

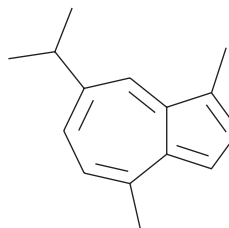
PRODUCT INFORMATION



Guaiazulene

Item No. 31506

CAS Registry No.: 489-84-9
Formal Name: 1,4-dimethyl-7-(1-methylethyl)-azulene
Synonym: NSC 4714
MF: C₁₅H₁₈
FW: 198.3
Purity: ≥95%
UV/Vis.: λ_{max}: 244, 284 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Guaiazulene is supplied as a solid. A stock solution may be made by dissolving the guaiazulene in the solvent of choice, which should be purged with an inert gas. Guaiazulene is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of guaiazulene in ethanol and DMSO is approximately 25 mg/ml and approximately 30 mg/ml in DMF.

Guaiazulene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, guaiazulene should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Guaiazulene 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

Guaiazulene is a sesquiterpene that has been found in *M. chamomilla* and has diverse biological activities.^{1,2,3} It inhibits lipid peroxidation in rat hepatic microsomes (IC₅₀ = 9.8 μM), as well as scavenges hydroxyl and 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in cell-free assays.¹ It inhibits LPS-induced nitric oxide production in RAW 264.7 cells (EC₅₀ = 10.1 μM) but is cytotoxic to RAW 264.7 cells at higher concentrations with cytotoxic concentration (CC₅₀) values of 29.8 and 30.8 μM in the presence and absence, respectively, of LPS.² Guaiazulene is cytotoxic to N2a neuroblastoma cells and primary rat neurons in a concentration-dependent manner.³ It inhibits decreases in hepatic glutathione (GSH) levels induced by paracetamol (acetaminophen; Item No. 10024) in rats when administered at a dose of 250 mg/kg.¹

References

1. Kourounakis, A.P., Rekka, E.A., and Kourounakis, P.N. Antioxidant activity of guaiazulene and protection against paracetamol hepatotoxicity in rats. *J. Pharm. Pharmacol.* **49**(9), 938-942 (1997).
2. Hashiba, K., Yokoyama, K., Wakabayashi, H., et al. Inhibition of LPS-stimulated NO production in mouse macrophage-like cells by azulenes. *Anticancer Res.* **24**(6), 3939-3944 (2004).
3. Togar, B., Turkez, H., Hacimuftuoglu, A., et al. Guaiazulene biochemical activity and cytotoxic and genotoxic effects on rat neuron and N2a neuroblastoma cells. *J. Intercult. Ethnopharmacol.* **4**(1), 29-33 (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, 11/16/2022

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