

Recombinant Murine EphrinB2-Fc

Ephrin-B2 (EPHB2) Fc Chimera is a fusion of EPHB2 to the C-terminal 6X histidine-tagged Fc region of IgG1. EPHB2 is a transmembrane protein-tyrosine kinase and member of the ephrin family. Implicated in developmental processes, particularly of the nervous system, as well as pattern formation and heart morphogenesis and angiogenesis through regulation of cell adhesion and migration.

Materials Provided

Catalog Number:	92-1089
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:	49600 g/mol
Source:	E. coli
Purity:	≥ 97% by SDS-PAGE
Endotoxin Level:	< 0.1 ng/µg (<1EU/µg).
Formulation:	Sterile filtered through a 0.2 micron filter. Lyophilized with no additives
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