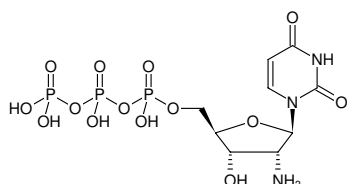


**2'NH₂-dUTP**

2'-Amino-dUTP

2'-Amino-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-242S	50 µl (100 mM)
NU-242L	5 x 50 µl (100 mM)

Structural formula of 2'NH₂-dUTP**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:** C₉H₁₆N₃O₁₄P₃ (free acid)**Molecular Weight:** 483.16 g/mol (free acid)**Exact Mass:** 482.98 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 262 nm, ε 10.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Govindan *et al.* (2012) P2Y receptor subtypes evoke different Ca²⁺ signals in cultured aortic smooth muscle cells. *Purinergic Signalling* **8**:763.Lauridsen *et al.* (2012) Enzymatic recognition of 2'-modified ribonucleoside 5'-triphosphates: towards the evolution of versatile aptamers. *ChemBioChem* **13**:19.Kanwar *et al.* (2011) Chimeric aptamers in cancer cell-targeted drug delivery. *Crit. Rev. Biochem. Mol. Biol.* **46** (6):459.Lin *et al.* (1994) Modified RNA sequence pools for in vitro selection. *Nucleic Acids Research* **22** (24):5229.