

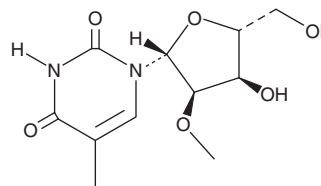
PRODUCT INFORMATION



2'-O-Methyl-5-methyluridine

Item No. 35357

CAS Registry No.: 55486-09-4
Formal Name: 5-methyl-2'-O-methyl-uridine
Synonyms: 2'-O-Methylribothymidine, 2'-O-Methylthymidine
MF: $C_{11}H_{16}N_2O_6$
FW: 272.3
Purity: $\geq 98\%$
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

2'-O-Methyl-5-methyluridine is supplied as a solid. A stock solution may be made by dissolving the 2'-O-methyl-5-methyluridine in the solvent of choice, which should be purged with an inert gas. 2'-O-Methyl-5-methyluridine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2'-O-methyl-5-methyluridine in these solvents is approximately 10 and 14 mg/ml, respectively. 2'-O-Methyl-5-methyluridine is also slightly soluble in ethanol.

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-methyl-5-methyluridine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 2'-O-methyl-5-methyluridine in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2'-O-Methyl-5-methyluridine is a modified nucleoside that has been found in tRNA isolated from rabbit liver.¹

Reference

1. Gross, H.J., Simsek, M., Raba, M., *et al.* 2'-O-methyl ribothymidine: A component of rabbit liver lysine transfer RNA. *Nucleic Acids Res.* **1(1)**, 35-43 (1974).

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, 12/12/2022

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