

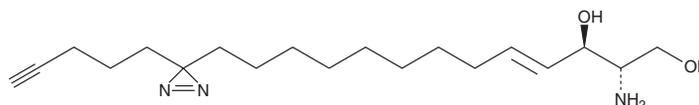
PRODUCT INFORMATION



Photoclick Sphingosine

Item No. 25364

CAS Registry No.: 1823021-14-2
Formal Name: (2S,3R,4E)-2-amino-13-[3-(4-pentyn-1-yl)-3H-diazirin-3-yl]-4-tridecene-1,3-diol
Synonyms: Click Tag™ Sphingosine, pacSph, pacSphingosine (d18:1)
MF: C₁₉H₃₃N₃O₂
FW: 335.5
Purity: ≥95%
UV/Vis.: λ_{max}: 349 nm
Supplied as: A 5 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

Photoclick sphingosine is a photoreactive probe for the study of sphingosine metabolism.¹ It contains a terminal clickable alkyne moiety and a photoactivatable diazirine group that allow its interaction with sphingosine binding proteins and photo-activated cross-linking of its interacting partners, respectively. Photoclick sphingosine has been used to monitor the metabolism of sphingosine and visualize protein-lipid interaction complexes in *S1PL*^{-/-} mouse embryonic fibroblasts (MEFs).

Reference

1. Haberkant, P., Stein, F., Höglinger, D., *et al.* Bifunctional sphingosine for cell-based analysis of protein-sphingolipid interactions. *ACS Chem. Biol.* **11**(1), 222-230 (2016).

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, 02/05/2026

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