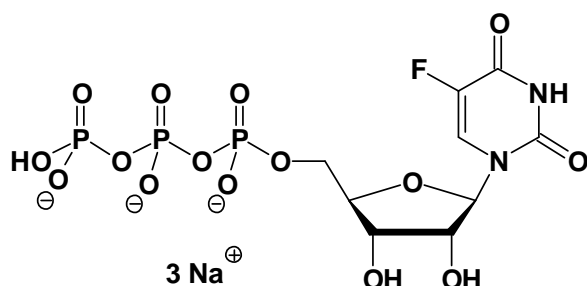


## 5F-UTP

### 5-Fluoro-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1115S	100 Units
NU-1115L	500 Units



**Cat. No.:** NU-1115

**Molecular Formula:** C<sub>9</sub>H<sub>11</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub>F (Anion)

**Molecular Weight:** 499.11 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

#### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

#### For research use only!

1 unit = 1 µl of a 10 mM solution

#### Selected References:

Au *et al.* (1982) Reversed-phase ion-pair high-performance liquid-chromatographic assay of 5-fluorouracil, 5'-deoxy-5-fluorouridine, their nucleosides, monophosphate, diphosphate, and triphosphate nucleotides with a mixture of quaternary ammonium-ions. *J. Chromatogr.* **228**:245.

Glazer *et al.* (1980) The effect of 5-fluorouridine 5'-triphosphate on RNA transcribed in isolated-nuclei invitro. *Mol. Pharmacol.* **17** (2):279.

Anukarahanonta *et al.* (1979) Enhancement of 5-fluorouridine action in hepatoma-cells by galactosamine-induced uridine triphosphate deficiency. *H-S Z. Physiol. Chem.* **360** (3):225.