

APP monoclonal antibody (1G6)

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

Manuals, SDS & CofA [View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Amyloid precursor protein, Amyloid A4 protein
Application	ELISA, IHC (PS), WB
Clone	1G6
Formulation	Liquid. In PBS. Does not contain any preservatives or carrier proteins.
Host	Mouse
Isotype	IgG2b
Purity Detail	Purified.
Species Reactivity	Human
Specificity	Epitope lies within amino acids 573-596 of amyloid precursor protein.
UniProt ID	P05067
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-fr@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