

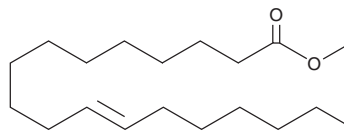
PRODUCT INFORMATION



trans-Vaccenic Acid methyl ester

Item No. 26731

CAS Registry No.: 6198-58-9
Formal Name: (11E)-11-octadecenoic acid methyl ester
Synonyms: C18:1 methyl ester, Methyl *trans*-11-Octadecenoate, Methyl *trans*-Vaccenate, *trans*-11-Octadecenoic Acid methyl ester, SFE 19:1
MF: C₁₉H₃₆O₂
FW: 296.5
Purity: ≥98%
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.

Laboratory Procedures

trans-Vaccenic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of *trans*-vaccenic acid methyl ester in these solvents is approximately 30 mg/ml.

Description

trans-Vaccenic acid methyl ester is an esterified form of *trans*-vaccenic acid (Item No. 15301). It has been used as a standard for the identification and quantification of fatty acids in Uruguayan Colonia cheese by GC-MS.¹

Reference

1. Hirigoyen, D., de los Santos, R., Calvo, M.F., *et al.* Chemical composition and seasonal changes in the fatty acid profile of Uruguayan "colonia" cheeses. *Grasas Y Aceites* **69**(2), e254 (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, 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