

Gastrin-releasing peptide

Gastrin-releasing peptide (GRP) is a 27 amino acid regulatory neuropeptide that is an agonist for gastrin-releasing peptide receptor and is implicated in both physiological and pathophysiological processes, including release of gastrin from G cells of the stomach and the regulation of gastric acid secretion and enteric motor function. It also has a role in the circadian system, regulating light signaling to the master circadian oscillator in the hypothalamus. Negative feedback regulator of fear, and may also mediate stress.

Materials Provided

Catalog Number:	92-1057
Quantity Supplied:	1 vial(s), 0.5 mg 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:	2859.4 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	0.700 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