

Fentanyl succinimidyl ester

Catalog Number: 50560

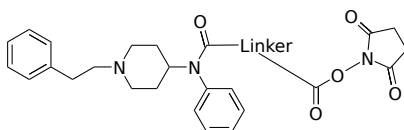
Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid off- white
Molecular Weight	491.59
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

Fentanyl succinimidyl ester can be conveniently used to make biological conjugates of fentanyl that are used to develop fentanyl assays and tests. It specifically reacts with aliphatic amines such as lysine residue of proteins under slightly basic conditions (pH 7.5-9.5).