

# Anti-N-Terminal A $\beta$ Mab 3A1 Preferentially Recognizes A $\beta$ Aggregates and Does Not Cross-React with APP

Jeffrey L. Frost<sup>1</sup>, James S. Foster<sup>2</sup>, Peggy Taylor<sup>1</sup>, Brian O'Nuallain<sup>1</sup>

<sup>1</sup>BioLegend, San Diego, CA 92121, and <sup>2</sup>University of Tennessee Graduate School of Medicine, Knoxville, TN 37920, USA

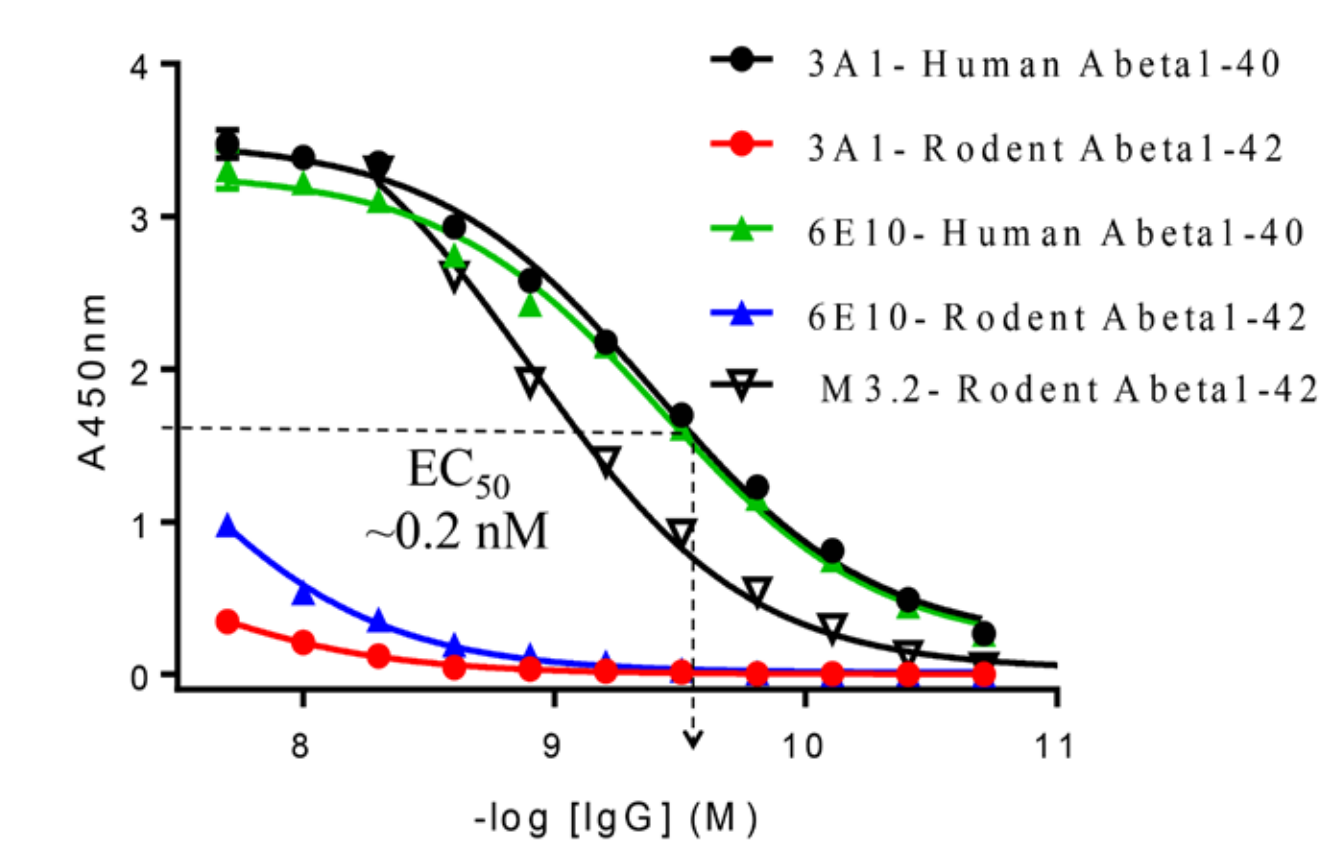
## Introduction

A $\beta$ 's N-terminus is highly immunogenic and antibodies that target this region are standard reagents for Alzheimer's disease (AD) research, and have been utilized as investigational AD therapeutics. A murine anti-N-terminal A $\beta$  monoclonal antibody (mAb), 3A1, was generated against dityrosine cross-linked A $\beta$ 1-40 protein species (CAPS), and its epitope mapped to the peptide's first 15 amino acids. 3A1 has demonstrated activity *in vivo* by decreasing plaque burden and increasing the levels of plasma A $\beta$  in an APPsw/PS1 $\Delta$ E9 transgenic mouse model of AD [Frost *et al.* (2015) *Neurobiol Aging*. 36(12): 3187]. To better understand the mAb's specificity for A $\beta$  we established *in vitro* the antibody's ability to recognize A $\beta$ 1-40 conformers (monomers, dimers, protofibrils), APP, and rodent A $\beta$ .

