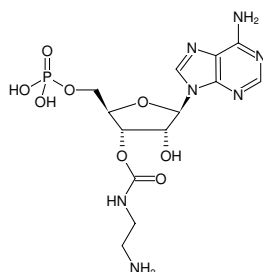
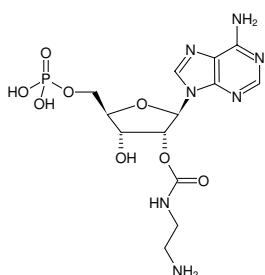


**EDA-AMP**

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-909-1	1 mg
NU-909-5	5 mg



Structural formula of EDA-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:** C₁₃H₂₀N₇O₈P (free acid)**Molecular Weight:** 433.11 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)