

## Calcium chloride dihydrate

Calcium chloride is a water soluble ionic compound common in labs. Uses include brine refrigeration, ice and dust control, water hardness, as a sequestrant and firming agent in food, for treatment of hydrofluoric acid burns and magnesium intoxication, and as a desiccator.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1220</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>	147.01 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 99%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	RT Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	1 M
<b>Reconstitution Volume:</b>	0.068 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)