

Norfentanyl-HRP Conjugate

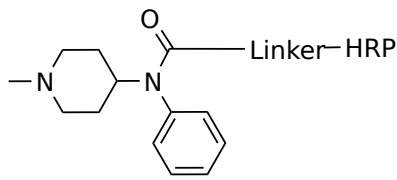
Catalog Number: 50631

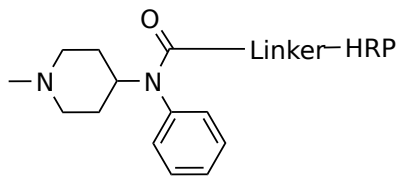
Unit Size: 1 mg

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

Norfentanyl-HRP is an excellent probe to develop assays and tests for detecting norfentanyl substances. AAT Bioquest offers the largest collection of HRP substrates that can be used with Norfentanyl-HRP for detecting norfentanyl substances. Norfentanyl is a monocarboxylic acid amide resulting from the formal condensation of the aryl amino group of 4-(N'-phenyl)piperidin-4-amine with propanoic acid. It is a major metabolite of fentanyl. It has a role as an opioid analgesic and a drug metabolite. It is an anilide, a member of piperidines and a monocarboxylic acid amide.