

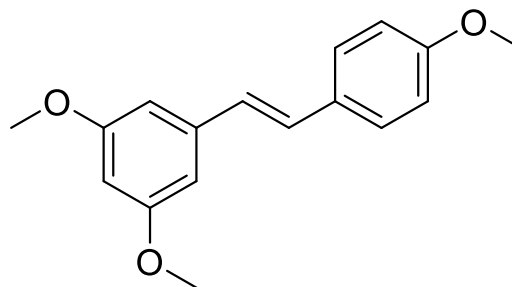
## Trismethoxyresveratrol

Catalog number: B-0022

Molecular Formula:  $C_{17}H_{18}O_3$

MW: 270.3

CAS: 22255-22-7



Alternate Name: 3,5,4'-Trimethoxystilbene

Solubility: Ethanol (~10 mg/mL), DMSO, DMF (~50 mg/mL), DMSO:PBS (1:2) (pH 7.2) - 0.25 mg/mL (DMSO solution diluted with buffer), sparingly soluble in water.

Storage and Handling: Store dry at -20°C. Stock solutions can be prepared in organic solvent and flushed with inert gas. Aqueous solutions should not be store longer than 1 day

Background: 3,5,4'-Trimethoxystilbene induces microtubule disassembly and tubulin depolymerization. it is also a powerful antiangiogenic agent and inhibits endothelial cell proliferation, morphogenesis, sprouting and collagen gel invasion.

References: M. Belleri *et al.* Antiangiogenic and vascular-targeting activity of the microtubule-destabilizing trans-resveratrol derivative 3,5,4'-trimethoxystilbene. *Mol. Pharmacol.* 2005 67 1451

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

---

Technical Data Sheet, Rev 2, 10-9-18 – For research use only. Not intended for diagnostic or therapeutic use.

