

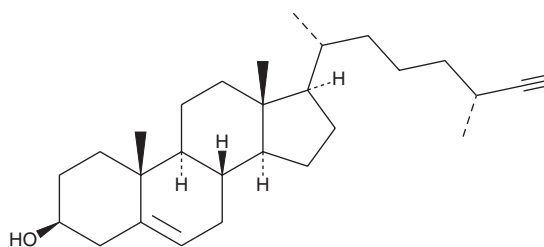
PRODUCT INFORMATION



27-alkyne Cholesterol

Item No. 24319

CAS Registry No.: 1527467-07-7
Formal Name: 20R-[(4R)-4-methyl-5-hexyn-1-yl]-pregn-5-en-3 β -ol
Synonym: Click Tag™ 27-alkyne Cholesterol
MF: C₂₈H₄₄O
FW: 396.6
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

27-alkyne Cholesterol is supplied as a crystalline solid. A stock solution may be made by dissolving the 27-alkyne cholesterol in the solvent of choice. 27-alkyne Cholesterol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 27-alkyne cholesterol in these solvents is approximately 20, 0.1, and 2 mg/ml, respectively.

27-alkyne Cholesterol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 27-alkyne cholesterol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 27-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

27-alkyne Cholesterol is an alkyne derivative of cholesterol. It has been used to track cholesterol metabolism and localization using click chemistry and various reporter azides in fixed or living cells.¹ It is a substrate for enzymes from a variety of species including bacteria, yeast, rat, and human.

Reference

1. Hofmann, K., Thiele, C., Schött, H.F., *et al.* A novel alkyne cholesterol to trace cellular cholesterol metabolism and localization. *J. Lipid Res.* **55(3)**, 583-591 (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, 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