

Recombinant Human Artemin

Artemin is a neurotrophic factor in the glial cell line-derived neurotrophic factor family of ligands which are a group of ligands within the TGF-beta superfamily of signaling molecules. This protein is a ligand for the RET receptor and uses GFR-alpha 3 as a co-receptor.

Materials Provided

Catalog Number:	92-1346
Quantity Supplied:	1 vial(s), 5 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:	24,200 g/mol
Source:	Recombinant
Purity:	≥ 98% by SDS-PAGE gel and HPLC analyses
Endotoxin Level:	< 0.1 ng/µg of protein
Formulation:	Lyophilized
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:	4 mM HCl

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