

Secretin, Human

Secretin is a 27 amino acid GI peptide hormone responsible for regulation of water homeostasis, secretions of the stomach and pancreas, and pH of the duodenum. May also play a role in osmoregulation of the hypothalamus, pituitary, and kidney.

Materials Provided

Catalog Number:	92-1181
Quantity Supplied:	1 vial(s), 250 µg per vial

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:	3039.41 g/mol
Source:	Synthetic
Purity:	≥ 97%
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:	0.822 mL
Reconstitution Solvent:	H ₂ O

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