

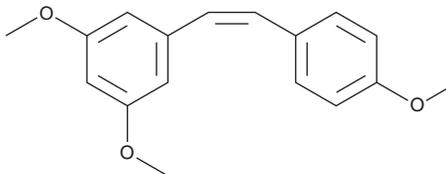
PRODUCT INFORMATION



cis-trimethoxy Resveratrol

Item No. 13199

CAS Registry No.: 94608-23-8
Formal Name: 1,3-dimethoxy-5-[(1Z)-2-(4-methoxyphenyl)ethenyl]-benzene
Synonym: *cis*-Trimethoxy Stilbene
MF: C₁₇H₁₈O₃
FW: 270.3
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 285 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

cis-trimethoxy Resveratrol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of *cis*-trimethoxy resveratrol in these solvents is approximately 50 mg/ml.

cis-trimethoxy Resveratrol is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Resveratrol is a potent antioxidant found in grapes and red wine that also has anti-proliferative, anti-neoplastic, and anti-angiogenic activities. In addition, resveratrol activates sirtuins and, in yeast, extends lifespan.^{1,2} *cis*-trimethoxy Resveratrol is a potent anti-mitotic drug that is 100-fold more active than resveratrol at inhibiting the growth of human colon cancer Caco-2 cells.³ It inhibits tubulin polymerization in a dose-dependent manner (IC₅₀ = 4 μM) and inhibits enzymes involved in the synthesis of the polyamines, putrescine, and spermidine.^{3,4} *trans*-trimethoxy Resveratrol has superior pharmacokinetic characteristics when compared with resveratrol, including greater plasma exposure, longer elimination half-life, and lower clearance.⁵

References

1. Borra, M.T., Smith, B.C., and Denu, J.M. *J. Biol. Chem.* **280**(17), 17187-17195 (2005).
2. Howitz, K.T., Bitterman, K.J., Cohen, H.Y., et al. *Nature* **425**, 191-196 (2003).
3. Schneider, Y., Chabert, P., Stutzmann, J., et al. *Int. J. Cancer* **107**, 189-196 (2003).
4. Cushman, M., Nagarathnam, D., Gopal, D., et al. *J. Med. Chem.* **34**, 2579-2588 (1991).
5. Lin, H.-S. and Ho, P.C. *J. Pharm. Biomed. Analysis* **49**, 387-392 (2009).

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, 02/26/2024

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