

Recombinant Human Angiopoietin-2

Angiopoietin-2 (ANGPT2) is a modulating antagonist of angiopoietin-1 (ANGPT1) and TIE2, binding to TIE2 and competing with ANGPT1. Can induce tyrosine phosphorylation of TIE2, and may facilitate endothelial cell apoptosis and vascular regression through loosening of cell-matrix contacts. May also contribute to angiogenic signaling via stimulation of endothelial cell migration and proliferation.

Materials Provided

Catalog Number:	92-1293
Quantity Supplied:	1 vial(s), 25 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:	56 kg/mol
Source:	Recombinant
Purity:	≥ 97%
Endotoxin Level:	<0.10EU/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:	0.25 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