

ApoE4 monoclonal antibody (4E4)

Apolipoproteins are plasma lipoproteins that function as transporters of lipids and cholesterol in the circulatory system. Human ApoE is an apolipoprotein that has three isoforms: ApoE2, ApoE3, and ApoE4.

The ApoE4 variant is the largest known genetic risk factor for late-onset Alzheimer's disease, and while the exact mechanism remains to be determined, it appears there may be an interaction with beta-amyloid. Alzheimer's disease is characterized by build-ups of aggregates of the peptide beta-amyloid. Apolipoprotein E enhances proteolytic break-down of this peptide, both within and between cells. The isoform ApoE4 is not as effective as the others at promoting these reactions, resulting in increased vulnerability to AD in individuals with that gene variation.

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

Ordering Information

[Order Online »](#)

ENZ-ABS295-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

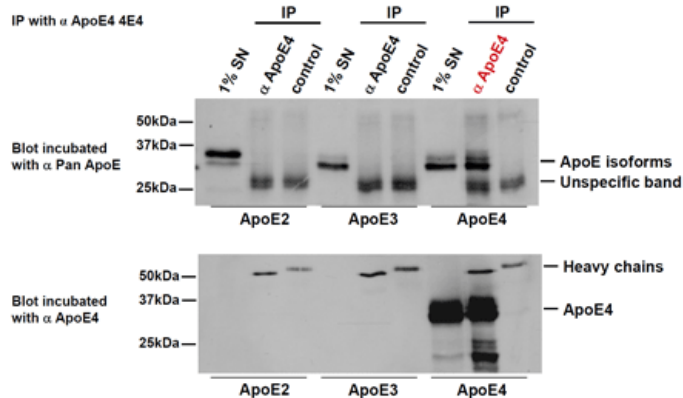


Figure 2. Immunoprecipitation. ApoE4, mAb (4E4) immunoprecipitates specifically ApoE4 but not ApoE2 or ApoE3.

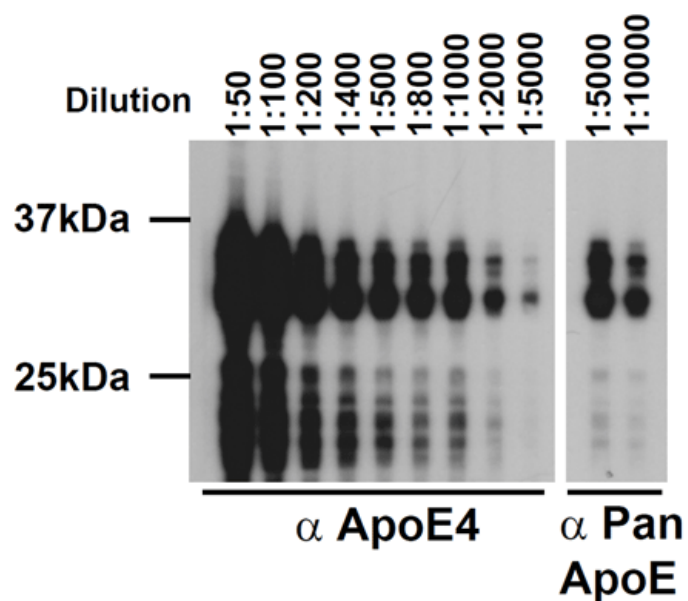


Figure 1. Western blot of ApoE4, mAb (4E4) (Prod. No. ENZ-ABS295). α ApoE4 4E4: Mouse monoclonal antibody 4E4 specific for ApoE4, cell culture supernatant.
 α Pan ApoE: Rabbit polyclonal serum recognizing all three ApoE proteins.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Apolipoprotein E4

Application ELISA, IP, WB

Clone 4E4

Crossreactivity Does not cross-react with ApoE2 or ApoE3.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to residues surrounding Arg110 of human ApoE4 protein.

Isotype IgG1κ

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Human

Specificity Recognizes endogenous levels of total ApoE4 protein.

UniProt ID P02649

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)