

PRODUCT INFORMATION



C-6 NBD

Item No. 18871

CAS Registry No.: 88235-25-0
Formal Name: 6-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-hexanoic acid

Synonyms: C₆-NBD, NBD Hexanoic Acid, NBD-X, 6-(7-Nitrobenzofurazan-4-ylamino)hexanoic Acid

MF: C₁₂H₁₄N₄O₅

FW: 294.3

Purity: ≥96%

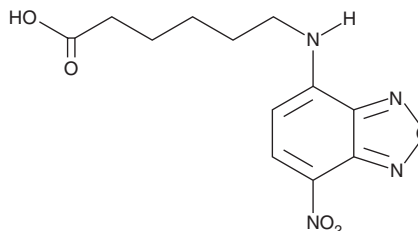
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: Protect from light and moisture

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



Laboratory Procedures

C-6 NBD is supplied as a solid. A stock solution may be made by dissolving the C-6 NBD in the solvent of choice, which should be purged with an inert gas. C-6 NBD is soluble in the organic solvent DMSO.

Description

C-6 NBD is a derivative of the environmentally sensitive fluorophore nitrobenzoxadiazole (NBD; absorbance 465 nm, emission 535 nm). It is modified with a hexanoic acid tail, which provides a reactive carboxylic acid site. C-6 NBD has been used to synthesize a variety of fluorescently tagged compounds for research purposes.¹⁻⁵

References

1. Lipsky, N.G. and Pagano, R.E. A vital stain for the golgi apparatus. *Science* **228**, 745-757 (1985).
2. Paul, P., Kamisaka, Y., Marks, D.L., *et al.* Purification and characterization of UDP-glucose: Ceramide glucosyltransferase from rat liver Golgi membranes. *J. Biol. Chem.* **271**, 2287-2293 (1996).
3. Zurzolo, C., van't Hof, W., van Meer, G., *et al.* VIP21/caveolin, glycosphingolipid clusters and the sorting of glycosylphosphatidylinositol-anchored proteins in epithelial cells. *EMBO J.* **13**, 42-53 (1994).
4. Engfeldt, T., Renberg, B., Brumer, J., *et al.* Chemical synthesis of triple-labelled three-helix bundle binding proteins for specific fluorescent detection of unlabelled protein. *ChemBioChem* **6(6)**, 1043-1050 (2005).
5. Escoffre, J.-M., Bellard, E., Faurie, C., *et al.* Membrane disorder and phospholipid scrambling in electroporabilized and viable cells. *Biochim. Biophys. Acta.* **1838(7)**, 1701-1709 (2014).

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, 11/02/2022

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