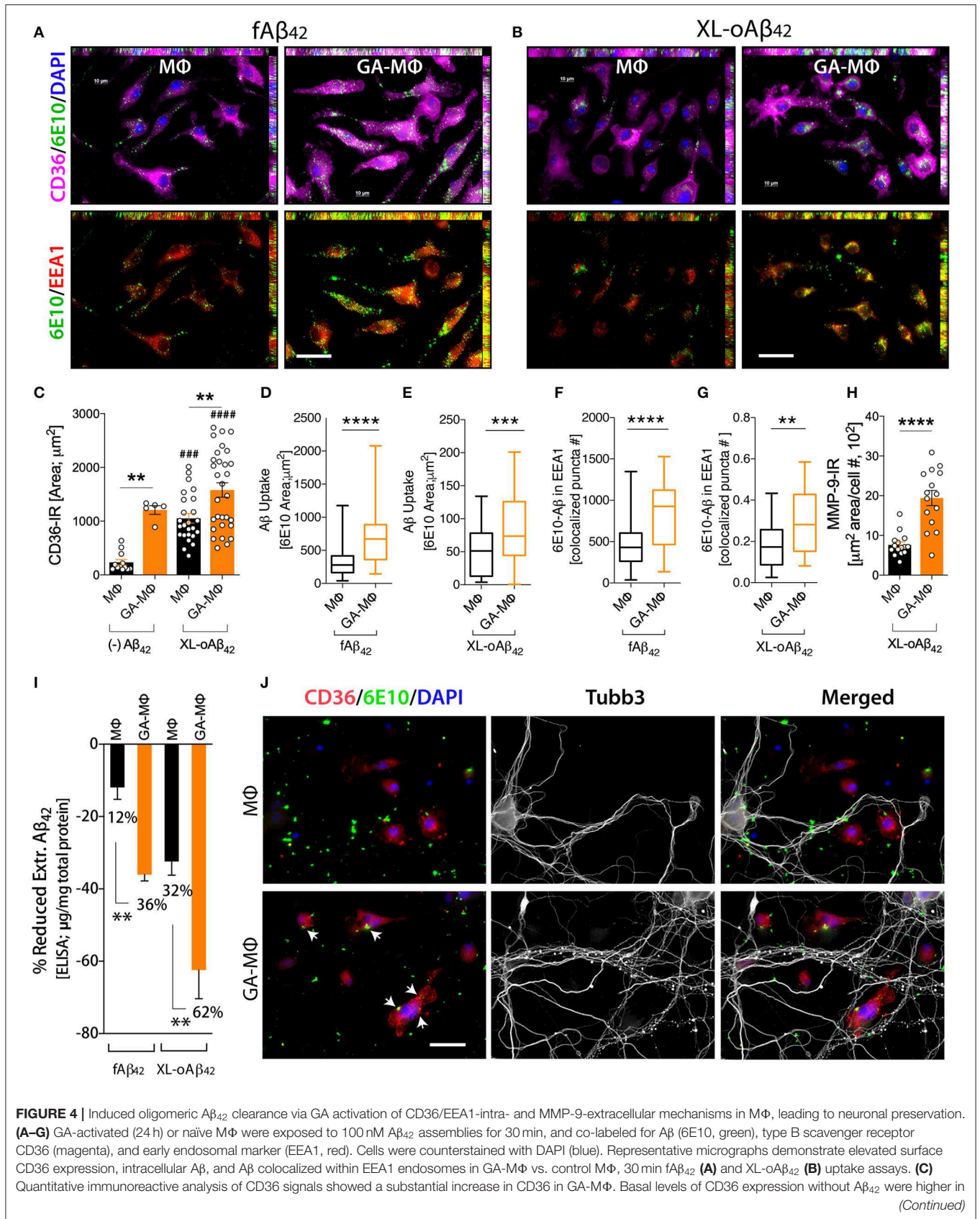


FIGURE 4 | GA-M Φ vs. control M Φ . This increase maintained its significance after exposure to XL-oA β_{42} . Regardless of GA effects, the presence of oligomeric A β_{42} substantially increased CD36 levels in both control M Φ and GA-M Φ . **(D,E)** Quantitative analysis of intracellular A β in M Φ following 30 min fA β_{42} **(D)** and XL-oA β_{42} **(E)**. **(F,G)** Quantitative analysis of fA β_{42} **(F)** and XL-oA β_{42} **(G)** within EEA1-positive vesicles in early endosomes in GA-M Φ vs. control M Φ . Co-labeled 6E10/EEA1-immunoreactive puncta number per M Φ is displayed. **(H)** Increased MMP-9 signal in GA-M Φ after exposure to XL-oA β_{42} . **(I)** ELISA quantification of extracellular A β_{42} after 24 h incubation with fA β_{42} or XL-oA β_{42} compared to 30 min basal peptide levels. **(J)** Representative microphotographs of co-cultured P1 neurons and M Φ exposed to XL-oA β_{42} for 24 h, stained with CD36, 6E10, Tubb3, and DAPI. P1 cortical neurons co-cultured with GA-M Φ vs. control M Φ exhibit increased neuritic outgrowth density with increased CD36-mediated A β uptake (arrows) and reduced extracellular A β_{42} . Data expressed as mean \pm s.e.m., with individual data point; $n = 48$ fields analyzed from 3 independent experiments; immunoreactive areas are normalized by cell number; ** $P < 0.01$, *** $P < 0.001$, **** $P < 0.0001$, GA-M Φ vs. M Φ with fA β_{42} or XL-oA β_{42} , and ### $P < 0.001$, #### $P < 0.0001$ for comparisons to M Φ without A β [(-) A β_{42}], by one-way ANOVA with Tukey's test or two-tailed student t -test. Scale bar = 20 μ m.

clearance, and moreover, MMP-9/extracellular degradation of synaptotoxic A β_{42} oligomers, ultimately protecting synapses, their dendritic structure, and neural function.

Severe Loss of Excitatory Pre- and Postsynaptic Terminals in the Entorhinal Cortex and Hippocampus of ADtg Mice

Given the crucial role of entorhinal cortex layers 2 and 3 in connecting brain cortical regions to the hippocampus, we sought to analyze pre- and postsynaptic VGLUT1 and PSD95 biomarkers in substructures of the entorhinal cortex (ENT) and hippocampus (HIPPO) of 10- and 13-month-old symptomatic ADtg vs. wild type (WT) mice (**Figures 5A–E**; **Supplementary Figures 5–8**). Due to lack of differences between the lateral (ML-L) and the medial (ML-M) blade molecular layers in pre- and postsynaptic areas across all experimental groups (**Supplementary Figure 5**), both regions were combined in our analyses (**Figures 5D,E**). Ten-month-old ADtg mice already exhibited significant pre- and postsynaptic losses in the molecular layers (ML) of the dentate gyrus (DG) and the stratum lacunosum-moleculare (SLM), but not in the stratum oriens (SO) nor in the stratum radiatum (SR) of the cornu ammonis 1 (CA1; **Supplementary Figure 6**). At 13 months, ADtg mice exhibited 50–62% presynaptic VGLUT1-immunoreactive (IR) area loss in all hippocampal regions as compared to WT mice (**Figure 5D**; $P < 0.05–0.001$). Similarly, analyses of postsynaptic PSD95-IR areas in 13-month-old ADtg mice vs. WT littermates indicated significant reductions by 28–42% in all hippocampal regions (**Figure 5E**; $P < 0.05$). In ENT layers 2 and 3, 68 and 53% of presynaptic VGLUT1-positive areas were lost (**Figure 5D**; $P < 0.001$), while 38 and 51% of postsynaptic PSD95 signals were reduced (**Figure 5E**; $P < 0.05$ and $P < 0.01$) in ADtg mice compared to WT mice, respectively. Overall, our data demonstrate a progressive and considerable loss of VGLUT1/PSD95 synaptic density, while the presynaptic VGLUT1 terminals appear more vulnerable to A β accumulation in the 13-month ADtg animals.

