

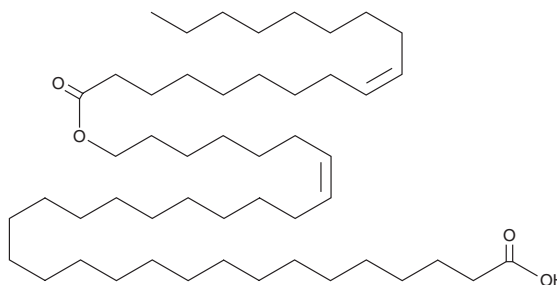
PRODUCT INFORMATION



FAHFA 18:1(9Z)/32-O-32:1(25Z)

Item No. 37879

Formal Name: (Z)-32-(oleoyloxy)dotriacont-25-enoic acid
Synonym: OAHFA 18:1(n-9, cis)/ω-O-32:1(n-7, cis)
MF: C₅₀H₉₄O₄
FW: 759.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

FAHFA 18:1(9Z)/32-O-32:1(25Z) is an (O-acyl)-ω-hydroxy fatty acid (OAHFA) containing the monounsaturated fatty acid oleic acid (Item Nos. 90260 | 24659) and the very long-chain ω-hydroxy fatty acid 25(Z)-dotriacontenoic acid.¹

Reference

1. Hancock, S.E., Ailuri, R., Marshall, D.L., *et al.* Mass spectrometry-directed structure elucidation and total synthesis of ultra-long chain (O-acyl)-ω-hydroxy fatty acids. *J. Lipid Res.* **59(8)**, 1510-1518 (2018).

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, 10/22/2025

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