

PRODUCT INFORMATION



Homobrassinolide

Item No. 36038

CAS Registry No.: 74174-44-0
Formal Name: (1R,3aS,3bS,6aS,8S,9R,10aR,10bS,12aS)-1-[(1S,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

Synonym: 28-Highbrassinolide

MF: C₂₉H₅₀O₆

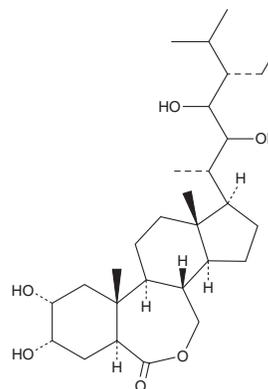
FW: 494.7

Purity: ≥85%

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

Homobrassinolide is supplied as a solid. A stock solution may be made by dissolving the homobrassinolide in the solvent of choice, which should be purged with an inert gas. Homobrassinolide is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of homobrassinolide in these solvents is approximately 5 and 1 mg/ml, respectively. Homobrassinolide is slightly soluble in ethanol.

Description

Homobrassinolide is a brassinosteroid.¹ Field application of homobrassinolide (0.12 g/ha) increases the number of fruits per plant, fruit length, and fruit breadth in tomato plants.

Reference

1. Sridhara, S., Ramesh, N., Gopakkali, P., *et al.* Application of homobrassinolide enhances growth, yield and quality of tomato. *Saudi J. Biol. Sci.* **28(8)**, 4800-4806 (2021).

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, 12/19/2022

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