Synaptic Rescue Following Enhanced Cerebral Recruitment of Macrophages via Immunomodulation Therapy in ADtg Mice

To better understand the role of peripheral Mo/M Φ in preservation of excitatory synapses within AD-relevant cortical and hippocampal regions, we analyzed both pre- and postsynaptic density in substructures of ENT and HIPPO

following weekly s.c. injections of GA immunization or monthly i.v. injections of CD115⁺ Mo^{BM} into the peripheral blood (schematic of *in vivo* experiment in **Figure 5C**). Rescue pre- and postsynaptic density was noted in ENT2/3 by GA immunization or blood enrichment with Mo^{BM} in ADtg mice (**Figures 5F,G**, $P < 0.05–0.001$; **Supplementary Figure 7**, $P < 0.05–0.01$). Similarly, GA led to rescue of pre-VGLUT1 and post-PSD95 synapses in HIPPO across various ML, SLM, SR, and SO subregions (**Figures 5F,G**, $P < 0.05–0.01$; **Supplementary Figure 8**). Notably, throughout the analyzed brain regions, both immunomodulation approaches rescued postsynaptic density (i.e., PSD95-IR area) in symptomatic ADtg mice to levels similar or higher than those observed in WT mice (**Figure 5G**; **Supplementary Figures 7B, 8B**). No cumulative effect was noted between a group of ADtg mice treated with both GA and Mo^{BM} (combined treatment) and groups of mice treated with GA or Mo^{BM} alone (data not shown). Taken together, our *in vitro* and *in vivo* data suggest that immunomodulation approaches that deliver macrophages to reduce pathogenic A β_{42} burden at synaptic clefts substantially preserve pre- and postsynaptic terminals.

To identify confounders that potentially predict excitatory synaptic loss and contribute to immunotherapy-based synaptic preservation, we carried out multiple Pearson's coefficient (r) correlation analyses in the 13-month-old mouse cohort (**Supplementary Figure 9**). Specifically, we explored possible relationships between synaptic integrity and parameters of disease stage (cognitive function in Barnes maze test, astrogliosis, and amyloidosis) and therapeutic response (cerebral M Φ recruitment and phenotype). A strong linear association was observed between cerebral astrogliosis (GFAP⁺ reactive astrocyte area) and impaired cognitive function, as measured by error count on day 7 (long-term memory retention) of Barnes maze test (**Supplementary Figure 9A**; $P = 0.0187$, Pearson's $r = 0.60$). Similar correlations were found between hippocampal and cortical astrogliosis and other parameters of cognitive deficit in the Barnes maze test (not shown), suggesting that astrogliosis is a strong predictor of cognitive decline. Comparable to these observations, tight correlations were detected between cortical astrogliosis and postsynaptic loss (**Supplementary Figures 9B–D**; $P = 0.028–0.007$ and Pearson's $r = -0.55$ to -0.61). These data show an association between astrogliosis and cognitive outcome.

In relation to A β -plaque pathology, both cortical and hippocampal regions further revealed close inverse associations with postsynaptic density (**Supplementary Figures 9E–H**; $P =$

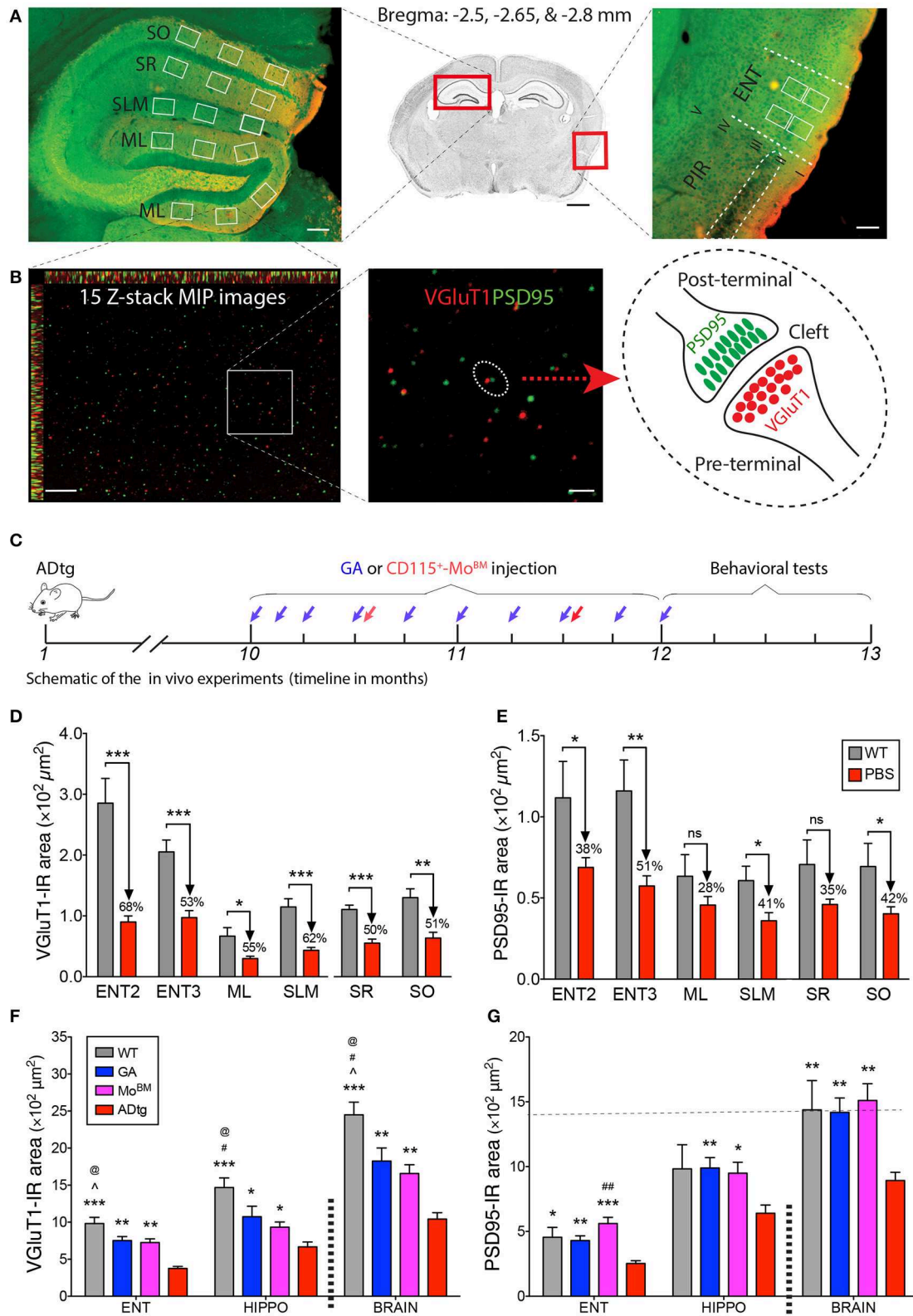


FIGURE 5 | Cortical and hippocampal synaptic rescue by CD115⁺-monocyte blood enrichment and GA immunization in ADtg mice. **(A,B)** VGluT1/PSD95 synaptic quantification scheme in entorhinal cortex (ENT) and hippocampus (HIPPO). **(A)** (Middle) Representative crystal violet stained image of a coronal brain section at (Continued)

FIGURE 5 | Bregma -2.65 mm. Scale bar = 1 mm. (Left) Representative microimage of HIPPO stained with the presynaptic (VGLuT1, red) and postsynaptic (PSD95, green) markers. Note that 3 fields are chosen in each of the lateral and medial blade molecular layers (ML) of dentate gyrus (DG), as well as in the stratum lacunosum-moleculare (SLM), stratum radiatum (SR), and stratum oriens (SO) of CA1. (Right) Representative microphotograph of ENT immunolabeled with the same pre- and postsynaptic markers. Note that ENT adjacent to piriform cortex (PIR) and 2 same fields are precisely selected from layers 2 and 3. Scale bar = $100\ \mu\text{m}$. **(B)** Representative microphotographs from the lateral blade ML immunostained with the same synaptic markers. Each high-magnification maximum intensity projection (MIP) image contains 15 Z-stack optical scanning (left). Scale bar = $10\ \mu\text{m}$. (Middle) Magnification for visualizing the proximity of pre- and postsynaptic signals. Scale bar = $2\ \mu\text{m}$. (Right) Schematic representation of pre- and post-terminal synapse. **(C)** Schematic of the *in vivo* experiments, where 10-month-old ADtg mice (all males) were injected with weekly s.c. GA immunization or monthly i.v. CD115⁺-Mo^{BM} injections for a duration of 2 months. Control groups were either naïve WT mice or monthly i.v. PBS injected ADtg mice. At the completion of the experiment, mice underwent behavioral testing, were euthanized, and brains were collected for analyses. **(D,E)** Quantification of presynaptic VGLuT1 **(D)** and postsynaptic PSD95 **(E)** areas in ENT2/3 and hippocampal substructures of 13-month-old ADtg mice vs. age-matched WT littermates. **(F,G)** Quantification of VGLuT1 **(F)** and PSD95 **(G)** immunoreactive areas in the entorhinal cortex, hippocampus, and whole brain of WT, GA-immunized, Mo^{BM}-injected, and PBS-control mice. Data expressed as mean \pm s.e.m.; $n = 6$ mice per group; * $P < 0.05$, ** $P < 0.01$, *** $P < 0.001$, and # $P < 0.001$, ## $P < 0.0001$, compared to ADtg control; @ $P < 0.05$ compared to Mo^{BM}, ^ $P < 0.05$ compared to GA, one-way ANOVA with Tukey's test.

