

Calcitonin, Human

Calcitonin is a polypeptide hormone produced in the thyroid by parafollicular cells (C-Cells). It reduces blood calcium levels in opposition to parathyroid hormone effects. It participates in both calcium and phosphorus metabolism, stimulates osteoblast activity, and inhibits intestinal calcium absorption, osteoclast activity, and renal tubular cell calcium absorption. Calcitonin levels in the blood are dictated by blood calcium levels.

Materials Provided

Catalog Number:	92-1054
Quantity Supplied:	1 vial(s), 1 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:	3417.9 g/mol
Source:	Synthetic
Purity:	≥ 95% 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:	500 µM
Reconstitution Volume:	0.585 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