

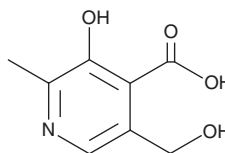
PRODUCT INFORMATION



4-Pyridoxic Acid

Item No. 33968

CAS Registry No.: 82-82-6
Formal Name: 3-hydroxy-5-(hydroxymethyl)-2-methyl-4-pyridinecarboxylic acid
Synonym: NSC 41814
MF: C₈H₉NO₄
FW: 183.2
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 323 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

4-Pyridoxic acid is supplied as a solid. A stock solution may be made by dissolving the 4-pyridoxic acid in the solvent of choice, which should be purged with an inert gas. 4-Pyridoxic acid is soluble in aqueous acid at a concentration of approximately 10 mg/ml.

Description

4-Pyridoxic acid is an inactive metabolite of pyridoxine (Item No. 20706).¹ It is formed from pyridoxine via pyridoxal 5'-phosphate and pyridoxal intermediates by pyridoxine kinase or pyridoxamine phosphate oxidase, as well as pyridoxine-5'-phosphate oxidase, aldehyde dehydrogenase, or aldehyde oxidase.²

References

1. Sharma, S.K. and Dakshinamurti, K. Determination of vitamin B₆ vitamers and pyridoxic acid in biological samples. *J. Chromatogr.* **578**(1), 45-51 (1992).
2. Kohlmeier, M. Water-soluble vitamins and nonnutrients. *Nutrient metabolism*. Kohlmeier, M., editor, 2nd edition, Elsevier (2015).

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, 11/17/2022

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