

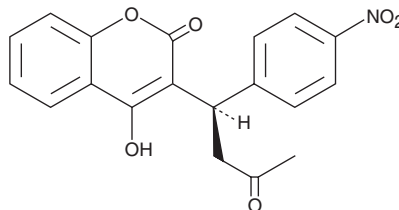
Product Information



(R)-(+)-Acenocoumarol

Catalog No. 9000336

CAS Registry No.: 66556-77-2
Formal Name: 4-hydroxy-3-[1-(4-nitrophenyl)-3-oxobutyl]-2H-1-benzopyran-2-one
Synonyms: (R)-(+)-Acenocoumarin, (R)-(+)-Nicoumalone
MF: C₁₉H₁₅NO₆
FW: 353.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that (R)-(+)-acenocoumarol be stored as supplied at -20°C. It should be stable for at least two years.

(R)-(+)-Acenocoumarol is supplied as a crystalline solid. A stock solution may be made by dissolving the (R)-(+)-acenocoumarol in an organic solvent purged with an inert gas. (R)-(+)-Acenocoumarol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (R)-(+)-acenocoumarol in these solvents is approximately 0.2, 10, and 20 mg/ml, respectively.

(R)-(+)-Acenocoumarol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (R)-(+)-acenocoumarol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (R)-(+)-Acenocoumarol has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Acenocoumarol is a short-lived oral anti-coagulant, which, like warfarin, functions by inhibiting vitamin K epoxide reductase. It has higher intrinsic anti-coagulant potency than warfarin and phenprocoumon, when evaluated *in vitro*. Acenocoumarol has a single chiral center that gives rise to two different enantiomeric forms. (R)-(+)-Acenocoumarol has a longer plasma elimination half-life (6.6 hours) and slower plasma clearance (1.9 L/hour), compared to the (S)-(-)-enantiomer (1.8 hours, 28.5 L/hour).¹ The R-enantiomer is rapidly absorbed from the gastrointestinal tract with essentially complete oral bioavailability, whereas (S)-(-)-acenocoumarol undergoes extensive first-pass metabolism.¹ Perhaps related to these pharmacokinetic characteristics, (R)-(+)-acenocoumarol is more potent *in vivo* as an anti-coagulant than the (S)-(-)-enantiomer. As the clearance of acenocoumarol is ~20-fold faster than that for warfarin, the plasma concentrations of acenocoumarol are substantially lower than those for warfarin in patients receiving long-term treatment.

Reference

1. Ufer, M. Comparative pharmacokinetics of vitamin K antagonists warfarin, phenprocoumon and acenocoumarol. *Clin. Pharmacokinet.* **44**(12), 1227-1246 (2005).

Related Products

(+)-Warfarin - Cat. No. 13526 • (-)-Warfarin - Cat. No. 13531 • (±)-Warfarin - Cat. No. 13566 • (S)-(-)-Acenocoumarol - Cat. No. 9000337 • Acenocoumarol - Cat. No. 10010569

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/01/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com