

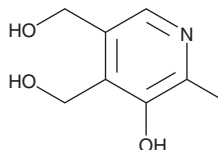
PRODUCT INFORMATION



Pyridoxine

Item No. 20706

CAS Registry No.: 65-23-6
Formal Name: 5-hydroxy-6-methyl-3,4-pyridinedimethanol
MF: $C_8H_{11}NO_3$
FW: 169.2
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 291 nm
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥ 4 years



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

Laboratory Procedures

Pyridoxine is supplied as a crystalline solid. A stock solution may be made by dissolving the pyridoxine in the solvent of choice, which should be purged with an inert gas. Pyridoxine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of pyridoxine in these solvents is approximately 10 and 0.25 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 pyridoxine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of pyridoxine in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pyridoxine is a 4-methanol form of vitamin B₆ that is converted to pyridoxal 5'-phosphate (Item No. 20352), the active form of vitamin B₆ and an important cofactor for metabolism.¹ Pyridoxine has been used in cell culture media as a precursor to pyridoxal 5'-phosphate.²

References

1. Vitamin B₆ (Pyridoxine; Pyridoxal 5'-Phosphate). *Alt. Med. Rev.* **6**(1), 87-92 (2001).
2. Block, G.D. Cell culture media for mammalian cells. *University of Pittsburgh US 7,022,520 B2* (2006).

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, 12/01/2022

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