

Isoproterenol

Isoproterenol is a synthetic small molecule non-selective beta-adrenergic agonist that acts as a sympathomimetic and is structurally similar to adrenaline. Activation of beta 2 andrenoceptor activates PKA and ERK downstream. Used as a medication for treatment of bradycardia, heart block , and asthma by relaxing blood vessels and airway.

Materials Provided

| | |
|---------------------------|---------------------------|
| Catalog Number: | 92-1119 |
| Quantity Supplied: | 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

| | |
|--------------------------|---|
| Molecular Weight: | 247.72 g/mol |
| Source: | Synthetic |
| Purity: | ≥ 98% by TLC |
| Endotoxin Level: | N/A |
| Formulation: | N/A |
| Storage: | ≤ -20°C Please avoid multiple freeze/thaw cycles. |

Reagent Preparation

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

| | |
|--------------------------------|----------------------|
| Stock Concentration: | 10 mM |
| Reconstitution Volume: | 4 mL |
| Reconstitution Solvent: | 0.3 mM Ascorbic Acid |

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com