

## IP1 Succinimidyl Ester

Catalog Number: 20340

Unit Size: 100 ug

### Product Details

---

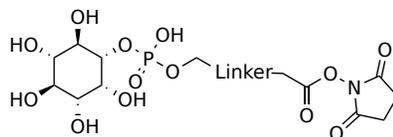
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid
Molecular Weight	471.35
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 amino compounds
- Might be used for detecting myo-inositol

IP1 Succinimidyl Ester is an excellent building block to develop probes for detecting 1D-myo-inositol 1-phosphate (IP1). It readily reacts with amino group-containing biomolecules (such as peptides, antibodies and amino-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 a myo-inositol. It is a conjugate acid of 1D-myo-inositol 1-phosphate.