

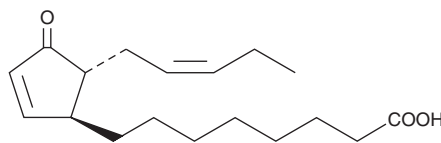
PRODUCT INFORMATION



13-*epi*-12-oxo Phytodienoic Acid

Item No. 10195

CAS Registry No.: 71606-07-0
Formal Name: 4-oxo-5R-(2Z)-2-pentenyl-2-cyclopentene-1S-octanoic acid
Synonym: 13-*epi*-12-oxo PDA
MF: C₁₈H₂₈O₃
FW: 292.4
Purity: ≥95%
UV/Vis.: λ_{max}: 220 nm
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

13-*epi*-12-oxo Phytodienoic acid (13-*epi*-12-oxo PDA) 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 13-*epi*-12-oxo PDA in these solvents is approximately 100, 10, and 20 mg/ml, respectively.

13-*epi*-12-oxo PDA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 13-*epi*-12-oxo PDA should be diluted with the aqueous buffer of choice. The solubility of 13-*epi*-12-oxo PDA in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

13-*epi*-12-oxo PDA is a lipoxygenase metabolite of α-linolenic acid in the leaves of green plants such as corn. ω-3 and ω-6 polyunsaturated fatty acids in plants are substrates for plant lipoxygenases.¹ 12-oxo PDA is one of the best studied end metabolites of this enzymatic pathway.² While the initial enzymatic product and major isomer of 12-oxo PDA contains side chains in the *cis* position, both being β to the ring, the upper side chain attached at C-13 can and frequently does isomerize when 12-oxo PDA is extracted, isolated or stored. 13-*epi*-12-oxo PDA is the product of this isomerization.

References

1. Crombie, L. and Mistry, K.M. Synthesis of 12-oxophytodienoic acid (12-oxo PDA) and the compounds of its enzymic degradation cascade in plants, OPC-8:0, -6:0, -4:0 and -2:0 (*epi*-jasmonic acid), as their methyl esters. *J. Chem. Soc. Perkin Trans. 1* **8**, 1981-1991 (1991).
2. Hamberg, M. and Gardner, H.W. Oxylipin pathway to jasmonates: Biochemistry and biological significance. *Biochim. Biophys. Acta* **1165**(1), 1-18 (1992).

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, 03/07/2024

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