

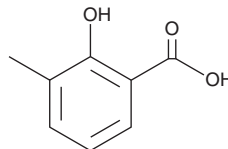
# PRODUCT INFORMATION



## 3-Methylsalicylic Acid

Item No. 36536

**CAS Registry No.:** 83-40-9  
**Formal Name:** 2-hydroxy-3-methyl-benzoic acid  
**Synonyms:** 3-methyl-2-Hydroxybenzoic Acid, NSC 1772, NSC 17561, NSC 50796, 3-methyl SA  
**MF:** C<sub>8</sub>H<sub>8</sub>O<sub>3</sub>  
**FW:** 152.1  
**Purity:** ≥98%  
**Supplied as:** A 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

3-Methylsalicylic acid is supplied as a solid. A stock solution may be made by dissolving the 3-methylsalicylic acid in the solvent of choice, which should be purged with an inert gas. 3-Methylsalicylic acid is soluble in the organic solvent DMSO.

### Description

3-Methylsalicylic acid is a derivative of salicylic acid. It has been used in the synthesis of salicylamides with *in vitro* anticancer activity.<sup>1</sup> 3-Methylsalicylic acid has also been used in the synthesis of pyrimethamine-salicylic acid salts.<sup>2</sup>

### References

1. Chae, H.-D., Cox, N., Capolicchio, S., *et al.* SAR optimization studies on modified salicylamides as a potential treatment for acute myeloid leukemia through inhibition of the CREB pathway. *Bioorg. Med. Chem. Lett.* **29(16)**, 2307-2315 (2019).
2. Ceborska, M., Kędra-Królik, K., Narodowicz, J., *et al.* Influence of hydroxyl group position and substitution pattern of hydroxybenzoic acid on the formation of molecular salts with the antifolate pyrimethamine. *Cryst. Growth Des.* **21(12)**, 6714-6726 (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, 10/04/2022

#### 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)