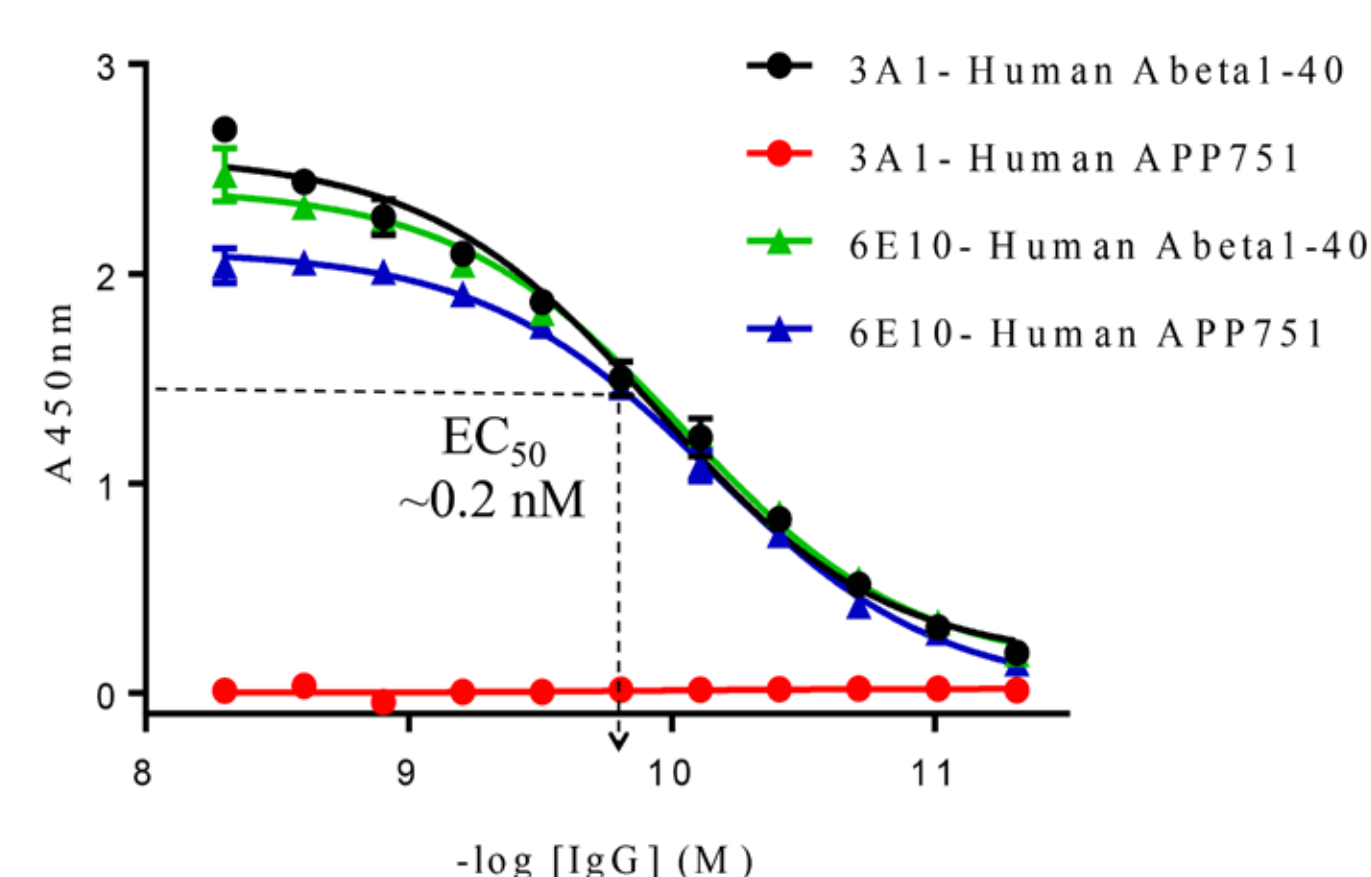
## Methods

3A1's avidity for A $\beta$  conformers, APP and rodent A $\beta$  were determined using ELISA, and Western blot assays. Anti-N-terminal A $\beta$  mAb, 6E10 [Kim *et al.* (1988) *Neurosci. Res. Comm.* 7:113], was used as a positive control. A $\beta$  conformers were generated as previously described [Welzel *et al.* (2012) *PLoS One* 7(11):e50317].

