

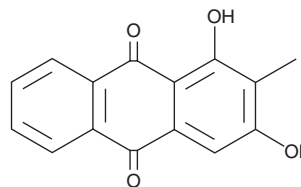
# PRODUCT INFORMATION



## Rubiadin

Item No. 11753

**CAS Registry No.:** 117-02-2  
**Formal Name:** 1,3-dihydroxy-2-methyl-9,10-anthracenedione  
**Synonym:** C.I. 75350  
**MF:** C<sub>15</sub>H<sub>10</sub>O<sub>4</sub>  
**FW:** 254.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 245, 279, 411 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Rubiadin is supplied as a crystalline solid. A stock solution may be made by dissolving the rubiadin in the solvent of choice, which should be purged with an inert gas. Rubiadin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of rubiadin in these solvents is approximately 0.5 mg/ml.

### Description

Rubiadin is an anthraquinone that is a minor component of certain *Rubia* species, as well as other plants. It demonstrates antioxidant activity.<sup>1</sup>

### Reference

1. Tripathi, Y.B., Sharma, M., and Manickam, M. Rubiadin, a new antioxidant from *Rubia cordifolia*. *Indian J. Biochem. Biophys.* **34(3)**, 302-306 (1997).

#### 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, 11/29/2022

#### 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