

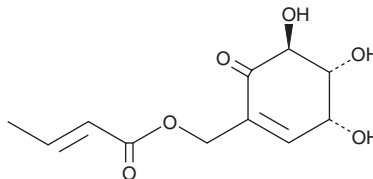
PRODUCT INFORMATION



6-*epi* COTC

Item No. 37746

CAS Registry No.: 959150-61-9
Formal Name: 2E-butenic acid, [(3R,4R,5S)-3,4,5-trihydroxy-6-oxo-1-cyclohexen-1-yl]methyl ester
MF: C₁₁H₁₄O₆
FW: 242.2
Purity: ≥90%
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

6-*epi* COTC is supplied as a solid. A stock solution may be made by dissolving the 6-*epi* COTC in the solvent of choice. 6-*epi* COTC is soluble in organic solvents such as DMSO, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

6-*epi* COTC is a derivative of COTC (Item No. 36408) with anticancer activity.¹ It is cytotoxic to A549 and H460 non-small cell lung cancer (NSCLC) cells (IC₅₀s = 170 and 158 μM, respectively).

Reference

1. Arthurs, C.L., Raftery, J., Whitby, H.L., et al. Arene *cis*-dihydrodiols: Useful precursors for the preparation of analogues of the anti-tumour agent, 2-crotonyloxymethyl-(4R,5R,6R)-4,5,6-trihydroxycyclohex-2-enone (COTC). *Bioorg. Med. Chem. Lett.* **17**(21), 5974-5977 (2007).

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, 02/14/2023

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