

## Neuropeptide S

Neuropeptide S (NPS) is a selective agonist of the G protein-coupled receptor NPSR that acts to suppress anxiety, fear, and appetite, and induce wakefulness and hyperactivity. Produced in the neurons of the amygdala, and possible uses include treatment for anxiety, sleep disorders, attention deficit hyperactivity disorder (ADHD), ADD, and asthma.

### Materials Provided

|                           |                          |
|---------------------------|--------------------------|
| <b>Catalog Number:</b>    | <b>92-1188</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> | 2187 g/mol                                        |
| <b>Source:</b>           | Synthetic                                         |
| <b>Purity:</b>           | ≥ 95% by HPLC                                     |
| <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>  | 0.457 mL              |
| <b>Reconstitution Solvent:</b> | Reconstitution Buffer |

***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)