

Presenilin 1 monoclonal antibody (APS 18)

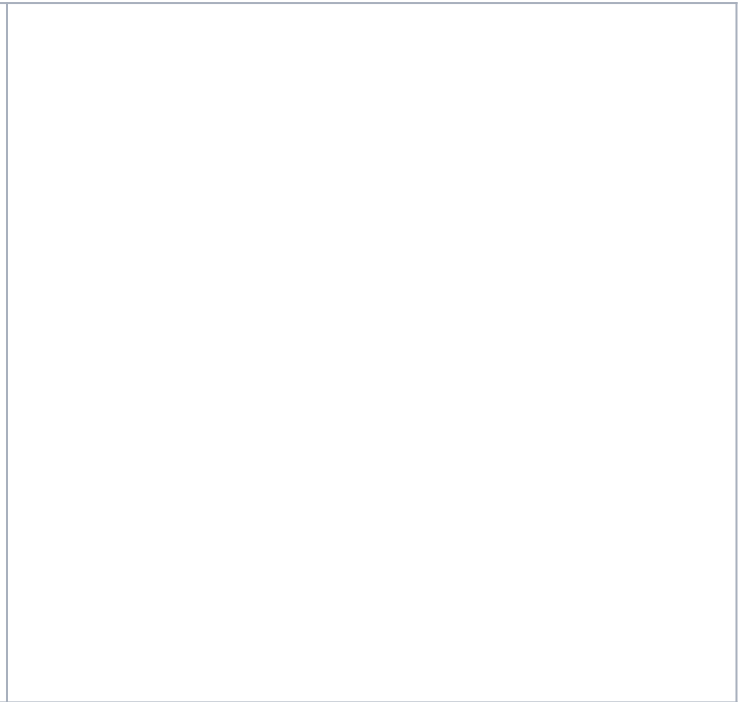
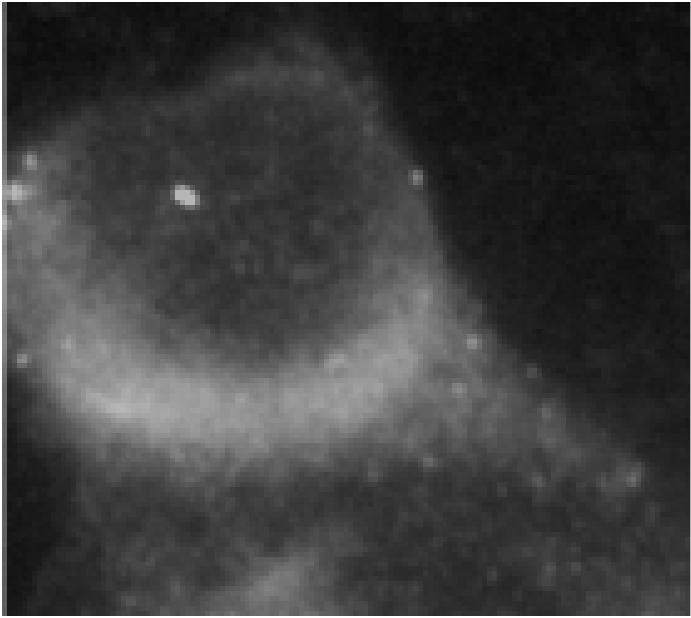
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling Avoid freeze/thaw cycles.

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 ~18kDa (C-terminal fragment) and ~46kDa (full-length) presenilin 1 by Western blot.

Clone APS 18

Crossreactivity Does not cross-react with presenilin 2.

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 313-334 (S313KYNAESTERESQDTVAENDDG334) of C-terminal human presenilin 1.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions ELISA (1µg/ml)Immunocytochemistry (6µg/ml)Western Blot (12µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Monkey, Mouse, Rat

UniProt ID P49768

Last modified: May 29, 2024



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)