

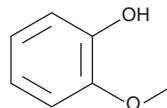
PRODUCT INFORMATION



Guaiacol

Item No. 70430

CAS Registry No.: 90-05-1
Formal Name: 2-methoxy phenol
MF: C₇H₈O₂
FW: 124.1
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 277 nm



Supplied as: A colorless liquid
Storage: Room temperature
Stability: ≥2 years

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

Laboratory Procedures

Guaiacol is supplied as a colorless liquid. A stock solution may be made by diluting the guaiacol in an organic solvent purged with an inert gas. Guaiacol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of guaiacol in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of guaiacol can be prepared by directly diluting the colorless liquid in aqueous buffers. The solubility of guaiacol in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Guaiacol is a phenolic natural product first isolated from Guaiac resin and the oxidation of lignin. Guaiacol is readily oxidized by the heme iron of peroxidases including the peroxidase of cyclooxygenase (COX) enzymes. It can therefore serve as a reducing co-substrate for COX reactions.¹ The color yield per peroxide at 470 nm is 6,650 (M peroxide reduced)⁻¹ cm⁻¹.²⁻⁴ Two moles of guaiacol are oxidized for each mole of hydroperoxide reduced by the peroxidase. The resulting guaiacol chromophore can be used for the colorimetric determination of hydroperoxidase activity.

References

1. Markey, C.M., Alward, A., Weller, P.E., *et al.* Quantitative studies of hydroperoxide reduction by prostaglandin H synthase. Reducing substrate specificity and the relationship of peroxidase to cyclooxygenase activities. *J. Biol. Chem.* **262**, 6266-6279 (1987).
2. Doerge, D.R., Divi, R.L., and Churchwell, M.I. Identification of the colored guaiacol oxidation product produced by peroxidases. *Anal. Biochem.* **250**, 10-17 (1997).
3. Kulmacz, R.J. and Lands, W.E.M. Quantitative similarities in the several actions of cyanide on prostaglandin H synthase. *Prostaglandins* **29**, 175-190 (1985).
4. Marquez, L.A. and Dunford, H.B. Mechanism of the oxidation of 3,5,3',5'-tetramethylbenzidine by myeloperoxidase determined by transient-and steady-state kinetics. *Biochemistry* **36**, 9349-9355 (1997).

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, 12/20/2022

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