## Mab 3A1 Recognized Human A $\beta$ and Did Not Bind to Rodent A $\beta$ or to Human APP



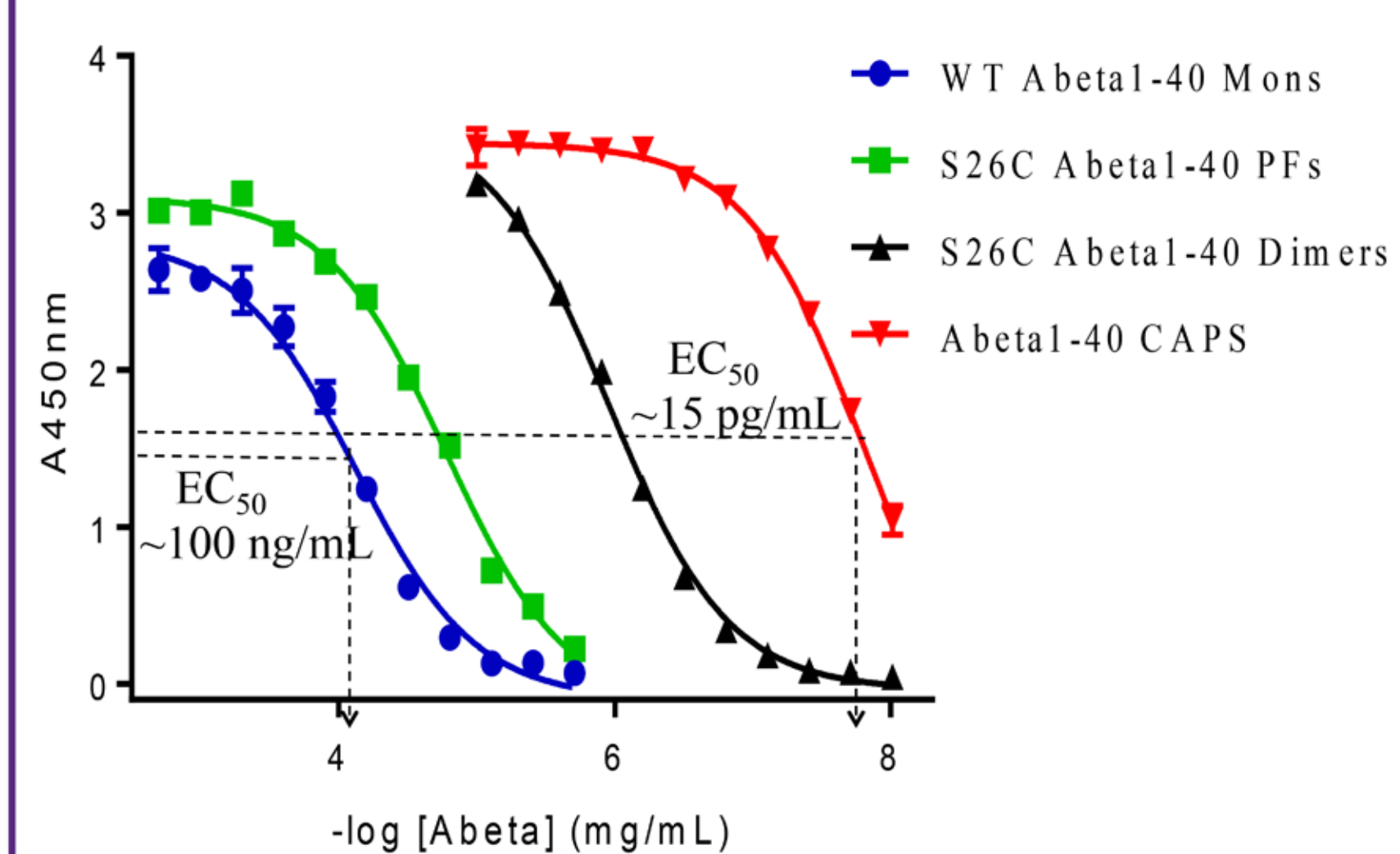
Both anti-N-terminal A $\beta$  mAbs, 3A1 and 6E10, bound to plate-immobilized human A $\beta$ 1-40 with EC<sub>50</sub> of ~0.2 nM. The mAbs had low to no binding to rodent A $\beta$ .



