

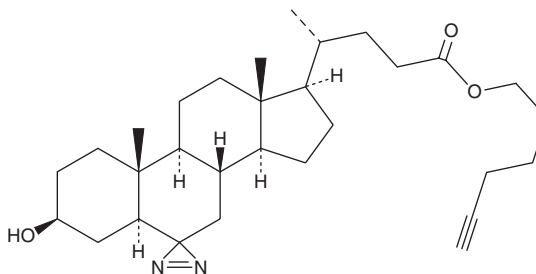
PRODUCT INFORMATION



Photoclick Cholesterol

Item No. 24528

CAS Registry No.: 1485490-47-8
Formal Name: (3' β ,5' α)-3'-hydroxy-spiro[3H-diazirine-3,6'-cholan]-24'-oic acid, 5-hexyn-1-yl ester
Synonym: Click Tag™ Cholesterol
MF: C₃₀H₄₆N₂O₃
FW: 482.7
Purity: ≥95%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: ≥1 year



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

Description

Photoclick cholesterol is a derivative of the sterol lipid cholesterol (Item No. 9003100) that contains a terminal clickable alkyne moiety and a photoactivatable diazirine group.¹ It has been used in quantitative proteomic screens for sterol-binding proteins (SBPs) in *S. cerevisiae*.

Reference

1. Chauhan, N., Sere, Y.Y., Sokol, A.M., *et al.* A PhotoClick cholesterol-based quantitative proteomics screen for cytoplasmic sterol-binding proteins in *Saccharomyces cerevisiae*. *Yeast* **37(1)**, 15-25 (2020).

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