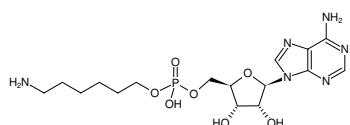


 **α -(6-Aminoethyl)-AMP** α -(6-Aminoethyl)-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-859-1	1 mg
NU-859-5	5 mg

Structural formula of α -(6-Aminoethyl)-AMP**For research use only!****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:** C₁₆H₂₇N₆O₇P (free acid)**Molecular Weight:** 446.40 g/mol (free acid)**Exact Mass:** 446.17 g/mol (free acid)**Purity:** \geq 95 % (HPLC)**Form:** solid**Color:** white to off-white