

Pancreatic polypeptide

Pancreatic polypeptide is a 95 amino acid polypeptide secreted by the pancreas that is cleaved into a 36 amino acid active hormone that regulates pancreatic secretion, and an icosapeptide of unknown function. It is an endogenous agonist of the NPY Y4 receptor ($K_i=0.056\text{nM}$), and may play an important part in gastrointestinal function.

Materials Provided

| | |
|-------------------------------|----------------------------|
| Catalog Number: | 92-1062 |
| 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: | 4181.8 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: | 300 μM |
| Reconstitution Volume: | 0.398 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