

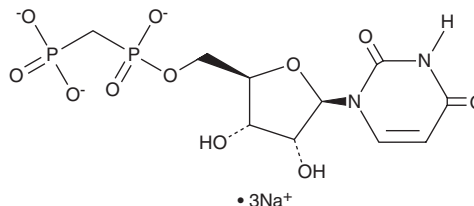
PRODUCT INFORMATION



Uridine 5'-methylenediphosphate (sodium salt)

Item No. 38691

Formal Name: 5'-[hydrogen P-(phosphonomethyl) phosphonate], uridine, trisodium salt
Synonyms: α,β -methylene UDP, UMP-CP
MF: $C_{10}H_{13}N_2O_{11}P_2 \cdot 3Na$
FW: 468.1
Purity: $\geq 95\%$
Supplied as: A solution in water
Storage: $-80^\circ C$
Stability: ≥ 2 years



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

Description

Uridine 5'-methylenediphosphate (UMP-CP) is an agonist of the purinergic P2Y₁₄ receptor and a hydrolytically stable derivative of UDP (Item No. 18137).^{1,2} It selectively inhibits cAMP production in cells expressing human P2Y₁₄ over those expressing human P2Y₆ (EC₅₀s = 11 and 339 nM, respectively).¹

References

1. Das, A., Ko, H., Burianek, L., *et al.* Human P2Y₁₄ receptor agonists: Truncation of the hexose moiety of uridine-5'-diphosphoglucose and its replacement with alkyl and aryl groups. *J. Med. Chem.* **53(1)**, 471-480 (2010).
2. Ko, H., Carter, R.L., Cosyn, L., *et al.* Synthesis and potency of novel uracil nucleotides and derivatives as P2Y₂ and P2Y₆ receptor agonists. *Bioorg. Med. Chem.* **16(12)**, 6319-6332 (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, 08/22/2023

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