Mab 3A1 did not bind to plate-immobilized human recombinant APP751 protein. Mab 6E10 bound similarly to A $\beta$ 1-40 and APP751 with EC<sub>50</sub>'s of ~0.2 nM.

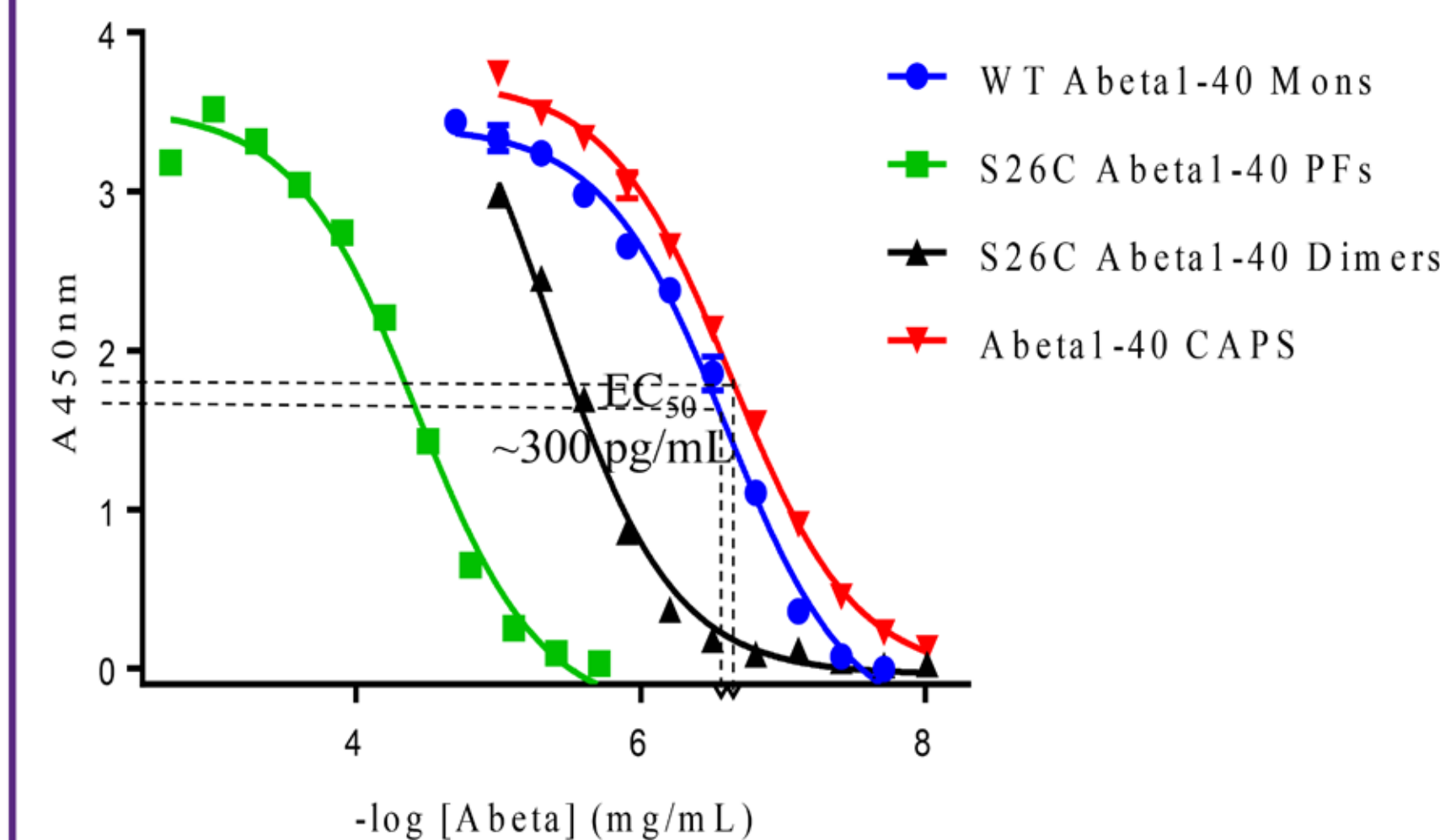
## Mab 3A1 Preferentially Captured Aggregated A $\beta$

