

Presenilin 1 monoclonal antibody (APS 11)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-554-C200	200µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

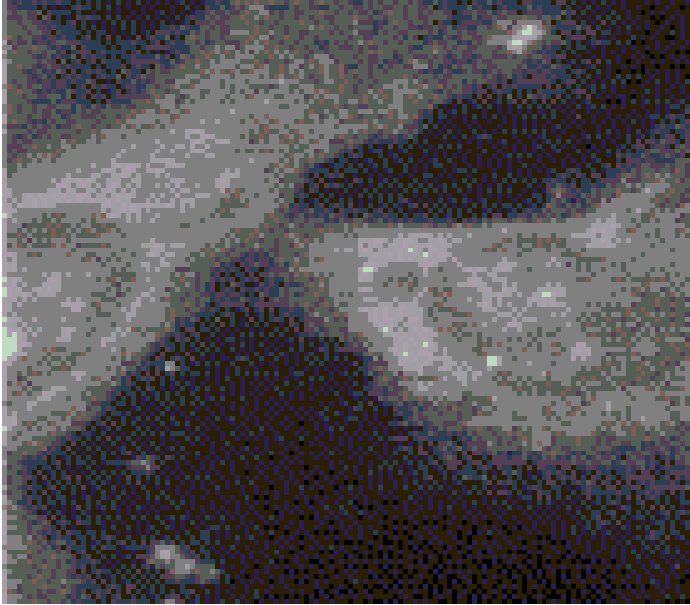


Figure: Immunocytochemistry on mouse fibroblast cell line stained with MAb to presenilin 1 (APS 11) (Prod. No. ALX-804-554).

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, IHC (PS), WB
Application Notes	Detects a band of ~28kDa by Western blot.
Clone	APS 11
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 21-34 (CH21LSNTVRSQNDNRE34) of N-terminal human presenilin 1.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	ELISA (4.5µg/ml)Immunocytochemistry (0.75µg/ml)Immunohistochemistry (paraffin sections, 0.75µg/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, Mouse, Primate, 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](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)