

Endothelin-1

Endothelin 1 is an endothelium derived vasoconstrictive peptide produced in vascular endothelial cells. It is one of the three isoforms of endothelin, and a potent agonist of the endothelin receptors type A and type B (ETA and ETB). Widely expressed in all tissues, and responsible for maintenance of vascular tone.

Materials Provided

Catalog Number:	92-1164
Quantity Supplied:	1 vial(s), 10 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:	2491.9 g/mol
Source:	Synthetic
Purity:	≥ 97% by HPLC
Endotoxin Level:	N/A
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:	30 µM
Reconstitution Volume:	0.133 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