

# PRODUCT INFORMATION

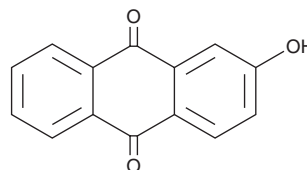


## 2-Hydroxyanthraquinone

Item No. 32906

**CAS Registry No.:** 605-32-3  
**Formal Name:** 2-hydroxy-9,10-anthracenedione  
**Synonyms:** 2-hydroxy-9,10-Anthraquinone,  
2-hATQ,  $\beta$ -Hydroxyanthraquinone,  
NSC 2595

**MF:** C<sub>14</sub>H<sub>8</sub>O<sub>3</sub>  
**FW:** 224.2  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 246, 272 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Plant/*Spermacoce latifolia*



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

### Laboratory Procedures

2-Hydroxyanthraquinone is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxyanthraquinone in the solvent of choice, which should be purged with an inert gas. 2-Hydroxyanthraquinone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2-hydroxyanthraquinone in these solvents is approximately 30 and 5 mg/ml, respectively.

### Description

2-Hydroxyanthraquinone is an anthraquinone that has been found in *Spermacoce latifolia* and has antibacterial and estrogenic activities.<sup>1,2</sup> It is active against *B. subtilis* and *B. cereus* (MICs = 1.9 and 62.5  $\mu$ g/ml, respectively).<sup>1</sup> 2-Hydroxyanthraquinone (19  $\mu$ M) induces estrogen receptor  $\alpha$  (ER $\alpha$ ) activation in a yeast two-hybrid assay.<sup>2</sup>

### References

1. Luo, Y., Shen, H.-Y., Shen, Q.-X., *et al.* A new anthraquinone and a new naphthoquinone from the whole plant of *Spermacoce latifolia*. *J. Asian Nat. Prod. Res.* **19(9)**, 869-876 (2017).
2. Kurihara, R., Shiraishi, F., Tanaka, N., *et al.* Presence and estrogenicity of anthracene derivatives in coastal Japanese waters. *Environ. Toxicol. Chem.* **24(8)**, 1984-1993 (2005).

#### 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, 10/03/2022

#### 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