

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 »](#)

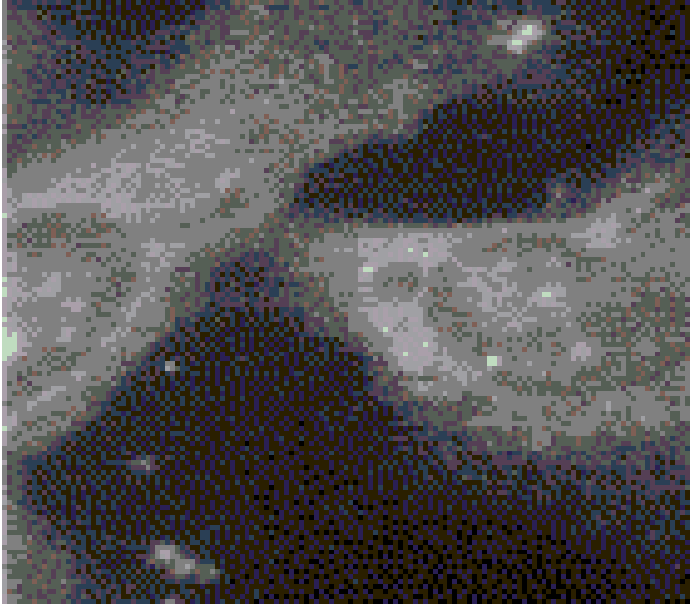


Figure: Immunocytochemistry on mouse fibroblast cell line stained with MAb to presenilin 1 (APS 11) (Prod. No. ALX-804-554).

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 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

ENZO LIFE SCIENCES

INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

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 1 72 44 01 555
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