

Cloprostenol

(+/-)-Cloprostenol is a synthetic water-soluble prostaglandin F_{2α} analog that can terminate the pregnancies of rats and hamsters 200 times more potently than PGF_{2α}. Potent agonist of the FP receptor (EC₅₀=0.84nM), optically active, and a potent luteolytic agent in rats and hamsters.

Materials Provided

Catalog Number:	92-1180
Quantity Supplied:	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

Molecular Weight:	446.9 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	5 mM
Reconstitution Volume:	0.447 mL
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