

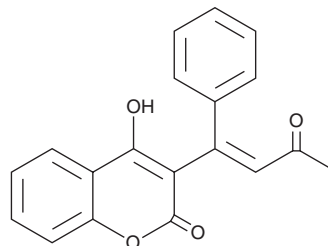
# PRODUCT INFORMATION



## Dehydro Warfarin

Item No. 33995

**CAS Registry No.:** 67588-18-5  
**Formal Name:** 4-hydroxy-3-(3-oxo-1-phenyl-1-buten-1-yl)-2H-1-benzopyran-2-one  
**Synonym:** NSC 289346  
**MF:** C<sub>19</sub>H<sub>14</sub>O<sub>4</sub>  
**FW:** 306.3  
**Purity:** ≥95%  
**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

Dehydro warfarin is supplied as a solid. A stock solution may be made by dissolving the dehydro warfarin in the solvent of choice, which should be purged with an inert gas. Dehydro warfarin is slightly soluble in methanol and chloroform.

### Description

Dehydro warfarin is a metabolite of (±)-warfarin (Item No. 13566).<sup>1</sup> It is formed from (±)-warfarin by rat liver microsomes.

### Reference

1. Kaminsky, L.S., Fasco, M.J., and Guengerich, F.P. Comparison of different forms of liver, kidney, and lung microsomal cytochrome P-450 by immunological inhibition of regio- and stereoselective metabolism of warfarin. *J. Biol. Chem.* **254**(19), 9657-9662 (1979).

#### 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, 11/17/2022

#### 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