

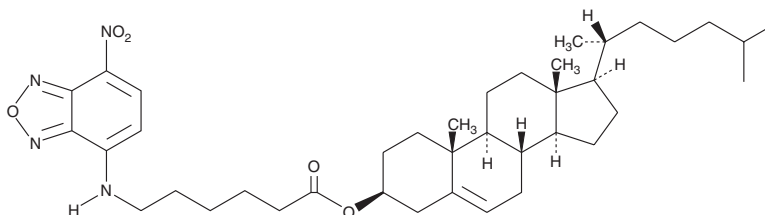
PRODUCT INFORMATION



3-hexanoyl-NBD Cholesterol

Item No. 13221

CAS Registry No.: 201731-19-3
Formal Name: 6-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-cholest-5-en-3-ol
Synonym: 3-C₆-NBD Cholesterol
MF: C₃₉H₅₈N₄O₅
FW: 662.9
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 332, 465 nm
Ex./Em. Max: 473/536 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

3-hexanoyl-NBD cholesterol is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-hexanoyl-NBD cholesterol in the solvent of choice, which should be purged with an inert gas. 3-hexanoyl-NBD cholesterol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-hexanoyl-NBD cholesterol is approximately 1 mg/ml in ethanol and DMF and approximately 0.25 mg/ml in DMSO.

Description

3-hexanoyl-NBD cholesterol is a fluorescently tagged cholesterol derivative with the hydrophilic NBD fluorophore attached to carbon 3, at the hydrophilic end of cholesterol, separated by a 6-carbon spacer. This design allows the cholesterol to properly orient in membrane bilayers while the fluorescent tag is presented outside of the bilayer. This should model the behavior of cholesterol in membranes better than the previously-used 25-NBD cholesterol, which positions NBD directly on the 25th carbon of cholesterol at the hydrophobic terminus. 3-hexanoyl-NBD cholesterol has excitation/emission maxima of 473/536 nm, respectively, in vesicles comprised of dioleoylphosphatidylcholine (DOPC), however, the maxima will vary depending on membrane composition.¹ NBD has excitation/emission maxima of 465/535 nm, respectively. Fluorescently tagged lipids have been used to study their interactions with proteins, their utilization by cells and liposomes, and for the development of assays for lipid metabolism.²⁻⁶

References

1. Ramirez, D.M.C., Ogilvie, W.W., and Johnston, L.J. NBD-cholesterol probes to track cholesterol distribution in model membranes. *Biochim. Biophys. Acta.* **1798(3)**, 558-568 (2010).
2. Kumagai, K., Yasuda, S., Okemoto, K., et al. CERT mediates intermembrane transfer of various molecular species of ceramides. *J. Biol. Chem.* **280(8)**, 6488-6495 (2005).
3. Luo, M., Jones, S.M., Peters-Golden, M., et al. Nuclear localization of 5-lipoxygenase as a determinant of leukotriene B₄ synthetic capacity. *Proc. Nat. Acad. Sci. USA* **100(21)**, 12165-12170 (2003).
4. Moreno, M.J., Estronca, L.M.B.B., and Vaz, W.L.C. Translocation of phospholipids and dithionite permeability in liquid-ordered and liquid-disordered membranes. *Biophys. J.* **91(3)**, 873-881 (2006).
5. Loidl, A., Claus, R., Deigner, H.P., et al. High-precision fluorescence assay for sphingomyelinase activity of isolated enzymes and cell lysates. *J. Lipid Res.* **43(5)**, 815-823 (2002).
6. Tani, M., Okino, N., Mitsutake, S., et al. Specific and sensitive assay for alkaline and neutral ceramidases involving C₁₂-NBD-ceramide. *J. Biochem.* **125(4)**, 746-749 (1999).

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, 02/18/2020

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