

Recombinant Human Leptin

Leptin is an adipose-derived cytokine hormone encoded by the ob (obese) gene. The actions of leptin involve reducing body weight, blood glucose levels, and food consumption as well as regulating energy balance, suppressing the appetite, increasing thermogenesis.

Materials Provided

Catalog Number:	92-1327
Quantity Supplied:	1 vial(s), 200 µg per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	16,000 g/mol
Source:	Recombinant
Purity:	>98%
Endotoxin Level:	<1.0 EU per 1 µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	1 mg/mL
Reconstitution Volume:	0.200 ml
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com