

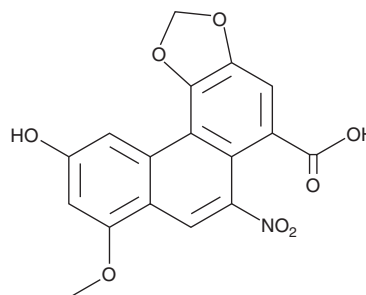
PRODUCT INFORMATION



Aristolochic Acid D

Item No. 35492

CAS Registry No.: 17413-38-6
Formal Name: 10-hydroxy-8-methoxy-6-nitro-phenanthro
[3,4-d]-1,3-dioxole-5-carboxylic acid
Synonym: Aristolochic Acid IVa
MF: C₁₇H₁₁NO₈
FW: 357.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Aristolochia manshuriensis*



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

Laboratory Procedures

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

Description

Aristolochic acid D is an alkaloid and a derivative of aristolochic acid (Item No. 21520) originally isolated from *Aristolochia indica*.¹

Reference

1. Kupchan, S.M. and Merianos, J.J. The isolation and structural elucidation of novel derivatives of aristolochic acid from *Aristolochia indica*. *J. Org. Chem.* **33**(10), 3735-3738 (1968).

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, 12/12/2022

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