

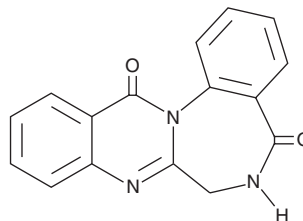
PRODUCT INFORMATION



Sclerotigenin

Item No. 23492

CAS Registry No.: 65641-84-1
Formal Name: 6,7-dihydro-quinazolino[3,2-a][1,4]benzodiazepine-5,13-dione
Synonym: (±)-Sclerotigenin
MF: C₁₆H₁₁N₃O₂
FW: 277.3
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

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

Description

Sclerotigenin is a natural diazepine originally isolated from *P. sclerotigenum* that acts as an insect growth regulator.¹ It reduces the growth rate of first instar *H. zea* larvae by 42% relative to controls when administered in the standard test diet at 200 ppm.

Reference

1. Joshi, B.K., Gloer, J.B., Wicklow, D.T., *et al.* Sclerotigenin: A new antiinsectan benzodiazepine from the sclerotia of *penicillium sclerotigenum*. *J. Nat. Prod.* **62**(4), 650-652 (1999).

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, 11/16/2022

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