

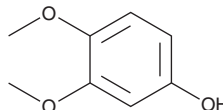
PRODUCT INFORMATION



3,4-Dimethoxyphenol

Item No. 45665

CAS Registry No.: 2033-89-8
Synonym: NSC 140927
MF: $C_8H_{10}O_3$
FW: 154.2
Purity: $\geq 95\%$
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

3,4-Dimethoxyphenol is supplied as a solid. A stock solution may be made by dissolving the 3,4-dimethoxyphenol in the solvent of choice, which should be purged with an inert gas. 3,4-Dimethoxyphenol is soluble (≥ 10 mg/ml) in ethanol and DMSO.

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 3,4-dimethoxyphenol can be prepared by directly dissolving the solid in aqueous buffers. 3,4-Dimethoxyphenol is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

3,4-Dimethoxyphenol is a phenol that has been found in *W. nobilis* and has diverse biological activities.¹⁻³ It scavenges DPPH (Item No. 14805) radicals in a cell-free assay and is active against *S. aureus*, *M. luteus*, and *C. albicans*.^{1,2}

References

1. Seal, A.N., Pratley, J.E., Haig, T.J., et al. Plants with phytotoxic potential: Wollemi pine (*Wollemia nobilis*). *Agric. Ecosyst. Environ.* **135**(1-2), 52-57 (2010).
2. Furtado, N.A., Said, S., Ito, I.Y., et al. The antimicrobial activity of *Aspergillus fumigatus* is enhanced by a pool of bacteria. *Microbiol. Res.* **157**(3), 207-211 (2002).
3. Liu, Q., Zhang, Y.-J., Yang, C.-R., et al. Phenolic antioxidants from green tea produced from *Camellia crassicolumna* Var. *multiplex*. *J. Agric. Food Chem.* **57**(2), 586-590 (2009).

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, 06/08/2026

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