

# PRODUCT INFORMATION **Thermo Scientific** Biotin DecaLabel DNA Labeling Kit

for \_\_\_\_\_reactions #\_\_\_ Lot: Expiry Date: \_\_\_ Store at -20°C

#### www.thermoscientific.com/onebio

#### COMPONENTS OF THE KIT

Component	<b>#K0651</b> 10 rxns	# <b>K0652</b> 30 rxns
Klenow Fragment, exo⁻, 5 u/µl	15 µl	30 µl
Decanucleotide in 5X Reaction Buffer	100 µl	300 µl
Biotin Labeling Mix 1 mM dGTP, 1 mM dATP, 1 mM dCTP, 0.65 mM dTTP, 0.35 mM Biotin-11-dUTP aqueous solution	50 µl	150 µl
Control Template λ DNA/HindIII fragments,10 ng/µI	50 µl	125 µl
<b>Biotin-labeled DNA</b> 50 μl of biotin-labeled DNA (λ DNA/HindIII), 5 ng/μl	50 µl	125 µl
Water, nuclease-free	1.25 ml	2x1.25 ml

# CERTIFICATE OF ANALYSIS

All components of the kit were tested in a control labeling reaction. Labeled DNA probe was used for a spot hybridization. 0.3-0.1 pg of homologous DNA was detected after 16 hours color development with a streptavidin conjugated to alkaline phosphatase which catalyzes a color reaction with NBT/BCIP.



### DESCRIPTION

Thermo Scientific Biotin DecaLabel DNA Labeling Kit is an advanced system for the efficient synthesis of biotin-labeled DNA probes, based on an improved random-primed labeling method originally developed by Feinberg and Vogelstein (1, 2). The primary improvement over the traditional randomprimed method involves the use of random decamers instead of hexamers to ensure more efficient annealing with DNA at 37°C. Klenow Fragment, exo- is also included in the kit; this genetically engineered enzyme has no detectable exonuclease activity. Therefore, the enzyme does not degrade the labeled probe during reaction, which results in a high labeling yield even with low amounts of template. As a result, DNA fragments of any length can be uniformly labeled. Biotin-labeled DNA is detected with the Biotin Chromogenic Detection Kit (#K0661) or with conventional biotinavidin or biotin-streptavidin detection systems.

### DNA LABELING with BIOTIN-11-dUTP

1. Add the following components into 1.5 ml microcentrifuge tube:

DNA template (100 ng - 1 µg)	10 µl
Decanucleotide in 5X Reaction Buffer	10 µl
Water, nuclease-free	to 44 µl

Vortex the tube and spin down in a microcentrifuge for 3-5 s.

Incubate the tube in a boiling water bath for 5-10 min and cool it on ice. Spin down quickly.

2. Add the following components in the same tube:

Biotin Labeling Mix	5 µl
Klenow fragment, exo- (5 u)	1 µl

Shake the tube and spin down in a microcentrifuge for 3-5 s. Incubate for 1 hour at 37°C. Incubation at 37°C for up to 20 hours increases the yield of labeled DNA.

- 3. Stop the reaction by the addition of 1  $\mu$ l of 0.5 M EDTA, pH 8.0.
- 4. The labeled DNA is used directly for hybridization or stored at -20°C. Removal of the unincorporated label is not necessary for most applications. If required, the unincorporated dNTP can be removed by chromatography on Sephadex<sup>®</sup> G-50 or by selective precipitation of DNA with ethanol in the presence of ammonium acetate (3).

## CONTROL LABELING REACTION

1. Add the following components into 1.5 ml microcentrifuge tube:

Control Template, 10 ng/µl	25 µl
Decanucleotide in 5X Reaction Buffer	10 µl
Water, nuclease-free	9 µl

Vortex the tube and spin down in a microcentrifuge for 3-5 s.

Incubate the tube in a boiling water bath for 5-10 min and cool it on ice. Spin down quickly.

2. Add the following components in the same tube:

Biotin Labeling Mix	5 µl
Klenow fragment, exo- (5 u)	1 µl

Shake the tube and spin down in a microcentrifuge for 3-5 s. Incubate for 1 hour at 37°C. Incubation at 37°C for up to 20 hours increases the yield of labeled DNA.

3. Stop the reaction by the addition of 1  $\mu$ l of 0.5 M EDTA, pH 8.0.

## References

- 1. Feinberg, A.P., Vogelstein, B., Biochem. 132, 6-13, 1983.
- 2. Feinberg, A.P., Vogelstein, B., Biochem. 137, 266-267, 1984.
- 3. Sambrook, J., Fritsch, E.F. and Maniatis, T., Molecular Cloning: A Laboratory Manual; Second Edition, Cold Spring Horbor laboratory, Cold Spring Harbor, N. Y., 1989.

#### PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively for research purposes and *in vitro* use only. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to <u>www.thermoscientific.com/onebio</u> for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Saphadex is a registered trademark of GE Healthcare Ltd. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.