

Leptin, Pegylated (mouse), (recombinant)

Ordering Information

[Order Online »](#)

RLT-500-047	1mg
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Biological Activity	Recombinant mouse pegylated leptin is capable of stimulating proliferation of BAF/3 cells stably transfected with the long form of human leptin receptor. Its in vitro activity is only slightly lower than the non-pegylated recombinant mouse leptin but in vivo it has profound weight reducing effect (as compared to the non-pegylated recombinant mouse leptin), resulting mainly from reduced food intake.
Endotoxin Content	< 0.1 ng/μg of protein (
Formulation	Lyophilized.
MW	35.6 KkDa
Purity	> 98.0% as determined by Gel filtration and SDS-PAGE gel.
Source	Produced in <i>E. coli</i> .
UniProt ID	P41160



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)