

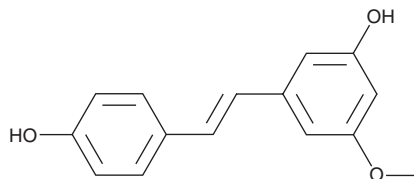
PRODUCT INFORMATION



Pinostilbene

Item No. 14044

CAS Registry No.: 42438-89-1
Formal Name: 3-[(1E)-2-(4-hydroxyphenyl)ethenyl]-5-methoxy-phenol
MF: $C_{15}H_{14}O_3$
FW: 242.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 218, 307 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

Pinostilbene is supplied as a crystalline solid. A stock solution may be made by dissolving the pinostilbene in the solvent of choice, which should be purged with an inert gas. Pinostilbene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pinostilbene in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Pinostilbene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pinostilbene should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Pinostilbene has a solubility of approximately 0.15 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pinostilbene is a stable, monomethoxylated resveratrol derivative with enhanced bioavailability compared to resveratrol.¹ At 0.1-10 μM pinostilbene reduces neurotoxicity in SH-SY5Y cells induced by the parkinsonian mimetic 6-hydroxydopamine and scavenges 2,2-diphenyl-1-picrylhydrazyl radicals with an IC_{50} value of 47.1 μM .²

References

1. Wilson, M.A., Rimando, A.M., and Wolkow, C.A. Methoxylation enhances stilbene bioactivity in *Caenorhabditis elegans*. *BMC Pharmacol.* **8**(15), (2008).
2. Chao, J., Li, H., Cheng, K.W., et al. Protective effects of pinostilbene, a resveratrol methylated derivative, against 6-hydroxydopamine-induced neurotoxicity in SH-SY5Y cells. *J. Nutr. Biochem.* **21**(6), 482-489 (2010).

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, 11/07/2022

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