

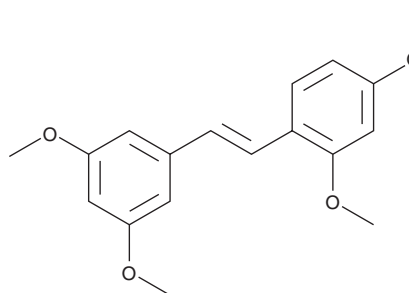
PRODUCT INFORMATION



TMS

Item No. 10038

CAS Registry No.: 24144-92-1
Formal Name: 1-[(1E)-2-(3,5-dimethoxyphenyl)ethenyl]-2,4-dimethoxy-benzene
Synonym: 2,3',4,5'-Tetramethoxystilbene
MF: C₁₈H₂₀O₄
FW: 300.4
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 302, 326 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

TMS is supplied as a crystalline solid. A stock solution may be made by dissolving the TMS in the solvent of choice, which should be purged with an inert gas. TMS is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TMS in these solvents is approximately 0.4, 20, and 30 mg/ml, respectively.

TMS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TMS should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TMS has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CYP1B1 is mainly an extrahepatic enzyme which oxidatively metabolizes both endogenous (steroids; eicosanoids) and exogenous xenobiotics such as polycyclic aromatic hydrocarbons. TMS is a potent and selective inhibitor of CYP1B1, with an IC₅₀ of 6 nM.¹ It is 50-fold selective for the inhibition of CYP1B1 versus CYP1A1, making it a useful tool to differentiate between various CYP450 families.¹ In cultured human colon cancer cells, TMS induces apoptosis and inhibits cell growth with an IC₅₀ of 0.8 µg/ml.²

References

1. Kim, S., Ko, H., Park, J.E., *et al.* Design, synthesis, and discovery of novel trans-stilbene analogues as potent and selective human cytochrome P450 1B1 inhibitors. *J. Med. Chem.* **45(1)**, 160-164 (2002).
2. Nam, K.A., Kim, S., Heo, Y.H., *et al.* Resveratrol analog, 3,5,2',4'-tetramethoxy-trans-stilbene, potentiates the inhibition of cell growth and induces apoptosis in human cancer cells. *Arch. Pharm. Res.* **24(5)**, 441-445 (2001).

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, 04/26/2024

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