

## Prostaglandin D2

Prostaglandin D2 is a naturally occurring eicosanoid product of mast cells that potently activates the prostaglandin D2 and CRTH2 receptors ( $K_i=21$  and  $2.4\text{nM}$ , respectively). Is released during allergic and asthmatic anaphylaxis, causing vasodilation, flushing, hypotension, inflammation, syncopal episodes, and immune cell recruitment.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1160</b>
<b>Quantity Supplied:</b>	1 vial(s), 1 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>	352.47 g/mol
<b>Source:</b>	Synthetic
<b>Purity:</b>	≥ 95%
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	≤ -20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	2.5 mM
<b>Reconstitution Volume:</b>	1.135 mL
<b>Reconstitution Solvent:</b>	Ethanol

***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)