

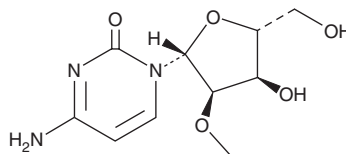
# PRODUCT INFORMATION



## 2'-O-Methylcytidine

Item No. 33411

CAS Registry No.: 2140-72-9  
Formal Name: 2'-O-methyl-cytidine  
MF: C<sub>10</sub>H<sub>15</sub>N<sub>3</sub>O<sub>5</sub>  
FW: 257.2  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 271 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

2'-O-Methylcytidine is supplied as a crystalline solid. A stock solution may be made by dissolving the 2'-O-methylcytidine in the solvent of choice, which should be purged with an inert gas. 2'-O-Methylcytidine is soluble in organic solvents such as DMSO at a concentration of approximately 0.2 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 2'-O-methylcytidine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2'-O-methylcytidine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2'-O-Methylcytidine is a nucleoside derivative and an inhibitor of hepatitis C virus (HCV) non-structural protein 5B (NS5B; IC<sub>50</sub> = 3.8 μM).<sup>1</sup> It inhibits HCV replication in a replicon assay (IC<sub>50</sub> = 21.2 μM).

### Reference

1. Migliaccio, G., Tomassini, J.E., Carroll, S.S., *et al.* Characterization of resistance to non-obligate chain-terminating ribonucleoside analogs that inhibit hepatitis C virus replication *in vitro*. *J. Bio. Chem.* **278**(49), 49164-49170 (2003).

#### 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, 11/16/2022

#### 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