

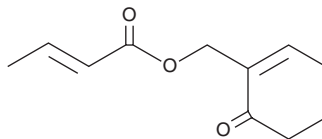
PRODUCT INFORMATION



COMC

Item No. 37736

CAS Registry No.: 106281-45-2
Formal Name: 2-butenic acid, (6-oxo-1-cyclohexen-1-yl) methyl ester
Synonym: 2-Crotonyloxymethyl-2-cyclohexenone
MF: C₁₁H₁₄O₃
FW: 194.2
Purity: ≥90%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

COMC is supplied as a neat oil. A stock solution may be made by dissolving the COMC in the solvent of choice, which should be purged with an inert gas. COMC is soluble in acetonitrile, chloroform, and DMSO.

Description

COMC is an anticancer agent and a derivative of the bacterial metabolite COTC (Item No. 36408).¹ It forms adducts with glutathione and nucleic acids.^{2,3} COMC inhibits the proliferation of B16 murine melanoma cells (IC₅₀ = 0.041 μM).¹

References

1. Joseph, E., Eiseman, J.L., Hamilton, D.S., *et al.* Molecular basis of the antitumor activities of 2-crotonyloxymethyl-2-cycloalkenones. *J. Med. Chem.* **46**(1), 194-196 (2003).
2. Joseph, E., Ganem, B., Eiseman, J.L., *et al.* Selective inhibition of MCF-7^{piGST} breast tumors using glutathione transferase-derived 2-methylene-cycloalkenones. *J. Med. Chem.* **48**(21), 6549-6552 (2005).
3. Zhang, Q., Ding, Z., Creighton, D.J., *et al.* Alkylation of nucleic acids by the antitumor agent COMC. *Org. Lett.* **4**(9), 1459-1462 (2002).

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, 01/13/2023

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