

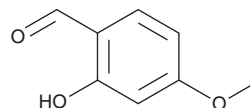
PRODUCT INFORMATION



2-Hydroxy-4-methoxybenzaldehyde

Item No. 30091

CAS Registry No.: 673-22-3
Synonyms: 4-Methoxysalicylaldehyde, NSC 155334
MF: C₈H₈O₃
FW: 152.1
Purity: ≥95%
Supplied as: A crystalline solid
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.

Laboratory Procedures

2-Hydroxy-4-methoxybenzaldehyde is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-hydroxy-4-methoxybenzaldehyde in the solvent of choice, which should be purged with an inert gas. 2-Hydroxy-4-methoxybenzaldehyde is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2-hydroxy-4-methoxybenzaldehyde in these solvents is approximately 30 mg/ml.

Description

2-Hydroxy-4-methoxybenzaldehyde is a phenol found in the root bark essential oil of *P. sepium* that has diverse biological activities, including antibacterial, antifungal, fungicidal, antioxidant, anti-virulence, and metal chelating properties.¹⁻³ It is active against Gram-positive (MICs = 100-200 µg/ml) and Gram-negative bacteria (MICs = 125-200 µg/ml), as well as the fungus *C. albicans* and the phytopathogenic fungus *M. oryzae* (MICs = 150 and 300 µg/ml, respectively). 2-Hydroxy-4-methoxybenzaldehyde scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH) radicals in a cell-free assay (IC₅₀ = 9.04 mg/ml) and acts as a metal chelator, inhibiting ferrozine-Fe²⁺ complex formation with an IC₅₀ value of 2.31 mg/ml. It reduces the production of the virulence factor staphyloxanthin in clinical isolates of methicillin-resistant *S. aureus* (MRSA) when used at a concentration of 200 µg/ml and decreases the expression of the MRSA virulence regulatory genes *saeS*, *geh*, *crtM*, and *sigB*.² 2-Hydroxy-4-methoxybenzaldehyde, when used in combination with the antibiotics amikacin (Item No. 15405), gentamicin, cefotaxime (Item No. 16040), vancomycin (Item No. 15327), or tetracycline (Item No. 14328), increases MRSA antibiotic sensitivity. It is also an intermediate in the synthesis of Schiff base-metal complexes with antiproliferative activity.³

References

1. Wang, J., Liu, H., Zhao, J., *et al.* Antimicrobial and antioxidant activities of the root bark essential oil of *Periploca sepium* and its main component 2-hydroxy-4-methoxybenzaldehyde. *Molecules* **15**(8), 5807-5817 (2010).
2. Kannappan, A., Srinivasan, R., Nivetha, A., *et al.* Anti-virulence potential of 2-hydroxy-4-methoxybenzaldehyde against methicillin-resistant *Staphylococcus aureus* and its clinical isolates. *Appl. Microbiol. Biotechnol.* **103**(16), 6747-6758 (2019).
3. Tyagi, P., Tyagi, M., Agrawal, S., *et al.* Synthesis, characterization of 1,2,4-triazole Schiff base derived 3d-metal complexes: Induces cytotoxicity in HepG2, MCF-7 cell line, BSA binding fluorescence and DFT study. *Spectrochim. Acta. A. Mol. Biomol. Spectrosc.* **171**, 246-257 (2017).

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, 05/13/2020

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