

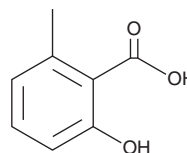
# PRODUCT INFORMATION



## 2-hydroxy-6-Methylbenzoic Acid

Item No. 19199

**CAS Registry No.:** 567-61-3  
**Formal Name:** 2-hydroxy-6-methyl-benzoic acid  
**Synonyms:** 2,6-Cresotic Acid, 6-Methylsalicylic Acid, 6-MSA, NSC 403256, 6-hydroxy-*o*-Toluic Acid  
**MF:** C<sub>8</sub>H<sub>8</sub>O<sub>3</sub>  
**FW:** 152.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 209, 242, 308 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

2-hydroxy-6-Methylbenzoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-hydroxy-6-methylbenzoic acid in the solvent of choice. 2-hydroxy-6-Methylbenzoic acid is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 2-hydroxy-6-methylbenzoic acid in these solvents is approximately 5 mg/ml.

2-hydroxy-6-Methylbenzoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-hydroxy-6-methylbenzoic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 2-hydroxy-6-Methylbenzoic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2-hydroxy-6-Methylbenzoic acid is a constituent of *G. anandria*, a well-known traditional Chinese medicinal herb, and also produced in various fungi, including *P. patulum*, where it is a precursor to the toxin patulin (Item No. 11346).<sup>1,2</sup> This polyketide is a key structural moiety for many different antibiotic and anticancer agents, including chlorothricin (Item No. 15501), maduropeptin, neocarzinostatin, and terreic acid.<sup>2,3</sup>

### References

1. He, F., Wang, M., Gao, M., *et al.* Chemical composition and biological activities of *Gerbera anandria*. *Molecules* **19**(4), 4046-4057 (2014).
2. Parascandolo, J.S., Havermann, J., Potter, H.K., *et al.* Insights into 6-methylsalicylic acid bio-assembly by using chemical probes. *Angew. Chem. Int. Ed. Engl.* **55**(10), 3463-3467 (2016).
3. Guo, C.-J., Sun, W.-W., Bruno, K.S., *et al.* Molecular genetic characterization of terreic acid pathway in *Aspergillus terreus*. *Org. Lett.* **16**(20), 5250-5253 (2014).

#### 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, 11/07/2022

#### 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](http://WWW.CAYMANCHEM.COM)