

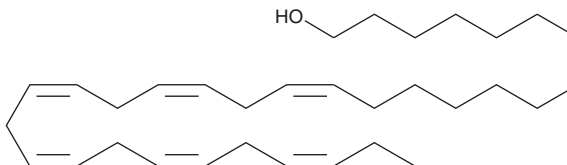
PRODUCT INFORMATION



all-cis-14,17,20,23,26,29-Dotriacontahexaenol

Item No. 10562

CAS Registry No.: 2649103-52-4
Formal Name: 14Z,17Z,20Z,23Z,26Z,29Z-dotriacontahexaen-1-ol
Synonyms: CAY10631, 14(Z),17(Z),20(Z),23(Z),26(Z),29(Z)-Dotriacontahexaenol, FOH 32:6
MF: $C_{32}H_{54}O$
FW: 454.8
Purity: $\geq 95\%$
Supplied as: A 1 mg/ml solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

all-cis-14,17,20,23,26,29-Dotriacontahexaenol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of all-cis-14,17,20,23,26,29-dotriacontahexaenol in these solvents is approximately 10 and 20 mg/ml, respectively.

Description

all-cis-14,17,20,23,26,29-Dotriacontahexaenol is a very long-chain fatty alcohol. It is an intermediate in the synthesis of the very long-chain PUFA all-cis-14,17,20,23,26,29-dotriacontahexaenoic acid.¹

Reference

1. Wade, A., Rallabandi, R., Lucas, S., *et al.* The synthesis of the very long chain polyunsaturated fatty acid (VLC-PUFA) 32:6 n-3. *Org. Biomol. Chem.* **19(25)**, 5563-5566 (2021).

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, 11/19/2025

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