

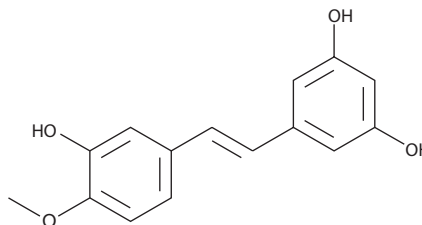
PRODUCT INFORMATION



Rhapontigenin

Item No. 13293

CAS Registry No.: 500-65-2
Formal Name: 5-[(1E)-2-(3-hydroxy-4-methoxyphenyl)ethenyl]-1,3-benzenediol
MF: C₁₅H₁₄O₄
FW: 258.3
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 326 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

Rhapontigenin is supplied as a crystalline solid. A stock solution may be made by dissolving the rhapontigenin in the solvent of choice, which should be purged with an inert gas. Rhapontigenin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of rhapontigenin in these solvents is approximately 30 mg/ml.

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

Description

Rhapontigenin is a natural analog of resveratrol with antioxidant and anti-cancer activity. It is a mechanism-based, selective inactivator of cytochrome P450 1A1 (IC₅₀ = 400 nM), an aryl hydrocarbon hydroxylase which activates polycyclic aromatic hydrocarbons that act as procarcinogens.¹ At higher concentrations, rhapontigenin inhibits the proliferation of HepG2 and HL-60R cancer cell lines (IC₅₀ = 48 μM).^{2,3}

References

1. Chun, Y.J., Ryu, S.Y., Jeong, T.C., *et al.* Mechanism-based inhibition of human cytochrome P450 1A1 by rhapontigenin. *Drug Metab. Dispos.* **29(4)**, 389-393 (2001).
2. Roupe, K.A., Helms, G.L., Halls, S.C., *et al.* Preparative enzymatic synthesis and HPLC analysis of rhapontigenin: Applications to metabolism, pharmacokinetics and anti-cancer studies. *J. Pharm. Pharm. Sci.* **8(3)**, 374-386 (2005).
3. Roberti, M., Pizzirani, D., Simoni, D., *et al.* Synthesis and biological evaluation of resveratrol and analogues as apoptosis-inducing agents. *J. Med. Chem.* **46(16)**, 3546-3554 (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, 09/30/2022

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