

# NBD-PE [N-(7-Nitrobenz-2-Oxa-1,3-Diazol-4-yl)-1,2-Dihexadecanoyl-sn-Glycero-3-Phosphoethanolamine, Triethylammonium Salt]

Catalog Number: 23305

Unit Size: 5 mg

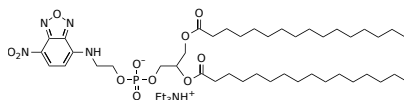
## Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

## Chemical Properties

Appearance	Solid brown
Molecular Weight	956.24
Soluble In	DMSO

Chemical Structure



## Spectral Properties

Excitation Wavelength	467 nm
Emission Wavelength	539 nm

## Applications

NBD-PE (N-(7-Nitrobenz-2-oxa-1,3-diazol-4-yl)-1,2-dihexadecanoyl-sn-glycero-3-phosphoethanolamine, triethylammonium salt) is a fluorescent phospholipid commonly used in cell biology research for studying membrane dynamics, including membrane fluidity and trafficking. Its ability to integrate seamlessly into cellular membranes and fluoresce upon excitation makes it a valuable tool for investigating the physical and functional properties of lipid bilayers.