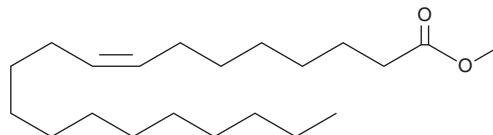


PRODUCT INFORMATION

8(Z)-Eicosenoic Acid methyl ester

Item No. 35428

CAS Registry No.: 69119-99-9
Formal Name: 8Z-eicosenoic acid, methyl ester
Synonyms: *cis*-8-Eicosenoic Acid methyl ester, SFE 21:1
MF: C₂₁H₄₀O₂
FW: 324.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

8(Z)-Eicosenoic acid methyl ester is an esterified form of 8(Z)-eicosenoic acid (Item No. 22579).¹ It has been used as a standard for the quantification of 8(Z)-eicosenoic acid by gas chromatography.

Reference

1. Stránský, K., Jursík, T., Vitek, A., *et al.* An improved method of characterizing fatty acids by equivalent chain length values. *J. High Resolut. Chrom.* **15(11)**, 730-740 (1992).

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, 02/13/2024

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