

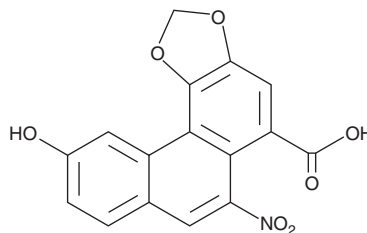
# PRODUCT INFORMATION



## Aristolochic Acid C

Item No. 28505

**CAS Registry No.:** 4849-90-5  
**Formal Name:** 10-hydroxy-6-nitro-phenanthro[3,4-d]-1,3-dioxole-5-carboxylic acid  
**Synonym:** Aristolochic Acid IIIa  
**MF:** C<sub>16</sub>H<sub>9</sub>NO<sub>7</sub>  
**FW:** 327.2  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/Aristolochiaceae



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

### Laboratory Procedures

Aristolochic acid C is supplied as a solid. A stock solution may be made by dissolving the aristolochic acid C in the solvent of choice, which should be purged with an inert gas. Aristolochic acid C is slightly soluble in methanol (warmed) and DMSO.

### Description

Aristolochic acid C is an alkaloid originally isolated from *Aristolochia argentina* and a derivative of aristolochic acid (Item No. 21520) that has anticancer activities.<sup>1-3</sup> It inhibits the growth of LLC-PK1, P388, HT-29, and HL-60 cancer cells (IC<sub>50</sub>s = 85, 8.8, 17.1, and 40 μM, respectively).<sup>1,2</sup>

### References

1. Cai, Y. and Cai, T.-G. Two new aristolochic acid derivatives from the roots of *Aristolochia fangchi* and their cytotoxicities. *Chem. Pharm. Bull. (Tokyo)* **58(8)**, 1093-1095 (2010).
2. Wu, T.-S., Chan, Y.-Y., and Leu, Y.-L. The constituents of the root and stem of *Aristolochia cucurbitifolia* Hayata and their biological activity. *Chem. Pharm. Bull. (Tokyo)* **48(7)**, 1006-1009 (2000).
3. Kumar, V., Poonam, Prasad, A.K., et al. Naturally occurring aristolactams, aristolochic acids and dioxoaporphines and their biological activities. *Nat. Prod. Rep.* **20(6)**, 565-583 (2003).

#### 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, 11/07/2022

#### 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