

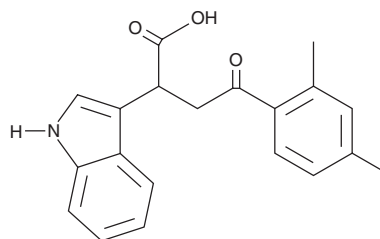
PRODUCT INFORMATION



Auxinole

Item No. 40334

CAS Registry No.: 86445-22-9
Formal Name: α -[2-(2,4-dimethylphenyl)-2-oxoethyl]-1H-indole-3-acetic acid
MF: C₂₀H₁₉NO₃
FW: 321.4
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Auxinole is supplied as a solid. A stock solution may be made by dissolving the auxinole in the solvent of choice, which should be purged with an inert gas. Auxinole is soluble in acetonitrile and DMSO.

Description

Auxinole is an antagonist of the auxin receptor transport inhibitor response 1 (TIR1).¹ It inhibits indole 3-acetic acid-induced root shoot growth in *Arabidopsis* seedlings when used at a concentration of 20 μ M. Auxinole (5 or 10 μ M) prevents indole 3-acetic acid-induced increases in root hair growth, root gravitropism, the number of lateral roots, and indole 3-acetic acid-induced decreases in primary root length in *Arabidopsis* seedlings. It reduces the number of programmed cell death-induced perforations in *A. madagascariensis* leaves when used at a concentration of 25 μ M.²

References

- Hayashi, K., Neve, J., Hirose, M., *et al.* Rational design of an auxin antagonist of the SCF(TIR1) auxin receptor complex. *ACS Chem. Biol.* **7(3)**, 590-598 (2012).
- Denbigh, G.L., Dauphinee, A.N., Fraser, M.S., *et al.* The role of auxin in developmentally regulated programmed cell death in lace plant. *Am. J. Bot.* **107(4)**, 577-586 (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, 03/08/2024

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