INSTRUCTIONS



North2South[®] Chemiluminescent Hybridization and Detection Kit

| 17097 | 1758.1 |
|--------|---|
| Number | Description |
| 17097 | North2South[®] Chemiluminescent Hybridization and Detection Kit, contains sufficient reagents for 1,000 cm ² of membrane |
| | Kit Contents: |
| | Stabilized Streptavidin-Horseradish Peroxidase Conjugate, 1.5 ml |
| | Chemiluminescent Substrate, stable for six months at room temperature or one year at 4°C |
| | Luminol/Enhancer Solution, 80 ml |
| | Stable Peroxide Solution, 80 ml |
| | Blocking Buffer, 500 ml |
| | 4X Wash Buffer, 500 ml |
| | Substrate Equilibration Buffer, 500 ml, store at room temperature or 4°C |
| | North2South [®] Hybridization Buffer, 125 ml |
| | North2South [®] Hybridization Stringency Wash Buffer (2X), 375 ml |

Storage: Upon receipt store kit at 4°C. Kit is shipped at ambient temperature.

Introduction

The North2South[®] Chemiluminescent Hybridization and Detection Kit is a complete system for chemiluminescent detection of nucleic acids in Northern and Southern blot applications. This system combines an enhanced luminol substrate for horseradish peroxidase (HRP) with optimized hybridization and blocking conditions that ensure consistent results with sensitivity equal to or exceeding ³²P. This robust system greatly reduces processing and film exposure time. Post-hybridization processing has been reduced from the standard 2.5 hours to 1 hour. Film exposure times range from 0.5 to 10 minutes with the substrate emitting light with relatively constant intensity for six hours, thus allowing for multiple exposures.

Additional Materials Required

- Biotinylated nucleic acid probe (can be produced using Product No. 17075)
- Positively charged nylon transfer membrane (Product No. 77016)
- X-ray film (Product No. 34090 or 34091) or a CCD imaging instrument

Procedure for Nucleic Acid Hybridization and Detection

A. Pre-Hybridization and Hybridization

Notes:

- Equilibrate kit buffers to room temperature before use. If there is a precipitate in any of the kit buffers, heat buffer in a 37°C water bath until precipitate disappears.
- Heat incubator to appropriate temperature (e.g., 55°C for DNA hybrids or 65°C for RNA:RNA hybrids).
- Quantify probe using a spectrophotometer.^{1, 2}



1. Equilibrate the North2South[®] Hybridization Buffer to room temperature (RT).

Note: This buffer has been specifically optimized for use with the North2South® Kit

- 2. Place blot in a container such as a 50 ml centrifuge tube and add sufficient Hybridization Buffer to completely cover the blot. Use at least 0.1 ml per cm² of membrane.
- 3. Seal the container and pre-hybridize the membrane with shaking or rotating for at least 30 minutes. For DNA hybrids incubate at 55°C; for RNA:RNA hybrids incubate at 65°C.
- 4. While pre-hybridizing, denature the biotinylated DNA probe. Heat DNA probe at 100°C for 10 minutes and place on ice for 5 minutes.
- 5. After pre-hybridization, add the denatured biotinylated probe. For biotinylated RNA probes, add 3-5 ng of probe per milliliter of hybridization buffer; for biotinylated DNA probes, add ~30 ng of probe per milliliter of hybridization buffer.
- 6. Incubate overnight with shaking or rotating at 55°C for DNA hybrids or 65°C for RNA:RNA hybrids.

B. Stringency Washes

1. On the next day, equilibrate the North2South[®] Hybridization Stringency Wash Buffer (2X) to RT. Once the Wash Buffer is fully in solution, add an equal volume of sterile ultrapure water. The resulting 1X buffer contains 2X SSC/0.1% SDS.

Note: If required use other wash conditions to increase or decrease stringency. In general, stringency increases with increasing temperature and decreasing ionic strength.

2. Wash blot three times for 15-20 minutes per wash with agitation. Use 0.2 ml of 1X Stringency Wash Buffer per cm² of membrane and perform washes at 55°C for DNA hybrids or 65°C for RNA:RNA hybrids.

C. Probe Detection

Note: Use clean forceps to handle only the corners of the membrane. For optimal results, rinse forceps with ethanol and allow them to dry between steps.

