

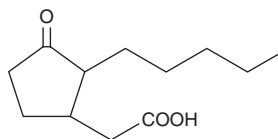
PRODUCT INFORMATION



Dihydrojasmonic Acid

Item No. 35362

CAS Registry No.: 3572-64-3
Formal Name: 3-oxo-2-pentyl-cyclopentaneacetic acid
Synonym: (±)-9,10-Dihydrojasmonic Acid
MF: $C_{12}H_{20}O_3$
FW: 212.3
Purity: $\geq 95\%$ (mixture of isomers)
Supplied as: A neat liquid
Storage: -20°C
Stability: ≥ 2 years
Item Origin: Synthetic



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

Laboratory Procedures

Dihydrojasmonic acid is supplied as a neat liquid. A stock solution may be made by dissolving the dihydrojasmonic acid in the solvent of choice, which should be purged with an inert gas. Dihydrojasmonic acid is soluble in chloroform.

Description

Dihydrojasmonic acid is a plant growth regulator.¹ It inhibits lettuce seed germination and reduces the radicle length of germinated seedlings when used at 10 mM.

Reference

1. Ravid, U., Ikan, R., and Sachs, R.M. Structures related to jasmonic acid and their effect on lettuce seedling growth. *J. Agric. Food Chem.* **23(5)**, 835-838 (1975).

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, 10/06/2022

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