

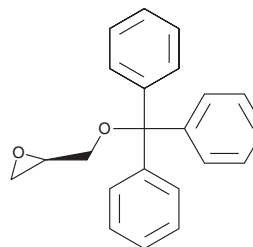
PRODUCT INFORMATION



(R)-(+)-Trityl glycidyl ether

Item No. 34131

CAS Registry No.: 65291-30-7
Formal Name: 2R-[(triphenylmethoxy)methyl]-oxirane
Synonym: (R)-Trityl Glycidol
MF: C₂₂H₂₀O₂
FW: 316.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(R)-(+)-Trityl glycidyl ether is supplied as a solid. A stock solution may be made by dissolving the (R)-(+)-trityl glycidyl ether in the solvent of choice, which should be purged with an inert gas. (R)-(+)-Trityl glycidyl ether is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of (R)-(+)-trityl glycidyl ether in DMSO and DMF is approximately 1 and 10 mg/ml, respectively.

Description

(R)-(+)-Trityl glycidyl ether is a precursor that has been used in the synthesis of glycerophospholipids, as well as compounds with antiviral and antimalarial activities.¹⁻³

References

1. Hendrickson, H.S. and Hendrickson, E.K. A facile asymmetric synthesis of glycerol phospholipids via trityl glycidol prepared by the asymmetric epoxidation of allyl alcohol. Thiolester and thioether analogs of phosphatidylcholine. *Chem. Phys. Lipids* **53(1)**, 115-120 (1990).
2. Baszczyński, O., Jansa, P., Dračinský, M., et al. Synthesis and antiviral activity of N⁹-[3-fluoro-2-(phosphonomethoxy)propyl] analogues derived from N⁶-substituted adenines and 2,6-diaminopurines. *Bioorg. Med. Chem.* **19(7)**, 2114-2124 (2011).
3. Sato, E., Morita, M., Ogawa, H., et al. Design, synthesis and anti-malarial activities of synthetic analogs of biselyngbyolide B, a Ca²⁺ pump inhibitor from marine cyanobacteria. *Bioorg. Med. Chem. Lett.* **28(3)**, 298-301 (2018).

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, 11/17/2022

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