

Recombinant Human Ephrin-A3

Ephrin-A3 is a membrane protein tyrosine kinase and member of the ephrin (EPH) family. Implicated in cell migration, repulsion, and adhesion during neuronal, vascular, and epithelial development. Potent agonist of the EPH receptors.

Materials Provided

Catalog Number:	92-1142
Quantity Supplied:	1 vial(s), 200 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:	47700 g/mol
Source:	Recombinant
Purity:	≥ 90% by SDS-PAGE
Endotoxin Level:	< 0.1 ng/µg (<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:	200 µg/mL
Reconstitution Volume:	1 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