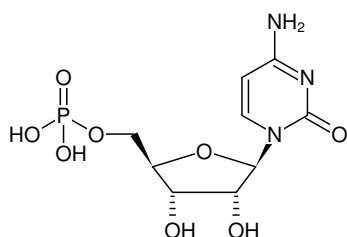


**CMP - Solid**

Cytidine-5'-monophosphate, Disodium salt

Cat. No.	Amount
NU-1032-5G	5 g
NU-1032-25G	25 g



Structural formula of CMP - 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:** C₉H₁₄N₃O₈P (free acid)**Molecular Weight:** 323.20 g/mol (free acid)**Exact Mass:** 323.05 g/mol (free acid)**CAS#:** 6757-06-8**EC number:** 229-819-3**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)