

## Dexamethasone

Dexamethasone is an anti-inflammatory and immunosuppressant corticosteroid medication. Its glucocorticoid effect is 25 times more potent than cortisol. Some uses of Dexamethasone include inflammation relief, arthritis, rheumatologic problems, skin diseases, allergies, asthma, chronic obstructive lung disease, and certain cancers.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1132</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>	392.46 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 97%
<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>	2.548 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)