

Isotonitazene HRP Conjugate

Catalog Number: 50642

Unit Size: 1 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 24 months upon receiving

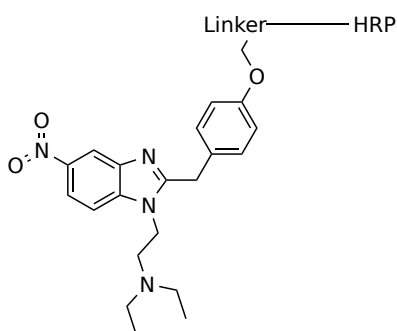
Chemical Properties

Appearance Solid

Molecular Weight N/A

Soluble In Water

Chemical Structure



Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

Applications

Isotonitazene-HRP 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.