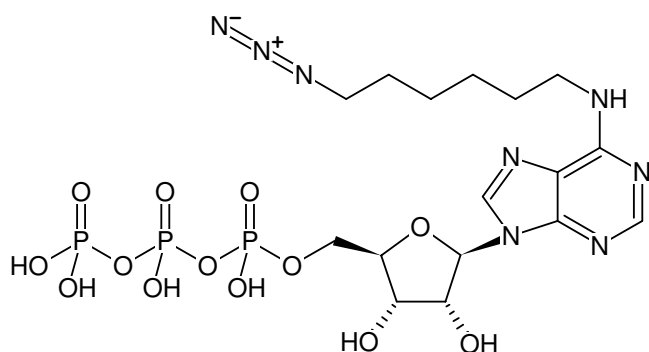




### N<sup>6</sup>-(6-Azido)hexyl-ATP

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

Cat. No.	Amount
CLK-NU-003-1	1 mg
CLK-NU-003-5	5 mg



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

**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>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 632.35 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**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)