





T4 dNMP (deoxy-Nucleotide Monophosphate) Kinase

bacteriophage T4, recombinant, E. coli

Cat. No.	Amount
PR-340	100 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 27 kDa

Accession number: AAD42414

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 100 mM Tris-HCl pH 7.5, 10 mM $\rm MgCl_2$ and 50 % glycerol)

Description:

NMP kinases catalyse the phosphorylation of nucleotide monophosphates resulting in the corresponding nucleotide diphosphates by using ATP as a phosphate donor. T4 dNMP kinase is the only member of the family of NMP kinases that recognizes three structurally dissimilar nucleotides: dGMP, dTMP and 5-hydroxymethyl-dCMP (5-OH-dCMP).

Activity Assay: 1 μg T4 dNMP kinase, 2 mM ATP and 2 mM dGMP in 100 mM Tris-HCl pH 7.5, 10 mM MgCl_2.

Activity:

One unit of enzyme catalyses the phosphorylation of 1 nmol dGMP to dGDP per minute at 37°C. Specific activity: 100 $u/\mu g$.

Selected References:

Teplyakov *et al.* (1996) Crystal structure of bacteriophage T4 deoxynucleotide kinase with its substrates dGMP and ATP. *EMBO J.* **15**:3487.

Brush *et al.* (1993) Chemical modification of bacteriophage T4 deoxynucleotide kinase. Evidence of a single catalytic region. *J. Biol. Chem.* **268**:1603.

Brush *et al.* (1990) Bacteriophage T4 deoxynucleotide kinase: gene cloning and enzyme purification. *J. Bacteriol.* **172**:2935.

