

## Neuromedin B

Neuromedin B (NMB) is a mammalian peptide related to bombesin that is present in the CNS and GI tract. Regulator of exocrine and endocrine secretions, cell growth, body temperature, smooth muscle contraction, and blood pressure and glucose levels. Potent agonist of the neuromedin B receptor (IC<sub>50</sub>=2nM).

### Materials Provided

<b>Catalog Number:</b>	<b>92-1233</b>
<b>Quantity Supplied:</b>	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

<b>Molecular Weight:</b>	1132.3 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 97.5%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	1 mM
<b>Reconstitution Volume:</b>	0.883 mL
<b>Reconstitution Solvent:</b>	H <sub>2</sub> O

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)