0.011–0.002 and Pearson's $r = -0.57$ to -0.70). Similar correlations were also observed for presynaptic areas (not shown). These data suggest that cerebral A β -plaque burden are important indicators of synaptic loss in this murine model. We then established the relationship between excitatory synaptic integrity and cognition, showing significant associations between hippocampal and cortical synaptic density and cognitive function (**Supplementary Figures 9I,J**; $P = 0.015$ – 0.009 and Pearson's $r = -0.45$ to -0.53). The latter cognitive performance was determined by escape latency times in the learning and memory reversal phase on day 9 of the Barnes maze test. These data suggest that synaptic integrity is tightly associated with cognitive function, and that both astrogliosis and amyloidosis adversely affect it.

We previously showed that effective immunomodulation approaches in ADtg mice, such as GA immunization or blood enrichment with Mo^{BM}, reduced A β -plaque pathology and preserved cognition (50, 52, 63, 70). These therapeutic effects were achieved via cerebral recruitment of Iba-1⁺/CD115⁺/CD45^{hi} Mo/M Φ that expressed high levels of osteopontin (OPN; encoded by the gene secreted phosphoprotein 1 or SPP1) (63, 70). OPN/SPP1 is an immunoregulatory cytokine shown to both promote synaptogenesis (93–95) and be crucial for M Φ -mediated A β ₄₂ clearance (70). Significant and strong associations were observed between cerebral and hippocampal Iba-1⁺/CD45^{hi} Mo/M Φ (Mo/M Φ cell count) and cognitive function, as measured by latency times and error counts on day 9 of the Barnes maze test (**Supplementary Figures 9K,L**; $P = 0.0472$ – 0.011 and Pearson's $r = -0.64$ – -0.76). We further found a significant and robust association between OPN and postsynaptic but not presynaptic density areas (**Supplementary Figures 9M,N**; $P = 0.0133$ and Pearson's $r = 0.69$). These analyses suggest that increased Mo/M Φ -mediated OPN/SPP1 presence in the brain is linked with synaptic integrity and cognitive preservation.

DISCUSSION

This study supports the conclusion that activated M Φ ^{BM} can reduce A β ₄₂ oligomers and protect synaptic integrity and neuronal structure. Initially, we determined the impact of defined and stabilized populations of soluble A β ₄₂ oligomers on

excitatory synapses, neuritic arborization, and neuronal function as compared to pure fibrils. We indicated that synaptotoxicity was not elicited by non-aggregating sA β ₄₂ peptides but was dependent on their conformation state; cortical neurons were far more vulnerable to A β ₄₂ oligomers than to fibrils. A β ₄₂ oligomers caused severe synaptic loss, neurite arborization retraction, and hyperactivity, as well as changes in two critical features of the extracellular waveform in cortical neurons. We further found that M Φ ^{BM} effectively protect against synaptic loss and neurite retraction caused by A β ₄₂ fibrils yet provide insufficient protection against oligomers. Importantly, M Φ activation with glatiramer acetate (GA, an FDA-approved drug) fully protected neurons from detrimental effects triggered by both oligomers and fibrils. We identified distinct M Φ -mediated mechanisms for clearance of oligomers as compared to fibrils leading to synaptic protection. While eradication of A β ₄₂ oligomers predominately involves increased extracellular/MMP-9 degradation, fibrils are primarily eliminated via CD36/EEA1⁺-endosomal intracellular proteolysis. Strikingly, both pathways are boosted in M Φ stimulated with GA. Finally, our *in vivo* studies corroborated the *in vitro* findings, demonstrating that loss of cortical and hippocampal excitatory synapses in transgenic mouse models of AD is reversed by increased cerebral recruitment of Mo-derived M Φ , achieved via GA immunomodulation or peripheral-blood enrichment of BM-derived monocytes. Overall, these findings should guide future therapeutic interventions for AD to boost M Φ -mediated clearance of synaptotoxic A β ₄₂ oligomers and preserve synapses and cognition.

Our data in mouse primary cortical neurons provide evidence to support the notion that A β ₄₂ oligomers pose a greater threat to the neuronal network than do fibrils. These results align with previous studies reporting that A β ₄₂ oligomers impair neuronal plasticity and activation of excitatory mGluR hippocampal neurons, as well as exhibit direct synaptotoxic effects leading to cognitive impairments (8, 13, 15, 23–29, 96). As compared to postsynaptic terminals, presynaptic terminals displayed greater vulnerability to A β ₄₂ assemblies in both *in vivo* and *in vitro* models. We found that compared to fibrils, stabilized A β ₄₂ oligomers caused greater synaptic loss and neuritic retraction along with hyperactivity and altered action potential shape in primary cortical neurons. Due to the massive reduction in synapses ($\sim 70\%$) observed in response to A β ₄₂ oligomers as early as 12 h after incubation, a corresponding

reduction in neuronal activity was anticipated. To the contrary, however, we found elevated levels of spontaneous activity after incubation with oligomers. These observations are in accordance with previous studies in murine models of AD that showed increased neuronal excitability near A β plaques or in A β -depositing mouse brains, using Ca²⁺ imaging and *in vivo* intracellular recording (97, 98). Combined with the modified action potential shape we documented, this suggests that oligomeric A β ₄₂ directly affects the biophysical (intrinsic) properties of neurons, possibly via altered ion channel densities, in addition to its effects on synapses. Because oligomeric A β ₄₂ species are highly metastable and exist in dynamically changing mixtures, the conclusions of previous reports on A β ₄₂ oligomers may have been confounded by the existence of heterogeneous A β populations. By using defined and stabilized A β ₄₂ oligomers, this study establishes the precise neuropathological consequences caused by oligomers vs. fibrils and supports therapeutic avenues that clear cerebral A β ₄₂ oligomers to prevent such neuronal and synaptic damage.

One powerful method to clear A β ₄₂ oligomers and protect synapses was revealed in this study. Cortical neurons that were severely impaired by A β ₄₂ oligomers showed partial recovery of pre- and postsynaptic density and neurite arborization when co-cultured with M Φ ^{BM}, and full recovery with GA-M Φ . While naïve M Φ were able to effectively protect against synaptic and neurite length loss caused by the fibrils, they exhibited poor capacity to prevent oligomeric-induced synaptotoxicity. Yet, GA-M Φ successfully eliminated both fibrils and oligomers, essentially protecting neuronal network integrity despite these adverse conditions. Postsynaptic terminals, which were less vulnerable to A β ₄₂ species, were fully rescued by the presence of GA-M Φ . However, it would be interesting to look at the alteration of inhibitory gephyrin immunosignal expression following the administration of GA or GA-M Φ . We discovered that the EEA1⁺-early endosomal/lysosomal pathway was far more involved in the M Φ -mediated engulfment of fibrils than of oligomers. In contrast, reduction of A β ₄₂ in the extracellular space of M Φ was substantially more efficient for oligomers than for fibrils, potentially through enhanced release and activity of proteolytic enzymes (i.e., MMP-9). These *in vitro* data corroborate our *in vivo* studies in ADtg mice that demonstrated enhanced MMP-9 expression by infiltrating macrophages surrounding cerebral A β plaques in GA-immunized animals (63, 70). Notably, GA enhanced both internalization and extracellular pathways in M Φ to break down synaptotoxic A β ₄₂ forms. Recent studies indicated that depleting microglia with colony-stimulating factor 1 receptor (CSF1R) signaling inhibitors could impair formation of the plaques in an AD model (99, 100). Microglial depletion was further found not only to lead to a complete rescue of deficits in proliferation, differentiation and survival of adult hippocampal neural progenitor cells (101), but also significantly reduce ApoE, which is a major constituent of amyloid plaques and promotes their aggregation and deposition (102). It would be interesting to look at changes of synapses, A β deposition and cognition when microglial depletion plus GA immunotherapy is applied to this ADtg model. Additional studies are needed to

further establish direct structure-function connection between these and other A β ₄₂ forms (e.g., dimers, trimers, etc.), synaptotoxicity, and macrophage-mediated synaptic protection. A limitation of this study is that it does not allow for conclusions regarding potential rescue of neuronal function by macrophages due to lack of electrophysiological studies in macrophages and neurons co-cultures. Yet, since this study clearly demonstrates neuroprotection by morphological (neurite length) and synaptic biomarker measurements, future studies should evaluate whether these are accompanied with functional protection.

