

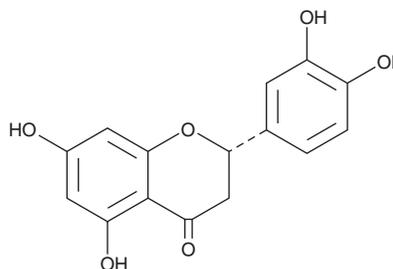
# PRODUCT INFORMATION



## 2S-Eriodictyol

Item No. 26803

**CAS Registry No.:** 552-58-9  
**Formal Name:** (2S)-2-(3,4-dihydroxyphenyl)-2,3-dihydro-5,7-dihydroxy-4H-1-benzopyran-4-one  
**Synonym:** (+)-Eriodictyol  
**MF:** C<sub>15</sub>H<sub>12</sub>O<sub>6</sub>  
**FW:** 288.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 289 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/Unidentified sp.



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

### Laboratory Procedures

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

2S-Eriodictyol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2S-eriodictyol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2S-Eriodictyol has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2S-Eriodictyol a flavanone that has been found in *Citrus* and has antioxidant activity.<sup>1</sup> It is protective against hydrogen peroxide-induced cell injury and apoptosis in EA.hy926 cells when used at a concentration of 20 μM.

### Reference

1. Li, H., Li, C., Shen, T., *et al.* R-eriodictyol and S-eriodictyol exhibited comparable effect against H<sub>2</sub>O<sub>2</sub>-induced oxidative stress in EA.hy926 cells. *Drug Discov. Ther.* **8(5)**, 218-224 (2014).

#### 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, 09/14/2023

#### 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