

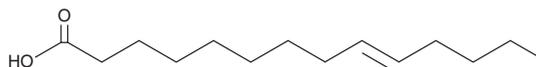
PRODUCT INFORMATION



Myristelaidic Acid

Item No. 22361

CAS Registry No.: 50286-30-1
Formal Name: (9E)-tetradecenoic acid
Synonyms: FA 14:1, *trans*-9-Tetradecenoic Acid
MF: C₁₄H₂₆O₂
FW: 226.4
Purity: ≥98%
UV/Vis.: λ_{max}: 201 nm
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

Myristelaidic acid is supplied as a solution in ethanol. A stock solution may be made by dissolving the myristelaidic acid in the solvent of choice, which should be purged with an inert gas. Myristelaidic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of myristelaidic acid in ethanol and DMF is approximately 15 mg/ml and approximately 12 mg/ml in DMSO.

Description

Myristelaidic acid is a 14-carbon monounsaturated ω-5 long-chain fatty acid. It is the *trans* isomer of myristoleic acid (Item No. 9002461). Myristelaidic acid is converted to myristelaidoyl CoA at a faster rate than myristoleic acid to myristoleoyl CoA.¹

Reference

1. Heuckeroth, R.O., Glaser, L., and Gordon, J.I. Heteroatom-substituted fatty acid analogs as substrates for N-myristoyltransferase: An approach for studying both the enzymology and function of protein acylation. *Proc. Natl. Acad. Sci. U.S.A.* **85(23)**, 8795-9799 (1988).

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, 01/25/2024

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