

PRODUCT INFORMATION



NBD-Phosphatidylethanolamine (ammonium salt)

Item No. 45094

CAS Registry No.: 477225-98-2

Synonym: NBD-PE

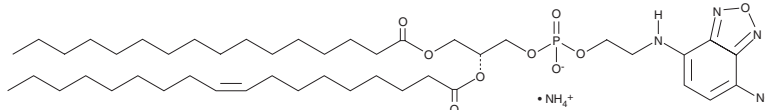
MF: $C_{45}H_{76}N_4O_{11}P \cdot NH_4$

FW: 898.1

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

NBD-Phosphatidylethanolamine is a phosphatidylethanolamine isolated from chicken egg that has been labeled with an N-linked fluorescent nitrobenzoxadiazole (NBD) group. NBD-phosphatidylethanolamines can be used in assays that monitor membrane fusion and aggregation in cells and to study the role of lipids in membrane fluidity.¹ NBD displays excitation/emission maxima of 460/538 nm, respectively. This product contains L- α -phosphatidylethanolamine molecular species with primarily C16:0, C18:0, C18:1, C18:2, and C20:4 fatty acyl chains.

Reference

1. Chattopadhyay, A. Chemistry and biology of N-(7-nitrobenz-2-oxa-1,3-diazol-4-yl)-labeled lipids: Fluorescent probes of biological and model membranes. *Chem. Phys. Lipids* **53**(1), 1-15 (1990).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/23/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM