

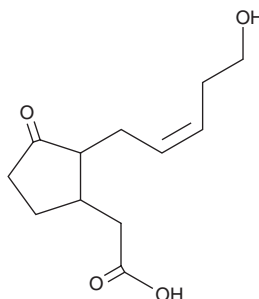
PRODUCT INFORMATION



(±)-Tuberonic Acid

Item No. 35750

CAS Registry No.: 221682-86-6
Formal Name: 2-[(Z)-5-hydroxy-2-penten-1-yl]-3-oxo-cyclopentaneacetic acid
Synonyms: 12-hydroxy Jasmonic Acid, 12-hydroxy *epi*-Jasmonic Acid
MF: C₁₂H₁₈O₄
FW: 226.3
Purity: ≥70%
Supplied as: A solution in methanol
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

(±)-Tuberonic acid is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. (±)-Tuberonic acid is slightly soluble in methanol.

Description

(±)-Tuberonic acid is a derivative of the plant growth regulator jasmonic acid (Item No. 88300) that has been found in *M. oleifera*.^{1,2}

References

1. Seto, Y., Hamada, S., Ito, H., *et al.* Tobacco salicylic acid glucosyltransferase is active toward tuberonic acid (12-hydroxyjasmonic acid) and is induced by mechanical wounding stress. *Biosci. Biotechnol. Biochem.* **75(12)**, 2316-2320 (2011).
2. Luetragoon, T., Sranujit, R.P., Noysang, C., *et al.* Bioactive compounds in *Moringa oleifera* Lam. leaves inhibit the pro-inflammatory mediators in Lipopolysaccharide-induced human monocyte-derived macrophages. *Molecules* **25(1)**, 191 (2020).

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, 01/03/2023

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