

## N-Cyclopentyladenosine

N-Cyclopentyladenosine (CPA) is a centrally active synthetic drug compound that acts as an anticonvulsant, as well as a cardiovascular and sleep regulator. Potent and selective agonist of the adenosine receptor A1, with activity for A2A and A3 also ( $K_i=2.3, 790, \text{ and } 43\text{nM}$ , respectively).

### Materials Provided

<b>Catalog Number:</b>	<b>92-1195</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 mg (re-packaged) 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>	335.4 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 98%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	2-8°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

<b>Stock Concentration:</b>	50 mM
<b>Reconstitution Volume:</b>	0.596 mL
<b>Reconstitution Solvent:</b>	DMSO

***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)