

Peptide YY, Human

Pancreatic peptide tyrosine tyrosine (peptide YY or PYY) is a 36 amino acid peptide released in the ileum and colon in response to feeding. It acts through neuropeptide Y (NPY) receptors to reduce appetite, inhibit gastric motility, and increase water and electrolyte absorption in the colon. It is a selective agonist for NPY Y2 over Y1 (0.11nM and 1050nM, respectively).

Materials Provided

Catalog Number:	92-1063
Quantity Supplied:	1 vial(s), 0.5 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:	4309.8 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:	200 µM
Reconstitution Volume:	0.580 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