

Isotonitazene BSA Conjugate

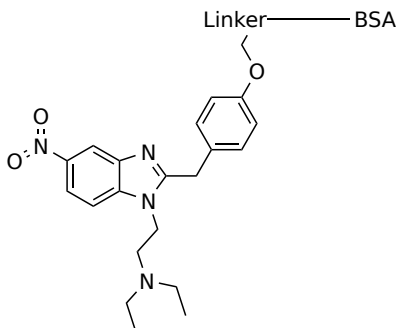
Catalog Number: 50640

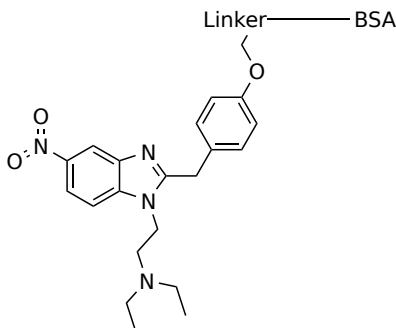
Unit Size: 1 mg

Product Details

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

Chemical Properties

Appearance	Solid white
Molecular Weight	N/A
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Isotonitazene-BSA conjugate is an excellent probe for developing ELISA 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.