

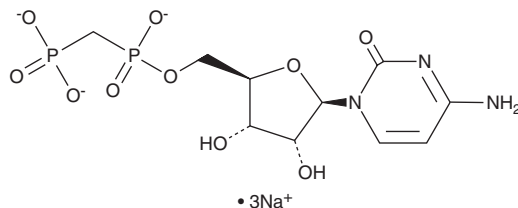
PRODUCT INFORMATION



Cytidine 5'-methylenediphosphate (sodium salt)

Item No. 38372

Formal Name: cytidine, 5'-[hydrogen P-(phosphonomethyl) phosphonate], trisodium salt
Synonyms: CBP, CMP-CP, Cytidine-5'-O-(α,β -methylene)diphosphate
MF: $C_{10}H_{14}N_3O_{10}P_2 \cdot 3Na$
FW: 467.2
Purity: $\geq 95\%$
Supplied as: A solution in water
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Description

Cytidine 5'-methylenediphosphate is an inhibitor of ecto-5'-nucleotidase, also known as CD73 ($K_i = 898$ nM), an enzyme that converts AMP to adenosine.¹

Reference

1. Junker, A., Renn, C., Dobelmann, C., *et al.* Structure-activity relationship of purine and pyrimidine nucleotides as ecto-5'-nucleotidase (CD73) inhibitors. *J. Med. Chem.* **62**(7), 3677-3695 (2019).

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, 04/13/2026

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