In this study, we demonstrated that progressive cortical and hippocampal pre- and postsynaptic loss tightly correlated with increased cerebral A β burden in the double-transgenic APP_{SWE}/PS1 Δ E9 murine model. The direct association between pre- and postsynaptic biomarkers and cognitive scores further emphasizes a potential link between A β ₄₂, synaptotoxicity, and cognitive decline. Moreover, the unexpected strong association between synaptic density and GFAP⁺ astrogliosis suggested that beyond direct A β toxicity, chronically reactive astrocytes not only fail to protect synapses, but also detrimentally impair them and lead to cognitive decline. Future investigations are warranted to better understand the impact of astrocytes and their phenotype on synaptic and cognitive decline in AD.

Here, we showed that both immunotherapies—GA immunization or blood-enrichment with CD115⁺ monocytes—restored presynaptic and fully rescued postsynaptic density in ADtg mice exhibiting significant synaptic loss. In our previous studies, these immunomodulation approaches have led to increased cerebral infiltration of monocyte-derived macrophages directly involved in A β clearance, leading to improved hippocampal-based memory and learning (52, 63). In this study, cognitive function, which directly correlated with synaptic integrity, also strongly associated with cerebral recruitment of (GFP⁺) Iba1⁺/CD45^{high} monocytes and macrophages. These data suggest that M Φ activity in the brain associated with decreased synaptotoxic A β levels can enhance synaptic protection and regeneration. One interesting question would be what happens if microglia as residential macrophages are treated with GA. Immunophenotypes of microglia expressed differently in dementia with AD pathology. For example, Iba1 was negatively associated with AD while levels of CD68, MSR-A, and CD64 expression were positively related to the presence of dementia (103). In a previous paper we determined a significant reduction of GFAP expression in the cortex of AD mice following GA or/and GA-M Φ immunotherapy (63), and found GFAP expression in this model was negatively associated with pre-synapse VGluT1. We now further demonstrated that the relationship of GFAP levels vs. post-synapse PSD95 in the cortex of AD mice exhibits the same behavior pattern as that of GFAP vs. VGluT1.

Cerebral clearance of A β by innate immune cells has been shown to take place via multiple mechanisms, including phagocytosis and enzymatic degradation (48, 49, 51, 54, 56, 57, 61, 104–106). Increasing evidence over the past decade has established the notion that peripheral M Φ contribute to the phagocytosis of cerebral A β plaque, an idea that is

further supported by reports of defective A β phagocytosis by M Φ isolated from AD patients (52, 53, 58, 60, 63, 104, 107). Several studies corroborating this idea have shown the involvement of scavenger receptors such as CD36 and SCARA-1 in facilitating fibrillar A β engulfment by innate immune cells. Rare variants in the *TREM2* gene were also associated with diminished M Φ phagocytosis and increased risk for AD (108–110). Indeed, parallel to the aforementioned investigations, this study's findings indicated elevated levels of surface scavenger receptors (e.g., CD36) in response to not only fibrillar but also oligomeric A β_{42} , with increased uptake by GA-M Φ . Importantly, the role of M Φ and, moreover, GA-activated M Φ in the clearance of non-fibrillar oligomeric A β forms has not been previously studied (111, 112). Enhanced colocalization of oligomers and fibrils in EEA1⁺-vesicles in GA-M Φ imply that these A β assemblies are redirected to lysosomal degradation rather than for perinuclear or endoplasmic reticulum pathologic accumulation (113). Therefore, this FDA-approved drug emerges again as a promising modality to safely stimulate phagocytic M Φ and improve ability to resist synaptotoxic A β forms associated with AD.

In the absence of A β , cortical neurons supplemented with GA-M Φ exhibited elevated levels of synaptic density and neuritic length that were higher than that of the baseline levels detected in cortical neurons alone. These results suggest that increased synaptic density in the presence of GA-M Φ is a result of both A β_{42} clearance and upregulation of neurotrophic factors stimulating synaptogenesis, such as insulin-like growth factor 1 (IGF-1), transforming growth factor beta 1 (TGF β 1), and/or OPN/SPP1 (50, 66, 70). In our previous study, we observed that OPN/SPP1 was markedly elevated in M Φ following GA treatment *in vitro* and *in vivo* (70). Here, a strong correlation (Pearson's $r = 0.7$) was found between cerebral OPN/SPP1 levels and postsynaptic density, which can explain increased synaptogenesis in the presence of M Φ^{BM} overexpressing OPN in the brains of immunized ADtg mice (70). In support of these findings, we reported that harnessing monocyte-derived macrophages by GA- or MOG-45D-immunization in ADtg mice enhanced cerebral expression of neurotrophic factors such as IGF-1, TGF β 1, and nuclear transcription factor early growth response 1 (Egr-1) (50, 58, 66), concomitant with recovered hippocampal neurogenesis. Hence, this immunomodulation approach holds great promise to positively affect neural regeneration as well as synaptic and cognitive preservation through combined M Φ activity of eradicating pathogenic A β_{42} forms and increasing neurotrophic support.

In conclusion, therapeutic approaches designed to prevent synaptotoxicity or even rescue synapses are becoming more prominent. Our extended data from both *in vitro* and *in vivo* studies may signify that GA-stimulated macrophages confer a potential synaptoprotective phenotype, at least in part through augmented capacity to eliminate highly toxic A β_{42} oligomers and induce synaptic preservation and regeneration. Hence, this study

encourages the development of a macrophage-based approach as therapeutic intervention for AD.

DATA AVAILABILITY STATEMENT

The raw data and protocols used to generate datasets for this study are presented in the article text, figures and tables. Supplementary data are also available online. Extended data can be made available upon reasonable request from the corresponding author to any qualified researcher.

ETHICS STATEMENT

The animal study was reviewed and approved by Cedars-Sinai Medical Center Institutional Animal Care and Use Committee (IACUC).

AUTHOR CONTRIBUTIONS

MK-H: study conception, design, study supervision, and final approval of the manuscript. SL: major experimental contributor, data acquisition and design. VG, YK, D-TF, JS, AR, and TT: experimental contributors with assistance from MK-H and SL. EH, JS, and DT: preparation of A β_{42} assemblies. DD: correlational analysis with guidance from MK-H. SL, VG, and MK-H: data analysis, interpretation, and presentation. MK-H, SL, EH, D-TF, TT, YK, UR, KB, and DT: discussion on intellectual content and manuscript editing. MK-H, SL, EH, VG, D-TF, UR, and DT: manuscript drafting, editing, and revisions.

FUNDING

This study was supported by the BrightFocus Foundation A2013328S00 grant (formerly AHAF; MK-H), The Coins for Alzheimer's Research Trust (C.A.R.T.; MK-H) Fund, NIH/NIA R01AG056478 and R01AG055865 (MK-H), The Cheryl and Haim Saban Foundation (MK-H), The Marciano Family Foundation (MK-H), National Natural Science Foundation of China 81671229, 81871032 (SL) and the National Center for Advancing Translational Sciences through CTSI (UL1TR000124; MK-H).

ACKNOWLEDGMENTS

We thank Mia Oviatt for assistance with editing this manuscript and Prof. Jevgenij Raskatov for his helpful comments. This paper is dedicated to Dr. Salomon Moni Hamaoui and Lillian Jones Black, both of whom died from Alzheimer's disease.

