

# Amyloid $\beta$ A4 (CT, 1-42) monoclonal antibody (8G7)

The  $\beta$ -amyloid peptide ( $\beta$  A4), proteolytically released from the amyloid precursor protein (APP), provides the principal component of senile plaques in Alzheimer's disease. Cleavage of APP by  $\alpha$ -secretase or alternatively by  $\beta$ -secretase leads to generation and extracellular release of soluble APP peptides, S-APP- $\alpha$  and S-APP- $\beta$ , respectively, and the retention of corresponding membrane-anchored C-terminal fragments, C83 and C99. Subsequent processing of C83 by  $\gamma$ -secretase yields P3 peptides. This is the major secretory pathway and is nonamyloidogenic. Alternatively, presenilin/ nicastrin-mediated  $\gamma$ -secretase processing of C99 releases the amyloid  $\beta$  proteins, amyloid- $\beta$  40 ( $A\beta$ 40) and amyloid- $\beta$  42 ( $A\beta$ 42), major components of amyloid plaques, and the cytotoxic C-terminal fragments,  $\gamma$ -CTF(50),  $\gamma$ -CTF(57) and  $\gamma$ -CTF(59).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

[View Online »](#)

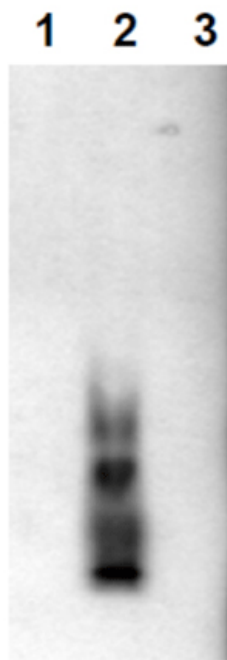
## Ordering Information

[Order Online »](#)

ADI-905-804-100	100 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of (1)  $\beta$  A4 (1-40), (2)  $\beta$  A4 (1-42), and (3)  $\beta$  A4 (1-43), probed with 2 $\mu$ g/ml of Amyloid  $\beta$  A4 (42) mAb (8G7).

## Handling & Storage

**Use/Stability** Stable at -80°C up to 1 year, at 4°C up to 3 months.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** ELISA, ICC, WB

**Application Notes** Detects a band of ~4kDa by Western blot.

**Clone** 8G7

**Crossreactivity** Does not cross react with  $\beta$  A4 (1-40).

**Formulation** Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.

**Host** Mouse

**Immunogen** Synthetic peptide corresponding to aa 1-42 of amyloid  $\beta$  A4.

**Isotype** IgG1

**Purity Detail** Thiophilic adsorption and size exclusion chromatography purified.

**Recommendation Dilutions/Conditions** ELISA (0.05 $\mu$ g/ml)Immunocytochemistry (0.1-1.0 $\mu$ g/ml)Western Blot (1 $\mu$ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Reconstitution** Reconstitute with 1ml water (15 minutes at room temperature).

**Species Reactivity** Human

## Technical Info / Product Notes

UniProt ID

Worry-free Guarantee

Last modified: May 29, 2024

## Application Note:

[Analysis of the Effect of Aggregated  \$\beta\$ -Amyloid on Cellular Signaling Pathways Critical for Memory in Alzheimer's disease](#)

P05067

This antibody is covered by our [Worry-Free Guarantee](#)

.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)