

PRODUCT INFORMATION



Jacaric Acid methyl ester

Item No. 27098

Formal Name: methyl (8Z,10E,12Z)-octadeca-8,10,12-trienoate

Synonyms: Methyl jacarate, Methyl 8(Z),10(E),12(Z)-Octadecatrienoate, SFE 19:3

MF: $C_{19}H_{32}O_2$

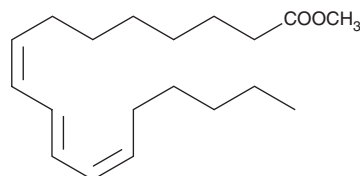
FW: 292.5

Purity: $\geq 96\%$

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

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

Description

Jacaric acid methyl ester is a methyl ester form of the conjugated PUFA jacaric acid (Item No. 16036).

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, 03/04/2024

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