

PRODUCT INFORMATION



22(S),23(S)-Homobrassinolide

Item No. 32884

CAS Registry No.: 80483-89-2
Formal Name: (1R,3aS,3bS,6aS,8S,9R,10aR,10bS,12aS)-1-
[[1S,2S,3S,4S)-4-ethyl-2,3-dihydroxy-1,5-
dimethylhexyl]hexadecahydro-8,9-dihydroxy-
10a,12a-dimethyl-6H-benz[c]indeno[5,4-e]
oxepin-6-one

Synonyms: (22S,23S)-28-Homobrassinolide,
Isohomobrassinolide

MF: C₂₉H₅₀O₆

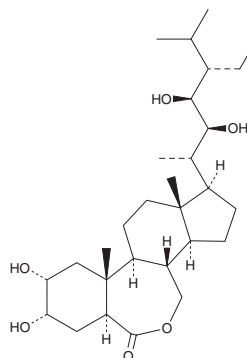
FW: 494.7

Purity: ≥90%

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

22(S),23(S)-Homobrassinolide is supplied as a solid. A stock solution may be made by dissolving the 22(S),23(S)-homobrassinolide in the solvent of choice, which should be purged with an inert gas. 22(S),23(S)-Homobrassinolide is soluble in the organic solvent dimethyl formamide at a concentration of approximately 5 mg/ml. 22(S),23(S)-Homobrassinolide is slightly soluble in DMSO.

Description

22(S),23(S)-Homobrassinolide is a synthetic brassinosteroid.^{1,2} It increases the rooting percentage in *P. abies* cuttings when applied topically at a concentration of 60 ppm.¹

References

1. Rönsch, H., Adam, G., Matschke, J., *et al.* Influence of (22S,23S)-homobrassinolide on rooting capacity and survival of adult Norway spruce cuttings. *Tree Physiol.* **12(1)**, 71-80 (1993).
2. McMorris, T.C., Patil, P.A., Chavez, R.G., *et al.* Synthesis and biological activity of 28-homobrassinolide and analogues. *Phytochemistry* **36(3)**, 585-589 (1994).

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