

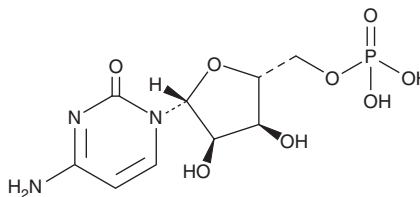
# PRODUCT INFORMATION



## Cytidine 5'-monophosphate

Item No. 35116

**CAS Registry No.:** 63-37-6  
**Formal Name:** 5'-cytidylic acid  
**Synonyms:** CMP, 5'-CMP, Cytidine 5'-phosphate  
**MF:** C<sub>9</sub>H<sub>14</sub>N<sub>3</sub>O<sub>8</sub>P  
**FW:** 323.2  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 276 nm  
**Supplied as:** A 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

Cytidine 5'-monophosphate (CMP) is supplied as a solid. Aqueous solutions of CMP can be prepared by directly dissolving the solid in aqueous buffers. The solubility of CMP in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

CMP is a pyrimidine nucleotide.<sup>1</sup> It is used in the synthesis of RNA and can be further phosphorylated to form cytidine 5'-diphosphate (CDP) and cytidine 5'-triphosphate (CTP). CMP is also an intermediate in the synthesis of phospholipids.

### Reference

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. *Biochemistry*. 5<sup>th</sup> ed., W.H. Freeman, New York (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, 10/05/2022

#### 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