

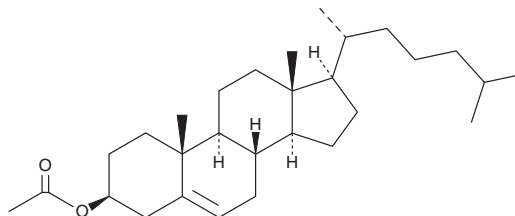
PRODUCT INFORMATION



Cholesteryl Acetate

Item No. 30816

CAS Registry No.: 604-35-3
Formal Name: cholest-5-en-3 β -ol, 3-acetate
Synonyms: 3 β -acetoxy-5-Cholestene,
3 β -hydroxy-5-Cholestene 3-acetate,
Cholesteryl Acetic Acid, NSC 8799
MF: C₂₉H₄₈O₂
FW: 428.7
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

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

Description

Cholesteryl acetate is a cholesterol ester.¹ It has been used as an internal standard for the quantification of cholesterol, as well as cholesterol derivatives, by GC or GLC.^{2,3} Cholesteryl acetate has also been derivatized from, and used for the quantification of, cholesterol by ESI-MS/MS.⁴ It has also been used in the manufacturing of liquid crystal displays (LCDs), wrist watches, and thermometers.⁵

References

1. Huang, Z., Kawi, S., and Chiew, Y.C. Solubility of cholesterol and its esters in supercritical carbon dioxide with and without cosolvents. *J. Supercrit. Fluids* **30(1)**, 25-39 (2004).
2. Jones Owen, V.M., Ho, F.K., Mazzuchin, A., et al. Cholesterol in amniotic fluid, determined by gas chromatography. *Clin. Chem.* **22(2)**, 224-226 (1976).
3. Gray, G.M. and Yardley, H.J. Lipid compositions of cells isolated from pig, human, and rat epidermis. *J. Lipid Res.* **16(6)**, 434-440 (1975).
4. Liebis, G., Binder, M., Schifferer, R., et al. High throughput quantification of cholesterol and cholesteryl ester by electrospray ionization tandem mass spectrometry (ESI-MS/MS). *Biochim. Biophys. Acta* **176(1)**, 121-128 (2006).
5. Vijayakumar, V.N., Rajasekaran, T.R., and Baskar, K. Influence of hydrogen bond on thermal and phase transitions of binary complex liquid crystals. *Russ. J. Phys. Chem. A* **91(13)**, 2578-2584 (2017).

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, 10/06/2022

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