

Isotonitazene NHS Ester

Catalog Number: 50635

Unit Size: 1 mg

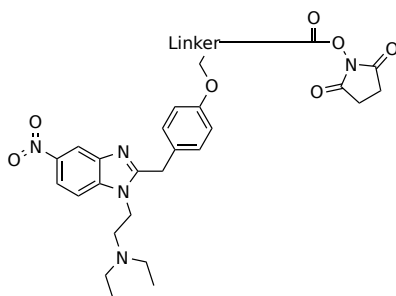
Product Details

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

Chemical Properties

Appearance	Solid brown
Molecular Weight	593.68
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Isotonitazene NHS ester is an excellent amino-reactive building block for developing assays and tests for detecting Isotonitazene. Isotonitazene is a potent synthetic opioid. This product is solely used for developing tools or analytical methods for detecting Isotonitazene.