

PRODUCT INFORMATION



cis-9,10-Methyleneoctadecanoic Acid methyl ester

Item No. 24825

CAS Registry No.: 3971-54-8

Formal Name: (1R,2S)-*rel*-2-octyl-cyclopropaneoctanoic acid, methyl ester

Synonyms: Dihydrosterculic Acid methyl ester,
Methyl *cis*-9,10-Methyleneoctadecanoate,
SFE 20:1

MF: C₂₀H₃₈O₂

FW: 310.5

Purity: ≥98%

Supplied as: A liquid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

cis-9,10-Methyleneoctadecanoic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the *cis*-9,10-methyleneoctadecanoic acid methyl ester in the solvent of choice. *cis*-9,10-Methyleneoctadecanoic acid methyl ester is soluble in organic solvents such as chloroform, ethanol, methanol, and hexane, which should be purged with an inert gas.

Description

cis-9,10-Methyleneoctadecanoic acid methyl ester is a cyclopropane fatty acid methyl ester.¹ It stimulates DNA synthesis in rat hepatocytes. *cis*-9,10-Methyleneoctadecanoic acid methyl ester has been used as a standard for the quantification of *cis*-9,10-methyleneoctadecanoic acid (Item No. 24824) in bacterial membranes by co-chromatography.²

References

1. Scarpelli, D.G. Experimental modification of the mitogenic effect of methyl sterculate on rat liver. *Virchows Arch. B Cell Path.* **16(2)**, 211-220 (1974).
2. Montgomerie, J.Z., Kalmanson, G.M., and Guze, L.B. Fatty acid composition of membranes of L-forms of *Streptococcus faecalis* and *Proteus mirabilis* cultured at different osmolalities. *Curr. Microbiol.* **6(5)**, 301-303 (1981).

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, 04/05/2024

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