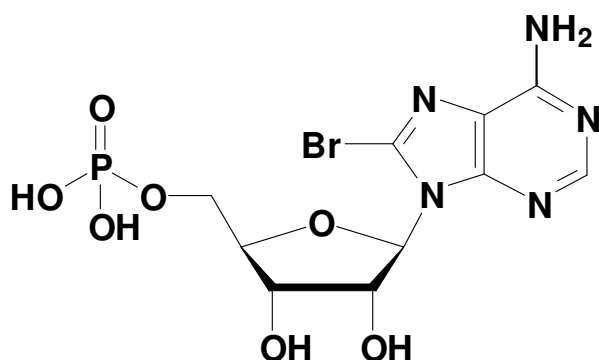


## 8Br-AMP, lyophilized

8-Bromo-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-156-5	5 mg
NU-156-25	25 mg



**Cat. No.:** NU-156

**Molecular Formula:**  $C_{10}H_{13}BrN_5O_7P$  (free acid)

**Molecular Weight:** 426.12 (free acid)

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

**Spectroscopic Properties:**  $\lambda_{max}$  264 nm;  
 $\epsilon$  17,000  $M^{-1}cm^{-1}$  Tris HCl pH 7.5

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at  $< -20^{\circ}C$ . 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:

Ke *et al.* (2007) Crystal structures of Phosphodiesterases and Implications on Substrate Specificity and Inhibitor Selectivity. *Curr. Topics Medicinal Chem.* **7 (4)**:391.

Lascu *et al.* (1979) Enzymatic Properties of 8-Bromoadenine Nucleotides. *Biochemistry* **18 (22)**:4818.