

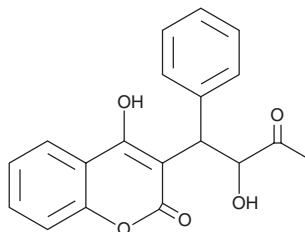
PRODUCT INFORMATION



10-hydroxy Warfarin

Item No. 33929

CAS Registry No.: 83219-99-2
Formal Name: 4-hydroxy-3-(2-hydroxy-3-oxo-1-phenylbutyl)-2H-1-benzopyran-2-one
MF: C₁₉H₁₆O₅
FW: 324.3
Purity: ≥95% (mixture of diastereomers)
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

10-hydroxy Warfarin is supplied as a solid. A stock solution may be made by dissolving the 10-hydroxy warfarin in the solvent of choice, which should be purged with an inert gas. 10-hydroxy Warfarin is slightly soluble in methanol and DMSO.

Description

10-hydroxy Warfarin is a metabolite of (R)-warfarin (Item No. 13526).¹ It is formed from (R)-warfarin by the cytochrome P450 (CYP) isoform CYP3A4. 10-hydroxy Warfarin is an inhibitor of CYP2C9 (IC₅₀ = 1.6 μM), the enzyme that converts (S)-warfarin (Item No. 13531) to its 6-hydroxy warfarin (Item No. 9000632) and 7-hydroxy warfarin (Item No. 9000672) metabolites.²

References

1. Kaminsky, L.S. and Zhang, Z.-Y. Human P450 metabolism of warfarin. *Pharmacol. Ther.* **73(1)**, 67-74 (1997).
2. Jones, D.R., Kim, S.-Y., Guderyon, M., *et al.* Hydroxywarfarin metabolites potently inhibit CYP2C9 metabolism of S-warfarin. *Chem. Res. Toxicol.* **23(5)**, 939-945 (2010).

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, 11/08/2022

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