

Leptin (human), (recombinant) (>95%)

Leptin (obese gene product) is a ligand for the OB receptor.

Ordering Information

[Order Online »](#)

BML-SE161-0001	1mg
BML-SE161-0005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name OB gene product

CAS 177404-21-6

Formulation Lyophilized.

MW 16 kDa

Purity ≥95% (HPLC & SDS-PAGE)

Purity Detail Highly purified and biologically active.

Source Produced in *E. coli*.

UniProt ID P41159

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