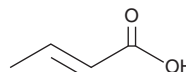


PRODUCT INFORMATION

Crotonic Acid

Item No. 43903

CAS Registry No.: 107-93-7
 Formal Name: 2E-butenoic acid
 Synonyms: *trans*-2-Butenoic acid, NSC 8751
 MF: C₄H₆O₂
 FW: 86.1
 Purity: ≥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

Crotonic acid is supplied as a solid. A stock solution may be made by dissolving the crotonic acid in the solvent of choice, which should be purged with an inert gas. Crotonic acid is soluble (≥10 mg/ml) in ethanol and DMSO.

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 crotonic acid can be prepared by directly dissolving the solid in aqueous buffers. Crotonic acid is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Crotonic acid is a short-chain unsaturated carboxylic acid that has been used as a building block in the synthesis of various small molecules and polymers.^{1,2}

References

- Chen, Z., Geng, D.-W., Yuan, T.-B., *et al.* Design and evaluation of novel N-substituent diphenylamine derivatives as tubulin colchicine binding site inhibitors. *Bioorg. Med. Chem. Lett.* **115**, 130031 (2025).
- Hu, X., Wang, Y., Zhang, L., *et al.* Design of a pH-sensitive magnetic composite hydrogel based on salean graft copolymer and Fe₃O₄@SiO₂ nanoparticles as drug carrier. *Int. J. Biol. Macromol.* **107(Pt B)**, 1811-1820 (2018).

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, 07/28/2025

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