

## Oxytocin

Oxytocin is a nine amino acid neurohypophysial hormone that acts as a neuromodulator in the brain. It is produced in the hypothalamus and secreted from the pituitary, and controls key parts of the reproductive system and brain, especially active during and after childbirth.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1060</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>	1007.2 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>	2.5 mM
<b>Reconstitution Volume:</b>	1.99 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)