

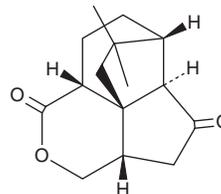
PRODUCT INFORMATION



Quadrone

Item No. 26688

CAS Registry No.: 66550-08-1
Formal Name: (3aS,5aR,6R,8aR,8bR)-octahydro-10,10-dimethyl-6,8b-ethano-8bH-cyclopenta[de]-2-benzopyran-1,4-dione
Synonym: (-)-Quadrone
MF: C₁₅H₂₀O₃
FW: 248.3
Purity: ≥98%
Supplied as: A residue
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Quadrone is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Quadrone is a sesquiterpene originally isolated from *A. terreus*.^{1,2} It is cytotoxic to human carcinoma KB cells *in vitro* (EC₅₀ = 1.3 µg/ml).¹

References

1. Ranieri, R.L. and Calton, G.J. Quadrone, a new antitumor agent from *Aspergillus terreus*. *Tetrahedron Lett.* **19(6)**, 499-502 (1978).
2. Calton, G.J., Ranieri, R.L., and Espenshade, M.A. Quadrone, a new antitumor substance produced by *Aspergillus terreus*. Production, isolation and properties. *J. Antibiot. (Tokyo)* **31(1)**, 38-42 (1978).

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, 08/08/2023

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