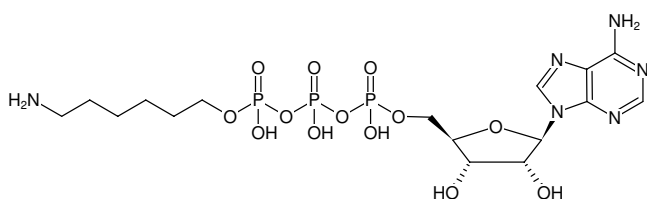




### $\gamma$ -(6-Aminoethyl)-ATP

$\gamma$ -(6-Aminoethyl)-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-833S	50 $\mu$ l (10 mM)
NU-833L	5 x 50 $\mu$ l (10 mM)



Structural formula of  $\gamma$ -(6-Aminoethyl)-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>29</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 606.36 g/mol (free acid)

**Exact Mass:** 606.10 g/mol (free acid)

**Purity:**  $\geq$  95 % (HPLC)

**Form:** clear aqueous solution

**Concentration:** 10 mM - 11 mM

**pH:** 7.5  $\pm$  0.5

**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  15.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)