

APLP2 (human), (recombinant) (His- tag)

APLP2 is a member of an evolutionarily conserved APP family. APLP2 functions as a novel G(0)-protein-coupled receptor for Müllerian inhibiting substance (MIS) in cell survival. APLP2 is required for proper cell cycle exit of neuronal progenitors, and has a distinct role in priming cortical progenitors for neuronal differentiation. APLP2, a closely related homolog to AbetaPP, share overlapping anticoagulant functions with regard to regulating thrombosis after cerebral vascular injury.

Ordering Information

[Order Online »](#)

ENZ-PRT332-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa M

116

66.2

45.0

35.0

25.0

18.4

14.4



As a result of glycosylation, the recombinant protein migrates as an approximately 107.6 kDa protein in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Amyloid beta precursor like protein 2

Formulation Lyophilized from sterile PBS, pH 7.4.

MW ~107.6 kDa (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

Source Produced in HEK293 cells. A DNA sequence encoding the human APLP2 (Met1-Ser692) was expressed with a polyhistidine tag at the C-terminus.

UniProt ID Q06481-1



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)