



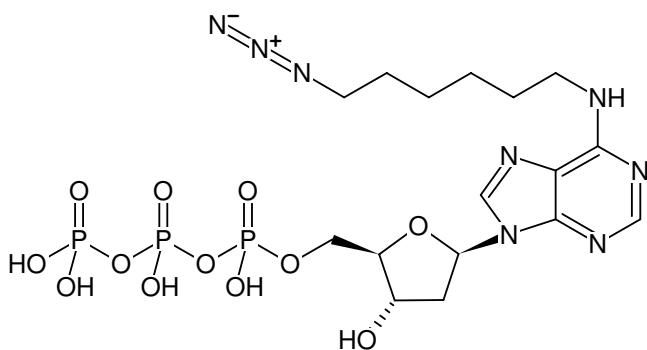
## N<sup>6</sup>-(6-Azido)hexyl-dATP

N<sup>6</sup>-(6-Azido)hexyl-2'-deoxy-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
CLK-NU-002-S	50 µl (10 mM)
CLK-NU-002-L	5 x 50 µl (10 mM)

### Selected References:

Ando *et al.* (2019) ELTA: Enzymatic Labeling of Terminal ADP-Ribose. *Mol. Cell.* **73** (4):845.



Structural formula of N<sup>6</sup>-(6-Azido)hexyl-dATP

### For research use only!

**Shipping:** shipped on blue ice

**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>16</sub>H<sub>27</sub>N<sub>8</sub>O<sub>12</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 616.35 g/mol (free acid)

**Exact Mass:** 616.10 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** sterile solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>max</sub> 266 nm, ε 16.2 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)