

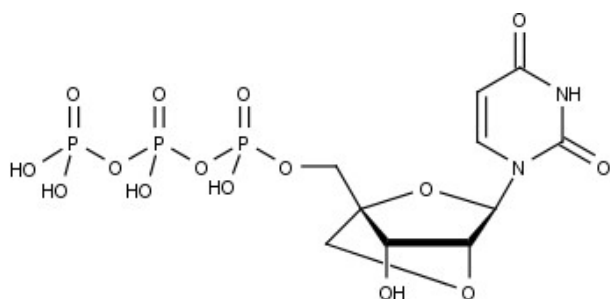
**LNA-UTP**

LNA-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-985	10 µl (100 mM)

Related Products:

HighFidelity LNA PCR Labeling Kit, #APP-101-LNA

Selected References:Veedu *et al.* (2010) Polymerase-directed synthesis of C5-ethynyl locked nucleic acids. *Bioorg. Med. Chem. Lett.* **20** (22):565.

Structural formula of LNA-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:** C₁₀H₁₅N₂O₁₅P₃ (free acid)**Molecular Weight:** 496.15 g/mol (free acid)**Exact Mass:** 495.97 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} 262 nm, ε 9.8 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)