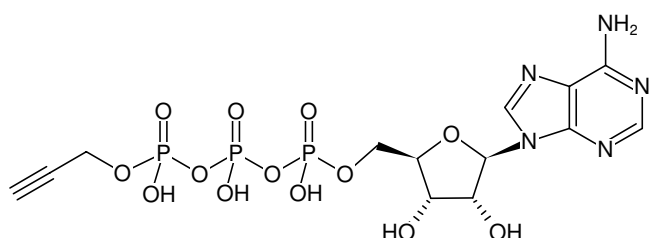


**γ-Propargyl-ATP**

Adenosine-5'-[γ-(propargyl)]triphosphate, Sodium salt

| Cat. No.  | Amount |
|-----------|--------|
| CLK-T10-1 | 1 mg   |



Structural formula of γ-Propargyl-ATP

**For research use only!****Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dry and under inert gas

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>13</sub>H<sub>18</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 545.23 g/mol (free acid)**Exact Mass:** 545.01 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)