

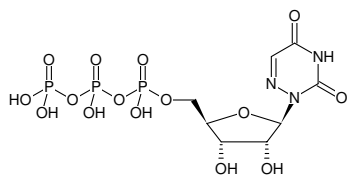
**6-Aza-UTP**

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 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>8</sub>H<sub>14</sub>N<sub>3</sub>O<sub>15</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 485.13 g/mol (free acid)**Exact Mass:** 484.96 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<sub>max</sub> 260 nm, ε 6.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)