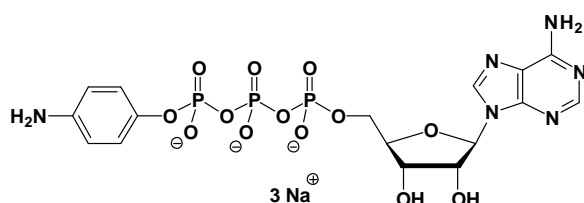


## $\gamma$ -Aminophenyl-ATP

Adenosine-5'-[ $\gamma$ -(4-aminophenyl)]triphosphate, Sodium salt

Cat. No.	Amount
NU-801S	3 mg
NU-801L	15 mg



**Cat. No.:** NU-801

**Molecular Formula:** C<sub>16</sub>H<sub>18</sub>N<sub>6</sub>O<sub>13</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 574.32 (Anion)

**Purity:** > 95%, lyophilized solid

### 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!**

### Selected References:

Haystead *et al.* (1993) Gamma-phosphate-linked ATP-Sepharose for the affinity purification of protein-kinases - rapid purification to homogeneity of skeletal-muscle mitogen-activated protein-kinase kinase. *Eur. J. Biochem.* **214** (2):459.

Trayer *et al.* (1974) Preparation of adenosine nucleotide derivatives suitable for affinity chromatography. *Biochem. J.* **139** (3):609.