

Presenilin 1 monoclonal antibody (APS 18)

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

Citations: 3

[View Online »](#)

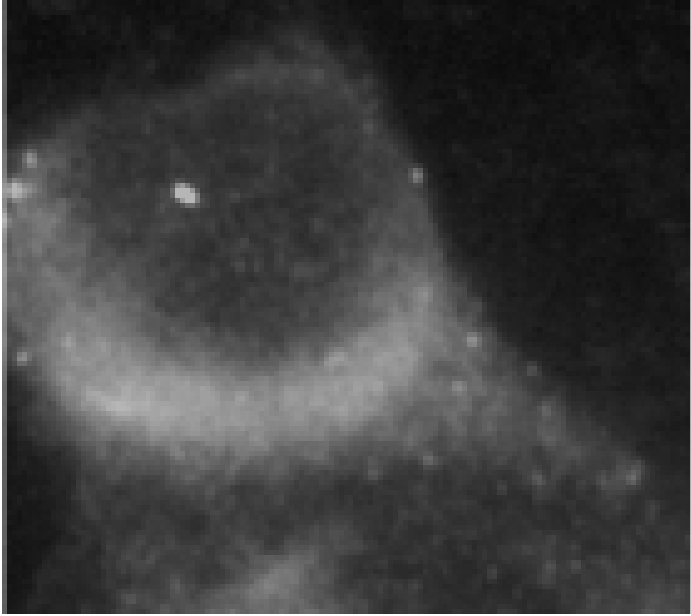
Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ALX-804-555-C200 | 200µg |
|------------------|-------|

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 |

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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