

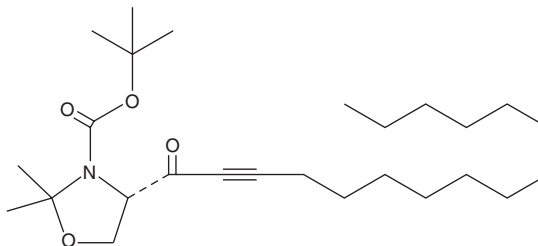
PRODUCT INFORMATION



CAY10621

Item No. 13371

CAS Registry No.: 120005-55-2
Formal Name: (4S)-2,2-dimethyl-4-(1-oxo-2-hexadecyn-1-yl)-3-oxazolidinecarboxylic acid, 1,1-dimethylethyl ester
Synonyms: SKI 5C, SPHK1 Inhibitor 5C
MF: C₂₆H₄₅NO₄
FW: 435.6
Purity: ≥98%
UV/Vis.: λ_{max}: 224 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

CAY10621 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of CAY10621 in these solvents is approximately 2 mg/ml.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10621 is an inhibitor of sphingosine kinase 1 (SPHK1; IC₅₀ = 3.3 μM).¹ It is selective for SPHK1 over SPHK2 at 10 μM and PKC at concentrations less than 100 μM. CAY10621 inhibits SPHK1 activity by 70% in U937 cells when used at a concentration of 5 μM.

Reference

1. Wong, L., Tan, S.L., Lam, Y., *et al.* Synthesis and evaluation of sphingosine analogues as inhibitors of sphingosine kinases. *J. Med Chem.* **52(12)**, 3618-3626 (2009).

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, 03/08/2024

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