

Isotonitazene KLH Conjugate

Catalog Number: 50641

Unit Size: 1 mg

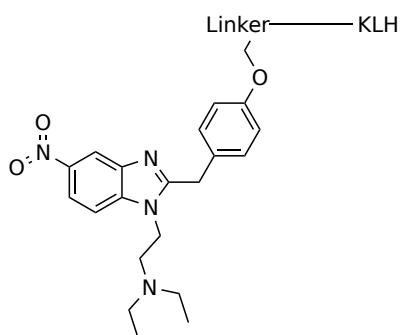
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	Water

Chemical Structure



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Isotonitazene-KLH conjugate is an excellent carrier protein for developing anti-Isotonitazene antibodies for detecting Isotonitazene. Isotonitazene is a potent synthetic opioid. This product is solely used for developing tools or analytical methods for detecting Isotonitazene.