

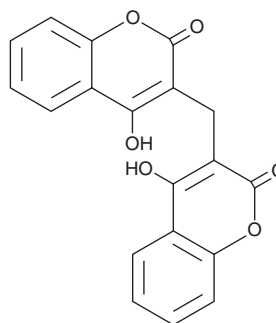
# PRODUCT INFORMATION



## Dicoumarol

Item No. 20764

**CAS Registry No.:** 66-76-2  
**Formal Name:** 3,3'-methylenebis[4-hydroxy-2H-1-benzopyran-2-one]  
**Synonyms:** NSC 17860, NSC 41834, NSC 221570  
**MF:** C<sub>19</sub>H<sub>12</sub>O<sub>6</sub>  
**FW:** 336.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 309 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

Dicoumarol is supplied as a crystalline solid. A stock solution may be made by dissolving the dicoumarol in the solvent of choice. Dicoumarol is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of dicoumarol in these solvents is approximately 2.5 and 1.25 mg/ml, respectively. Dicoumarol is also slightly soluble in ethanol.

Dicoumarol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dicoumarol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dicoumarol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Dicoumarol is a competitive inhibitor of NADH:quinone oxidoreductase (NQO1) with IC<sub>50</sub> values of 2.6 and 404 nM in the absence and presence of 2 μM BSA, respectively.<sup>1</sup> It has antiproliferative activity against MIA PaCa-2 pancreas and HCT116 colon carcinoma cells (IC<sub>50</sub>s = 52 and 19 μM, respectively, after a 96 hour incubation). Dicoumarol inhibits stress-activated protein kinase (SAPK) in HEK293 cells (IC<sub>50</sub> = 19-33 μM) at a point upstream of MEKK1 and downstream of TNF receptor-associated factor 2 (TRAF2), and it inhibits TNF-α and LPS-induced NF-κB activation in HeLa cells.<sup>2</sup> It also has antiproliferative activity against FL5.12 lymphocytic and MCF-7 breast carcinoma cells (100 μM) by suppressing JNK activation.<sup>3</sup>

### References

1. Nolan, K.A., Doncaster, J.R., Dunstan, M.S., *et al.* Synthesis and biological evaluation of coumarin-based inhibitors of NAD(P)H: Quinone oxidoreductase-1 (NQO1). *J. Med. Chem.* **52**(22), 7142-7156 (2009).
2. Cross, J.V., Deak, J.C., Rich, E.A., *et al.* Quinone reductase inhibitors block SAPK/JNK and NFκB pathways and potentiate apoptosis. *J. Biol. Chem.* **274**(44), 31150-31154 (1999).
3. Krause, D., Lyons, A., Fennelly, C., *et al.* Transient activation of Jun N-terminal kinases and protection from apoptosis by the insulin-like growth factor I receptor can be suppressed by dicoumarol. *J. Biol. Chem.* **276**(22), 19244-19252 (2001).

#### 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, 11/04/2022

#### 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