

Growth Hormone Releasing Factor (1-29)

Growth Hormone Releasing Factor (GRF or GHRF) is a preproprotein produced in the hypothalamus that acts on the adenohypophyse to stimulate growth hormone secretion from the pituitary gland. Antagonists of these receptors can inhibit tumor growth. GRF levels are regulated by somatostatin, growth hormone, and insulin-like growth factor 1.

Materials Provided

Catalog Number:	92-1058
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:	3357.9 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:	1 mM
Reconstitution Volume:	0.149 mL
Reconstitution Solvent:	H ₂ O

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