

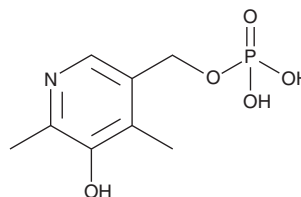
PRODUCT INFORMATION



4-Deoxyribose Phosphate

Item No. 29973

CAS Registry No.: 883-84-1
Formal Name: 5-hydroxy-4,6-dimethyl-3-pyridinemethanol, 3-(dihydrogen phosphate)
Synonyms: 4-Deoxyribose Phosphate, DOP-P, DOP Phosphate, NSC 29870
MF: C₈H₁₂NO₅P
FW: 233.2
Purity: ≥90%
UV/Vis.: λ_{max}: 283 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

4-Deoxyribose phosphate is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-deoxyribose phosphate in water. We do not recommend storing the aqueous solution for more than one day.

Description

4-Deoxyribose phosphate is a vitamin B₆ analog and an inhibitor of sphingosine-1-phosphate (S1P) lyase.¹ It induces accumulation of [³H]sphingosine-1-phosphate ([³H]dhS1P), an S1P lyase substrate, in IT-79MTNC3 cells that endogenously express high levels of S1P lyase (EC₅₀s = <1 and >100 μM in the absence and presence of vitamin B₆, respectively).

Reference

- Ohtoyo, M., Tamura, M., Machinaga, N., *et al.* Scintillation proximity assay to detect the changes in cellular dihydrosphingosine 1-phosphate levels. *Lipids* **51(10)**, 1207-1216 (2016).

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, 06/12/2020

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