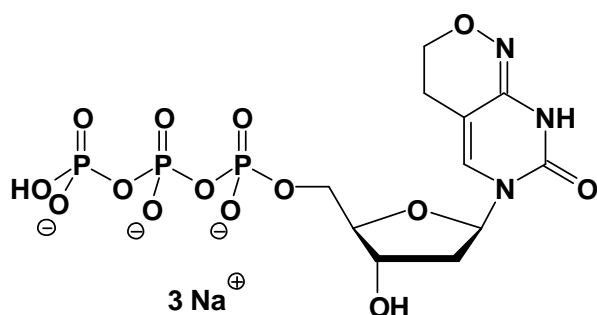


dPTP

6H,8H-3,4-Dihydro-pyrimido[4,5-c][1,2]oxazin-7-one-8-
 β -D-2'-deoxy-ribofuranosid-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1119S	30 Units
NU-1119L	150 Units



Cat. No.: NU-1119

Molecular Formula: C₁₁H₁₅N₃O₁₄P₃ (Anion)

Molecular Weight: 506.17 (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 μ l of a 10 mM solution

Selected References:

Cummins *et al.* (2001) Synthesis and study of the fluorescein conjugate of the nucleotide dPTP. *Nucleos. Nucleot. Nucl.* **20** (4-7):1049.

Zaccolo *et al.* (1996) An approach to random mutagenesis of DNA using mixtures of Triphosphate derivatives of Nucleoside analogues. *J. Mol. Biol.* **255** (4):589.

Zaccolo, M. *et al.* (1996) *J. Mol. Biol.* **255**:589

Lethal mutagenesis:
 Loeb, L. *et al.* (1999) *Proc. Natl. Acad. Sci. USA* **96**:1492