

Dopamine

Dopamine hydrochloride is an endogenous neurotransmitter with sympathomimetic properties that is a potent agonist of the dopamine receptors D1-5. Increases heartbeat and blood flow to the body's tissues, also playing a role in reward-motivated behavior. Addictive drugs tend to increase dopamine activity, and multiple nervous system diseases are associated with its dysfunction, including Parkinson's and schizophrenia. Precursor in noradrenalin and adrenalin biosynthesis.

Materials Provided

Catalog Number:	92-1161
Quantity Supplied:	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

Molecular Weight:	189.64 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	2-8°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 mM
Reconstitution Volume:	527 µL with 10 mg vial
Reconstitution Solvent:	H ₂ O

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