

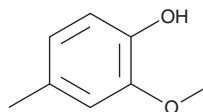
PRODUCT INFORMATION



2-Methoxy-4-methylphenol

Item No. 36533

CAS Registry No.: 93-51-6
Formal Name: 2-methoxy-4-methyl-phenol
Synonyms: *p*-Creosol, 4-methyl Guaiacol, *p*-Methylguaiacol, NSC 4969
MF: C₈H₁₀O₂
FW: 138.2
Purity: ≥98%
Supplied as: A neat liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

2-Methoxy-4-methylphenol is supplied as a neat liquid. A stock solution may be made by dissolving the 2-methoxy-4-methylphenol in water. 2-Methoxy-4-methylphenol is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

2-Methoxy-4-methylphenol is a phenolic pyrolysis product of plant biomass.^{1,2} It is a volatile component of black olives and roasted coffee beans.^{3,4}

References

1. Galletti, G.C., Ward, R.S., and Pelter, A. Analysis of pyrolysis products of lignans using capillary gas chromatography and ion trap detection. *J. Anal. Appl. Pyrolysis* **21(3)**, 281-292 (1991).
2. Theapparat, C.Y., Khongthong, S., Rodjan, P., *et al.* Physicochemical properties and in vitro antioxidant activities of pyroligneous acid prepared from brushwood biomass waste of Mangosteen, Durian, Rambutan, and Langsat. *J. For. Res.* **30(3)**, 1139-1148 (2018).
3. Sánchez, R., Martín-Tornero, E., Lozano, J., *et al.* Electronic nose application for the discrimination of sterilization treatments applied to Californian-style black olive varieties. *J. Sci. Food Agric.* **102(6)**, 2232-2241 (2022).
4. Chindapan, N., Puangngoen, C., and Devahastin, S. Profiles of volatile compounds and sensory characteristics of Robusta coffee beans roasted by hot air and superheated steam. *Int. J. Food Sci.* **56(8)**, 3814-3825 (2021).

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, 03/28/2022

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