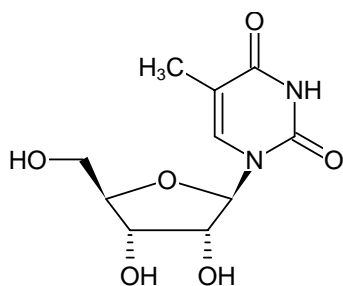


**5-Methyl-uridine**

Cat. No.	Amount
N-1057	100 mg



Structural formula of 5-Methyl-uridine

For research use only!**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 24 months after date of delivery**Molecular Formula:** C₁₀H₁₄N₂O₆**Molecular Weight:** 258.23 g/mol**CAS#:** 1463-10-1**EC number:** 215-973-9**Purity:** > 98 %**Form:** solid**Selected References:**Bullinger *et al.* (2007) Metabolic signature of breast cancer cell line MCF-7: profiling of modified nucleosides via LC-IT MS coupling. *BMC Biochem* **29**:25.Glenn *et al.* (1971) Analysis of 5-Methyluridine Function in the Transfer RNA of *Escherichia coli*. *Cancer Research* **31**:706.