

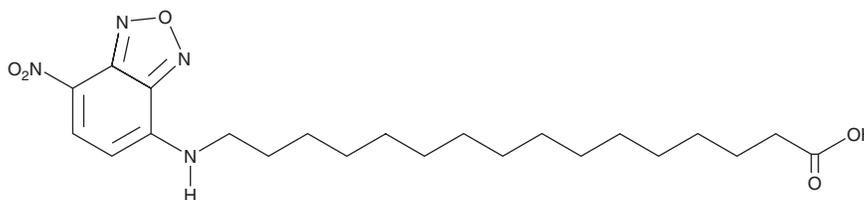
PRODUCT INFORMATION



NBD-Palmitic Acid

Item No. 44455

CAS Registry No.: 2260669-56-3
Formal Name: 16-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-hexadecanoic acid
Synonyms: NBD-C16:0, NBD-Cetylic Acid, NBD-FA 16:0, NBD-Hexadecanoic Acid
MF: C₂₂H₃₄N₄O₅
FW: 434.5
Purity: ≥98%
Ex./Em. Max.: 460/538 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

NBD-palmitic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the NBD-palmitic acid in the solvent of choice, which should be purged with an inert gas. NBD-palmitic acid is soluble (≥10 mg/ml) in DMSO.

Description

NBD-palmitic acid is a biologically active derivative of palmitic acid (Item Nos. 10006627 | 43561) tagged with a fluorescent nitrobenzoxadiazole (NBD) group at carbon 16. It has been used in the study of lipid uptake and lipid-binding proteins.¹⁻³ NBD displays excitation/emission maxima of 460/538 nm, respectively.

References

1. Das, S., Manna, A., Majumdar, O., *et al.* M-O-M mediated denaturation resistant P2 tetramer on the infected erythrocyte surface of malaria parasite imports serum fatty acids. *iScience* **27(5)**, 109760 (2024).
2. Ha, H.T., Liu, S., Nguyen, X.T., *et al.* Lack of SPNS1 results in accumulation of lysolipids and lysosomal storage disease in mouse models. *JCI Insight* **9(8)**, e175462 (2024).
3. Osawa, T., Kotani, T., Kawaoka, T., *et al.* Atg2 mediates direct lipid transfer between membranes for autophagosome formation. *Nat. Struct. Mol. Biol.* **26(4)**, 281-288 (2019).

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, 12/02/2025

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