

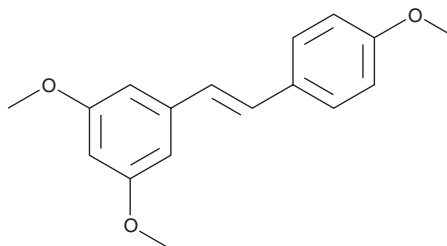
Product Information



trans-trimethoxy Resveratrol

Item No. 10188

CAS Registry No.: 22255-22-7
Formal Name: 1,3-dimethoxy-5-[(1E)-2-(4-methoxyphenyl)ethenyl]-benzene
Synonym: (E)-5-[2-(4-hydroxyphenyl)ethenyl]-1,3-benzene diol
MF: C₁₇H₁₈O₃
FW: 270.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 218, 306, 319 nm



Laboratory Procedures

For long term storage, we suggest that *trans*-trimethoxy resveratrol be stored as supplied at -20°C. It should be stable for at least two years.

trans-trimethoxy Resveratrol is supplied as a crystalline solid. A stock solution may be made by dissolving the *trans*-trimethoxy resveratrol in an organic solvent purged with an inert gas. *trans*-trimethoxy Resveratrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *trans*-trimethoxy resveratrol in these solvents is approximately 10 mg/ml in ethanol and 50 mg/ml in DMSO and DMF.

trans-trimethoxy Resveratrol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, *trans*-trimethoxy resveratrol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. *trans*-trimethoxy Resveratrol has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. 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 Resveratrol is a potent phenolic antioxidant found in grapes and red wine that also has antiproliferative and anti-inflammatory activity.⁴ When the three phenolic hydroxyl groups of resveratrol are converted to methyl ethers, the inhibition of cell growth and pro-apoptotic activities of resveratrol are enhanced.⁵

References

- Frémont, L., Belguendouz, L., and Delpal, S. Antioxidant activity of resveratrol and alcohol-free wine polyphenols related to LDL oxidation and polyunsaturated fatty acids. *Life Sci.* **64**, 2511-2521 (1999).
- Johnson, J.L. and Maddipati, K.R. Paradoxical effects of resveratrol on the two prostaglandin H synthases. *Prostaglandins and Other Lipid Mediators* **56**, 131-143 (1998).
- Miller, N.J. and Rice-Evans, C. Antioxidant activity of resveratrol in red wine. *Clin. Chem.* **41**, 1789 (1995).
- Rotondo, S., Rajtar, G., Manarini, S., *et al.* Effect of *trans*-resveratrol, a natural polyphenolic compound, on human polymorphonuclear leukocyte function. *Brit. J. Pharmacol.* **123**, 1691-1699 (1998).
- Pettit, G.R., Grealish, M.P., Jung, M.K., *et al.* Antineoplastic agents. 465. Structural modification of resveratrol: Sodium resverastatin phosphate. *J. Med. Chem.* **45**, 2534-2542 (2002).

Related Products

TMS - Item No. 10038 • *trans*-trimethoxy Resveratrol-d₄ - Item No. 13129 • *trans*-Resveratrol-d₄ - Item No. 13130 • Resveratrol - Item No. 70675 • 3,4',5'-Trimethoxybenzophenone - Item No. 10004185 • *cis*-Resveratrol - Item No. 10004235 • CAY10512 - Item No. 10009536 • CAY10591 - Item No. 10009797 • SIRT1 Direct Fluorescent Screening Assay Kit - Item No. 10010401 • SIRT1 (human recombinant) - Item No. 10011190 • SIRT3 (human recombinant) - Item No. 10011194 • SIRT3 Direct Fluorescent Screening Assay Kit - Item No. 10011566

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, 03/08/2011

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