

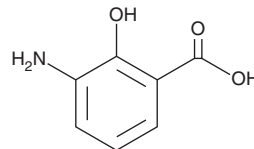
# PRODUCT INFORMATION



## 3-Aminosalicylic Acid

Item No. 36543

**CAS Registry No.:** 570-23-0  
**Formal Name:** 3-amino-2-hydroxy-benzoic acid  
**Synonyms:** 3-ASA, NSC 285111  
**MF:** C<sub>7</sub>H<sub>7</sub>NO<sub>3</sub>  
**FW:** 153.1  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224 nm  
**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-Aminosalicylic acid is supplied as a solid. A stock solution may be made by dissolving the 3-aminosalicylic acid in the solvent of choice, which should be purged with an inert gas. 3-Aminosalicylic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 3-aminosalicylic acid in these solvents is approximately 1 and 2 mg/ml, respectively.

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-aminosalicylic acid can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 3-aminosalicylic acid in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

3-Aminosalicylic acid is a salicylic acid derivative and potential impurity in commercial preparations of 5-aminosalicylic acid.<sup>1,2</sup> It protects against apoptosis, as well as reverses reductions in mitochondrial complex I, also known as NADH dehydrogenase, activity and increases in reactive oxygen species (ROS) production induced by manganese in SK-N-MC cells when used at a concentration of 1 μM.<sup>1</sup>

### References

1. Yoon, H., Lee, G.-H., Kim, D.-S., *et al.* The effects of 3, 4 or 5 amino salicylic acids on manganese-induced neuronal death: ER stress and mitochondrial complexes. *Toxicol. In Vitro* **25(7)**, 1259-1268 (2011).
2. Gotti, R., Pomponio, R., Bertucci, C., *et al.* Determination of 5-aminosalicylic acid related impurities by micellar electrokinetic chromatography with an ion-pair reagent. *J. Chromatogr. A* **916(1-2)**, 175-183 (2001).

#### 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, 11/11/2022

#### 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