

TIP-39

Tuberoinfundibular peptide of 39 residues (TIP 39) is a neuropeptide protein related to parathyroid hormone that is a potent and selective agonist of parathyroid hormone 2 receptor (PTH2R), resulting in adenylyl cyclase activation and elevation of intracellular calcium levels. Also plays a role in protein kinase C beta activation, beta-arrestin recruitment, PTH2R internalization, cell proliferation inhibition, and possibly spermatogenesis and nociceptive circuit activation. Has little influence on PTH1R.

Materials Provided

Catalog Number:	92-1071
Quantity Supplied:	1 vial(s), 1 mg 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:	4505.3 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:	0.444 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