

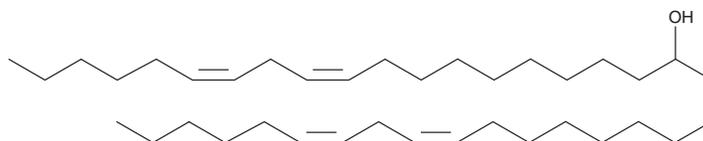
PRODUCT INFORMATION



DLin-MeOH

Item No. 41432

CAS Registry No.: 1169768-28-8
Formal Name: 6Z,9Z,28Z,31Z-heptatriacontatetraen-19-ol
MF: $C_{37}H_{68}O$
FW: 528.9
Purity: $\geq 98\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 1 year



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

Laboratory Procedures

DLin-MeOH is supplied as a solution in ethanol. To change the solvent, simply evaporate the DLin-MeOH under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO purged with an inert gas can be used. DLin-MeOH is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DLin-MeOH is a synthetic intermediate in the synthesis of the ionizable cationic amino lipid DLin-MC3-DMA (Item No. 34364).¹

Reference

1. Jayaraman, M., Ansell, S.M., Mui, B.L., *et al.* Maximizing the potency of siRNA lipid nanoparticles for hepatic gene silencing in vivo. *Angew. Chem. Int. Ed. Engl.* **51(34)**, 8529-8533 (2012).

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, 06/26/2024

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