INSTRUCTIONS

North2SouthTM Biotin Random Prime Labeling Kit

| 17075 | | 0810.4 |
|--------|--|--------|
| Number | Description | |
| 17075 | North2South Biotin Random Prime Labeling Kit, sufficient reagents for 10 biotin-labeling reactions, each with 100ng DNA. | |
| | Kit Contents: | |
| | Heptanucleotide mix (5X), 100µL | |
| | dNTP mix (5X), 1mM dATP, 1mM dTTP, 1mM dGTP, 1mM dCTP, 80µL | |
| | Reaction Buffer (10X), 50µL | |
| | DNA Polymerase, Klenow fragment (3'-5' exo-), 5 units/µL, 10µL | |
| | Control DNA, 250ng/µL, 5µL | |
| | EDTA, 500mM, pH 8.0, 1mL | |
| | Nuclease-free water, 1mL | |
| | Ammonium acetate (NH ₄ OAc), 5M, 1mL | |
| | Biotin-11-dUTP, 20µL | |
| | Storage: Upon receipt store all components at -20°C in a non-defrosting freezer. Product is sh | ipped |

Introduction

The Thermo ScientificTM North2SouthTM Biotin Random Prime Labeling Kit provides the components and procedure to successfully prepare biotin-labeled DNA probes for use in Southern blotting and other DNA hybridization methods involving streptavidin detection systems. The random prime labeling method is based on the procedure of Feinberg and Vogelstein,^{1,2} wherein random heptanucleotides containing all possible sequences anneal to a denatured DNA template and act as primers for complementary strand synthesis by DNA Polymerase (Klenow fragment, 3'-5' exo⁻). Inclusion of biotinylated nucleotides in the reaction mixture ensures that the newly synthesized DNA strands are labeled with biotin. This protocol yields biotinlabeled DNA of high activity for use as probes in experiments using the North2South Chemiluminescent Hybridization and Detection Kit (Product No. 17097).

Additional Materials Required

- Heating block
- 1X TE Buffer (10mM Tris•HCl, 1mM EDTA, pH 8.0)

with dry ice.

- Ethanol
- Linearized template DNA, i.e., the DNA to be labeled





Procedure for Biotin Random Prime Labeling

A. Prepare Complete 5X dNTP mix

Before first use of the kit, prepare a complete 5X dNTP mix by adding entire contents of the Biotin-11-dUTP (20μ L) to the tube of 5X dNTP mix (80μ L). This 100 μ L complete stock solution will henceforth be referred to simply as '5X dNTP mix'. After first use, store the prepared 5X dNTP mix at -20°C with the other kit components. In order to estimate your experimental probe labeling efficiency the control DNA will need to be labeled. Once labeled, this probe can be stored at -20°C for use in future probe labeling assessments.

B. Perform Labeling Reaction

- 1. In a microcentrifuge tube, dilute ~100ng of linear DNA for labeling (template DNA) to a final volume of 24μ L in nuclease-free water. For the positive control, dilute 0.5 μ L of control DNA (125ng) to 24μ L.
- 2. Add 10µL of heptanucleotide mix and denature the DNA template by boiling the tube for 5 minutes.
- 3. Quickly freeze the denatured DNA solution by placing the tube in a dry ice/ethanol bath for 5 minutes.
- 4. Thaw DNA solution at 4°C and briefly centrifuge to collect liquid in the bottom of the tube. Place the tube on ice.
- 5. In the order stated, add the following components to the prepared sample solution on ice:
 - 10µL of 5X dNTP mix
 - 5µL of 10X Reaction Buffer
 - 1µL Klenow fragment
 - (Final volume = 50μ L)
- 6. Mix contents by flicking or briefly vortexing the tube. Centrifuge briefly to collect liquid in the bottom of the tube.
- 7. Incubate labeling reaction for 60 minutes at 37°C.
- 8. Inactivate the enzyme by adding 2μ L of 500mM EDTA, pH 8.0.

C. Ethanol Precipitation to Remove Unincorporated Nucleotides

Note: As an alternative to the ethanol precipitation method described here, size exclusion chromatography with an appropriate gel filtration column may be used to remove unincorporated nucleotides.

- 1. Adjust the reaction to contain 0.5M ammonium acetate. For example, add 5μ L of 5M NH₄OAc to a 50μ L labeling reaction and mix well.
- 2. Add 2 volumes of 100% ethanol and mix well. For example, add 110μ L to the 55 μ L volume from Step 1.
- 3. Chill tube at -20°C for 15 minutes.
- 4. Centrifuge tube at 4°C for 30-60 minutes at maximum speed (> $10,000 \times g$).
- 5. Carefully aspirate to remove (and discard) the supernatant.
- 6. Wash the DNA pellet once by adding ice-cold 70% ethanol and centrifuging for 30-60 minutes at maximum speed. Carefully remove and discard the supernatant.
- 7. Dissolve the DNA pellet in 100μL of 1X TE or nuclease-free water and store at -20°C. For long-term storage, prepare single-use aliquots and store at -70°C.

