

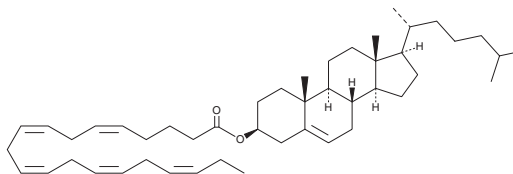
PRODUCT INFORMATION



Cholesteryl Eicosapentaenoate

Item No. 22719

CAS Registry No.: 74892-97-0
Formal Name: cholest-5-en-3 β -ol,
3-[(5Z,8Z,11Z,14Z,17Z)-
5,8,11,14,17-eicosapentaenoate]
Synonym: 20:5 Cholesteryl ester
MF: C₄₇H₇₄O₂
FW: 671.1
Purity: \geq 95%
Supplied as: A solution in chloroform
Storage: -20°C
Stability: \geq 2 years



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

Description

Cholesteryl eicosapentaenoate is a cholesterol ester.¹ It induces cytotoxicity in human monocyte-macrophages in a concentration-dependent manner at concentrations ranging from 88 to 880 μ M. Cholesteryl eicosapentaenoate levels are elevated 1.89-fold in the serum of pediatric patients with extrahepatic biliary atresia and decreased in the serum and plasma of patients with Alzheimer's disease and abetalipoproteinemia, respectively, compared with healthy individuals.²⁻⁴

References

1. Hardwick, S.J., Carpenter, K.L., Law, N.S., *et al.* Toxicity of polyunsaturated fatty acid esters for human monocyte-macrophages: The anomalous behaviour of cholesteryl linolenate. *Free Radic. Res.* **26(4)**, 351-362 (1997).
2. Robberecht, E., Koletzko, B., and Christophe, A. Several mechanisms contribute to the abnormal fatty acid composition of serum phospholipids and cholesterol esters in cholestatic children with extrahepatic biliary atresia. *Prostaglandins Leukot. Essent. Fatty Acids* **56(3)**, 199-204 (1997).
3. Tully, A.M., Roche, H.M., Doyle, R., *et al.* Low serum cholesteryl ester-docosahexaenoic acid levels in Alzheimer's disease: A case-control study. *Br. J. Nutr.* **89(4)**, 483-489 (2003).
4. Subbaiah, P.V., Banerji, B., Gregg, R.E., *et al.* Molecular species of cholesteryl esters formed in abetalipoproteinemia: Effect of apoprotein B-containing lipoproteins. *J. Lipid Res.* **31(5)**, 927-932 (1990).

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, 10/24/2022

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