

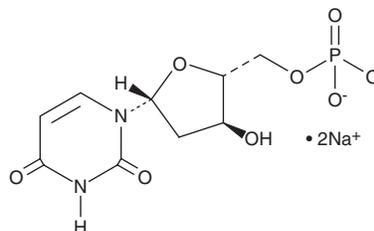
# PRODUCT INFORMATION



## Deoxyuridine 5'-monophosphate (sodium salt)

Item No. 34314

**CAS Registry No.:** 42155-08-8  
**Formal Name:** 2'-deoxy-5'-uridylic acid, disodium salt  
**Synonym:** dUMP  
**MF:**  $C_9H_{11}N_2O_8P \cdot 2Na$   
**FW:** 352.1  
**Purity:**  $\geq 95\%$   
**UV/Vis.:**  $\lambda_{max}$ : 263 nm  
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Deoxyuridine 5'-monophosphate (dUMP) (sodium salt) is supplied as a solid. Aqueous solutions of dUMP (sodium salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of dUMP (sodium salt) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

dUMP is a nucleotide intermediate in the biosynthesis of 5'-thymidylic acid (dTMP; Item No. 33549).<sup>1</sup>

### Reference

1. Berg, J.M., Tymoczko, J.L., and Stryer, L. Nucleotide Biosynthesis. *Biochemistry*. 5<sup>th</sup> edition, 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, 10/18/2022

#### 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