

Recombinant Human FGF-19

Fibroblast growth factor-19 (FGF-19 or FGF-J) protein functions as a hormone, regulating bile acid synthesis, with effects on glucose and lipid metabolism. FGF-19 also plays a role in prenatal development and postnatal growth and regeneration of a variety of tissues, by promoting cellular proliferation and differentiation.

Materials Provided

Catalog Number:	92-1301
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:	21800 g/mol
Source:	Recombinant
Purity:	≥ 95%
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:	0.05 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