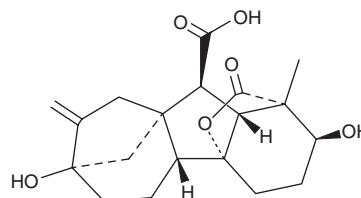


PRODUCT INFORMATION



Gibberellin A₁ Item No. 34291

CAS Registry No.: 545-97-1
Formal Name: (1 α ,2 β ,4 α ,4 β ,10 β)-2,4a,7-trihydroxy-1-methyl-8-methylene-gibbane-1,10-dicarboxylic acid, 1,4a-lactone
Synonym: GA₁
MF: C₁₉H₂₄O₆
FW: 348.4
Purity: $\geq 90\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

Gibberellin A₁ is supplied as a solid. A stock solution may be made by dissolving the gibberellin A₁ in the solvent of choice, which should be purged with an inert gas. Gibberellin A₁ is soluble in methanol, DMSO, and a 1:1 solution of acetone:water. We do not recommend storing the aqueous solution for more than one day.

Description

Gibberellin A₁ is a plant hormone that has been found in *Phaseolus multiflorus* and has growth stimulatory activities.¹⁻³ It reverses inhibition of spinach stem elongation induced by the plant growth retardants AMO-1618 or BX-112 (prohexadione; Item No. 33922).² Gibberellin A₁ increases hypocotyl, but not radicle, growth of cucumber seedlings when used at concentrations of 30 and 100 μM .³

References

1. MacMillan, J., Seaton, J.C., and Suter, P.J. Plant hormones—I: Isolation of gibberellin A₁ and gibberellin A₅ from *Phaseolus multiflorus*. *Tetrahedron* **11**(1-2), 60-66 (1960).
2. Zeevaart, J.A., Gage, D.A., and Talon, M. Gibberellin A₁ is required for stem elongation in spinach. *Proc. Natl. Acad. Sci. USA* **90**(15), 7401-7405 (1993).
3. Halevy, A.H. and Cathey, H.M. Effects of structure and concentration of gibberellins on the growth of cucumber seedlings. *Bot. Gaz.* **122**(1), 63-67 (1960).

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, 12/13/2022

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