

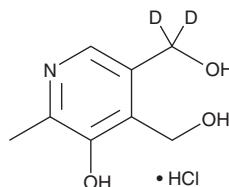
PRODUCT INFORMATION



Pyridoxine-d₂ (hydrochloride)

Item No. 45847

CAS Registry No.: 5027-82-7
Formal Name: 5-hydroxy-6-methyl-3,4-pyridinedimethan- α^3, α^3 -d₂-ol, monohydrochloride
MF: C₈H₉D₂NO₃ • HCl
FW: 207.7
Chemical Purity: ≥98% (pyridoxine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₂); ≤1% d₀
Supplied as: A 1 mg/ml solution in methanol
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

Pyridoxine-d₂ (hydrochloride) is intended for use as an internal standard for the quantification of pyridoxine (Item No. 20706) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

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 and pyridoxal 5'-phosphate). *Altern. Med. Rev.* **6**(1), 87-92 (2001).
2. Block, G.D. Cell culture media for mammalian cells. *University of Pittsburgh*. **US7,022,520B2** (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, 06/23/2026

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