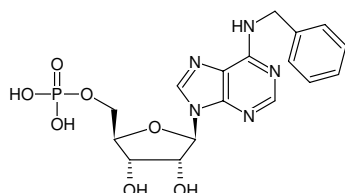


**N⁶-Benzyl-AMP**N⁶-Benzyl-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-1195S	300 µl (10 mM)
NU-1195L	5 x 300 µl (10 mM)

Selected References:

Eguchi *et al.* (1977) Synthesis of AMP analogs and their use for studies on the allosteric site of rabbit muscle glycogen phosphorylase b. *J. Biochem.* **81** (5):1401.

Structural formula of N⁶-Benzyl-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:** C₁₇H₂₀N₅O₇P (free acid)**Molecular Weight:** 437.35 g/mol (free acid)**Exact Mass:** 437.11 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 269 nm, ε 20.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)