

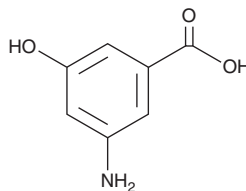
PRODUCT INFORMATION



3-Amino-5-hydroxybenzoic Acid

Item No. 35831

CAS Registry No.: 76045-71-1
Formal Name: 3-amino-5-hydroxy-benzoic acid
Synonym: AHBA
MF: C₇H₇NO₃
FW: 153.1
Purity: ≥98%
UV/Vis.: λ_{max}: 227 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-Amino-5-hydroxybenzoic acid is supplied as a solid. A stock solution may be made by dissolving the 3-amino-5-hydroxybenzoic acid in the solvent of choice, which should be purged with an inert gas. 3-Amino-5-hydroxybenzoic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 3-amino-5-hydroxybenzoic acid in these solvents is approximately 3 and 1 mg/ml, respectively.

Description

3-Amino-5-hydroxybenzoic acid is a biosynthetic precursor of the mC₇N unit of ansamycin antibiotics, including ansatrienin A (Item No. 21995) and geldanamycin (Item No. 13355), as well as the antitumor antibiotic mitomycin C (Item No. 11435).^{1,2}

References

1. Kim, C.-G., Kirschning, A., Bergon, P., *et al.* Biosynthesis of 3-amino-5-hydroxybenzoic acid, the precursor of mC₇N units in ansamycin antibiotics. *J. Am. Chem. Soc.* **118**(32), 7486-7491 (1996).
2. Mao, Y., Varoglu, M., and Sherman, D.H. Molecular characterization and analysis of the biosynthetic gene cluster for the antitumor antibiotic mitomycin C from *Streptomyces lavendulae* NRRL 2564. *Chem. Biol.* **6**(4), 251-263 (1999).

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, 12/15/2022

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