

(Tyr[SO₃]²⁷)CCK-8

(Tyr[SO₃]²⁷)CCK-8 (cholecystokinin) is an 8 amino acid GI peptide hormone and member of the CCK family that stimulates digestion of the fat and protein in the GI system. Research suggests it plays a role in opioid drug tolerance and hunger suppression. Agonist of the CCK A receptor.

Materials Provided

Catalog Number:	92-1152
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:	1143.27 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:	250 µM
Reconstitution Volume:	0.875 mL
Reconstitution Solvent:	0.05M NH ₄ OH

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