

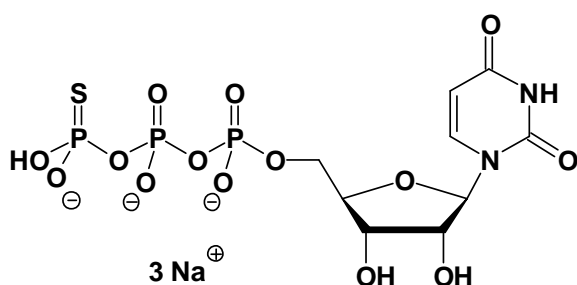
## UTP $\gamma$ S

Uridine-5'-( $\gamma$ -thio)-triphosphate, Sodium salt

Cat. No.	Amount
NU-416S	20 Units
NU-416L	100 Units

### Selected References:

Read *et al.* (2005) UTP-dependent turnover of *Trypanosoma brucei* mitochondrial mRNA requires UTP polymerization and involves the RET1 TUTase. *RNA* **11**:763.



**Cat. No.:** NU-416

**Molecular Formula:** C<sub>9</sub>H<sub>12</sub>N<sub>2</sub>O<sub>14</sub>P<sub>3</sub>S (Anion)

**Molecular Weight:** 497.18 (Anion)

**Purity:** 80 %, clear aqueous solution, pH 7.5  
contains ca. 20 % UDP, which UTP $\gamma$ S  
decomposes to at room temperature

### 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 3 months after date of delivery.

### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution