

IP1 Maleimide

Catalog Number: 20341

Unit Size: 100 ug

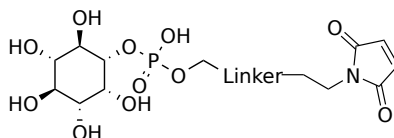
Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	454.32
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

- An excellent building block to develop IP1 probes
- Readily reacts with thiol compounds
- Might be used for detecting myo-inositol

IP1 Maleimide is an excellent building block to develop probes for detecting 1D-myo-inositol 1-phosphate (IP1). It readily reacts with thiol-containing biomolecules (such as peptides, antibodies and thiol-modified oligos) under extremely mild conditions. IP1 is an inositol having myo- configuration substituted at position 1 by a phosphate group. It has a role as a human metabolite, an Escherichia coli metabolite and a mouse metabolite. IP1 is functionally related to myo-inositol. It is a conjugate acid of 1D-myo-inositol 1-phosphate.