

Kisspeptin-10/Metastin

Kisspeptin-10 is a protein and potent endogenous ligand for the kisspeptin receptor, GPR54 ($K_i=2.33\text{nM}$). The kisspeptin-GPR54 signaling assists in gonadotropin-releasing hormone secretion at puberty. It is the smallest fully bioactive fragment of kisspeptin, and was originally identified as a metastasis suppressor for melanoma and breast cancer.

Materials Provided

Catalog Number:	92-1121
Quantity Supplied:	1 vial(s), 1 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:	1302.44 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:	0.768 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