

Thermo Scientific Phosphorylation of DNA with T4 Polynucleotide Kinase

This protocol is for the Thermo Scientific™ Phosphorylation™ of DNA with T4 Polynucleotide Kinase.

1. Prepare the following reaction mixture:

Linear ds DNA or Oligonucleotide	1-20 pmol of 5'-termini or 10-50 pmol
10x reaction buffer A for T4 Polynucleotide Kinase	2 µL
ATP, 10 mM*	2 µL
T4 Polynucleotide Kinase	1 µL (10 U)
Water, nuclease-free	to 20 µL
Total volume	20 µL

2. *Prepare 10 mM ATP solution by combining 10 µL of 100 mM ATP solution and 90 µL of Water, nuclease-free.
3. Mix thoroughly, spin briefly and incubate at 37 °C for 20 minutes.
4. Heat at 75 °C for 10 minutes.

Note

Visit [REviewer™](#) for molar calculations.

thermoscientific.com/onebio

© 2013 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Europe

Customer Service
cs.molbio.eu@thermofisher.com

Technical Support
ts.molbio.eu@thermofisher.com

Tel 00800 222 00 888
Fax 00800 222 00 889

United States

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 235 9880
Fax 800 292 6088

Canada

Customer Service
cs.molbio@thermofisher.com

Technical Support
ts.molbio@thermofisher.com

Tel 800 340 9026
Fax 800 472 8322

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific