

## (-)-Epinephrine (+)-bitartrate salt

(-)-Epinephrine (+)-bitartrate is the sympathomimetic adrenal medulla hormone and agonist for the adrenergic receptors alpha and beta. Stimulates the heart, including systemic vasoconstriction, and causes GI relaxation and bronchial dilation. Used for asthma and cardiac failure, as well as to delay absorption of local anesthetics.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1284</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>	333.29 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>	10 mM
<b>Reconstitution Volume:</b>	3 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)