

DNA 3'-end Labeling by Fill-in of 5'-overhangs with Klenow Fragment* or Bsm DNA Polymerase, Large Fragment

1. Prepare the following reaction mixture:

Linear DNA (aqueous solution)	0.1-4 µg
10X reaction buffer for Klenow Fragment or 10X Bsm buffer	2 µl
[alpha-32P]-dNTP, ~15-30 TBq/mmol (400-800 Ci/mmol) or [alpha-32P]-dNTP, ~110 TBq/mmol (3000 Ci/mmol)	0.74 MBq (20 µCi) 2.96 MBq (80 µCi)
3 dNTP Mix, 2 mM each (without a labeled dNTP)	2.5 µl (0.25 mM final concentration)
Klenow Fragment or Klenow Fragment, exo- or Bsm DNA Polymerase, Large Fragment	0.1 µl (1 u) 0.2 µl (1 u) 0.125 µl (1 u)
Water, nuclease-free	to 20 µl
Total volume	20 µl

2. Incubate at 37°C for 15 min.
3. Stop the reaction by heating at 75°C for 10 min.

Note

* This protocol is suitable labeling of the following Fermentas DNA markers, composed of DNA fragments with 5'-overhangs:

Lambda DNA EcoRI Marker, #SM028

Lambda DNA HindIII Marker, #SM0101

Lambda DNA EcoRI/HindIII Marker, #SM0191

Lambda DNA Eco91I Marker, #SM0111

phiX174 DNA HinfI Marker, #SM0261

- The modified version of this protocol can be used for non-radioactive labeling of DNA markers. Substitute a part of dTTP nucleotide with a modified nucleotide (e.g. Biotin-11-dUTP or Fluorescein-12-dUTP) at a molar ratio of 1:2.
- For estimation of pmol of DNA ends, see [REviewer™](#).

thermoscientific.com

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**North America
Technical Services:**
ts.molbio@thermofisher.com

Customer Services:
cs.molbio@thermofisher.com
Tel 800 235 9880
Fax 800 292 6088

**Europe and Asia
Technical Services:**
ts.molbio.eu@thermofisher.com

Customer Services:
cs.molbio.eu@thermofisher.com

Thermo
SCIENTIFIC