### 3A1 Capture: HRP-4G8 Detection



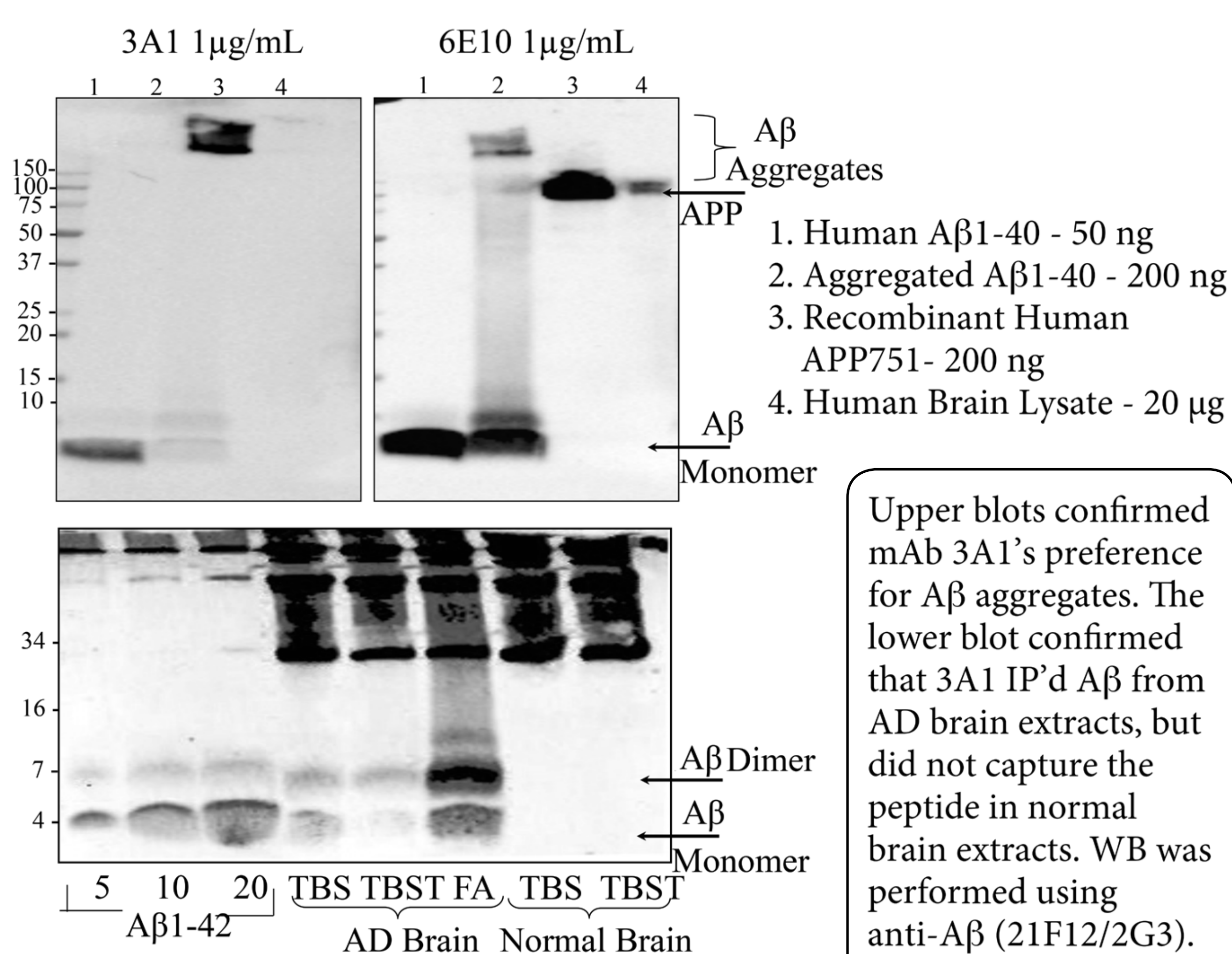
Mab 3A1 preferentially recognized A $\beta$  aggregates compared with the monomeric peptide, with ~700-fold stronger binding to dityrosine cross-linked A $\beta$ 1-40

