

## Orexin A

Orexin A is a neuropeptide hormone released by the hypothalamus that excites neurons and affects hunger and wakefulness by modulating the dopamine, norepinephrine, histamine, and acetylcholine systems. Lack of orexin A tends to result in narcolepsy. Potent orexin receptor agonist ( $K_i=20$  and  $38\text{nM}$  for OX1 and OX2, respectively).

### Materials Provided

<b>Catalog Number:</b>	<b>92-1123</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>	3561.1 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	$\geq 97\%$
<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>	300 $\mu\text{M}$
<b>Reconstitution Volume:</b>	0.936 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)