SUPPLEMENTARY MATERIAL

The Supplementary Material for this article can be found online at: <https://www.frontiersin.org/articles/10.3389/fimmu.2020.00049/full#supplementary-material>

REFERENCES

1. James BD, Leurgans SE, Hebert LE, Scherr PA, Yaffe K, Bennett DA. Contribution of Alzheimer disease to mortality in the United States. *Neurology*. (2014) 82:1045–50. doi: 10.1212/WNL.0000000000000240
2. Alzheimer's Association. 2018 Alzheimer's disease facts and figures. *Alzheimer's Dementia*. (2018) 14:367–429. doi: 10.1016/j.jalz.2018.02.001
3. Butterfield DA. Phosphoproteomics of Alzheimer disease brain: insights into altered brain protein regulation of critical neuronal functions and their contributions to subsequent cognitive loss. *Biochim Biophys Acta Mol Basis Dis*. (2019) 1865:2031–9. doi: 10.1016/j.bbadis.2018.08.035
4. Cardozo PL, De Lima IBQ, Maciel EMA, Silva NC, Dobransky T, Ribeiro FM. Synaptic Elimination in Neurological Disorders. *Curr Neuropharmacol*. (2019) 17:1071–95. doi: 10.2174/1570159X17666190603170511
5. Selkoe DJ. Alzheimer's disease results from the cerebral accumulation and cytotoxicity of amyloid beta-protein. *J Alzheimers Dis*. (2001) 3:75–80. doi: 10.3233/JAD-2001-3111
6. Hardy J, Selkoe DJ. The amyloid hypothesis of Alzheimer's disease: progress and problems on the road to therapeutics. *Science*. (2002) 297:353–6. doi: 10.1126/science.1072994
7. Selkoe DJ. Soluble oligomers of the amyloid beta-protein impair synaptic plasticity and behavior. *Behav Brain Res*. (2008) 192:106–13. doi: 10.1016/j.bbr.2008.02.016
8. Shankar GM, Li S, Mehta TH, Garcia-Munoz A, Shepardson NE, Smith I, et al. Amyloid-beta protein dimers isolated directly from Alzheimer's brains impair synaptic plasticity and memory. *Nat Med*. (2008) 14:837–42. doi: 10.1038/nm1782
9. Koronyo-Hamaoui M, Koronyo Y, Ljubimov AV, Miller CA, Ko MK, Black KL, et al. Identification of amyloid plaques in retinas from Alzheimer's patients and noninvasive *in vivo* optical imaging of retinal plaques in a mouse model. *Neuroimage*. (2011) 54(Suppl. 1):S204–17. doi: 10.1016/j.neuroimage.2010.06.020
10. La Morgia C, Ross-Cisneros FN, Koronyo Y, Hannibal J, Gallassi R, Cantalupo G, et al. Melanopsin retinal ganglion cell loss in Alzheimer disease. *Ann Neurol*. (2016) 79:90–109. doi: 10.1002/ana.24548
11. Koronyo YB, Biggs D, Barron E, Boyer DS, Pearlman JA, Au WJ. Retinal amyloid pathology and proof-of-concept imaging trial in Alzheimer's disease. *JCI Insight*. (2017) 2:e93621. doi: 10.1172/jci.insight.93621
12. Ferreira ST, Lourenco MV, Oliveira MM, De Felice FG. Soluble amyloid-beta oligomers as synaptotoxins leading to cognitive impairment in Alzheimer's disease. *Front Cell Neurosci*. (2015) 9:191. doi: 10.3389/fncel.2015.00191
13. Shankar GM, Bloodgood BL, Townsend M, Walsh DM, Selkoe DJ, Sabatini BL. Natural oligomers of the Alzheimer amyloid-beta protein induce reversible synapse loss by modulating an NMDA-type glutamate receptor-dependent signaling pathway. *J Neurosci*. (2007) 27:2866–75. doi: 10.1523/JNEUROSCI.4970-06.2007
14. Walsh DM, Selkoe DJ. A beta oligomers - a decade of discovery. *J Neurochem*. (2007) 101:1172–84. doi: 10.1111/j.1471-4159.2006.04426.x
15. Li S, Hong S, Shepardson NE, Walsh DM, Shankar GM, Selkoe D. Soluble oligomers of amyloid beta protein facilitate hippocampal long-term depression by disrupting neuronal glutamate uptake. *Neuron*. (2009) 62:788–801. doi: 10.1016/j.neuron.2009.05.012
16. Hayden EY, Conovaloff JL, Mason A, Bitan G, Teplow DB. Preparation of pure populations of covalently stabilized amyloid beta-protein oligomers of specific sizes. *Anal Biochem*. (2017) 518:78–85. doi: 10.1016/j.ab.2016.10.026
17. Li S, Jin M, Liu L, Dang Y, Ostaszewski BL, Selkoe DJ. Decoding the synaptic dysfunction of bioactive human AD brain soluble Abeta to inspire novel therapeutic avenues for Alzheimer's disease. *Acta Neuropathol Commun*. (2018) 6:121. doi: 10.1186/s40478-018-0626-x
18. Brinkmalm G, Hong W, Wang Z, Liu W, O'malley TT, Sun X, et al. Identification of neurotoxic cross-linked amyloid-beta dimers in the Alzheimer's brain. *Brain*. (2019) 142:1441–57. doi: 10.1093/brain/awz066
19. Raskatov JA. What is the "Relevant" Amyloid beta42 concentration? *Chembiochem*. (2019) 20:1725–6. doi: 10.1002/cbic.201900097
20. Yasumoto T, Takamura Y, Tsuji M, Watanabe-Nakayama T, Imamura K, Inoue H, et al. High molecular weight amyloid beta1-42 oligomers induce neurotoxicity via plasma membrane damage. *FASEB J*. (2019) 33:9220–34. doi: 10.1096/fj.201900604R
21. Shankar GM, Walsh DM. Alzheimer's disease: synaptic dysfunction and Abeta. *Mol Neurodegener*. (2009) 4:48. doi: 10.1186/1750-1326-4-48
22. Yu L, PetyukVA, Tasaki S, Boyle PA, Gaiteri C, Schneider JA, et al. Association of cortical beta-amyloid protein in the absence of insoluble deposits with Alzheimer disease. *JAMA Neurol*. (2019) 76:818–26. doi: 10.1001/jamaneurol.2019.0834
23. Lacor PN, Buniel MC, Furlow PW, Clemente AS, Velasco PT, Wood M, et al. Abeta oligomer-induced aberrations in synapse composition, shape, and density provide a molecular basis for loss of connectivity in Alzheimer's disease. *J Neurosci*. (2007) 27:796–807. doi: 10.1523/JNEUROSCI.3501-06.2007
24. Bitner RS, Bunnelle WH, Anderson DJ, Briggs CA, Buccafusco J, Curzon P, et al. Broad-spectrum efficacy across cognitive domains by alpha7 nicotinic acetylcholine receptor agonism correlates with activation of ERK1/2 and CREB phosphorylation pathways. *J Neurosci*. (2007) 27:10578–87. doi: 10.1523/JNEUROSCI.2444-07.2007
25. Vossel KA, Zhang K, Brodbeck J, Daub AC, Sharma P, Finkbeiner S, et al. Tau reduction prevents Abeta-induced defects in axonal transport. *Science*. (2010) 330:198. doi: 10.1126/science.1194653
26. Tang Y, Scott DA, Das U, Edland SD, Radomski K, Koo EH, et al. Early and selective impairments in axonal transport kinetics of synaptic cargoes induced by soluble amyloid beta-protein oligomers. *Traffic*. (2012) 13:681–93. doi: 10.1111/j.1600-0854.2012.01340.x
27. Walsh DM, Klyubin I, Fadeeva JV, Cullen WK, Anwyl R, Wolfe MS, et al. Naturally secreted oligomers of amyloid beta protein potently inhibit hippocampal long-term potentiation *in vivo*. *Nature*. (2002) 416:535–9. doi: 10.1038/416535a
28. Tu S, Okamoto S, Lipton SA, Xu H. Oligomeric Abeta-induced synaptic dysfunction in Alzheimer's disease. *Mol Neurodegener*. (2014) 9:48. doi: 10.1186/1750-1326-9-48
29. Umeda T, Ramser EM, Yamashita M, Nakajima K, Mori H, Silverman MA, et al. Intracellular amyloid beta oligomers impair organelle transport and induce dendritic spine loss in primary neurons. *Acta Neuropathol Commun*. (2015) 3:51. doi: 10.1186/s40478-015-0230-2
30. Selkoe DJ. Developing preventive therapies for chronic diseases: lessons learned from Alzheimer's disease. *Nutr Rev*. (2007) 65:S239–43. doi: 10.1111/j.1753-4887.2007.tb00370.x
31. Lee EB, Leng LZ, Zhang B, Kwong L, Trojanowski JQ, Abel T, et al. Targeting amyloid-beta peptide (Abeta) oligomers by passive immunization with a conformation-selective monoclonal antibody improves learning and memory in Abeta precursor protein (APP) transgenic mice. *J Biol Chem*. (2006) 281:4292–9. doi: 10.1074/jbc.M511018200
32. Klyubin I, Betts V, Welzel AT, Blennow K, Zetterberg H, Wallin A, et al. Amyloid beta protein dimer-containing human CSF disrupts synaptic plasticity: prevention by systemic passive immunization. *J Neurosci*. (2008) 28:4231–7. doi: 10.1523/JNEUROSCI.5161-07.2008
33. Nitsch RM, Hock C. Targeting beta-amyloid pathology in Alzheimer's disease with Abeta immunotherapy. *Neurotherapeutics*. (2008) 5:415–20. doi: 10.1016/j.nurt.2008.05.013
34. Cho JE, Kim JR. Recent approaches targeting beta-amyloid for therapeutic intervention of Alzheimer's disease. *Recent Pat CNS Drug Discov*. (2011) 6:222–33. doi: 10.2174/157488911796958002
35. Zuroff L, Daley D, Black KL, Koronyo-Hamaoui M. Clearance of cerebral Abeta in Alzheimer's disease: reassessing the role of microglia and monocytes. *Cell Mol Life Sci*. (2017) 74:2167–201. doi: 10.1007/s00018-017-2463-7
36. Guedes JR, Lao T, Cardoso AL, El Khoury J. Roles of microglial and monocyte chemokines and their receptors in regulating Alzheimer's disease-associated amyloid- β and tau pathologies. *Front Neurol*. (2018) 9:549. doi: 10.3389/fneur.2018.00549
37. Hickman SE, Allison EK, El Khoury J. Microglial dysfunction and defective beta-amyloid clearance pathways in aging Alzheimer's disease mice. *J Neurosci*. (2008) 28:8354–60. doi: 10.1523/JNEUROSCI.0616-08.2008

