

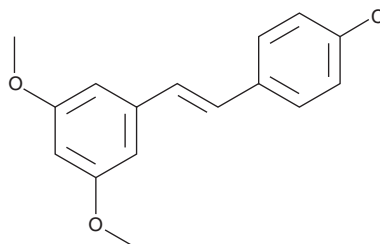
PRODUCT INFORMATION



PDM 11

Item No. 10006341

CAS Registry No.: 1032508-03-4
Formal Name: (E)-5-[2-(4-chlorophenyl)ethenyl]-1,3-dimethoxyphenyl
MF: C₁₆H₁₅ClO₂
FW: 274.7
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 303 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

PDM 11 is supplied as a crystalline solid. A stock solution may be made by dissolving the PDM 11 in the solvent of choice. PDM 11 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of PDM 11 in these solvents is approximately 2, 20, and 30 mg/ml, respectively.

PDM 11 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PDM 11 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PDM 11 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PDM 11 is a derivative of the antioxidant *trans*-resveratrol that is inactive *in vitro* in assays of resveratrol activity.¹ It does not activate quinone reductase 1, inhibit quinone reductase 2, or affect nitric oxide production or quenching of free radicals. It does not interact with estrogen receptors or affect the activity of COX-1 and COX-2. PDM 11 does not affect proliferation of K562, HT-29, and HepG2 cells.²

References

1. Kondratyuk, T.P., Park, E.J., Marler, L.E., *et al.* Resveratrol derivatives as promising chemopreventive agents with improved potency and selectivity. *Mol. Nutr. Food Res.* **55**(8), 1249-1265 (2011).
2. Chen, Y., Hu, F., Gao, Y., *et al.* Design, synthesis, and evaluation of methoxylated resveratrol derivatives as potential antitumor agents. *Res. Chem. Intermed.* **41**(5), 2725-2738 (2015).

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, 11/09/2023

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