Procedures for Estimating Labeling Efficiency

Labeling efficiency can be determined by dot blotting using a hand spotting method. The following dot blot protocol was developed for use with the detection reagents in our **Chemiluminescent Nucleic Acid Detection Module (Product No. 89880).**

Additional Materials Required

- Positively-charged nylon membrane (e.g., Biodyne™ B Membrane, Product No. 77016)
- TE Buffer: 10mM Tris•HCl, 1mM EDTA, pH 8.0
- Microcentrifuge tubes and microcentrifuge



- UV lamp or crosslinking device equipped with 254nm bulbs or 312nm transilluminator
- Chemiluminescent Nucleic Acid Detection Module (Product No. 89880)
- X-ray film or CCD camera

D. Dot Blot by Hand Spotting

- 1. Hydrate/equilibrate a positively-charged nylon membrane in TE Buffer for at least 10 minutes.
- 2. Pipette 20µL TE Buffer into a micro centrifuge tube labeled #1 and 10µL into micro centrifuge tubes labeled #2-7 for the control reaction and each experimental sample.
- 3. Prepare dilutions of the labeled DNA from both the control DNA and your experimental probe with TE. Dilute the labeled samples 10-fold in TE Buffer by adding 2μ L of the stated samples into the #1 tubes and serially diluting the sample serially into the other tubes using 10μ L per dilution.
- 4. Place the equilibrated membrane onto a clean, dry paper towel. Allow excess buffer to absorb into the membrane, but do not allow the membrane dry out.
- 5. Spot 2µL of samples and standards onto the hydrated membrane. Allow the samples to absorb into the membrane.

E. UV Crosslink the Dot Blot

- 1. Immediately UV crosslink the membrane by one of the following three methods:
- **Option 1**: Crosslink at 120mJ/cm² using a commercial UV-light crosslinking instrument equipped with 254nm bulbs (45-60 second exposure using the auto crosslink function).
- **Option 2**: Crosslink at a distance of approximately 0.5cm from the membrane for 5-10 minutes with a hand-held UV lamp equipped with 254nm bulbs.
- **Option 3**: Crosslink for 10-15 minutes with the membrane face down on a transilluminator equipped with 312nm bulbs.
- 2. Continue with the detection and analysis immediately or store the membrane dry at room temperature until the detection protocol can be performed.

F. Detection and Analysis

Detect the spotted standards and samples using the procedure and reagents for the North2South Chemiluminescent Hybridization and Detection Kit (Product No. 17097) or the Chemiluminescent Nucleic Acid Detection Module (Product No. 89880). To determine the labeling efficiency, compare spot intensities of the sample lanes to those of the labeled control DNA.

- If using X-ray film to document results, compare the row of sample spots to the row of labeled control DNA spots to estimate relative efficiency. Densitometry of the film can also be performed.
- If using a cooled CCD camera to document results, spot densitometry can be performed (choose a dilution that is within the linear exposure range) assess the labeling efficiency of the unknown samples.

Compare the spot intensities of experimental labeled probes to the labeled Control DNA. The intensity of your experimental probe should be approximately 50% of the labeled control DNA to yield satisfactory hybridization results. If the probe signal is weaker than this it will be difficult to detect signals from low abundance species. The amount of probe required for downstream applications must be empirically determined.

Troubleshooting

A poor quality starting DNA template is the most likely reason for poor probe synthesis labeling when compared to that of the purified control DNA supplied in the kit. In some cases, using less than the recommended starting DNA template is better because it reduces the concentration of inhibitors in the reaction. Alternatively, the DNA template may require further purification before use in the labeling reaction.



Related Thermo Scientific Products

| North2South Chemiluminescent Hybridization and Detection Kit |
|--|
| Chemiluminescent Nucleic Acid Detection Module (this module includes the same components as Product No. 17097, except the hybridization and stringency wash buffers.) |
| LightShift TM Chemiluminescent EMSA Kit |
| North2South Chemiluminescent Substrate for HRP, 100mL |
| Biodyne B Nylon Membranes, 8 × 12cm, 25 sheets, positively charged membrane with 0.45µm pore size |
| |

Cited References

- 1. Feinberg, A.P. and Vogelstein, B. (1983). Anal Biochem 132:6.
- 2. Feinberg, A.P. and Vogelstein, B. (1983). Anal Biochem 137:266.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at <u>www.thermoscientific.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2013 Thermo Fisher Scientific Inc. All rights reserved. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.