Glowing products for science

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

Call us: 800-304-5357

lam	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 November 23, 2025 at 03:04:15 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/