- 1. Decant the Stringency Wash Buffer and add sufficient Blocking Buffer to generously cover the membrane. Use at least 0.25 ml/cm² of membrane. Incubate with shaking or rotating for 15 minutes at RT.
- 2. Determine the amount of Streptavidin-HRP to add to obtain a 1:300 final dilution in the tube with the blot but do not add it yet. Decant a portion of the Blocking Buffer solution from the tube containing the membrane into a separate tube.
- 3. Add the Streptavidin-HRP conjugate from step 2 to the separated buffer. Add the buffer/Streptavidin-HRP mix to the tube containing the membrane and incubate for 15 minutes at RT with agitation.

Note: Performing steps 2 and 3 as described prevents the undiluted conjugate from coming in direct contact with the membrane, which would produce undesirable results.

- 4. Dilute Wash Buffer (4X) to 1X with sterile ultrapure water. Wash the membrane four times for 5 minutes each with 1X Wash Buffer at RT with agitation. Place membrane into a clean wash container for the next step.
- 5. Add Substrate Equilibration Buffer to container with the blot. Use 0.25 ml/cm² of membrane. Incubate blot for 5 minutes at RT with agitation.

D. Substrate Development

1. Prepare the Substrate Working Solution by mixing equal volumes of the Luminol/Enhancer Solution and Stable Peroxide Solution. Prepare enough solution to completely cover the membrane (i.e., ~0.1 ml/cm²).

Note: The Working Solution is stable for 6 hours at room temperature. Exposure to the sun or any other intense light can harm the Working Solution. For best results keep the Working Solution in an amber bottle and avoid prolonged exposure to any intense light. Short-term exposure to typical laboratory lighting will not harm the Working Solution.

- 2. Place the moist membrane on a tray or a piece of plastic wrap and cover with the Substrate Working Solution. Incubate for 5 minutes at RT. Make sure the membrane is fully covered with substrate.
- 3. Drain substrate from the membrane surface and transfer the moist membrane to a sheet protector or wrap in clear plastic wrap. Remove any trapped air bubbles or wrinkles within the plastic wrap and blot any substrate that may have leaked at the edges.



- 4. Expose blot to film for 1 minute. Adjust the exposure time as needed to obtain the desired signal.
- 5. Develop the film according to the manufacturer's instructions.

Note: For images that have been overexposed, or that show high background and/or speckling, use the Erase-It[®] Background Eliminator Kit (Product No. 21065).

Note: For best results, use a new membrane for each hybridization procedure. If stripping and reuse of the membrane is required, test the stripping protocol by incubating the stripped membrane in Substrate Working Solution, placing it in a sheet protector and exposing it to film for at least 30 minutes. If no bands are visible upon developing the film, rinse the membrane with Wash Buffer and re-probe.

Troubleshooting

| Problem | Possible Cause | Solution |
|-------------------------|---|--|
| High background | Probe is too concentrated | Reduce probe concentration during hybridization |
| | Too much HRP conjugate used | Dilute stabilized conjugate stock \geq 1:300 in Blocking Buffer |
| | Undiluted probe contacted the membrane | Thoroughly mix the probe in Hybridization Buffer before allowing it to contact the membrane |
| | Wrong blocking or hybridization buffer was used | Use the buffers supplied with the kit |
| | Wrong type of membrane was used | Use Biodyne [®] B Nylon Membrane (Product No. 77016) |
| Speckling background | Precipitate in the HRP conjugate | Filter conjugate through a 0.2 µm filter before use, or centrifuge 1 minute at maximum speed, or use Erase-It [®] Background Eliminator (Product No. 21065) |
| Clear (ghost) bands | Too much HRP | Dilute the HRP conjugate an additional 25- to 50-fold |
| Nonspecific bands | Degradation of target nucleic acid | Prevent nuclease contamination by using nuclease-free reagents or nuclease inhibitors |
| | Too much target nucleic acid present | Decrease the target concentration |
| | Too much probe was used | Reduce probe concentration during hybridization |
| | Stringency washes not sufficient | Increase stringency by decreasing the SSC concentration in the wash solution to 1X or 0.1X and/or by increasing temperature |
| Weak or no signal | Insufficient film exposure | Expose to film for 30 minutes or longer |
| | HRP conjugate concentration incorrect | Use a 1:300 dilution of a stabilized conjugate stock |
| | Low amount of target nucleic acid present | Increase probe concentration during hybridization and/or increase film exposure time |
| | DNA probe not completely denatured | Heat probe to 100°C (boil) for 10 minutes and chill on ice for 5 minutes |
| | Poor membrane transfer | Check transfer (blotting) protocol |
| | Inactive HRP conjugate | Test HRP by spotting 1 μ l of the HRP conjugate onto a small piece of membrane, wash and detect with the substrate |
| | Probe is degraded or not labeled | Test probe by spotting it onto membrane and detecting with Streptavidin-HRP as outlined in protocol |
| | Probe does not hybridize to target | Use Hybridization and other buffers supplied with the kit Reduce the hybridization temperature |



