

I-BOP

I-BOP is a small molecule that induces platelet aggregation and vascular smooth muscle contraction. It is a potent TP receptor agonist, with an EC50 of 0.34nM.

Materials Provided

Catalog Number:	92-1038
Quantity Supplied:	1 vial(s), 100 µg 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:	512.4 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	1.95 mM in 100 µL ethanol
Reconstitution Volume:	Follow Handling Instructions
Reconstitution Solvent:	N/A

Handling Instructions:

To change the solvent, simply evaporate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of I-BOP is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers.

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