

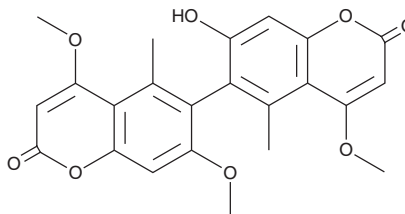
PRODUCT INFORMATION



Isokotanin B

Item No. 25082

CAS Registry No.: 154160-09-5
Formal Name: (+)-7-hydroxy-4,4',7'-trimethoxy-5,5'-dimethyl-[6,6'-bi-2H-1-benzopyran]-2,2'-dione
MF: $C_{23}H_{20}O_8$
FW: 424.4
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 215, 326 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungi/*Aspergillus* sp.



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

Laboratory Procedures

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

Description

Isokotanin B is a bicoumarin fungal metabolite originally isolated from *A. alliaceus*.¹ Dietary administration of isokotanin B (100 ppm) reduces *C. hemipterus* larvae feeding by 21%.

Reference

1. Laakso, J.A., Narske, E.D., Gloer, J.B., et al. Isokotanins A-C: New bicoumarins from the sclerotia of *Aspergillus alliaceus*. *J. Nat. Prod.* **57**(1), 128-133 (1994).

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, 12/20/2022

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