

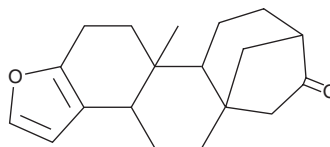
PRODUCT INFORMATION



16-Oxocafestol

Item No. 39940

CAS Registry No.: 108664-98-8
Formal Name: (3bS,5aS,8R,10aR,10bS)-3b,4,5,8,9,10,10a,10b,11,12-decahydro-10b-methyl-5a,8-methano-5aH-cyclohepta[5,6]naphtho[2,1-b]furan-7(6H)-one
MF: C₁₉H₂₄O₂
FW: 284.4
Purity: ≥95%
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

16-Oxocafestol is supplied as a solid. A stock solution may be made by dissolving the 16-oxocafestol in the solvent of choice, which should be purged with an inert gas. 16-Oxocafestol is slightly soluble in chloroform and methanol.

Description

16-Oxocafestol is a derivative of the diterpene cafestol (Item No. 13999).¹ It increases cytosolic glutathione S-transferase (GST) activity and acid-soluble sulfhydryl levels in mouse liver and small bowel mucosa when administered at a dose of 10 μmol/animal once per day for three days.

Reference

1. Lam, L.K., Sparnins, V.L., and Wattenberg, L.W. Effects of derivatives of kahweol and cafestol on the activity of glutathione S-transferase in mice. *J. Med. Chem.* **30(8)**, 1399-1403 (1987).

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, 02/13/2024

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