

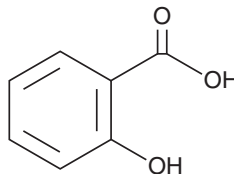
# PRODUCT INFORMATION



## Salicylic Acid

Item No. 32991

**CAS Registry No.:** 69-72-7  
**Formal Name:** 2-hydroxy-benzoic acid  
**Synonyms:** 2-Hydroxybenzoic Acid, NSC 180  
**MF:** C<sub>7</sub>H<sub>6</sub>O<sub>3</sub>  
**FW:** 138.1  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Salicylic acid (Item No. 32991) is an analytical reference standard categorized as a COX inhibitor metabolite.<sup>1</sup> Salicylic acid is a metabolite of aspirin. It has been found as an adulterant in seized heroin samples.<sup>2</sup> It has also been found as an adulterant in milk used to increase shelf life.<sup>3</sup> This product is intended for research and forensic applications.

### References

1. Klessig, D.F. Newly identified targets of aspirin and its primary metabolite, salicylic acid. *DNA Cell Biol.* **35(4)**, 163-166 (2016).
2. Lurie, I.S., Driscoll, S.E., Cathapermal, S.S., *et al.* Determination of heroin and basic impurities for drug profiling by ultra-high-pressure liquid chromatography. *Forensic Sci. Int.* **231(1-3)**, 300-305 (2013).
3. Nascimento, C.F., Santos, P.M., Pereira-Filho, E.R., *et al.* Recent advances on determination of milk adulterants. *Food Chem.* **221**, 1232-1244 (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, 07/05/2023

#### 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