

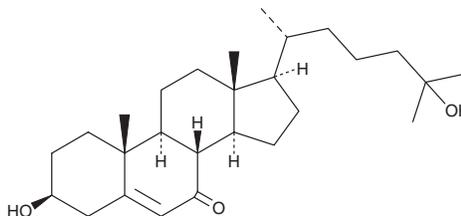
PRODUCT INFORMATION



7-keto-25-hydroxy Cholesterol

Item No. 25973

CAS Registry No.: 64907-23-9
Formal Name: 3 β ,25-dihydroxy-cholest-5-en-7-one
MF: C₂₇H₄₄O₃
FW: 416.6
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

7-keto-25-hydroxy Cholesterol is supplied as a solid. A stock solution may be made by dissolving the 7-keto-25-hydroxy cholesterol in the solvent of choice, which should be purged with an inert gas. 7-keto-25-hydroxy Cholesterol is soluble in organic solvents such as chloroform and methanol.

Description

7-keto-25-hydroxy Cholesterol is an oxysterol and a proposed metabolite of 7-keto cholesterol (Item No. 16339) and 7-dehydro cholesterol (Item No. 14612).¹ It binds to and activates Smoothed (Smo) at the conserved extracellular cysteine-rich domain (CRD) in a reporter assay. It also increases EGF high affinity binding site numbers in normal rat kidney (NRK) cells.²

References

1. Myers, B.R., Sever, N., Chong, Y.C., *et al.* Hedgehog pathway modulation by multiple lipid binding sites on the smoothed effector of signal response. *Dev. Cell.* **26(4)**, 346-357 (2013).
2. Astruc, M.E. and Lahoua, Z. Potentiation by cholesterol and vitamin D₃ oxygenated derivatives of arachidonic acid release and prostaglandin E₂ synthesis induced by the epidermal growth factor in NRK 49F cells: The role of protein kinase C. *Cell Signal.* **6(7)**, 763-775 (1994).

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, 12/20/2022

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