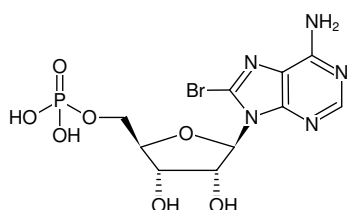


**8-Bromo-AMP**

(8Br-AMP)

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

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



Structural formula of 8-Bromo-AMP

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₁₀H₁₃BrN₅O₇P (free acid)**Molecular Weight:** 426.12 g/mol (free acid)**Exact Mass:** 424.97 g/mol (free acid)**CAS#:** 23567-96-6**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 264 nm, ε 17.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**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.