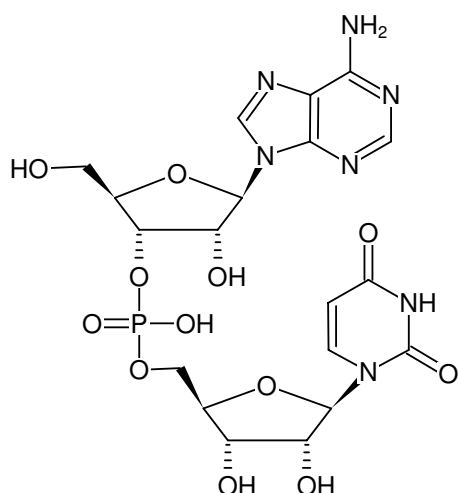


**ApU**

RNA Dinucleotide (5'→3'), Sodium salt

Cat. No.	Amount
NU-917S	10 µl (100 mM)
NU-917L	5 x 10 µl (100 mM)



Structural formula of ApU

For research use only!**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:** C₁₉H₂₄N₇O₁₂P (free acid)**Molecular Weight:** 573.41 g/mol (free acid)**Exact Mass:** 573.12 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm, ε 25.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)