

RFRP-3 (Human)

RFRP-3 is an inhibitory hormone that is a homolog of the gonadotropin-inhibitory hormone (GnIH) and blocks release of gonadotropin from GnRH in the pituitary. Inhibits cAMP accumulation and progesterone production from FSH, LH, and forskolin in granulosa-lutein cells. Agonist of the G protein-coupled receptor 147 (GPR147 or NPFF1, IC₅₀=0.7nM).

Materials Provided

Catalog Number:	92-1198
Quantity Supplied:	1 vial(s), 500 µg 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:	969.14 g/mol
Source:	Synthetic
Purity:	≥ 97% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°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:	0.516 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