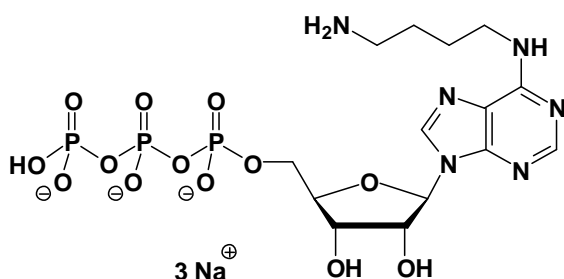


## N<sup>6</sup>-(4-Amino)butyl-ATP

N<sup>6</sup>-(4-Amino)butyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-804S	50 Units
NU-804L	250 Units



**Cat. No.:** NU-804

**Molecular Formula:** C<sub>14</sub>H<sub>22</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 575.28 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

### For research use only!

1 unit = 1 μl of a 10 mM solution