

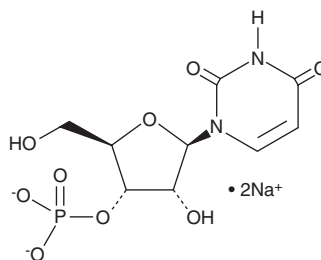
PRODUCT INFORMATION



Uridine-3'-monophosphate (sodium salt)

Item No. 16774

CAS Registry No.: 35170-03-7
Formal Name: 3'-uridylic acid, disodium salt
Synonyms: 3'-UMP, 3'-Uridylic Acid
MF: C₉H₁₁N₂O₉P • 2Na
FW: 368.1
Purity: ≥98%
UV/Vis.: λ_{max}: 262 nm
Supplied as: A crystalline 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

Uridine-3'-monophosphate is supplied as a crystalline solid. Aqueous solutions of uridine-3'-monophosphate can be prepared by directly dissolving the solid in aqueous buffers. The solubility of uridine-3'-monophosphate in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Uridine-3'-monophosphate is a nucleotide used in the biochemical synthesis of ribothymidine-3'-phosphate.¹ It has been used as a bioavailable source of uridine in research studies designed to enhance learning and memory.^{2,3}

References

1. Bowler, F.R., Chan, C.K.W., Duffy, C.D., *et al.* Prebiotically plausible oligoribonucleotide ligation facilitated by chemoselective acetylation. *Nat. Chem.* **5(5)**, 383-389 (2013).
2. Bronk, J.R. and Hastewell, J.G. The transport and metabolism of the uridine mononucleotides by rat jejunum *in vitro*. *J. Physiol.* **408**, 129-135 (1989).
3. Holguin, S., Martinez, J., Chow, C., *et al.* Dietary uridine enhances the improvement in learning and memory produced by administering DHA to gerbils. *FEBS J.* **22(11)**, 3938-3946 (2008).

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, 11/30/2022

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