

NBD-X acid *CAS 88235-25-0*

Catalog Number: 828

Unit Size: 1 g

Product Details

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

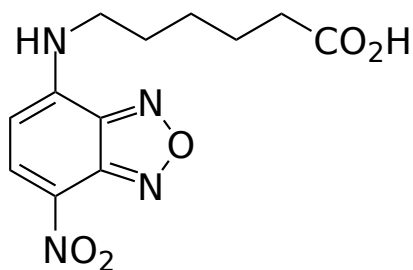
Chemical Properties

Appearance Solid yellow

Molecular Weight 294.27

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.