

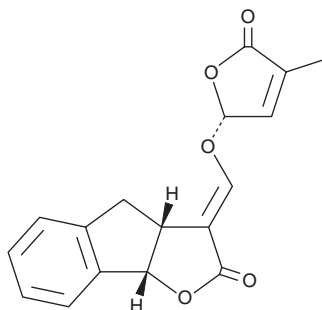
PRODUCT INFORMATION



GR24

Item No. 13210

CAS Registry No.: 76974-79-3
Formal Name: (3E,3aR,8bS)-rel-3-[[[(2R)-2,5-dihydro-4-methyl-5-oxo-2-furanyl]oxy]methylene]-3,3a,4,8b-tetrahydro-2H-indeno[1,2-b]furan-2-one
MF: C₁₇H₁₄O₅
FW: 298.3
Purity: ≥95%
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

GR24 is supplied as a solid. A stock solution may be made by dissolving the GR24 in the solvent of choice, which should be purged with an inert gas. GR24 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

GR24 is a synthetic strigolactone and plant growth regulator.^{1,2} It stimulates the germination of *Striga* and *Orobanch*e parasitic weed seeds (EC₅₀s = 5-50 nM).¹ GR24 increases root-hair length and decreases the number of lateral roots in *Arabidopsis*.²

References

1. Wigchert, S.C., Kuiper, E., Boelhouwer, G.J., *et al.* Dose-response of seeds of the parasitic weeds *Striga* and *Orobanch*e toward the synthetic germination stimulants GR 24 and Nijmegen 1. *J. Agric. Food Chem.* **47**(4), 1705-1710 (1999).
2. Kapulnik, Y., Delaux, P.M., Resnick, N., *et al.* Strigolactones affect lateral root formation and root-hair elongation in *Arabidopsis*. *Planta* **233**(1), 209-216 (2011).

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, 10/24/2022

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