

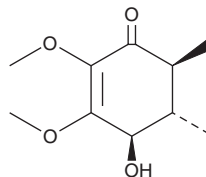
PRODUCT INFORMATION



Glioroseinol

Item No. 31679

Formal Name: 4S-hydroxy-2,3-dimethoxy-5S,6S-dimethylcyclohex-2-en-1-one
MF: C₁₀H₁₆O₄
FW: 200.2
Purity: ≥95%
Supplied as: A residue
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Gliocladium* sp.



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

Laboratory Procedures

Glioroseinol is supplied as a residue. A stock solution may be made by dissolving the glioroseinol in the solvent of choice, which should be purged with an inert gas. Glioroseinol is soluble in methanol and DMSO.

Description

Glioroseinol is a fungal metabolite and derivative of gliorosein (Item No. 24921) that has been found in *Gliocladium*.

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, 08/11/2023

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