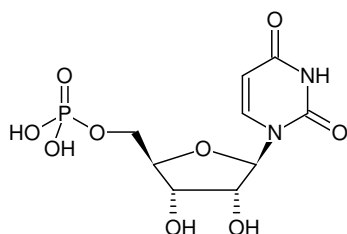


**UMP - Solid**

Uridine-5'-monophosphate, Disodium salt

Cat. No.	Amount
NU-1033-5G	5 g
NU-1033-25G	25 g



Structural formula of UMP - Solid

**For research use only!****Shipping:** shipped on blue ice**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>13</sub>N<sub>2</sub>O<sub>9</sub>P (free acid)**Molecular Weight:** 324.18 g/mol (free acid)**Exact Mass:** 324.04 g/mol (free acid)**CAS#:** 3387-36-8**EC number:** 222-211-9**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)