38. Tejera D, Heneka MT. Microglia in Alzheimer's disease: the good, the bad and the ugly. *Curr Alzheimer Res.* (2016) 13:370–80. doi: 10.2174/1567205013666151116125012
39. Hong S, Beja-Glasser VF, Nfonoyim BM, Frouin A, Li S, Ramakrishnan S, et al. Complement and microglia mediate early synapse loss in Alzheimer mouse models. *Science.* (2016) 352:712–6. doi: 10.1126/science.ad8373
40. Colonna M, Butovsky O. Microglia function in the central nervous system during health and neurodegeneration. *Annu Rev Immunol.* (2017) 35:441–68. doi: 10.1146/annurev-immunol-051116-052358
41. Keren-Shaul H, Spinrad A, Weiner A, Matcovitch-Natan O, Dvir-Szternfeld R, Ulland TK, et al. A Unique microglia type associated with restricting development of Alzheimer's Disease. *Cell.* (2017) 169:1276–90 e1217. doi: 10.1016/j.cell.2017.05.018
42. Krasemann S, Madore C, Cialic R, Baufeld C, Calcagno N, El Fatimy R, et al. The TREM2-APOE pathway drives the transcriptional phenotype of dysfunctional microglia in neurodegenerative diseases. *Immunity.* (2017) 47:566–81 e569. doi: 10.1016/j.immuni.2017.08.008
43. Song WM, Colonna M. The identity and function of microglia in neurodegeneration. *Nat Immunol.* (2018) 19:1048–58. doi: 10.1038/s41590-018-0212-1
44. Wendeln AC, Degenhardt K, Kaurani L, Gertig M, Ulas T, Jain G, et al. Innate immune memory in the brain shapes neurological disease hallmarks. *Nature.* (2018) 556:332–8. doi: 10.1038/s41586-018-0023-4
45. Ziegler-Waldkirch S, D'errico P, Sauer JF, Erny D, Savanthrapadian S, Loreth D, et al. Seed-induced Abeta deposition is modulated by microglia under environmental enrichment in a mouse model of Alzheimer's disease. *Embo J.* (2018) 37:167–82. doi: 10.15252/embj.2017.97021
46. Wang S, Colonna M. Microglia in Alzheimer's disease: a target for immunotherapy. *J Leukoc Biol.* (2019) 106:219–27. doi: 10.1002/JLB.MR0818-319R
47. Frautschy SA, Cole GM, Baird A. Phagocytosis and deposition of vascular beta-amyloid in rat brains injected with Alzheimer beta-amyloid. *Am J Pathol.* (1992) 140:1389–99.
48. Paresce DM, Ghosh RN, Maxfield FR. Microglial cells internalize aggregates of the Alzheimer's disease amyloid beta-protein via a scavenger receptor. *Neuron.* (1996) 17:553–65. doi: 10.1016/S0896-6273(00)80187-7
49. Wyss-Coray T, Lin C, Yan F, Yu GQ, Rohde M, Mconlogue L, et al. TGF-beta1 promotes microglial amyloid-beta clearance and reduces plaque burden in transgenic mice. *Nat Med.* (2001) 7:612–8. doi: 10.1038/87945
50. Butovsky O, Koronyo-Hamaoui M, Kunis G, Ophir E, Landa G, Cohen H, et al. Glatiramer acetate fights against Alzheimer's disease by inducing dendritic-like microglia expressing insulin-like growth factor 1. *Proc Natl Acad Sci USA.* (2006) 103:11784–9. doi: 10.1073/pnas.0604681103
51. Simard AR, Rivest S. Neuroprotective properties of the innate immune system and bone marrow stem cells in Alzheimer's disease. *Mol Psychiatry.* (2006) 11:327–35. doi: 10.1038/sj.mp.4001809
52. Butovsky O, Kunis G, Koronyo-Hamaoui M, Schwartz M. Selective ablation of bone marrow-derived dendritic cells increases amyloid plaques in a mouse Alzheimer's disease model. *Eur J Neurosci.* (2007) 26:413–6. doi: 10.1111/j.1460-9568.2007.05652.x
53. El Khoury J, Toft M, Hickman SE, Means TK, Terada K, Geula C, et al. Ccr2 deficiency impairs microglial accumulation and accelerates progression of Alzheimer-like disease. *Nat Med.* (2007) 13:432–8. doi: 10.1038/nm1555
54. Takata K, Kitamura Y, Yanagisawa D, Morikawa S, Morita M, Inubushi T, et al. Microglial transplantation increases amyloid-beta clearance in Alzheimer model rats. *FEBS Lett.* (2007) 581:475–8. doi: 10.1016/j.febslet.2007.01.009
55. Majumdar A, Chung H, Dolios G, Wang R, Asamoah N, Lobel P, et al. Degradation of fibrillar forms of Alzheimer's amyloid beta-peptide by macrophages. *Neurobiol Aging.* (2008) 29:707–15. doi: 10.1016/j.neurobiolaging.2006.12.001
56. Richard KL, Filali M, Prefontaine P, Rivest S. Toll-like receptor 2 acts as a natural innate immune receptor to clear amyloid beta 1-42 and delay the cognitive decline in a mouse model of Alzheimer's disease. *J Neurosci.* (2008) 28:5784–93. doi: 10.1523/JNEUROSCI.1146-08.2008
57. Bates KA, Verdile G, Li QX, Ames D, Hudson P, Masters CL, et al. Clearance mechanisms of Alzheimer's amyloid-beta peptide: implications for therapeutic design and diagnostic tests. *Mol Psychiatry.* (2009) 14:469–86. doi: 10.1038/mp.2008.96
58. Koronyo-Hamaoui M, Ko MK, Koronyo Y, Azoulay D, Seksenyan A, Kunis G, et al. Attenuation of AD-like neuropathology by harnessing peripheral immune cells: local elevation of IL-10 and MMP-9. *J Neurochem.* (2009) 111:1409–24. doi: 10.1111/j.1471-4159.2009.06402.x
59. Hickman SE, El Khoury J. Mechanisms of mononuclear phagocyte recruitment in Alzheimer's disease. *CNS Neurol Disord Drug Targets.* (2010) 9:168–73. doi: 10.2174/187152710791011982
60. Lebson L, Nash K, Kamath S, Herber D, Carty N, Lee DC, et al. Trafficking CD11b-positive blood cells deliver therapeutic genes to the brain of amyloid-depositing transgenic mice. *J Neurosci.* (2010) 30:9651–8. doi: 10.1523/JNEUROSCI.0329-10.2010
61. Lai AY, McLaurin J. Clearance of amyloid-beta peptides by microglia and macrophages: the issue of what, when and where. *Future Neurol.* (2012) 7:165–76. doi: 10.2217/fnl.12.6
62. Bernstein KE, Koronyo Y, Salumbides BC, Sheyn J, Pelissier L, Lopes DH, et al. Angiotensin-converting enzyme overexpression in myelomonocytes prevents Alzheimer's-like cognitive decline. *J Clin Invest.* (2014) 124:1000–12. doi: 10.1172/JCI66541
63. Koronyo Y, Salumbides BC, Sheyn J, Pelissier L, Li S, Ljubimov V, et al. Therapeutic effects of glatiramer acetate and grafted CD115(+) monocytes in a mouse model of Alzheimer's disease. *Brain.* (2015) 138:2399–422. doi: 10.1093/brain/awv150
64. He H, Altomare D, Ozer U, Xu H, Creek K, Chen H, et al. Cancer cell-selective killing polymer/copper combination. *Biomater Sci.* (2016) 4:115–20. doi: 10.1039/C5BM00325C
65. Frenkel D, Maron R, Burt DS, Weiner HL. Nasal vaccination with a proteosome-based adjuvant and glatiramer acetate clears beta-amyloid in a mouse model of Alzheimer disease. *J Clin Invest.* (2005) 115:2423–33. doi: 10.1172/JCI23241
66. Bakalash S, Pham M, Koronyo Y, Salumbides BC, Kramerov A, Seidenberg H, et al. Egr1 expression is induced following glatiramer acetate immunotherapy in rodent models of glaucoma and Alzheimer's disease. *Invest Ophthalmol Vis Sci.* (2011) 52:9033–46. doi: 10.1167/iiov.11-7498
67. Koronyo Y, Salumbides BC, Black KL, Koronyo-Hamaoui M. Alzheimer's disease in the retina: imaging retinal abeta plaques for early diagnosis and therapy assessment. *Neurodegener Dis.* (2012) 10:285–93. doi: 10.1159/000335154
68. Baruch K, Rosenzweig N, Kertser A, Deczkowska A, Sharif AM, Spinrad A, et al. Breaking immune tolerance by targeting Foxp3(+) regulatory T cells mitigates Alzheimer's disease pathology. *Nat Commun.* (2015) 6:7967. doi: 10.1038/ncomms8967
69. Kunis G, Baruch K, Miller O, Schwartz M. Immunization with a myelin-derived antigen activates the brain's choroid plexus for recruitment of immunoregulatory cells to the CNS and attenuates disease progression in a mouse model of ALS. *J Neurosci.* (2015) 35:6381–93. doi: 10.1523/JNEUROSCI.3644-14.2015
70. Rentsendorj A, Sheyn J, Fuchs DT, Daley D, Salumbides BC, Schubloom HE, et al. A novel role for osteopontin in macrophage-mediated amyloid-beta clearance in Alzheimer's models. *Brain Behav Immun.* (2018) 67:163–80. doi: 10.1016/j.bbi.2017.08.019
71. Khan UA, Liu L, Provenzano FA, Berman DE, Profaci CP, Sloan R, et al. Molecular drivers and cortical spread of lateral entorhinal cortex dysfunction in preclinical Alzheimer's disease. *Nat Neurosci.* (2014) 17:304–11. doi: 10.1038/nn.3606
72. Crisuolo C, Fontebasso V, Middei S, Stazi M, Ammassari-Teule M, Yan SS, et al. Entorhinal Cortex dysfunction can be rescued by inhibition of microglial RAGE in an Alzheimer's disease mouse model. *Sci Rep.* (2017) 7:42370. doi: 10.1038/srep42370
73. Frontinan-Rubio J, Sancho-Bielsa FJ, Peinado JR, Laferla FM, Gimenez-Llort L, Duran-Prado M, et al. Sex-dependent co-occurrence of hypoxia and beta-amyloid plaques in hippocampus and entorhinal cortex is reversed by long-term treatment with ubiquinol and ascorbic acid in the 3xTg-AD

