

Omentin 1 (human), (recombinant)

Has no effect on basal glucose uptake but enhances insulin-stimulated glucose uptake in adipocytes. Increases AKT phosphorylation in the absence and presence of insulin. May play a role in the defense system against microorganisms. May specifically recognize carbohydrate chains of pathogens and bacterial components containing galactofuranosyl residues, in a calcium-dependent manner. May be involved in iron metabolism.

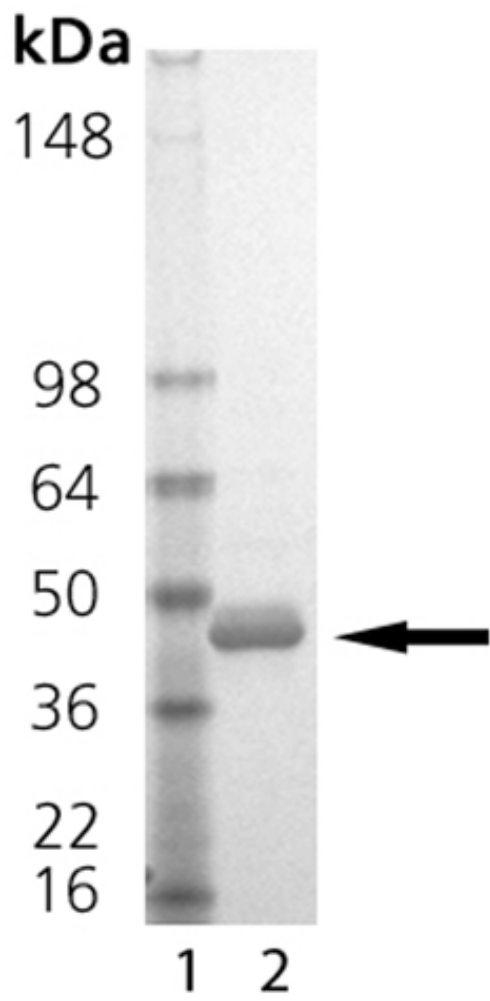
Omentin 1 (also named intelectin-1, endothelial lectin HL-1 and intestinal lactoferrin receptor) has been identified as a visceral fat (or omental fat) depot-specific adipokine. Omentin 1 is a secreted protein of ~38kDa containing a fibrinogen-related domain. Omentin 1 is highly expressed in the omental 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 1 triggers AKT signalling in the absence of insulin.

Citations: 5

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 1 μ g Omentin-1.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
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).
Formulation	Lyophilized. Contains PBS.
MW	~40kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100μl sterile water. PBS containing at least 0.1% BSA should be used for further dilutions.
Source	Produced in HEK 293 cells. Human omentin 1 (intelectin-1) (aa 17-313) 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	Q8WWA0



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: **May 29, 2024**
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com