

**5-Vinyl-UTP**

Sodium salt

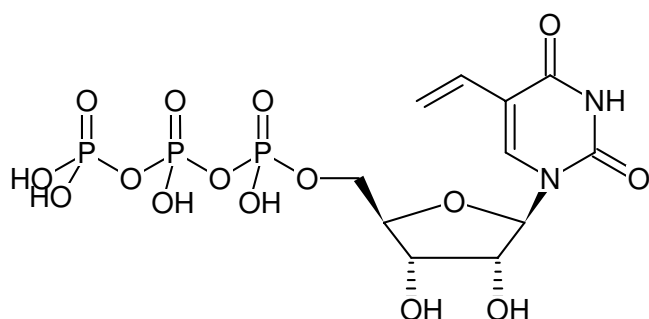
Cat. No.	Amount
CLK-069	50 µl (10 mM)

Applications:

Incorporation into RNA/cRNA by
- *in vitro* Transcription with T7 RNA polymerase ^[1]

Selected References:

[1] George *et al.* (2017) Vinyluridine as a Versatile Chemoselective Handle for the Posttranscriptional Chemical Functionalization of RNA. *Bioconjugate Chem.*:DOI: 10.1021/acs.bioconjchem.7b00169.



Structural formula of 5-Vinyl-UTP

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:** C₁₁H₁₇N₂O₁₅P₃ (free acid)**Molecular Weight:** 510.18 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** sterile solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 292 nm, ε 9.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)