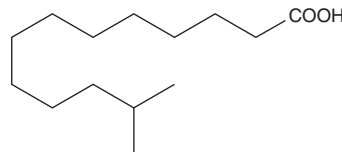


PRODUCT INFORMATION

12-methyl Tridecanoic Acid

Item No. 36461

CAS Registry No.: 2724-57-4
 Synonyms: FA 14:0, *iso*-14:0, *iso*-C14:0, 12-MTA
 MF: $C_{14}H_{28}O_2$
 FW: 228.4
 Purity: $\geq 98\%$
 Supplied as: A solid
 Storage: $-20^{\circ}C$
 Stability: ≥ 4 years



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

Laboratory Procedures

12-methyl Tridecanoic acid is supplied as a solid. A stock solution may be made by dissolving the 12-methyl tridecanoic acid in the solvent of choice, which should be purged with an inert gas. 12-methyl Tridecanoic acid is soluble in the organic solvent DMSO.

Description

12-methyl Tridecanoic acid is a methylated fatty acid that has been found in bovine milk.¹ Topical ocular application of 12-methyl tridecanoic acid (200 μM) decreases angiogenesis and corneal opacity in mouse models of alkali- or *P. aeruginosa*-induced eye injury.²

References

1. Ran-Ressler, R.R., Sim, D., O'Donnell-Megaro, A.M., *et al.* Branched chain fatty acid content of United States retail cow's milk and implications for dietary intake. *Lipids* **46**(7), 569-576 (2011).
2. Cole, N., Hume, E.B.H., Jalbert, I., *et al.* Effects of topical administration of 12-methyl tetradecanoic acid (12-MTA) on the development of corneal angiogenesis. *Angiogenesis* **10**(1), 47-54 (2007).

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