

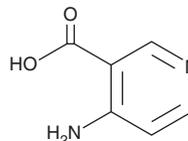
PRODUCT INFORMATION



4-amino Nicotinic Acid

Item No. 20733

CAS Registry No.: 7418-65-7
Formal Name: 4-amino-3-pyridinecarboxylic acid
Synonyms: 4-Aminopyridine-3-Carboxylic Acid, NSC 266130
MF: C₆H₆N₂O₂
FW: 138.1
Purity: ≥98%
UV/Vis.: λ_{max}: 261 nm
Supplied as: A crystalline 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

4-amino Nicotinic acid is supplied as a crystalline solid. Aqueous solutions of 4-amino nicotinic acid can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 4-amino nicotinic acid in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-amino Nicotinic acid is a synthetic intermediate that has been used in various syntheses.^{1,2}

References

1. Heber, D., Ravens, U., and Schulze, T. Synthesis and positive inotropic activity of several 5-aminopyrido[2,3-d]pyrimidines. Part 5: Compounds with positive inotropic activity. *Die. Pharmazie.* **48(7)**, 509-513 (1993).
2. San Feliciano, A., Caballero, E., Puebla, P., *et al.* Synthesis and pharmacological activities of some pyrido[2,1-b]oxazines. *Eur. J. Med. Chem.* **27(5)**, 527-535 (1992).

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