

Ghrelin, Human

Ghrelin is a neuropeptide GI hormone that regulates hunger and plays a role in energy use and distribution. It is secreted in the GI tract when the stomach is empty and stimulates gastric acid secretion and GI motility to prepare the body for food consumption. Acts in opposition to the "satiety hormone", leptin, and is an agonist for the same receptor in the brain, called the growth hormone secretagogue type 1A receptor, or GHSR.

Materials Provided

Catalog Number:	92-1345
Quantity Supplied:	1 vial(s), 0.5 mg per vial

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:	3370.9 g/mol
Source:	Synthetic
Purity:	96.2% (HPLC)
Endotoxin Level:	N/A
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:	250 μM
Reconstitution Volume:	0.594 mL
Reconstitution Solvent:	5% acetic acid

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