- mouse model of Alzheimer's disease. *Mol Cell Neurosci.* (2018) 92:67–81. doi: 10.1016/j.mcn.2018.06.005
74. Marzi SJ, Leung SK, Ribarska T, Hannon E, Smith AR, Pishva E, et al. A histone acetylo-me-wide association study of Alzheimer's disease identifies disease-associated H3K27ac differences in the entorhinal cortex. *Nat Neurosci.* (2018) 21:1618–27. doi: 10.1038/s41593-018-0253-7
 75. Nakazono T, Jun H, Blurton-Jones M, Green KN, Igarashi KM. Gamma oscillations in the entorhinal-hippocampal circuit underlying memory and dementia. *Neurosci Res.* (2018) 129:40–6. doi: 10.1016/j.neures.2018.02.002
 76. Bitan G, Kirkitadze MD, Lomakin A, Vollers SS, Benedek GB, Teplow DB. Amyloid beta -protein (Abeta) assembly: Abeta 40 and Abeta 42 oligomerize through distinct pathways. *Proc Natl Acad Sci USA.* (2003) 100:330–5. doi: 10.1073/pnas.222681699
 77. Rahimi F, Maiti P, Bitan G. Photo-induced cross-linking of unmodified proteins (PICUP) applied to amyloidogenic peptides. *J Vis Exp.* (2009) 12:1071. doi: 10.3791/1071
 78. Beaudoin GM 3rd, Lee SH, Singh D, Yuan Y, Ng YG, Reichardt LF, Arikath J. Culturing pyramidal neurons from the early postnatal mouse hippocampus and cortex. *Nat Protoc.* (2012) 7:1741–54. doi: 10.1038/nprot.2012.099
 79. Li S, Nie EH, Yin Y, Benowitz LI, Tung S, Vinters HV, et al. GDF10 is a signal for axonal sprouting and functional recovery after stroke. *Nat Neurosci.* (2015) 18:1737–45. doi: 10.1038/nn.4146
 80. Paxinos G, Franklin KB. *The Mouse Brain in Stereotaxic Coordinates.* 2nd ed. San Diego, CA: Academic Press (2001).
 81. Eroglu C, Allen NJ, Susman MW, O'Rourke NA, Park CY, Ozkan E, et al. Gabapentin receptor alpha2delta-1 is a neuronal thrombospondin receptor responsible for excitatory CNS synaptogenesis. *Cell.* (2009) 139:380–92. doi: 10.1016/j.cell.2009.09.025
 82. Ippolito DM, Eroglu C. Quantifying synapses: an immunocytochemistry-based assay to quantify synapse number. *J Vis Exp.* (2010) 19:2270. doi: 10.3791/2270
 83. Pool M, Thiemann J, Bar-Or A, Fournier AE. NeuriteTracer: a novel ImageJ plugin for automated quantification of neurite outgrowth. *J Neurosci Methods.* (2008) 168:134–9. doi: 10.1016/j.jneumeth.2007.08.029
 84. Bitan G, Teplow DB. Preparation of aggregate-free, low molecular weight amyloid-beta for assembly and toxicity assays. *Methods Mol Biol.* (2005) 299:3–9. doi: 10.1385/1-59259-874-9:003
 85. Bitan G, Lomakin A, Teplow DB. Amyloid beta-protein oligomerization: pre-nucleation interactions revealed by photo-induced cross-linking of unmodified proteins. *J Biol Chem.* (2001) 276:35176–84. doi: 10.1074/jbc.M102223200
 86. Bitan G, Teplow DB. Rapid photochemical cross-linking—a new tool for studies of metastable, amyloidogenic protein assemblies. *Acc Chem Res.* (2004) 37:357–64. doi: 10.1021/ar000214l
 87. Bitan G. Structural study of metastable amyloidogenic protein oligomers by photo-induced cross-linking of unmodified proteins. *Methods Enzymol.* (2006) 413:217–36. doi: 10.1016/S0076-6879(06)13012-8
 88. Kyte J, Doolittle RF. A simple method for displaying the hydropathic character of a protein. *J Mol Biol.* (1982) 157:105–32. doi: 10.1016/0022-2836(82)90515-0
 89. Borreguero JM, Urbanc B, Lazo ND, Buldyrev SV, Teplow DB, Stanley HE. Folding events in the 21-30 region of amyloid beta-protein (Abeta) studied in silico. *Proc Natl Acad Sci USA.* (2005) 102:6015–20. doi: 10.1073/pnas.0502006102
 90. Yamin G, Coppola G, Teplow DB. Design, characterization, and use of a novel amyloid beta-protein control for assembly, neurotoxicity, and gene expression studies. *Biochemistry.* (2016) 55:5049–60. doi: 10.1021/acs.biochem.6b00579
 91. Fradinger EA, Maji S, Lazo ND, Teplow DB. Studying amyloid β -protein assembly. In: Xia W, Xu H, editors. *Amyloid Precursor Protein, A Practical Approach.* Boca Raton, FL: CRC Press (2005). p. 83–110.
 92. Finkelstein A, Kunis G, Seksenyan A, Ronen A, Berkutzki T, Azoulay D, et al. Abnormal changes in NKT cells, the IGF-1 axis, and liver pathology in an animal model of ALS. *PLoS ONE.* (2011) 6:e22374. doi: 10.1371/journal.pone.0022374
 93. Lund SA, Giachelli CM, Scatena M. The role of osteopontin in inflammatory processes. *J Cell Commun Signal.* (2009) 3:311–22. doi: 10.1007/s12079-009-0068-0
 94. Rittling SR. Osteopontin in macrophage function. *Expert Rev Mol Med.* (2011) 13:e15. doi: 10.1017/S1462399411001839
 95. Chan JL, Reeves TM, Phillips LL. Osteopontin expression in acute immune response mediates hippocampal synaptogenesis and adaptive outcome following cortical brain injury. *Exp Neurol.* (2014) 261:757–71. doi: 10.1016/j.expneurol.2014.08.015
 96. Zhao J, Li A, Rajsombath M, Dang Y, Selkoe DJ, Li S. Soluble abeta oligomers impair dipolar heterodendritic plasticity by activation of mGluR in the hippocampal CA1 region. *iScience.* (2018) 6:138–50. doi: 10.1016/j.isci.2018.07.018
 97. Busche MA, Eichhoff G, Adelsberger H, Abramowski D, Wiederhold KH, Haass C, et al. Clusters of hyperactive neurons near amyloid plaques in a mouse model of Alzheimer's disease. *Science.* (2008) 321:1686–9. doi: 10.1126/science.1162844
 98. Kellner V, Menkes-Caspi N, Beker S, Stern EA. Amyloid-beta alters ongoing neuronal activity and excitability in the frontal cortex. *Neurobiol Aging.* (2014) 35:1982–91. doi: 10.1016/j.neurobiolaging.2014.04.001
 99. Spangenberg EE, Green KN. Inflammation in Alzheimer's disease: lessons learned from microglia-depletion models. *Brain Behav Immun.* (2017) 61:1–11. doi: 10.1016/j.bbi.2016.07.003
 100. Spangenberg E, Severson PL, Hohsfield LA, Crapser J, Zhang J, Burton EA, et al. Sustained microglial depletion with CSF1R inhibitor impairs parenchymal plaque development in an Alzheimer's disease model. *Nat Commun.* (2019) 10:3758. doi: 10.1038/s41467-019-11674-z
 101. Ortega-Martinez S, Palla N, Zhang X, Lipman E, Sisodia SS. Deficits in enrichment-dependent neurogenesis and enhanced anxiety behaviors mediated by expression of Alzheimer's disease-linked Ps1 variants are rescued by microglial depletion. *J Neurosci.* (2019) 39:6766–80. doi: 10.1523/JNEUROSCI.0884-19.2019
 102. Parhizkar S, Arzberger T, Brendel M, Kleinberger G, Deussing M, Focke C, et al. Loss of TREM2 function increases amyloid seeding but reduces plaque-associated ApoE. *Nat Neurosci.* (2019) 22:191–204. doi: 10.1038/s41593-018-0296-9
 103. Minett T, Classey J, Matthews FE, Fahrenhold M, Taga M, Brayne C, et al. Microglial immunophenotype in dementia with Alzheimer's pathology. *J Neuroinflammation.* (2016) 13:135. doi: 10.1186/s12974-016-0601-z
 104. Simard AR, Soulet D, Gowing G, Julien JP, Rivest S. Bone marrow-derived microglia play a critical role in restricting senile plaque formation in Alzheimer's disease. *Neuron.* (2006) 49:489–502. doi: 10.1016/j.neuron.2006.01.022
 105. Yamada K, Hashimoto T, Yabuki C, Nagae Y, Tachikawa M, Strickland DK, et al. The low density lipoprotein receptor-related protein 1 mediates uptake of amyloid beta peptides in an *in vitro* model of the blood-brain barrier cells. *J Biol Chem.* (2008) 283:34554–62. doi: 10.1074/jbc.M801487200
 106. Zlokovic BV, Deane R, Sagare AP, Bell RD, Winkler EA. Low-density lipoprotein receptor-related protein-1: a serial clearance homeostatic mechanism controlling Alzheimer's amyloid beta-peptide elimination from the brain. *J Neurochem.* (2010) 115:1077–89. doi: 10.1111/j.1471-4159.2010.07002.x
 107. Malm TM, Koistinaho M, Parepalo M, Vatanen T, Ooka A, Karlsson S, et al. Bone-marrow-derived cells contribute to the recruitment of microglial cells in response to beta-amyloid deposition in APP/PS1 double transgenic Alzheimer mice. *Neurobiol Dis.* (2005) 18:134–42. doi: 10.1016/j.nbd.2004.09.009
 108. Guerreiro R, Wojtas A, Bras J, Carrasquillo M, Rogava E, Majounie E, et al. TREM2 variants in Alzheimer's disease. *N Engl J Med.* (2013) 368:117–27. doi: 10.1056/NEJMoa1211851
 109. Jin SC, Benitez BA, Karch CM, Cooper B, Skorupa T, Carrell D, et al. Coding variants in TREM2 increase risk for Alzheimer's disease. *Hum Mol Genet.* (2014) 23:5838–46. doi: 10.1093/hmg/ddu277
 110. Finelli D, Rollinson S, Harris J, Jones M, Richardson A, Gerhard A, et al. TREM2 analysis and increased risk of Alzheimer's disease. *Neurobiol Aging.* (2015) 36:546. e549–13. doi: 10.1016/j.neurobiolaging.2014.08.001

