

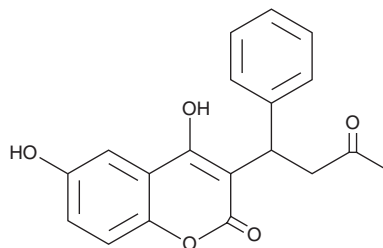
# PRODUCT INFORMATION



## 6-hydroxy Warfarin

Item No. 9000632

**CAS Registry No.:** 17834-02-5  
**Formal Name:** 4,6-dihydroxy-3-(3-oxo-1-phenylbutyl)-2H-1-benzopyran-2-one  
**MF:** C<sub>19</sub>H<sub>16</sub>O<sub>5</sub>  
**FW:** 324.3  
**Purity:** ≥98%  
**Supplied as:** A 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

6-hydroxy Warfarin is supplied as a solid. A stock solution may be made by dissolving the 6-hydroxy warfarin in the solvent of choice, which should be purged with an inert gas. 6-hydroxy Warfarin is soluble in organic solvents such as DMSO and methanol.

### Description

6-hydroxy Warfarin is a metabolite of (+)-warfarin (Item No. 13526), which is a weaker vitamin K antagonist than (-)-warfarin (Item No. 13531).<sup>1-3</sup> Both isomers of warfarin are metabolized by the cytochrome P450 isoform 2C9 (CYP2C9); however, stereoselective metabolism also occurs, with CYP1A1 and CYP1A2 primarily responsible for metabolism of (+)-warfarin to 6-hydroxy warfarin.<sup>1,2</sup>

### References

1. Zhang, Z., Fasco, M.J., Huang, Z., *et al.* Human cytochromes P4501A1 and P4501A2: R-warfarin metabolism as a probe. *Drug. Metab. Dispos.* **23(12)**, 1339-1346 (1995).
2. Kaminsky, L.S., and Zhang, Z.-Y. Human P450 metabolism of warfarin. *Pharmacol. Ther.* **73(1)**, 67-74 (1997).
3. Hirsh, J., Fuster, V., Ansell, J., *et al.* American Heart Association/American College of Cardiology Foundation guide to warfarin therapy. *J. Am. Coll. Cardiol.* **41(9)**, 1633-1652 (2003).

#### 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, 12/13/2022

#### 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