

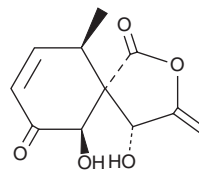
PRODUCT INFORMATION



Massarigenin C

Item No. 29772

CAS Registry No.: 496926-08-0
Formal Name: (4R,5R,6R,10R)-*rel*-(-)-4,6-dihydroxy-10-methyl-3-methylene-2-oxaspiro[4.5]dec-8-ene-1,7-dione
MF: C₁₁H₁₂O₅
FW: 224.2
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

Massarigenin C is supplied as a solid. A stock solution may be made by dissolving the massarigenin C in the solvent of choice, which should be purged with an inert gas. Massarigenin C is soluble in ethanol, methanol, DMSO, and dichloromethane.

Description

Massarigenin C is a fungal metabolite that has been found in *M. flavorosea* and has enzyme inhibitory activities.^{1,2} Massarigenin C inhibits neuraminidase *in vitro* (IC₅₀ = 4.15 μM).² It is also an inhibitor of yeast α-glucosidase (IC₅₀ = 1.25 mM).¹ It reduces the postprandial peak in blood glucose levels in an oral sucrose tolerance test in normo- and hyperglycemic mice when administered at doses of 3.2, 10, and 31.6 mg/kg.

References

1. Verastegui-Omaña, B., Rebollar-Ramos, D., Pérez-Vásquez, A., *et al.* α-Glucosidase inhibitors from *Malbranchea flavorose*. *J. Nat. Prod.* **80**(1), 190-195 (2017).
2. Zhang, G.F., Han, W.B., Cui, J.T., *et al.* Neuraminidase inhibitory polyketides from the marine-derived fungus *Phoma herbaru*. *Planta Med.* **78**(1), 76-78 (2012).

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, 08/16/2023

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