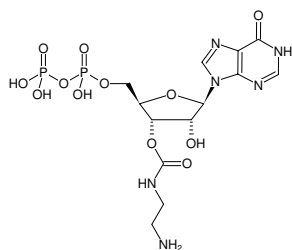
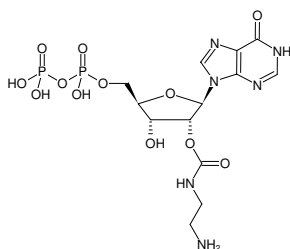


**EDA-IDP**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Inosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-837S	50 µl (10 mM)
NU-837L	5 x 50 µl (10 mM)



Structural formula of EDA-IDP

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:** C₁₃H₂₀N₆O₁₂P₂ (free acid)**Molecular Weight:** 514.28 g/mol (free acid)**Exact Mass:** 514.06 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