

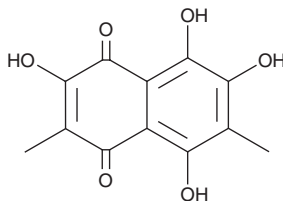
PRODUCT INFORMATION



Aureoquinone

Item No. 45482

CAS Registry No.: 321658-21-3
Formal Name: 2,5,7,8-tetrahydroxy-3,6-dimethyl-1,4-naphthalenedione
MF: $C_{12}H_{10}O_6$
FW: 250.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*Talaromyces* sp.



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

Laboratory Procedures

Aureoquinone is supplied as a solid. A stock solution may be made by dissolving the aureoquinone in the solvent of choice, which should be purged with an inert gas. Aureoquinone is slightly soluble (0.1-1 mg/ml) in methanol.

Description

Aureoquinone is a fungal metabolite originally isolated from *Aureobasidium* with protease inhibitory and antibacterial activity.¹ It inhibits trypsin, papain, thermolysin, collagenase, and zinc protease (IC_{50} s = 11.4, 14.5, 17.8, 7.1, and 8.7 $\mu\text{g/ml}$, respectively). Aureoquinone is active against *B. subtilis* (MIC = 120 $\mu\text{g/ml}$).

Reference

1. Berg, A., Görls, H., Dörfelt, H., *et al.* Aureoquinone, a new protease inhibitor from *Aureobasidium* sp. *J. Antibiot. (Tokyo)* **53**(11), 1293-1295.

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, 04/02/2026

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