

Non-radioactive Random-primed DNA Labeling with Klenow Fragment, exo-

This protocol is for Non-radioactive Random-primed DNA Labeling with Klenow Fragment, exo-.

1. Prepare the following reaction mixture:

DNA template	10 μ L (100 ng – 1 μ g)
10x reaction buffer for Klenow Fragment, exo-	5 μ L
6.0 A₂₆₀ units/mL (100 μM) Random Hexamer Primer	12.5 μ L
Water, nuclease-free	to 39 μ L
Total volume	39 μL

2. Incubate the mixture in a boiling water bath for 5-10 minutes and then chill on ice.
3. Add:

3 dNTP Mix, 1 mM each (without the dTTP)	5 μ L (0.1 mM final concentration)
dTTP	3.25 μ L (0.065 mM final concentration)
Biotin-11-dUTP*, 1 mM	1.75 μ L
Thermo Scientific Klenow Fragment, exo- (Cat #EP0421, #EP0422)	1 μ L (5 U)
Water, nuclease-free	to 20 μ L
Total volume	50 μL

* Fluorescein-12-dUTP, DIG-dUTP or Aminoallyl-dUTP can also be used with the same protocol.

4. Incubate the reaction mixture at 37 °C for 1 hour. Add 1 μ L 0.5 M EDTA, pH 8.0 to stop the reaction.
5. Remove 1 μ L of the reaction mixture and determine the percentage of label incorporated.
6. Optionally, purify by using Sephadex G-50 or Bio-Gel P-60 or size-cutoff appropriate Thermo Scientific GeneJET nucleic acid purification kit.

thermoscientific.com/onebio

© 2012 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