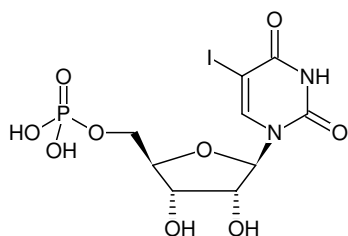


**5-Iodo-UMP**

(5I-UMP)

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

Cat. No.	Amount
NU-868S	50 µl (10 mM)
NU-868L	5 x 50 µl (10 mM)



Structural formula of 5-Iodo-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:** C<sub>9</sub>H<sub>12</sub>N<sub>2</sub>O<sub>9</sub>PI (free acid)**Molecular Weight:** 450.08 g/mol (free acid)**Exact Mass:** 449.93 g/mol (free acid)**CAS#:** 34198-43-1**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<sub>max</sub> 287 nm, ε 7.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)