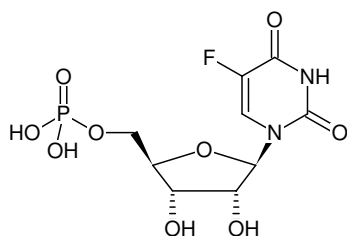


**5-Fluoro-UMP**

(5F-UMP)

5-Fluoro-uridine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-146S	50 µl (10 mM)
NU-146L	5 x 50 µl (10 mM)



Structural formula of 5-Fluoro-UMP

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

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

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₉H₁₂N₂O₉PF (free acid)**Molecular Weight:** 342.17 g/mol (free acid)**Exact Mass:** 342.03 g/mol (free acid)**CAS#:** 796-66-7**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 267 nm, ε 10.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**

Labesse *et al.* (2011) Structural and functional characterization of the Mycobacterium tuberculosis uridine monophosphate kinase: insights into the allosteric regulation. *Nucleic Acids Res.* **39** (8):3458.

Gilles *et al.* (2007) Regulatory Mechanisms Differ in UMP Kinases from Gram-negative and Gram-positive Bacteria. *J. Biol. Chem.* **282** (10):7242.