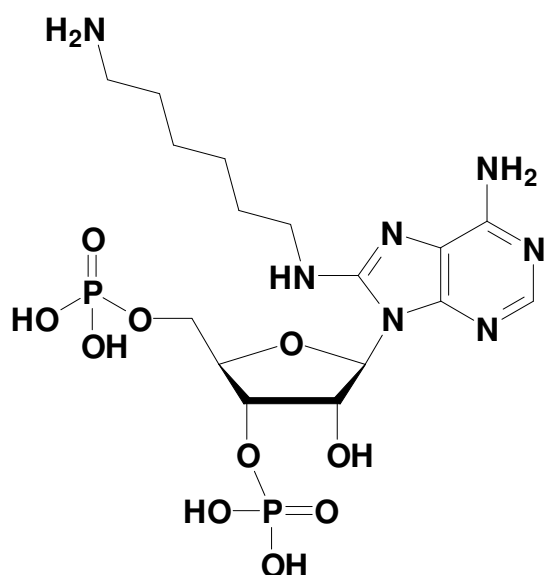


## 8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate

8-[(6-Amino)hexyl]-amino-adenosine-3'/5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-811S	30 Units
NU-811L	150 Units



**Cat. No.:** NU-811

**Molecular Formula:**  $C_{16}H_{29}N_7O_{10}P_2$  (free acid)

**Molecular Weight:** 541.39 (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  $\mu$ l of a 10 mM solution