

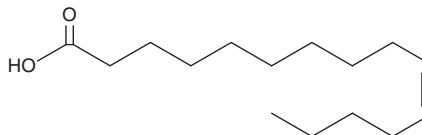
PRODUCT INFORMATION



10(Z)-Pentadecenoic Acid

Item No. 22362

CAS Registry No.: 84743-29-3
Formal Name: 10(Z)-pentadecenoic acid
Synonyms: FA 15:1, *cis*-10-Pentadecenoic acid
MF: C₁₅H₂₈O₂
FW: 240.4
Purity: ≥98%
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

10(Z)-Pentadecenoic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 10(Z)-pentadecenoic acid in these solvents is approximately 10 and 25 mg/ml, respectively.

10(Z)-Pentadecenoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 10(Z)-pentadecenoic acid should be diluted with the aqueous buffer of choice. 10(Z)-Pentadecenoic acid has a solubility of 0.25 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

10(Z)-Pentadecenoic acid is a 15-carbon, long-chain monounsaturated fatty acid. 10(Z)-Pentadecenoic acid inhibits IFN-γ-induced production of kynurenine in Thp-1 cells by 16%.¹

Reference

1. Costabile, M., Bassal, N.K., Gerber, J.P., *et al.* Inhibition of indoleamine 2,3-dioxygenase activity by fatty acids and prostaglandins: A structure function analysis. *Prostaglandins Leukot. Essent. Fatty Acids* **122**, 7-15 (2017).

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