

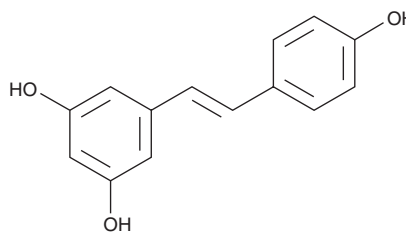
Product Information



trans-Resveratrol

Catalog 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%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 218, 235, 307, 321 nm



Laboratory Procedures

For long term storage, we suggest that *trans*-Resveratrol be stored as supplied at -20°C. It should be stable for at least one year.

trans-Resveratrol is supplied as a crystalline solid. A stock solution may be made by dissolving the Resveratrol in an organic solvent purged with an inert gas. *trans*-Resveratrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of *trans*-Resveratrol in these solvents is approximately 65 mg/ml.

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 compound in aqueous buffers. The solubility of *trans*-Resveratrol in PBS (pH 7.2) is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Phenolic compounds, particularly flavonoids, from plant sources have long been observed to have antioxidant activity with potential benefits for human health.¹⁻³ *trans*-Resveratrol is a potent phenolic antioxidant found in grapes and red wine that also has antiproliferative and anti-inflammatory activity.⁴ *trans*-Resveratrol is also a selective inhibitor of cyclooxygenase-1 (COX-1).⁵ It inhibits COX and peroxidase activities of COX-1 with ED₅₀ values of 15 and 3.7 µM, respectively; with essentially no inhibition of the COX activity of COX-2. Resveratrol also activates sirtuins⁶ and, in *C. elegans*, extends lifespan.⁷

References

1. Frémont, L., Belguendouz, L., and Delpal, S. *Life Sci.* **64**, 2511-2521 (1999).
2. Johnson, J.L. and Maddipati, K.R. *Prostaglandins and Other Lipid Mediators* **56**, 131-143 (1998).
3. Miller, N.J. and Rice-Evans, C. *Clin. Chem.* **41**, 1789 (1995).
4. Rotondo, S., Rajtar, G., Manarini, S., et al. *Br. J. Pharmacol.* **123**, 1691-1699 (1998).
5. Jang, M., Cai, L., Udeani, G.O., et al. *Science* **275**, 218-220 (1997).
6. Borra, M.T., Smith, B.C., and Denu, J.M. *J. Biol. Chem.* **280**(17), 17187-17195 (2005).
7. Howitz, K.T., Bitterman, K.J., Cohen, H.Y., et al. *Nature* **425**, 191-196 (2003).

Related Products

COX-PAK - Cat. No. 10186 • Trismethoxy Resveratrol - Cat. No. 10188 • Trismethoxy Resveratrol-d₄ - Cat. No. 13129 • *trans*-Resveratrol-d₄ - Cat. No. 13130 • Astaxanthin - Cat. No. 70685 • Carnosol - Cat. No. 89800 • Luteolin - Cat. No. 10004161 • 3,4',5'-Trismethoxybenzophenone - Cat. No. 10004185 • *cis*-Resveratrol - Cat. No. 10004235 • PDM 11 - Cat. No. 10006341 • PDM 2 - Cat. No. 10006342 • CAY10465 - Cat. No. 10006546 • Combrestatin A4 - Cat. No. 10007412 • CAY10503 - Cat. No. 10008872 • CAY10512 - Cat. No. 10009536 • CAY10591 - Cat. No. 10009797 • SIRT1 Direct Fluorescent Screening Assay Kit - Cat. No. 10010401 • SIRT1 FRET-Based Screening Assay Kit - Cat. No. 10010991 • SIRT3 Direct Fluorescent Screening Assay Kit - Cat. No. 10011566 • SIRT1 (human recombinant) - Cat. No. 10011190 • SIRT3 (human recombinant) - Cat. No. 10011194

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

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 under separate cover to the MSDS supervisor at 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, 02/18/2009