

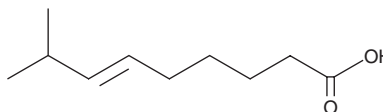
PRODUCT INFORMATION



(E)-8-Methyl-6-nonenoic Acid

Item No. 16609

CAS Registry No.: 59320-77-3
Formal Name: (6E)-8-methyl-6-nonenoic acid
Synonyms: trans-8-Methylnon-6-enoic Acid,
8-Methylnon-trans-6-enoic Acid
MF: C₁₀H₁₈O₂
FW: 170.3
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

(E)-8-Methyl-6-nonenoic acid is supplied as a neat oil. A stock solution may be made by dissolving the neat oil in an organic solvent purged with an inert gas. (E)-8-Methyl-6-nonenoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of (E)-8-methyl-6-nonenoic acid in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of (E)-8-methyl-6-nonenoic acid can be prepared by directly dissolving the neat oil in aqueous buffers. The solubility of (E)-8-methyl-6-nonenoic acid in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(E)-8-Methyl-6-nonenoic acid is a potential thermal degradant of capsaicin (Item No. 92350) with possible mutagenic effects.^{1,2}

References

1. Gannett, P.M., Nagel, D.L., Reilly, P.J., *et al.* The capsaicinoids: Their separation, synthesis, and mutagenicity. *J. Org. Chem.* **53(5)**, 1064-1071 (1988).
2. Hoffman, P.G., Lego, M.C., and Galetto, W.G. Separation and quantitation of red pepper major heat principles by reverse-phase high-pressure liquid chromatography. *J. Agric. Food Chem.* **31(6)**, 1326-1330 (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, 11/16/2022

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