

RAP (human), (recombinant)

RAP functions in ER maturation of lipoprotein receptor family members by acting as an antagonist. It blocks binding of PCSK9 to lipoprotein receptors and of all known ligands (including MMP-9 and apoE) to lipoprotein receptor-related protein 1 (LRP-1). Modulates amyloid deposition in mouse models of Alzheimer's disease. Reduces endocytosis of prion protein (PrP^C) by neuronal cells.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE552-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	LDL receptor associated protein, LRPAP1, LDL receptor-related protein-associated protein 1
Application Notes	Useful tool for binding studies, and testing <i>in vivo</i> and in cell culture.
Formulation	Liquid. In 0.0067M Phosphate, pH 7.4 containing 0.15M sodium chloride.
MW	~39kDa (SDS-PAGE)
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	P30533

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-
fr@enzolifesciences.com](mailto:info-fr@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)