

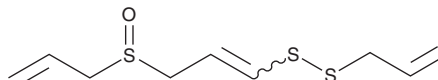
PRODUCT INFORMATION



(E/Z)-Ajoene

Item No. 34881

CAS Registry No.: 92285-01-3
Formal Name: 2-propen-1-yl [3-(2-propen-1-ylsulfanyl)-1-propen-1-yl] disulfide
MF: C₉H₁₄OS₃
FW: 234.4
Purity: ≥95%
Supplied as: A solution in ethyl acetate
Storage: -20°C
Stability: ≥2 years
Item Origin: Synthetic



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

Description

(E/Z)-Ajoene is a disulfide that has been found in *A. sativum* and has antitumor and antiplatelet activities.^{1,2} It is cytotoxic to a variety of cancer cells, including B16-BL6 murine melanoma, and human HT-29 colon adenocarcinoma, MDA-MB-231 mammary, and PANC-1 pancreatic adenocarcinoma cells (IC₅₀s = 18, 19, 7, and 38 μM, respectively).¹ (E/Z)-Ajoene (25 mg/kg) reduces tumor growth and decreases the number of lung metastases in B16-BL6 murine models of melanoma. It also prevents thrombus formation in heparinized animals in a porcine model of severe vascular damage induced by intimal stripping and tunica media exposure when administered at a dose of 25 mg/kg.²

References

1. Taylor, P., Noriega, R., Farah, C., *et al.* Ajoene inhibits both primary tumor growth and metastasis of B16/BL6 melanoma cells in C57BL/6 mice. *Cancer Lett.* **239**(2), 298-304 (2006).
2. Apitz-Castro, R., Badimon, J.J., and Badimon, L. A garlic derivative, ajoene, inhibits platelet deposition on severely damaged vessel wall in an *in vivo* porcine experimental model. *Thromb. Res.* **75**(3), 243-249 (1994).

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, 02/15/2024

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