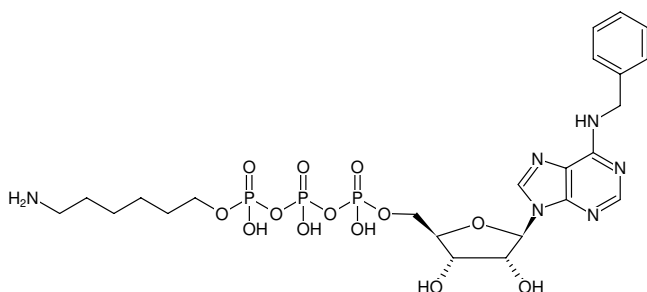


 **γ -(6-Aminoethyl)-N6-Benzyl-ATP** γ -(6-Aminoethyl)-N⁶-benzyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-911-1	1 mg

Structural formula of γ -(6-Aminoethyl)-N⁶-Benzyl-ATP**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:** C₂₃H₃₅N₆O₁₃P₃ (free acid)**Molecular Weight:** 696.48 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{\max} 269 nm, ϵ 20.5 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)