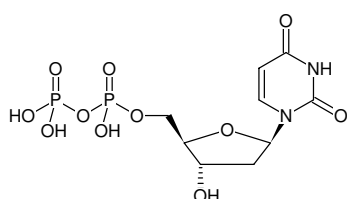


**dUDP**

2'-Deoxyuridine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-901S	150 µl (10 mM)
NU-901L	5 x 150 µl (10 mM)



Structural formula of dUDP

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:** C₉H₁₄N₂O₁₁P₂ (free acid)**Molecular Weight:** 388.16 g/mol (free acid)**Exact Mass:** 388.01 g/mol (free acid)**CAS#:** 4208-67-7**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_{max} 262 nm, ε 10.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Selected References:**Tóth *et al.* (2007) Kinetic Mechanism of Human dUTPase, an Essential Nucleotide Pyrophosphatase Enzyme. *J. Biol. Chem.* **282** (46):33572.Hidalgo-Zarco *et al.* (2001) Kinetic properties and inhibition of the dimeric dUTPase-dUDPase from *Leishmania major*. *Protein Sci.* **10** (7): 1426.Persson *et al.* (2001) Crystallization and preliminary crystallographic analysis of deoxyuridine 5'-triphosphate nucleotidohydrolyase from *Bacillus subtilis*. *Acta Cryst. D* **57**:876.Prasad *et al.* (2000) Structures of feline immunodeficiency virus dUTP pyrophosphatase and its nucleotide complexes in three crystal forms. *Acta Cryst. D* **56**:1100.Camacho *et al.* (2000) Properties of *Leishmania major* dUTP nucleotide-hydrolysing enzyme in kinetoplastids. *Biochem. J.* **346**:163.Dauter *et al.* (1999) Crystal structure of dUTPase from equine infectious anaemia virus, active site metal binding in a substrate analogue complex. *J. Mol. Biol.* **285** (2):655.Dauter *et al.* (1998) The refined structure of dUTPase from *Escherichia coli*. *Acta Cryst. D* **54**:735.Vertessy *et al.* (1998) The complete triphosphate moiety of non-hydrolyzable substrate analogues is required for a conformational shift of the flexible C-terminus in *E. coli* dUTP pyrophosphatase. *FEBS Lett.* **421** (1):83.Shao *et al.* (1997) Characterization and mutational studies of equine infectious anemia virus dUTPase. *BBA-Protein Struct. M* **1339** (2):181.Persson *et al.* (1997) Crystallization and preliminary X-ray crystallographic studies of dUTPase from equine infectious anemia virus. *Protein Peptide Lett.* **4** (2):145.Larsson *et al.* (1996) Kinetic characterization of dUTPase from *Escherichia coli*. *J. Biol. Chem.* **271** (39):24010.Larsson *et al.* (1996) Crystal structure of the *Escherichia coli* dUTPase in complex with a substrate analogue (dUDP). *Nat. Struct. Biol.* **3** (6):532.