

INSL 5, Murine

Insulin-like peptide 5 (INSL5) is a member of the insulin/relaxin superfamily that is a potent agonist for the G protein-coupled receptor RXFP4. Possible roles include gut contractility, thymic development and regulation, food intake, and glucose homeostasis.

Materials Provided

Catalog Number:	92-1228
Quantity Supplied:	1 vial(s), 20 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:	5117.04 g/mol
Source:	Synthetic
Purity:	≥ 90% 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:	25 µM
Reconstitution Volume:	0.156 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