### 6E10 Capture: HRP-4G8 Detection



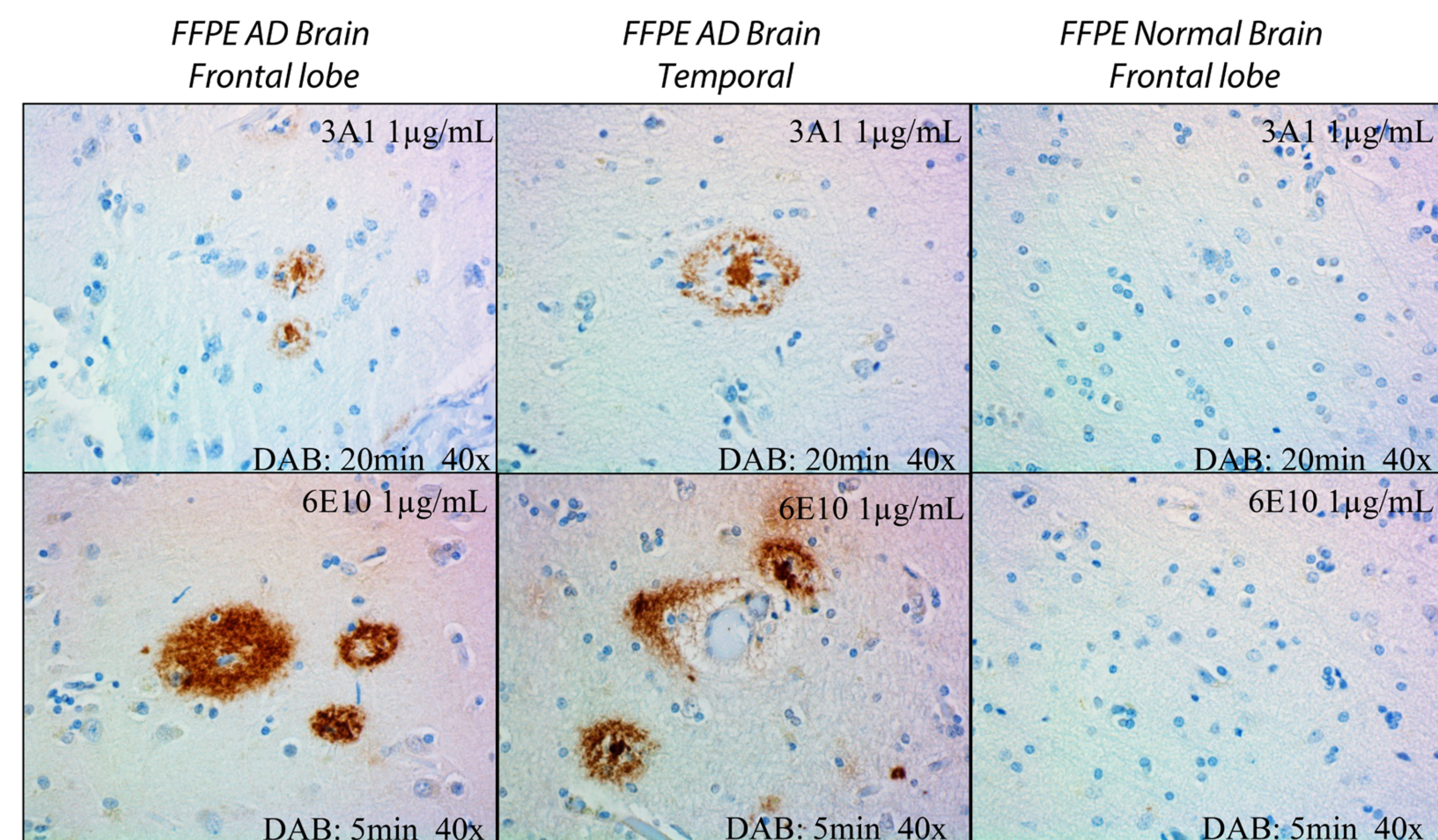
Mab 6E10 similarly bound to A $\beta$  monomers and CAPS, but bound significantly weaker to S26C A $\beta$  dimers and proto-fibrils (PFs).

## Mab 3A1 Preferentially Detected Aggregated A $\beta$ and Did Not Recognize APP



Upper blots confirmed mAb 3A1's preference for A $\beta$  aggregates. The lower blot confirmed that 3A1 IP'd A $\beta$  from AD brain extracts, but did not capture the peptide in normal brain extracts. WB was performed using anti-A $\beta$  (21F12/2G3).

## 3A1 Stained A $\beta$ Plaques in AD Brain Tissues



IHC was performed using 88% formic acid antigen retrieval, and by incubating the primary antibodies overnight at 4°C.

## Summary & Conclusions

Mab 3A1 is a novel anti-N-terminal A $\beta$  antibody that specifically detected human A $\beta$  conformers in ELISA, IP/WB, and IHC applications.

The antibody demonstrated up to a ~700-fold preference for aggregated compared with monomeric A $\beta$  in Capture/Sandwich ELISA. In contrast, another anti-N-terminal A $\beta$  antibody, 6E10, did not preferably capture A $\beta$  aggregates.

3A1's utility as an antibody research tool for AD was further demonstrated by its ability to avidly bind to human A $\beta$  conformers without appreciably cross-reacting with human APP or rodent A $\beta$ .

## Acknowledgement

We would like to thank the University of Tennessee Graduate School of Medicine for their contributions to this poster.

