

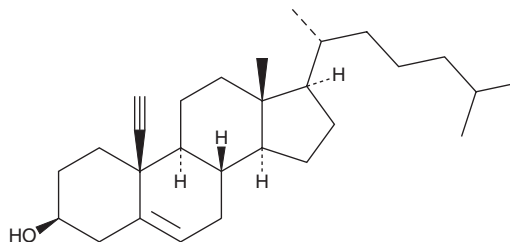
PRODUCT INFORMATION



19-alkyne Cholesterol

Item No. 16446

CAS Registry No.: 2058319-71-2
Formal Name: 10-ethynyl-19-norcholest-5-en-3 β -ol
Synonyms: 19-ethynyl Cholesterol,
Click Tag™ 19-alkyne Cholesterol
MF: C₂₈H₄₄O
FW: 396.7
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

19-alkyne Cholesterol is supplied as a crystalline solid. A stock solution may be made by dissolving the 19-alkyne cholesterol in the solvent of choice, which should be purged with an inert gas. 19-alkyne Cholesterol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 19-alkyne cholesterol in these solvents is approximately 20, 0.1, and 2 mg/ml, respectively.

19-alkyne Cholesterol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 19-alkyne Cholesterol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 19-alkyne Cholesterol has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

19-alkyne Cholesterol is an alkyne derivative of cholesterol (Item No. 9003100).¹ It has been used to track cholesterol localization and subcellular distribution in NIH3T3 cells and to study the role of cholesterol in hedgehog (Hh) signal transduction in Shh-LIGHT2 cells using click chemistry and various reporter azides.

Reference

1. Jao, C.Y., Nedelcu, D., Lopez, L.V., *et al.* Bioorthogonal probes for imaging sterols in cells. *ChemBiochem*. **16**(4), 611-617 (2015).

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, 08/24/2023

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