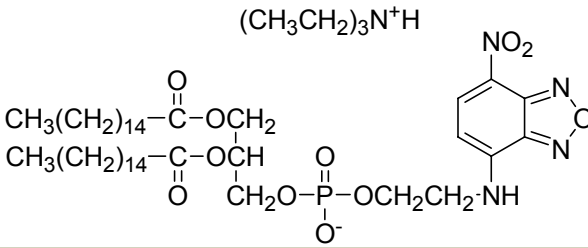


## N-(7-nitrobenz-2-oxa-1,3-diazol-4-yl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt

### Instruction Manual

Catalog Number	PK-CA707-60025
Description	NBD-PE in combination with Rhodamine-DHPE (Cat.No. PK-CA707-60026) or Texas Red™-DHPE (Cat.No. PK-CA707-60027) has been used to study membrane fusion via fluorescence resonance energy transfer (FRET). NBD-PE has also been used in photobleaching recovery measurement.
Quantity	10 mg
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (in MeOH) = 463/536 nm Extinction coefficient = 14,000
Molecular Structure	$(CH_3CH_2)_3N^+H$ 
Molecular Weight / Molecular Formula	956.24 Da; C <sub>49</sub> H <sub>90</sub> N <sub>5</sub> O <sub>11</sub> P
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Orange solid; soluble in chloroform.
Storage & Stability	Store at -20°C. Protect from light.
Applications	See Description
References	<ol style="list-style-type: none"> <li>1) Biochemistry 20, 4093(1981)</li> <li>2) Meth. Enzymol. 221, 239(1993)</li> <li>3) Meth. Enzymol. 171, 850(1989)</li> <li>4) Prog. Lipid Res. 33, 203(1994)</li> <li>5) J. Cell Biol. 122, 1253 (1993)</li> </ol>
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.