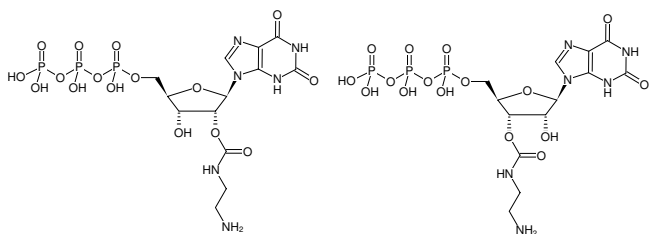


**EDA-XTP**

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

Cat. No.	Amount
NU-838S	100 µl (10 mM)
NU-838L	5 x 100 µl (10 mM)



Structural formula of EDA-XTP

**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:** C<sub>13</sub>H<sub>21</sub>N<sub>6</sub>O<sub>16</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 610.26 g/mol (free acid)**Exact Mass:** 610.02 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**Spectroscopic Properties:** λ<sub>max</sub> 276 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)