

Neuromedin U-25, Human

Neuromedin U-25 is a 25 amino acid neuropeptide in humans that has functions in smooth muscle contraction, blood pressure regulation, pain perception, appetite, bone growth, and hormone release. Activates the Neuromedin receptors. Multiple isoforms exist in different species, with as few as 8 amino acids in some mammals.

Materials Provided

Catalog Number:	92-1174
Quantity Supplied:	1 vial(s), 500 µg 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:	3080.41 g/mol
Source:	Synthetic
Purity:	≥ 97%
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:	0.325 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