

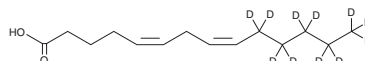
PRODUCT INFORMATION



5(Z),8(Z)-Tetradecadienoic Acid-d₁₁

Item No. 39711

Formal Name: (5Z,8Z)-tetradeca-5,8-dienoic-10,10,11,11,12,12,13,13,14,14,14-d₁₁ acid
Synonyms: C14:2(5Z,8Z)-d₁₁, C14:2 n-6-d₁₁, FA 14:2-d₁₁, Goshuyic Acid-d₁₁
MF: C₁₄H₁₃D₁₁O₂
FW: 235.4
Chemical Purity: ≥95% (5(Z),8(Z)-Tetradecadienoic Acid)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀
Supplied as: A 250 µg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Description

5(Z),8(Z)-Tetradecadienoic acid-d₁₁ is intended for use as an internal standard for the quantification of 5(Z),8(Z)-tetradecadienoic acid by GC- or LC-MS. 5(Z),8(Z)-Tetradecadienoic acid is an ω-6 polyunsaturated fatty acid (PUFA).

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, 04/06/2026

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