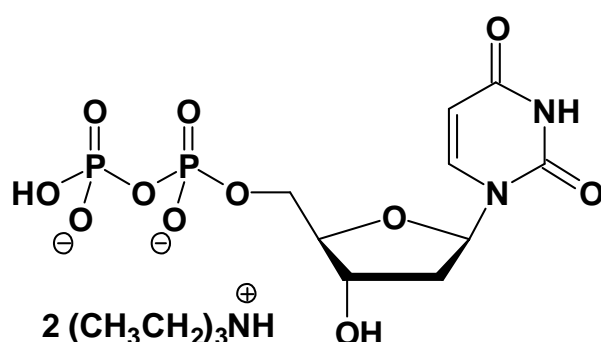


dUDP

2'-Deoxy-uridine-5'-diphosphate, Triethylammonium salt

Cat. No.	Amount
NU-901S	150 Units
NU-901L	750 Units



Cat. No.: NU-901

Molecular Formula: C₉H₁₂N₂O₁₁P₂ (Anion)

Molecular Weight: 386.14 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μl of a 10 mM solution

Selected References:

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 nucleotidohydrolase 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 nucleotidohydrolase, a distinct 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.