

# Strip Northern and Southern blots

TR0028.1

#### Introduction

Following exposure of a Northern or Southern blot (either by classical isotopic methods or Thermo Scientific North2South Reagents), researchers often wish to strip off the first probe and detecting reagents so that they may reprobe for the same or different target. This Tech Tip describes several commonly used conditions for stripping probe from a Northern or Southern blot while maintaining the nucleic acid target on the membrane (Table 1).

Be aware that any stripping procedure has the potential to denature sensitive nucleic acids, rendering them unrecognizable by the probe. Should this occur, reprobing will not be possible, and a new blot will have to be prepared for each probing experiment. Furthermore, any physical defects of a membrane will be exaggerated when stripped and reprobed.

Table 1. Possible stripping conditions for Northern and Southern blots.

Stripping Solution	Strength	DNA	RNA	Temperature & Time	
0.5% SDS	Mild	Yes	Yes	60°C for 60 minutes	
5 mM Tris•HCl, pH 8.0, 2 mM EDTA, 0.1X Denhardt's Reagent	Mild	Yes	Yes	65°C for 2 hours	
50% Formamide, 2X SSPE	Moderate	Yes	Yes	65°C for 1 hour	
0.4 M NaOH, 0.1% (w/v) SDS*	Moderate	Yes	No	45°C for 30 minutes, then room temperature for 2 x 10 minutes	
Boiling 0.1% SDS	Harsh	Yes	Yes	Pour on 100°C solution and let cool to room temperature; repeat 2 times	
0.1X SSPE or SSC/0.5% SDS	Harsh	Yes	Yes	Pour on 100°C solution and let cool to room temperature; repeat 2 times	

\*Blots must be neutralized in TE after stripping.

## **Tips for Stripping and Reprobing Blots**

- Wash blots with target side up
- Handle blots only at corners
- Keep blots wet between hybridization and stripping
- Keep blots wet after stripping
- Store blots in 0.1% SDS or 1X TE at 4°C
- Stripped blots may be dried and stored at room temperature for long-term storage. However, any probe remaining on the blot will become permanently bound once the blot is dried.

#### **Test for Stripping Efficiency**

For biotinylated probes detected using North2South<sup>®</sup> Chemiluminescent Hybridization and Detection Kit (Product No. 17097), test for stripping efficiency by reapplying streptavidin-HRP and detecting with the substrate as indicated in the protocol. If no signal is detected, then stripping was complete. If signal is detected, try a stronger stripping method.

### **Buffer Formulations**

50 X Denhardt's Reagent	<u>20 X SSC</u>	<u>20 X SSPE</u>	<u>TE, pH 7.4</u>
5 g Ficoll <sup>®</sup> Compound	800 ml H2O	800 ml H <sub>2</sub> O	10 mM Tris•HCI (pH 7.4)
5 g Polyvinylpyrrolidone	175.3 g NaCl	175.3 g NaCl	1 mM EDTA (pH 8.0)
5 g BSA	88.2 g sodium citrate	27.6 g NaH <sub>2</sub> PO <sub>4</sub> •H <sub>2</sub> O	
Adjust volume to 500 ml	Adjust pH to 7.0 with NaOH	Adjust pH to 7.4 with NaOH	
	Adjust volume to 1 L	Adjust volume to 1 L	



#### **Related Thermo Scientific Products**

77016 Biodyne<sup>®</sup> B Nylon Membranes, 25/pkg.
17097 North2South Chemiluminescent Hybridization and Detection Kit, complete set of hybridization buffers, streptavidin-HRP and substrate for use and detection of biotinylated probes
89880 Chemiluminescent Nucleic Acid Detection Module, streptavidin-HRP and substrate for detecting biotinylated probes
20148 LightShift<sup>®</sup> Chemiluminescent EMSA Kit, a complete kit for performing gel shift assays using biotin end-labeled DNA targets

Current versions of product instructions are available at www.thermo.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

Ficoll is a registered trademark of GE Healthcare.

© 2008 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.