

VIP, Human

Vasoactive intestinal peptide (VIP) is a 28 amino acid neuropeptide hormone that stimulates myocardial contractility, vasodilation and glycogenolysis, lowers arterial blood pressure, and relaxes the smooth muscles of the trachea, stomach, and gall bladder.

Materials Provided

Catalog Number:	92-1074
Quantity Supplied:	1 vial(s), 0.5 mg 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:	3552.9 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	500 µM
Reconstitution Volume:	0.280 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