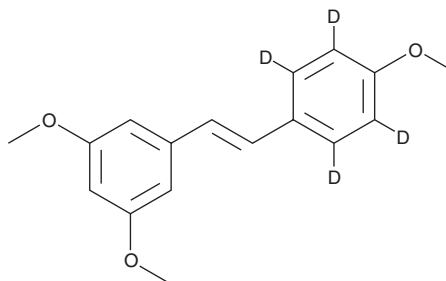


# PRODUCT INFORMATION



## *trans*-trimethoxy Resveratrol- $d_4$ Item No. 13129

**CAS Registry No.:** 1089051-64-8  
**Formal Name:** 3-[(1E)-2-(3,5-dimethoxyphenyl)ethenyl]-6-methoxy-benzene-1,2,4,5- $d_4$   
**Synonyms:** (E)-5-[2-(4-hydroxyphenyl)ethenyl]-1,3-benzene diol- $d_4$ , TMS- $d_4$ , *trans*-3,5,4'-Trimethoxystilbene- $d_4$   
**MF:**  $C_{17}H_{14}D_4O_3$   
**FW:** 274.4  
**Chemical Purity:**  $\geq 98\%$  (*trans*-trimethoxy Resveratrol)  
**Deuterium Incorporation:**  $\geq 99\%$  deuterated forms ( $d_1$ - $d_4$ );  $\leq 1\%$   $d_0$   
**UV/Vis.:**  $\lambda_{max}$ : 218, 305, 320 nm  
**Supplied as:** A solution in methyl acetate  
**Storage:**  $-20^\circ\text{C}$   
**Stability:**  $\geq 2$  years  
**Item Origin:** Plant/*Viola cuspidata*



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

### Laboratory Procedures

*trans*-trimethoxy Resveratrol- $d_4$  is intended for use as an internal standard for the quantification of *trans*-trimethoxy resveratrol (Item No. 10188) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

*trans*-trimethoxy Resveratrol- $d_4$  is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of *trans*-trimethoxy resveratrol- $d_4$  in ethanol is approximately 10 mg/ml and approximately 50 mg/ml in DMSO and DMF.

### Description

*trans*-trimethoxy Resveratrol is a polyketide synthase-derived stilbene originally isolated from *Viola cuspidata* that has diverse biological activities.<sup>1-4</sup> It is cytotoxic to several cancer cell lines, including PC3, KB, HT-29, SW480, and HL-60 cells ( $IC_{50}$ s = 3.6, 10.2, 16.1, 54, and 2.5  $\mu\text{M}$ , respectively).<sup>2</sup> *trans*-trimethoxy Resveratrol (15  $\mu\text{M}$ ) inhibits TNF- $\alpha$ -induced activation of NF- $\kappa\text{B}$  in HEK293T cells in a reporter assay.<sup>3</sup> It inhibits angiogenesis in zebrafish embryos when used at a concentration of 0.1  $\mu\text{M}$ .<sup>4</sup>

### References

1. Austin, M.B. and Noel, J.P. The chalcone synthase superfamily of type III polyketide synthases. *Nat. Prod. Rep.* **20**(1), 79-110 (2003).
2. Aldawsari, F.S. and Velázquez-Martínez, C.A. 3,4',5-*trans*-Trimethoxystilbene; a natural analogue of resveratrol with enhanced anticancer potency. *Invest. New Drugs* **33**(3), 775-786 (2015).
3. Heynekamp, J.J., Weber, W.M., Hunsaker, L.A., et al. Substituted *trans*-stilbenes, including analogues of the natural product resveratrol, inhibit the human tumor necrosis factor alpha-induced activation of transcription factor nuclear factor kappaB. *J. Med. Chem.* **49**(24), 7182-7189 (2006).
4. Belleri, M., Ribatti, D., Nicoli, S., et al. Antiangiogenic and vascular-targeting activity of the microtubule-destabilizing *trans*-resveratrol derivative 3,5,4'-trimethoxystilbene. *Mol. Pharmacol.* **67**(5), 1451-1459 (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, 03/08/2024

#### 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