## 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 insulinstimulated 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<sup>®</sup>-tag.

**Technical Info / Product Notes** FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID Q499T8



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