

## [Lys8,9]-Neurotensin (8-13)

[Lys8,9]-Neurotensin (8-13) is the smallest active fragment of Neurotensin, a 13 amino acid neuropeptide of the CNS that is implicated in luteinizing hormone (LH) and dopamine regulation, and prolactin release. Effects of Neurotensin activity include vasodilation, analgesia, hypothermia, and increased locomotion. This [Lys8,9]-analog of neurotensin enhances amiloride-sensitive Na<sup>+</sup>-dependent transmembrane H<sup>+</sup> flux.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1059</b>
<b>Quantity Supplied:</b>	1 vial(s), 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

<b>Molecular Weight:</b>	761 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 95%
<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>	10 mM
<b>Reconstitution Volume:</b>	0.675 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)