

NBD-PE

NBD-PE (N-(7-Nitrobenz-2-oxa-1,3-diazol-4-yl)-1,2-dihexadecanoyl-snglycero-3-triethylammonium salt) in combination with Rhodamine-DHPE (#60026) or Texas Red®-DHPE (#60027) has been used to study membrane fusion via fluorescence resonance energy transfer (FRET) (1-3). NBD-PE has also been used in photobleaching recovery measurement.



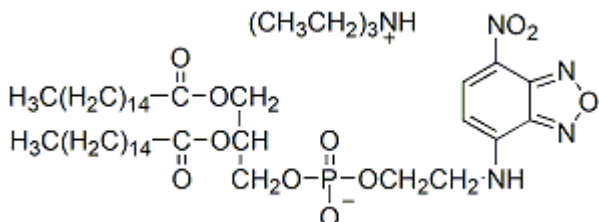
Product attributes

CAS number	178119-00-1
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Green
Excitation/Emission	463/536 nm
Conjugation	NBD

Product Description

NBD-PE (N-(7-Nitrobenz-2-oxa-1,3-diazol-4-yl)-1,2-dihexadecanoyl-snglycero-3-phosphoethanolamine, triethylammonium salt) in combination with Rhodamine-DHPE (#60026) or Texas Red®-DHPE (#60027) has been used to study membrane fusion via fluorescence resonance energy transfer (FRET) (1-3). NBD-PE has also been used in photobleaching recovery measurement.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 463/536 nm
- Orange solid soluble in chloroform
- Store at -20 °C and protect from light
- C₄₉H₉₀N₅O₁₁P
- MW: 956.24
- [178119-00-1]



Texas Red is a registered trademark of Thermo Fisher Scientific.

References

1. Biochemistry 20, 4093 (1981).
2. Meth Enzymol 221, 239 (1993).
3. Meth Enzymol 171, 850 (1989).
4. Prog Lipid Res 33, 203(1994).
5. J Cell Biol 122, 1253 (1993).

This datasheet was generated on June 20, 2026 at 11:09:18 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nbd-pe-n-7-nitrobenz-2-oxa-1,3-diazol-4-yl-1,2-dihexadecanoyl-snglycero-3-phosphoethanolamine-triethylammonium-salt/>