

PRODUCT INFORMATION



trans-Petroselinic Acid methyl ester

Item No. 36670

CAS Registry No.: 14620-36-1

Formal Name: 6E-octadecenoic acid, methyl ester

Synonyms: Δ^6 -*trans*-Octadecenoic Acid methyl ester,
trans-6-Octadecenoic Acid methyl ester,
Petroselaicid Acid methyl ester, SFE 19:1

MF: $C_{19}H_{36}O_2$

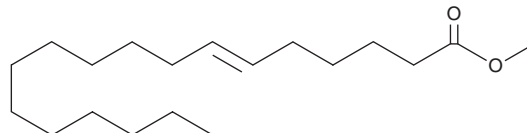
FW: 296.5

Purity: $\geq 95\%$

Supplied as: A neat 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

trans-Petroselinic acid methyl ester is supplied as a neat liquid. A stock solution may be made by dissolving the *trans*-petroselinic acid methyl ester in the solvent of choice, which should be purged with an inert gas. *trans*-Petroselinic acid methyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *trans*-petroselinic acid methyl ester in DMSO is approximately 14 mg/ml and approximately 33 mg/ml in ethanol and DMF.

Description

trans-Petroselinic acid methyl ester is a long-chain fatty acid methyl ester and esterified form of *trans*-petroselinic acid (Item No. 20026).

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, 02/14/2024

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