

## NECA

NECA (5'-N-ethylcarboxamidoadenosine) is a bioactive adenosine analog that is a 22,900 times more potent vasodilator than adenosine and inhibitor of platelet aggregation. Potent agonist of the adenosine receptors A1, A2A, A3, and A2B ( $K_i=14, 20,$  and  $6.2\text{nM}$  for A1, A2A, and A3 respectively, and  $EC_{50}=2.4\text{uM}$  for A2B).

### Materials Provided

<b>Catalog Number:</b>	<b>92-1191</b>
<b>Quantity Supplied:</b>	1 vial(s), 10 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>	326.32 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>	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>	40 mM
<b>Reconstitution Volume:</b>	0.765 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)