

PRODUCT INFORMATION



12-Tridecenoic Acid

Item No. 37601

CAS Registry No.: 6006-06-0

Synonyms: FA 13:1,
 ω -Tridecenoic acid

MF: $C_{13}H_{24}O_2$

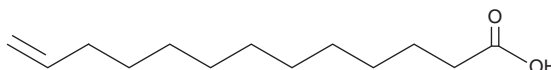
FW: 212.3

Purity: $\geq 95\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

12-Tridecenoic acid is supplied as a solid. A stock solution may be made by dissolving the 12-tridecenoic acid in the solvent of choice, which should be purged with an inert gas. 12-Tridecenoic acid is soluble in the organic solvent ethanol at a concentration of approximately 5 mg/ml.

Description

12-Tridecenoic acid is a monounsaturated fatty acid. It is active against *S. aureus* (MIC = 1.17 mM).¹ 12-Tridecenoic acid (5, 10, and 25 mg/kg) inhibits gastric juice secretion in a rat model of gastric ulcers induced by pylorus ligation.²

References

1. Zhang, H., Zhang, L., Peng, L.-j., *et al.* Quantitative structure-activity relationships of antimicrobial fatty acids and derivatives against *Staphylococcus aureus*. *J. Zhejiang Univ. Sci. B* **13**(2), 83-93 (2012).
2. Mimura, T., Kohda, I., Maeda, K., *et al.* Inhibitory effects of unsaturated fatty acids of *trans*-2- $C_{10:1}$ to *trans*-2- $C_{16:1}$ several $C_{18:n}$ and $C_{20:n}$ on gastric secretion and experimental ulceration in rats. *J. Pharmacobiodyn.* **6**(8), 527-538 (1983).

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, 02/20/2024

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