

# T4 Polynucleotide Kinase, 3'-phosphatase free

From phage T4 am N81 pse T1 infected *Escherichia coli* BB  
ATP: 5'-dephosphopolynucleotide 5'-phosphotransferase

**Cat. No. 10 709 557 001** 200 units

**Cat. No. 10 838 292 001** 1,000 units

 **Version 19**

Content version: February 2020

Store at -15 to -25°C

## Product overview

<b>Storage buffer</b>	50 mM Tris-HCl, 1 mM dithiothreitol, 0.1 mM EDTA, 1 μM ATP, 50% glycerol (v/v), pH approx. 7.5 (+25°C).
<b>Volume activity</b>	10 × 10 <sup>3</sup> U/ml. T4 Polynucleotide kinase (PNK), 3'-phosphatase free, is assayed acc. to (1). One unit is the enzyme activity which catalyzes the incorporation of 1 nmol [ <sup>32</sup> P] into acid-precipitable products within 30 min at +37°C.
<b>Specific activity</b>	≥ 40 × 10 <sup>3</sup> U/mg acc. to (1, 2).
<b>Source</b>	T4 PNK, 3'-phosphatase free, is prepared from phage T4 am N81 pse T1-infected <i>E. coli</i> BB cells.
<b>Purity</b>	≥ 2 μg of PNK, 3'-phosphatase free, migrate as a single band in SDS polyacrylamide gel electrophoresis according to (3).
<b>Supplied phosphorylation buffer</b>	Phosphorylation buffer for direct phosphorylation, 10 × conc.: 500 mM Tris-HCl, 100 mM MgCl <sub>2</sub> , 1 mM EDTA, 50 mM dithiothreitol, 1 mM spermidine, pH 8.2 (at +25°C).
<b>Properties</b>	This phage derivative lacks 3'-phosphatase activity. The enzyme catalyzes the transfer of the terminal phosphate group of ATP to the 5'-hydroxylated terminus of DNA or RNA (1). It also exchanges 5' terminal phosphate groups (4).
<b>Application</b>	T4 PNK, 3'-phosphatase free, is of interest for phosphorylating both 5'- and 3'-termini of RNA. The added 3'-phosphate prevents cyclization or self addition of the RNA. Another use of T4 PNK, 3-phosphatase free, is the 5' [ <sup>32</sup> P]-terminal labeling of 3'-CMP to give 5' [ <sup>32</sup> P]pCp. This substrate is commonly used for 3' end labeling of RNA with T4 RNA ligase (5). <b>Note:</b> The 3'-phosphatase activity of the wild-type kinase is not observed under optimal incubation conditions with the mutant T4 PNK.
<b>Stability</b>	The undiluted enzyme is stable at -15 to -25°C until the expiration date printed on the label.

## Protocols and required material

<b>Calculation of pmol ends</b>	Use the length and the concentration of your input DNA to calculate pmol ends. The conversion formula is described in our "Lab FAQs" described in Lab FAQs ( <a href="http://www.roche.com/labfaqs">http://www.roche.com/labfaqs</a> ) under commonly used formulas.
<b>Additional reagents required</b>	<ul style="list-style-type: none"> <li>20 pmol [<sup>32</sup>P] ATP, ≥10 μM (aqueous solution)</li> <li>dephosphorylated DNA-fragment solution (20 pmol 5'-OH-termini).</li> </ul>

## Protocol

Standard assay for 5'-end-labeling of free 5'-OH-termini.

Step	Action		
1	Pipet the following components to a reaction vial on ice:		
		<b>Reagent</b>	<b>Volume</b>
		Dephosphorylated DNA-fragment (20 pmol 5'-OH-termini)	variable
		20 pmol [ <sup>32</sup> P] ATP, (aqueous solution), ≥10 μM	variable
	Phosphorylation buffer, 10 × conc.	2 μl	
	<ul style="list-style-type: none"> <li>Add double dist. water to a final volume of 19 μl.</li> </ul>		
2	<ul style="list-style-type: none"> <li>Add 10 U T4 PNK 3'-phosphatase free.</li> <li>Mix well.</li> </ul>		
3	Incubate 30 min at +37°C.		
4	Stop the reaction by cooling in an ice bath.		

**Determination of labeling efficiency** The yield of the labeling reaction can be determined by trichloroacetic acid precipitation.

## Quality control

For lot-specific certificates of analysis, see section, **Contact and Support**.

---

**Changes to previous version** Update of the chapter Quality Control.

---

## References

---

- 1 Richardson, C. C. (1968) *Proc. Natl. Acad. Sci. USA* **54**, 158.
- 2 Bradford, M. (1976) *Anal. Biochem.* **72**, 248.
- 3 Laemmli, U. K. (1970) *Nature* **227**, 680.
- 4 Chaconas, G. & v. d. Sande, J. H. (1980). *Methods Enzymol.* **65**, 680.
- 5 Otsuka, E. *et al.* (1981), *Proc. Natl. Acad. Sci. USA* **78**, 5439.
- 6 Cameron, V. & Uhlenbeck, O. C. (1977) *Biochemistry* **16**, 5120.
- 7 Cameron, V. *et al.* (1978) *Nucleic Acids Res.* **5**, 825.
- 8 Soltis, D. A. & Uhlenbeck, O. C. (1982) *J. Biol. Chem.* **257**, 11332–11339.
- 9 Soltis, D. A. & Uhlenbeck, O. C. (1982) *J. Biol. Chem.* **257**, 11340–11345.
- 10 Vance J.R. & Wilson T.E. (2001) Uncoupling of 3'-Phosphatase and 5'-Kinase Functions in Budding Yeast. *J. Biol. Chem.* **276**, 15073-15081.
- 11 Vance J.R. & Wilson T.E. (2001) Repair of DNA Strand Breaks by the Overlapping Functions of Lesion-Specific and Non-Lesion-Specific DNA 3' Phosphatases. *Molecular and Cellular Biol.* **21**, 7191-7198.

---

**Regulatory Disclaimer** For life science research only. Not for use in diagnostic procedures.

---

**Trademarks** All third party product names and trademarks are the property of their respective owners.

---

**Disclaimer of License** For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

---

---

## Contact and Support

To ask questions, solve problems, suggest enhancements and report new applications, please visit our **Online Technical Support Site**.

To call, write, fax, or email us, visit [sigma-aldrich.com](http://sigma-aldrich.com), and select your home country. Country-specific contact information will be displayed.

---



Roche Diagnostics GmbH  
Sandhofer Strasse 116  
68305 Mannheim  
Germany