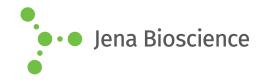
# **DATA SHEET**

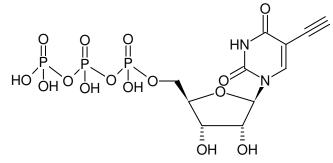




## ■ 5-Ethynyl-UTP (5-EUTP)

5-Ethynyl-uridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-T08-S	5 μl (100 mM)
CLK-T08-L	5 x 5 μl (100 mM)
CLK-T08-XL	50 μl (100 mM)



Structural formula of 5-Ethynyl-UTP (5-EUTP)

#### 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>11</sub>H<sub>15</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub> (free acid) **Molecular Weight:** 508.16 g/mol (free acid)

Exact Mass: 507.97 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:  $\lambda_{max}$  288 nm,  $\epsilon$  12.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl

pH 7.5)

#### **Applications:**

Incorporation into RNA by T7 RNA polymerase-mediated in vitro transcription.

The resulting alkyne-functionalized RNA can subsequently be processed via Cu(I)-catalyzed (azide-alkyne) click chemistry that offers the choice

- to introduce a Biotin group for subsequent purification tasks (via Azides of Biotin)
- to introduce fluorescent group for subsequent microscopic imaging (via Azides of fluorescent dyes)
- to crosslink the RNA to azide-functionalized biomolecules e.g.proteins

Presolski et al.<sup>[1]</sup> and Hong et al.<sup>[2]</sup> provide a general protocol for Cu(I)-catalyzed click chemistry reactions that may be used as a starting point for the set up and optimization of individual assays.

### **Related Products:**

Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-MI004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

#### Selected References:

[1] Presolski et al. (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. Current Protocols in Chemical Biology 3:153.

[2] Hong et al. (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. Angew. Chem. Int. Ed. **48**:9879.