

## Nociceptin

Nociceptin (Orphanin FQ) is a 17 amino acid opioid-related neuropeptide that is a potent anti-analgesic and is widely distributed in the CNS. Endogenous ligand for the nociceptin receptor (NOP,  $pK_i=8.6$ ), but has no activity at the opioid receptors  $\mu$ ,  $\delta$ , or  $\kappa$  ( $pK_i<6$ ). Plays a role in the regulation of reward and motivation pathways for substance abuse.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1182</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 mg per vial
<b>Reconstitution Buffer:</b>	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

<b>Molecular Weight:</b>	1808 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	$\geq 99\%$
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	$\leq -20^\circ\text{C}$ 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>	0.553 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)