

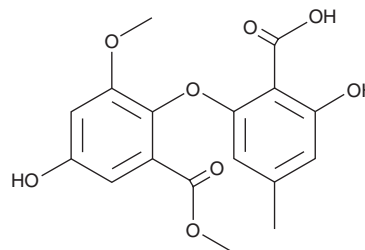
# PRODUCT INFORMATION



## Asterric Acid

Item No. 70630

**CAS Registry No.:** 577-64-0  
**Formal Name:** 2-(2-carboxy-3-hydroxy-5-methylphenoxy)-5-hydroxy-3-methoxy-benzoic acid, 1-methyl ester  
**Synonyms:** Dimethylsoic Acid, RES-1214-1, TAN 1415A  
**MF:** C<sub>17</sub>H<sub>16</sub>O<sub>8</sub>  
**FW:** 348.3  
**Purity:** ≥99%  
**UV/Vis.:** λ<sub>max</sub>: 214, 252, 315 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Aspergillus terreus*



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

### Laboratory Procedures

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

### Description

Asterric acid is an antibiotic fungal metabolite that completely inhibits the binding of the potent vasoconstrictor endothelin (ET)-1 to the ET<sub>A</sub> receptor in A10 cells at 0.1 μM.<sup>1</sup> Asterric acid derivatives have also been shown to inhibit VEGF-induced tube formation of human umbilical vein endothelial cells at 3-10 μM, which suggests its usefulness as an antiangiogenic agent.<sup>2</sup>

### References

1. Ohashi, H., Akiyama, H., Nishikori, K., *et al.* Asterric acid, a new endothelin binding inhibitor. *J. Antibiot.* **45**, 1684-1685 (1992).
2. Lee, H.J., Lee, J.H., Hwang, B.Y., *et al.* Fungal metabolites, asterric acid derivatives inhibit vascular endothelial growth factor (VEGF)-induced tube formation of HUVECs. *J. Antibiot. (Tokyo)* **55(6)**, 552-556 (2002).

#### 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, 08/25/2023

#### 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