

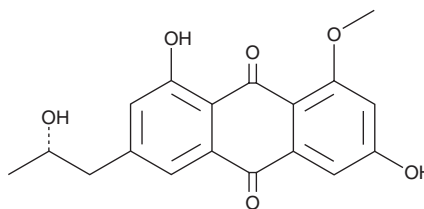
# PRODUCT INFORMATION



## Penipurdin A

Item No. 31657

CAS Registry No.: 1821668-16-9  
Formal Name: 1,6-dihydroxy-3-[(2S)-2-hydroxypropyl]-  
8-methoxy-9,10-anthracenedione  
MF:  $C_{18}H_{16}O_6$   
FW: 328.3  
Purity:  $\geq 95\%$   
Supplied as: A residue  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Fungus/*Penicillium* sp.



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

### Laboratory Procedures

Penipurdin A is supplied as a residue. A stock solution may be made by dissolving the penipurdin A in the solvent of choice, which should be purged with an inert gas. Penipurdin A is soluble in organic solvents such as methanol and DMSO.

### Description

Penipurdin A is an anthraquinone fungal metabolite originally isolated from *P. purpurogenum* that has antibacterial activity.<sup>1,2</sup> It is active against *M. tuberculosis* with a MIC value of 25  $\mu\text{g/ml}$ .<sup>2</sup>

### References

1. Xue, J., Fu, Y., Wu, P., *et al.* Two new anthraquinones from the soil fungus *Penicillium purpurogenum* SC0070. *J. Antibiot. (Tokyo)* **68**(9), 598-599 (2015).
2. Daengrot, C., Rukachaisirikul, V., Tadpetch, K., *et al.* Penicillanthone and penicillidic acids A-C from the soil-derived fungus *Penicillium aculeatum* PSU-RSPG105. *RSC Adv.* **6**(46), 39700-39709 (2016).

#### 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, 08/10/2023

#### 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