

Prostaglandin E2

Prostaglandin E2 is a naturally occurring product of the arachidonic acid/cyclooxygenase pathway that influences many biological processes, including inflammation, immunity, relaxation of smooth muscle, fertility, parturition, and gastric mucosal integrity. Potent agonist of the prostaglandin receptors EP1, EP2, EP3, and EP4 (KDs=1-10nM). It has been marketed as a medicine for labor induction by multiple laboratories, including Forest Laboratories, Ferring Pharmaceuticals, and Pfizer.

Materials Provided

Catalog Number:	92-1154
Quantity Supplied:	1 vial(s), 1 mg per vial
Reconstitution Buffer:	1 mL

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:	352.5 g/mol
Source:	Synthetic
Purity:	≥ 99%
Endotoxin Level:	N/A
Formulation:	
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:	0.284 mL
Reconstitution Solvent:	Ethanol

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