

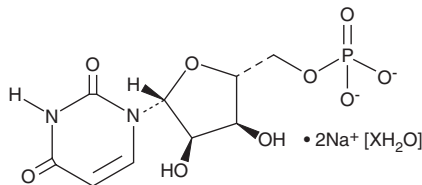
PRODUCT INFORMATION



Uridine-5'-monophosphate (sodium salt hydrate)

Item No. 34525

CAS Registry No.: 681435-27-8
Formal Name: 5'-uridylic acid, disodium salt, hydrate
Synonyms: UMP, 5'-UMP, 5'-Uridylic Acid
MF: $C_9H_{11}N_2O_9P \cdot 2Na [XH_2O]$
FW: 368.1
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 263 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Uridine-5'-monophosphate (UMP) (sodium salt hydrate) is supplied as a solid. Aqueous solutions of UMP (sodium salt hydrate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of UMP (sodium salt hydrate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

UMP is a ribonucleotide.¹ It is formed via decarboxylation of orotidine-5'-monophosphate (5'-OMP) by OMP decarboxylase. UMP is further phosphorylated by UMP-CMP kinase to form UDP (Item No. 18137) and UTP (Item No. 9003530) during nucleic acid biosynthesis. Formulations containing UMP have been used as dietary supplements.

Reference

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. Nucleotide Biosynthesis. *Biochemistry*. 5th ed., W.H. Freeman (2002).

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, 10/17/2022

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