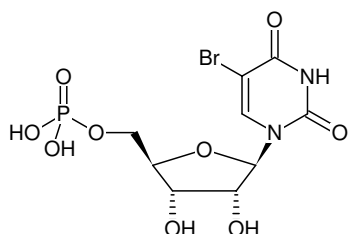


**5-Bromo-UMP**

(5Br-UMP)

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

Cat. No.	Amount
NU-135S	100 µl (10 mM)
NU-135L	5 x 100 µl (10 mM)



Structural formula of 5-Bromo-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:** C₉H₁₂BrN₂O₉P (free acid)**Molecular Weight:** 403.08 g/mol (free acid)**Exact Mass:** 401.95 g/mol (free acid)**CAS#:** 2149-79-3**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} 278 nm, ε 9.7 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.

Kwon *et al.* (1999) Bipartite modular structure of intrinsic, RNA hairpin-independent termination signal for phage RNA polymerases. *J. Biol. Chem.* **274** (41):29149.

Tourneux *et al.* (1998) Substitution of an alanine residue for glycine 146 in TMP kinase from *Escherichia coli* is responsible for bacterial hypersensitivity to bromodeoxyuridine. *J. Bacteriol.* **180** (16):4291.