

PTH(1-34)

Parathyroid hormone (PTH) is a polypeptide hormone secreted by the parathyroid that enhances bone calcium mobilization, enhances osteoblast activity (anabolic), increases blood calcium levels and calcium:phosphate ratio in the kidney, and promotes absorption of calcium by the intestines. The 1-34 amino acid fragment also induces BMP gene transcription. Agonist of the parathyroid hormone receptors PTH1 and PTH2.

Materials Provided

Catalog Number:	92-1082
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:	4117.74 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:	0.243 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