

## 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.



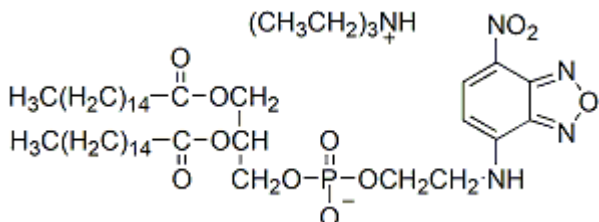
### Product attributes

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
- $C_{49}H_{90}N_5O_{11}P$
- 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.
5. J Cell Biol 122, 1253 (1993).

This datasheet was generated on January 19, 2026 at 02:12:21 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/nbd-pe-n-7-nitrobenz-2-oxa-13-diazol-4-yl-12-dihexadecanoyl-snglycero-3-phosphoethanolamine-triethylammonium-salt/>