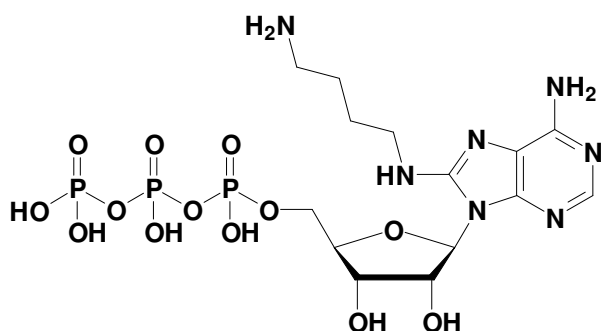


8-[(4-Amino)butyl]-amino-ATP

8-[(4-Amino)butyl]-amino-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-806S	50 Units
NU-806L	250 Units



Cat. No.: NU-806

Molecular Formula: $C_{14}H_{26}N_7O_{13}P_3$ (free acid)

Molecular Weight: 593.32 (free acid)

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 μ l of a 10 mM solution