

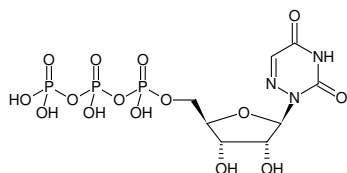
**6-Aza-UTP**

6-Azauridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-896S	10 µl (100 mM)
NU-896L	5 x 10 µl (100 mM)

Selected References:

Albulescu *et al.* (2014) An in vitro assay to study chikungunya virus RNA synthesis and the mode of action of inhibitors *J. Gen. Virol.* **95** (12):2683.



Structural formula of 6-Aza-UTP

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:** C₈H₁₄N₃O₁₅P₃ (free acid)**Molecular Weight:** 485.13 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm; ε 6.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)