

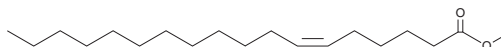
PRODUCT INFORMATION



cis-Petroselinic Acid methyl ester

Item No. 45681

CAS Registry No.: 2777-58-4
Formal Name: 6*Z*-octadecenoic acid, methyl ester
Synonyms: C18:1 methyl ester,
Methyl Octadecenoate,
cis-6-Octadecenoic
Acid methyl ester, SFE 19:1
MF: C₁₉H₃₆O₂
FW: 296.5
Purity: ≥98%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

cis-Petroselinic acid methyl ester is supplied as a liquid. A stock solution may be made by dissolving the *cis*-petroselinic acid methyl ester in the solvent of choice, which should be purged with an inert gas. *cis*-petroselinic acid methyl ester is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

cis-Petroselinic acid methyl ester is a long-chain fatty acid methyl ester and esterified form of *cis*-petroselinic acid (Item No. 20024). It has been found in *F. indica*.¹

Reference

1. Sulieman, A.M.E., Alanaizy, E., Alanaizy, N.A., *et al.* Unveiling chemical, antioxidant and antibacterial properties of *Fagonia indica* grown in the Hail Mountains, Saudi Arabia. *Plants (Basel)* **12**(6), 1354 (2023).

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, 07/06/2026

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