Related Thermo Scientific Products

| 37549North2South® Hybridization Buffer, 125 ml37555North2South® Hybridization Stringency Wash Buffer (2X), 375 ml17075North2South® Biotin Random Prime DNA Labeling Kit77016Biodyne® Membrane, 8 × 12 cm, 25/package34090CL-XPosure™ Film, 5 × 7 in, 100 sheets/package21065Erase-It® Background Eliminator Kit, for reducing background from exposed X-ray film | 89880 | Chemiluminescent Nucleic Acid Detection Module |
|--|-------|--|
| 17075North2South [®] Biotin Random Prime DNA Labeling Kit77016Biodyne [®] Membrane, 8 × 12 cm, 25/package34090CL-XPosure™ Film, 5 × 7 in, 100 sheets/package21065Erase-It [®] Background Eliminator Kit, for reducing background from exposed X-ray film | 37549 | North2South [®] Hybridization Buffer, 125 ml |
| 77016 Biodyne [®] Membrane, 8 × 12 cm, 25/package 34090 CL-XPosure™ Film, 5 × 7 in, 100 sheets/package 21065 Erase-It [®] Background Eliminator Kit, for reducing background from exposed X-ray film | 37555 | North2South [®] Hybridization Stringency Wash Buffer (2X), 375 ml |
| 34090CL-XPosure™ Film, 5 × 7 in, 100 sheets/package21065Erase-It® Background Eliminator Kit, for reducing background from exposed X-ray film | 17075 | North2South [®] Biotin Random Prime DNA Labeling Kit |
| 21065 Erase-It [®] Background Eliminator Kit, for reducing background from exposed X-ray film | 77016 | Biodyne[®] Membrane, 8 × 12 cm, 25/package |
| | 34090 | CL-XPosureTM Film, 5×7 in, 100 sheets/package |
| | 21065 | Erase-It [®] Background Eliminator Kit, for reducing background from exposed X-ray film |
| 17295 North2South [~] Chemiluminescent Substrate for HRP | 17295 | North2South [®] Chemiluminescent Substrate for HRP |

Additional Information

Please visit the web site for additional information on the product including the following items:

- Tech Tip protocol: Strip Northern and Southern blots
- Tech Tip resource: Northern/Southern blotting technical guide

Cited References

- 1. Sambrook, J., Fritsch, E.F. and Maniatis, T. (1989). Molecular Cloning: A Laboratory Manual, 2nd ed. Cold Spring Harbor Laboratory Press.
- 2. Ausubel, F.M., et al. (1995). Current Protocols in Molecular Biology. John Wiley & Sons, Inc.

Product References

- Adilakshmi, T. and Laine, R.O. (2002). Ribosomal protein S25 mRNA partners with MTF-1 and La to provide a p53-mediated mechanism for survival or death. J. Biol. Chem. 277(6):4147-51.
- Bach, S., *et al.* (2002). Gene expression profiling in AGS cells stimulated with *Heliobacter pylori* isogenic strains (*cagA* positive or *cagA* negative). *Infection and Immunity* **70**(2):988-92.
- Borst, E-M., *et al.* (1999). Cloning of the human cytomegalovirus (HCMV) genome as an infectious bacterial artificial chromosome in *Esherichia coli*: a new approach for construction of HCMV mutants. *J. Virology*. **73(10):**8320-9.
- Haertel-Wiesmann, M., *et al.* (2000). Regulation of cyclooxygenase-2 and periostin by Wnt-3 in mouse mammary epithelial cells. *J. Biol. Chem.* **275(41)**:32046-51.

Biodyne® is a registered trademark of Pall BioSupport Corporation.

This product ("Product") is 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") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. 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 the original purchaser of the Product ("Buyer").

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 replacement of or refund for the non-conforming Product(s).

There is no obligation to replace 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.

Current versions of product instructions are available at <u>www.thermo.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2010 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.