

NBD-X, succinimidyl ester *CAS 145195-58-0*

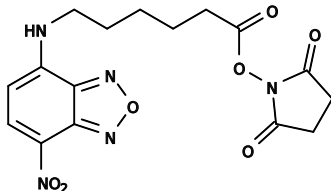
Catalog Number: 829

Unit Size: 100 mg

Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	391.33
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	467 nm
Emission Wavelength	539 nm

Applications

Compared to NBD chloride and fluoride, NBD-X acid gives better yield for labeling biopolymers. The fluorescence spectra of NBD derivatives are highly environment-sensitive, the fluorescence intensity decreases significantly in aqueous solutions.