

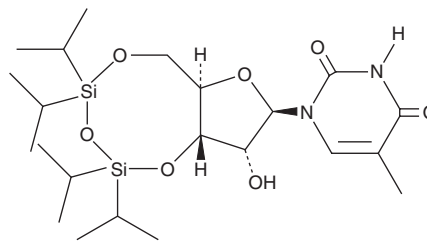
# PRODUCT INFORMATION



## 5',3'-TIPS 5-Methyluridine

Item No. 44624

**CAS Registry No.:** 130983-87-8  
**Formal Name:** 5-methyl-3',5'-O-[1,1,3,3-tetrakis(1-methylethyl)-1,3-disiloxanediyl]-uridine  
**MF:** C<sub>22</sub>H<sub>40</sub>N<sub>2</sub>O<sub>7</sub>Si<sub>2</sub>  
**FW:** 500.7  
**Purity:** ≥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

5',3'-TIPS 5-methyluridine is supplied as a solid. A stock solution may be made by dissolving the 5',3'-TIPS 5-methyluridine in the solvent of choice, which should be purged with an inert gas. 5',3'-TIPS 5-methyluridine is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

### Description

5',3'-TIPS 5-methyluridine is a modified nucleoside and derivative of 5-methyluridine (Item No. 27986). It inhibits the growth of U373 MG and Hs 683 glioma cells (GI<sub>50</sub>s = 32 and 29 μM, respectively).<sup>1</sup>

### Reference

1. Panayides, J.-L., Mathieu, V., Banuls, L.M., *et al.* Synthesis and in vitro growth inhibitory activity of novel silyl- and trityl-modified nucleosides. *Bioorg. Med. Chem.* **24(12)**, 2716-2724 (2016).

#### 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, 03/03/2026

#### 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