

Omentin (rat), (recombinant)

Omentin (also named intelectin-1, endothelial lectin HL-1 and intestinal lactoferrin receptor) has been identified as a visceral fat (or omental fat) depot-specific adipokine. Omentin is a secreted protein of ~38 kDa containing a fibrinogen-related domain. Omentin is highly expressed in the omental fat, and much less in intestine, lung and heart. Addition of recombinant omentin *in vitro* did not affect basal, but enhanced insulin-stimulated glucose uptake in subcutaneous as well as in omental human adipocytes. Omentin triggers AKT signalling in the absence of insulin.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-132-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Intelectin-1, Endothelial lectin HL-1, Intestinal lactoferrin receptor, Galactofuranose-binding lectin
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Associates of CAPE COD Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~38-40kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile water.
Source	Produced in HEK 293 cells. Rat omentin (aa 20-313) is fused at the N-terminus to a FLAG [®] -tag.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	Q499T8



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