

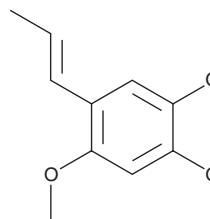
PRODUCT INFORMATION



α -Asarone

Item No. 11681

CAS Registry No.: 2883-98-9
Formal Name: 1,2,4-trimethoxy-5-(1E)-1-propen-1-yl-benzene
Synonym: *trans*-Asarone
MF: C₁₂H₁₆O₃
FW: 208.3
Purity: \geq 98%
UV/Vis.: λ_{\max} : 212, 258, 313 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Acorus calamus*



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

Laboratory Procedures

α -Asarone is supplied as a crystalline solid. A stock solution may be made by dissolving the α -asarone in the solvent of choice, which should be purged with an inert gas. α -Asarone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of α -asarone in these solvents is approximately 14, 11, and 16 mg/ml, respectively.

α -Asarone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, α -asarone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. α -Asarone 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

α -Asarone is a component of certain essential oils used in traditional Chinese herbal medicine. Although carcinogenic, α -asarone has been shown to be neuroprotective and radioprotective in mice when given at 50 mg/kg.¹⁻³ It also lowers serum cholesterol in rats, specifically reducing LDL-cholesterol levels and cholesterol saturation index, by directly inhibiting HMG-CoA reductase (IC₅₀ = 3 mM).^{4,5} α -Asarone also directly interacts with and inhibits certain isoforms of cytochrome P450 (IC₅₀ = 55.2 and 65.2 μ g/ml for CYP2D6 and CYP3A4, respectively).⁶

References

1. Limón, I.D., Mendieta, L., Díaz, A., *et al. Neurosci. Lett.* **453(2)**, 98-103 (2009).
2. Sandeep, D. and Nair, C.K. *Mutat. Res.* **722(1)**, 62-68 (2011).
3. Zuo, H.L., Yang, F.Q., Zhang, X.M., *et al. J. Anal. Methods Chem.* **2012**, 1-5 (2012).
4. Rodríguez-Páez, L., Juárez-Sánchez, M., Antúnez-Solís, J., *et al. Phytomedicine* **10(5)**, 397-404 (2003).
5. Antunez-Solis, J., Hernández-Derramadero, F., Aquino-Vega, M., *et al. J. Enzyme Inhib. Med. Chem.* **24(3)**, 903-909 (2009).
6. Pandit, S., Mukherjee, P.K., Ponnusankar, S., *et al. Fitoterapia* **82(3)**, 369-374 (2011).

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, 06/18/2024

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