

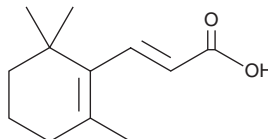
PRODUCT INFORMATION



β-Ionone Acid

Item No. 9004321

CAS Registry No.: 14393-45-4
Formal Name: (2E)-3-(2,6,6-trimethyl-1-cyclohexen-1-yl)-2-propenoic acid
MF: C₁₂H₁₈O₂
FW: 194.3
Purity: ≥98%
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

β-Ionone acid is supplied as a solid. A stock solution may be made by dissolving the β-ionone acid in the solvent of choice, which should be purged with an inert gas. β-Ionone acid is soluble in organic solvents such as acetonitrile, chloroform, and DMSO.

Description

β-Ionone acid is the acid form of the β-carotenoid derivative β-ionone.¹

Reference

1. Ansari, M. and Emami, S. β-Ionone and its analogs as promising anticancer agents. *Eur. J. Med. Chem.* **123**, 141-154 (2016).

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, 10/22/2025

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