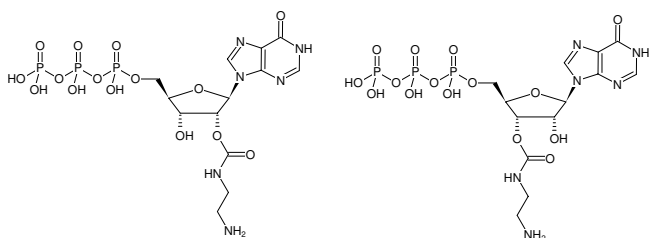


**EDA-ITP**

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

Cat. No.	Amount
NU-836S	100 µl (10 mM)
NU-836L	5 x 100 µl (10 mM)



Structural formula of EDA-ITP

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:** C₁₃H₂₁N₆O₁₅P₃ (free acid)**Molecular Weight:** 594.26 g/mol (free acid)**Exact Mass:** 594.03 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