





# 2-Ethynyl-ATP (2-EATP)

2-Ethynyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-NU-004S	100 μl (10 mM)
CLK-NU-004L	5 x 100 μl (10 mM)



Structural formula of 2-Ethynyl-ATP (2-EATP)

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>12</sub>H<sub>16</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub>

Molecular Weight: 531.20 g/mol

Exact Mass: 531.00 g/mol

**Purity:** ≥ 90 % (HPLC), contains approx. 6 % 2-Ethynyl-ADP

Form: solution in 100 mM Tris-HCl

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

**pH:** 7.5 ±0.5

Spectroscopic Properties:  $\lambda_{max}$  265 nm,  $\epsilon$  10.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

## **Applications:**

in vitro polyadenylation of RNA<sup>[1]</sup>

## **Description**:

2-Ethynyl-labeled adenosine triphosphate (2-EATP) is suitable for *in vitro* polyadenylation of RNA with recombinant poly(A) polymerase<sup>[1]</sup>.

The resulting Alkyne-functionalized RNA can subsequently be processed via Cu(I)-catalyzed Azide-Alkyne click chemistry (CUAAC) 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>[2]</sup> and Hong *et al.*<sup>[3]</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:**

5-Ethynyl-adenosine (5-EA), #CLK-N005 Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-M1004 Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010 Sodium Ascorbate (Na-Ascorbate), #CLK-M1005

#### Selected References:

[1] Curanovic *et al.* (2013) Global profiling of stimulus-induced polyadenylation in cells using a poly (A) trap. *Nat. Chem. Biol.* **9**:671.

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

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

