

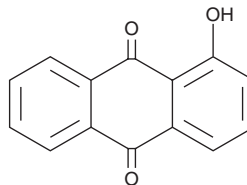
PRODUCT INFORMATION



1-Hydroxyanthraquinone

Item No. 32905

CAS Registry No.: 129-43-1
Formal Name: 1-hydroxy-9,10-anthracenedione
Synonym: NSC 8640
MF: C₁₄H₈O₃
FW: 224.2
Purity: ≥95%
UV/Vis.: λ_{max}: 217, 250 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

1-Hydroxyanthraquinone is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-hydroxyanthraquinone in the solvent of choice, which should be purged with an inert gas. 1-Hydroxyanthraquinone is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

1-Hydroxyanthraquinone is an anthraquinone that has been found in *Morinda officinalis* and has genotoxic and carcinogenic activities.^{1,2} 1-Hydroxyanthraquinone (5 and 500 μM) increases the percentage of isolated rat hepatocytes undergoing unscheduled DNA synthesis, indicating genotoxic activity.¹ Dietary administration of 1-hydroxyanthraquinone (1%) induces formation of cecum and colon adenocarcinomas, as well as hepatocellular carcinomas, in rats.

References

1. Kawai, K., Mori, H., Sugie, S., *et al.* Genotoxicity in the hepatocyte/DNA repair test and toxicity to liver mitochondria of 1-hydroxyanthraquinone and several dihydroxyanthraquinones. *Cell Biol. Toxicol.* **2(4)**, 457-467 (1986).
2. Mori, H., Yoshimi, N., Iwata, H., *et al.* Carcinogenicity of naturally occurring 1-hydroxyanthraquinone in rats: Induction of large bowel, liver and stomach neoplasms. *Carcinogenesis* **11(5)**, 799-802 (1990).

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, 11/16/2022

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