

Recombinant Human FGF-18

Fibroblast growth factor-18 (FGF-18) is a heparin-binding growth factor protein that possess broad mitogenic and cell survival activities, and is involved in a variety of biological processes, including embryonic development (prenatal development and postnatal growth), cell growth, morphogenesis, tissue repair, tumor growth, and invasion.

Materials Provided

Catalog Number:	92-1318
Quantity Supplied:	1 vial(s), 50 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:	20100 g/mol
Source:	Recombinant
Purity:	>95%
Endotoxin Level:	<0.10 EU
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:	500 µg/mL
Reconstitution Volume:	0.100 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