

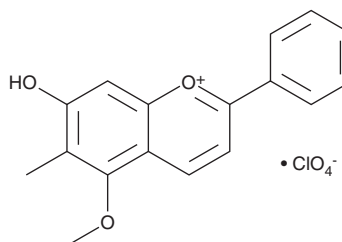
# PRODUCT INFORMATION



## Dracorhodin Perchlorate

Item No. 27294

**CAS Registry No.:** 125536-25-6  
**Formal Name:** 7-hydroxy-5-methoxy-6-methyl-2-phenyl-1-benzopyrylium, monopерchlorate  
**MF:** C<sub>17</sub>H<sub>15</sub>O<sub>3</sub> • ClO<sub>4</sub>  
**FW:** 366.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 272, 309, 445 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

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

Dracorhodin perchlorate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dracorhodin perchlorate should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dracorhodin perchlorate has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Dracorhodin perchlorate is a pro-apoptotic compound and a synthetic analog of dracorhodin, an anthocyanin that has been found in *D. draco* fruit.<sup>1</sup> Dracorhodin perchlorate is cytotoxic to U937, A375-S2, and MCF-7 cells in a concentration-dependent manner.<sup>1</sup> It is also cytotoxic to HeLa cells, an effect that can be reversed by Ac-YVAD-CMK (Item No. 10014), Z-DEVD-FMK (Item No. 14414), Z-IETD-FMK, and Z-LEHD-FMK, which are inhibitors of caspase-1, -3, -8, and -9, respectively.

### References

1. Xia, M., Wang, D., Wang, M., *et al.* Dracorhodin perchlorate induces apoptosis via activation of caspases and generation of reactive oxygen species. *J. Pharmacol. Sci.* **95(2)**, 273-283 (2004).

#### 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, 04/06/2023

#### 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