

Recombinant Human FGF-21

Fibroblast growth factor 21 (FGF-21) is a protein involved in many mitogenic and cell survival activities and biological processes, such as embryonic development, cell growth, morphogenesis, tissue repair, and tumor growth and invasion. Agonist of the glucose transporter SLC2A1/GLUT1 activity, stimulating glucose uptake in adipocytes.

Materials Provided

Catalog Number:	92-1231
Quantity Supplied:	1 vial(s), 5 µg 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:	19500 g/mol
Source:	Recombinant
Purity:	≥ 90% by SDS-PAGE
Endotoxin Level:	< 1EU/µg
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:	100 µg/mL
Reconstitution Volume:	0.050 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