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 insulinstimulated 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 »

Ordering Information

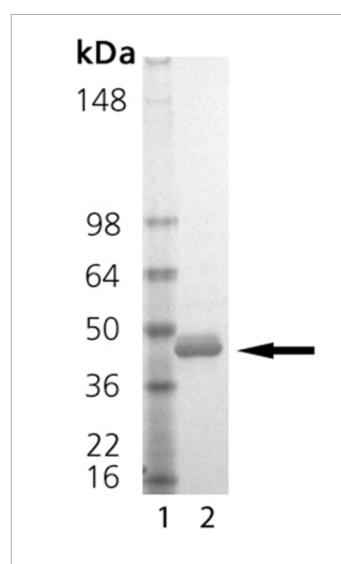
Order Online »

ALX-522-108-C010

10µg

Manuals, SDS & CofA

View Online »



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: $1\mu g$ 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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

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

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