

# Polynucleotide Kinase

From *Escherichia coli* E. D. pFDX

ATP: 5'-dephosphopolynucleotide 5'-phosphotransferase, EC 2.7.1.78

**Cat. No. 10 174 645 001** 200 U (10 U/  $\mu$ l)

**Cat. No. 10 633 542 001** 1,000 U (10 U/  $\mu$ l)

 **Version 19**

Content version: February 2020

Store at  $-15$  to  $-25^{\circ}\text{C}$

## Product description

**Phosphorylation buffer 10x conc.** 10x Incubation buffer for direct phosphorylation, supplied with the enzyme:

500 mM Tris-HCl, 100 mM MgCl<sub>2</sub>, 1 mM EDTA, 50 mM dithioerythritol, 1 mM spermidine, pH 8.2 (at  $+25^{\circ}\text{C}$ ).

**Storage buffer** 50 mM Tris-HCl, 1 mM dithioerythritol, 0.1 mM EDTA, 0.1  $\mu$ M ATP, 50% glycerol (v/v), pH approx. 7.5 ( $+25^{\circ}\text{C}$ ).

**Specific activity**  $> 40 \times 10^3$  U/mg acc. to (1, 2).

**Unit definition** Polynucleotide Kinase (PNK) 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  $+37^{\circ}\text{C}$ .

**Properties** Polynucleotide Kinase catalyzes the transfer of the terminal phosphate group of ATP to the 5'-hydroxylated terminus of DNA or RNA (1). The enzyme also catalyzes the exchange of 5'-terminal phosphate groups (4).

**Application** Labeling of 5'-hydroxyl ends of DNA or RNA with [<sup>32</sup>P]-ATP by direct phosphorylation or by an exchange reaction.  
Alternatively, polynucleotide kinase can be used for removal of 3'-phosphates from RNA or DNA at low pH conditions without altering the structure of the 5'-terminus.

**Storage/ stability** The undiluted enzyme is stable at  $-15$  to  $-25^{\circ}\text{C}$  through the expiration date printed on the label.

## Standard assay

**Protocol** Procedure for 5'-end-labeling of free 5'-OH-termini.

Step	Action										
1	<ul style="list-style-type: none"><li>Add to the dephosphorylated DNA-fragment solution (20 pmol 5'-OH-termini) on ice:<table border="1"><thead><tr><th>Reagent</th><th>Vol.</th></tr></thead><tbody><tr><td>Phosphorylation buffer, 10x conc.</td><td>2 <math>\mu</math>l</td></tr><tr><td>20 pmol [<sup>32</sup>P] ATP, 10 <math>\mu</math>mol (aqueous solution)</td><td>x <math>\mu</math>l</td></tr><tr><td>Add double-distilled water to a final volume of 19 <math>\mu</math>l.</td><td>x <math>\mu</math>l</td></tr><tr><td>10 U Polynucleotide Kinase</td><td>1 <math>\mu</math>l</td></tr></tbody></table></li><li>Mix and centrifuge briefly.</li><li>Incubate for 30 min at <math>+37^{\circ}\text{C}</math>.</li></ul>	Reagent	Vol.	Phosphorylation buffer, 10x conc.	2 $\mu$ l	20 pmol [ <sup>32</sup> P] ATP, 10 $\mu$ mol (aqueous solution)	x $\mu$ l	Add double-distilled water to a final volume of 19 $\mu$ l.	x $\mu$ l	10 U Polynucleotide Kinase	1 $\mu$ l
Reagent	Vol.										
Phosphorylation buffer, 10x conc.	2 $\mu$ l										
20 pmol [ <sup>32</sup> P] ATP, 10 $\mu$ mol (aqueous solution)	x $\mu$ l										
Add double-distilled water to a final volume of 19 $\mu$ l.	x $\mu$ l										
10 U Polynucleotide Kinase	1 $\mu$ l										
2	Stop the reaction by cooling in an ice bath.										

**Labeling yield** 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**.

## References

- Richardson, C. C. (1968) Proc. Natl. Acad. Sci. USA **54**, 158.
- Bradford, M. (1976) Anal. Biochem. **72**, 248.
- Laemmli, U. K. (1970) Nature **227**, 680.
- Chaconas, G. & v. d. Sande, J. H. (1980), Methods Enzymol. **65**, 680.

## Regulatory Disclaimer

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

## Changes to previous version

Update of the chapter Quality Control.

## 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.