111. Prokop S, Miller KR, Drost N, Handrick S, Mathur V, Luo J, et al. Impact of peripheral myeloid cells on amyloid-beta pathology in Alzheimer's disease-like mice. *J Exp Med.* (2015) 212:1811–8. doi: 10.1084/jem.20150479
112. Varvel NH, Grathwohl SA, Degenhardt K, Resch C, Bosch A, Jucker M, et al. Replacement of brain-resident myeloid cells does not alter cerebral amyloid-beta deposition in mouse models of Alzheimer's disease. *J Exp Med.* (2015) 212:1803–9. doi: 10.1084/jem.20150478
113. Chu T, Tran T, Yang F, Beech W, Cole GM, Frautschy SA. Effect of chloroquine and leupeptin on intracellular accumulation of amyloid-beta (A beta) 1-42 peptide in a murine N9 microglial cell line. *FEBS Lett.* (1998) 436:439–44. doi: 10.1016/S0014-5793(98)01161-2

Conflict of Interest: The authors declare that the research was conducted in the absence of any commercial or financial relationships that could be construed as a potential conflict of interest.

Copyright © 2020 Li, Hayden, Garcia, Fuchs, Sheyn, Daley, Rentsendorj, Torbati, Black, Rutishauser, Teplow, Koronyo and Koronyo-Hamaoui. This is an open-access article distributed under the terms of the Creative Commons Attribution License (CC BY). The use, distribution or reproduction in other forums is permitted, provided the original author(s) and the copyright owner(s) are credited and that the original publication in this journal is cited, in accordance with accepted academic practice. No use, distribution or reproduction is permitted which does not comply with these terms.