

PRODUCT INFORMATION



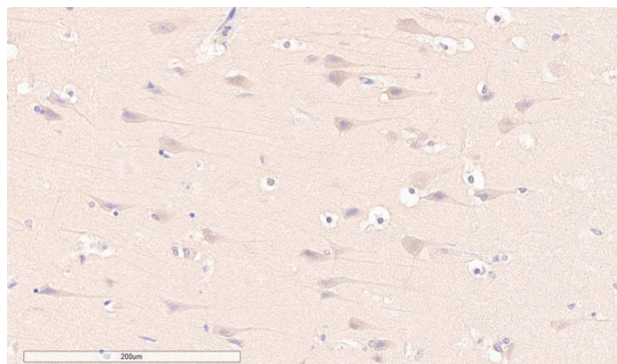
APP (C99 Fragment) Monoclonal Antibody (Clone 8G4)

Item No. 28636

Overview and Properties

Contents:	This vial contains 100 µg of protein G-purified monoclonal antibody.
Synonyms:	ABPP, Alzheimer's Disease Amyloid Protein, Amyloid-β A4 Protein, Amyloid Precursor Protein, APPI, Cerebral Vascular Amyloid Peptide, β C-terminal Fragment, βCTF, CVAP, PN-II, PreA4, Protease Nexin-II
Immunogen:	Peptide corresponding to the C99 fragment of human APP
Cross Reactivity:	(+) APP, C99 fragment, C47 fragment
Species Reactivity:	(+) Human; other species not tested
Uniprot No.:	P05067
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Clone:	8G4
Host:	Mouse
Isotype:	IgG2a
Applications:	ELISA and Immunohistochemistry (IHC); the recommended starting dilution is 1:1,000 for ELISA and 1:200 for IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human Alzheimer's disease brain tissue after heat-induced antigen retrieval in pH 6.0 citrate buffer. After incubation with APP (C99 Fragment) Monoclonal Antibody (Clone 8G4) at a dilution of 1:200, slides were incubated with a biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/07/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Amyloid precursor protein (APP) is a type I transmembrane protein that has a central role in the pathogenesis of Alzheimer's disease, as well as additional roles in brain development, neuronal plasticity, and memory.¹ APP is cleaved by β -secretase (BACE) in neuronal endosomes during amyloidogenic processing of APP, generating the C-terminal C99 fragment, which is localized to the endoplasmic membrane.^{2,3} C99 is further cleaved by γ -secretase, liberating the APP intracellular domain (AICD) and generating amyloid- β (A β) peptides of various lengths, including A β 40 (Item No. 21617) and A β 42 (Item No. 20574), which are hallmarks of Alzheimer's disease. APP can also be cleaved by α -secretase during non-amyloidogenic processing of APP, which occurs at the neuronal plasma membrane and generates the neuroprotective soluble APP fragment sAPP α , as well as a variety of other fragments, including the C47 fragment.^{1,4} Transgenic mice expressing mutant forms of APP exhibit extracellular A β deposits in the brain, as well as cognitive dysfunction, and are widely used models of Alzheimer's disease.⁵ Intraneuronal C99 levels are increased in the transgenic APP^{E693Q} mouse model of Alzheimer's disease, as well as postmortem frontal cortex from patients with sporadic Alzheimer's disease.^{3,6} Cayman's APP (C99 Fragment) Monoclonal Antibody (Clone 8G4) can be used for ELISA and immunohistochemistry (IHC) applications. The antibody recognizes the C-terminal region corresponding to the C99 fragment to detect intact APP, as well as the C47 APP fragment.

References

1. Müller, U.C., Deller, T., and Korte, M. Not just amyloid: Physiological functions of the amyloid precursor protein family. *Nat. Rev. Neurosci.* **18**(5), 281-298 (2017).
2. Zhao, J., Liu, X., Xia, W., et al. Targeting amyloidogenic processing of APP in Alzheimer's disease. *Front. Mol. Neurosci.* **13**, 137 (2020).
3. Lauritzen, I., Pardossi-Piquard, R., Bourgeois, A., et al. Does intraneuronal accumulation of carboxyl-terminal fragments of the amyloid precursor protein trigger early neurotoxicity in Alzheimer's disease? *Curr. Alzheimer Res.* **16**(5), 453-457 (2019).
4. Schrader-Fischer, G., Staufienbiel, M., and Paganetti, P.A. Insertion of lysosomal targeting sequences to the amyloid precursor protein reduces secretion of β A4. *J. Neurochem.* **68**(4), 1571-1580 (1997).
5. Sasaguri, H., Nilsson, P., Hashimoto, S., et al. APP mouse models for Alzheimer's disease preclinical studies. *EMBO J.* **36**(17), 2473-2487 (2017).
6. Pulina, M.V., Hopkins, M., Haroutunian, V., et al. C99 selectively accumulates in vulnerable neurons in Alzheimer's disease. *Alzheimers Dement.* **16**(2), 273-282 (2020).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM