

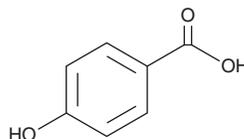
PRODUCT INFORMATION



4-Hydroxybenzoic Acid

Item No. 27468

CAS Registry No.: 99-96-7
Synonyms: *p*-Hydroxybenzoic Acid,
4-Salicyclic Acid, *p*-Salicyclic Acid
MF: C₇H₆O₃
FW: 138.1
Purity: ≥98%
UV/Vis.: λ_{max}: 256 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

4-Hydroxybenzoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-hydroxybenzoic acid in the solvent of choice, which should be purged with an inert gas. 4-Hydroxybenzoic acid is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of 4-hydroxybenzoic acid in these solvents is approximately 5 mg/ml.

4-Hydroxybenzoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-hydroxybenzoic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 4-Hydroxybenzoic 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

4-Hydroxybenzoic acid is a phenolic acid that has been found in a variety of fruit peels and leaves.^{1,2}

References

1. Kim, I., Ku, K.-H., Jeong, M.-C., *et al.* A comparison of the chemical composition and antioxidant activity of several new early- to mid-season apple cultivars for a warmer climate with traditional cultivars. *J. Sci. Food Agric.* (99)10, 4712-4724 (2019).
2. Zheng, Y., Huang, W., Yoo, J.-G., *et al.* Antibacterial compounds from *Siraitia grosvenorii* leaves. *Nat. Prod. Res.* 25(9), 890-897 (2011).

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, 11/28/2022

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