

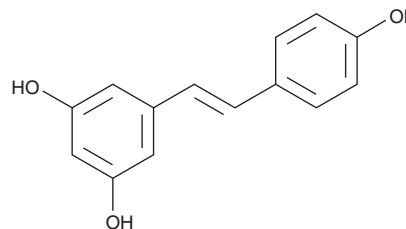
PRODUCT INFORMATION



trans-Resveratrol

Item No. 70675

CAS Registry No.: 501-36-0
Formal Name: 5-[(1E)-2-(4-hydroxyphenyl)ethenyl]-1,3-benzenediol
Synonym: (E)-Resveratrol
MF: C₁₄H₁₂O₃
FW: 228.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 307, 321 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Special Conditions: Light and pH sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

trans-Resveratrol is supplied as a crystalline solid. A stock solution may be made by dissolving the *trans*-resveratrol in the solvent of choice, which should be purged with an inert gas. *trans*-Resveratrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *trans*-resveratrol in ethanol and DMSO is approximately 50 mg/ml and approximately 100 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of *trans*-resveratrol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of *trans*-resveratrol in PBS, pH 7.2, is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

trans-Resveratrol is a polyphenol that has been found in grapes and has diverse biological activities.^{1,2} It inhibits the cyclooxygenase and hydroperoxidase activities of COX-1 (EC₅₀s = 15 and 3.7 μM, respectively) but not COX-2 (EC₅₀s = >100 μM and 85 μM, respectively).¹ *trans*-Resveratrol (200 μM) also activates sirtuin 1 (SIRT1), as well as inhibits a variety of targets including ERK1, JNK1, Src, PKCα, aromatase/CYP19, and DNA polymerases α and δ (IC₅₀s = 37, 50, 20, <10, 25, 3.3, and 5 μM, respectively) *in vitro*.^{2,3} It inhibits free radical formation induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in HL-60 cells and reduces the tumor incidence and number of tumors in a two-stage mouse model of skin cancer induced by TPA and 7,12-dimethyl-benz[*a*]anthracene (DMBA). *trans*-Resveratrol (3 and 8 mg/kg) inhibits carrageenan-induced paw edema in mice. Intravaginal administration of *trans*-resveratrol (12.5% v/v) inhibits herpes simplex virus 1 (HSV-1) and HSV-2 replication and delays the development of extrvaginal disease in mouse models of vaginal HSV infection.⁴ It also prolongs lifespan in model organisms including *C. elegans*, *D. melanogaster*, and mice.²

References

1. Jang, M., Cai, L., Udeani, G.O., *et al.* *Science* **275**(5297), 218-220 (1997).
2. Piroola, L. and Fröjdö, S. *IUBMB Life* **60**(5), 323-332 (2008).
3. Borra, M.T., Smith, B.C., and Denu, J.M. *J. Biol. Chem.* **280**(17), 17187-17195 (2005).
4. Docherty, J.J., Fu, M.M., Hah, J.M., *et al.* *Antiviral Res.* **67**(3), 155-162 (2005).

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, 11/11/2020

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