

# **Trans-Blot<sup>®</sup> Electrophoretic Transfer Cell**

## **Instruction Manual**

**Catalog Numbers  
170-3910, 170-3925,  
170-3926, 170-3939,  
170-3945, 170-3946,  
170-3950, 170-3951  
170-3953 and 170-3954**



## **Note**

### **1. Assembly & Disassembly**

To insure best performance from the Trans-Blot electrophoretic transfer cell, become fully acquainted with these operating instructions before using the cell to transfer samples. Bio-Rad recommends that you first read these instructions carefully. Then assemble and disassemble the cell completely. After these preliminary steps, you should be ready to transfer a sample.

### **2. Wash Cell Before Use**

Bio-Rad also recommends that all Trans-Blot electrophoretic transfer cell components and accessories be cleaned with a suitable laboratory cleaner (such as Bio-Rad Cleaning Concentrate, catalog number 161-0722) and rinsed thoroughly with distilled water, before use.

**Model** \_\_\_\_\_

**Catalog Number** \_\_\_\_\_

**Date of Delivery** \_\_\_\_\_

**Warranty Period** \_\_\_\_\_

**Serial Number** \_\_\_\_\_

**Invoice Number** \_\_\_\_\_

**Purchase Order Number** \_\_\_\_\_

## **Warranty**

Bio-Rad Laboratories warrants the Trans-Blot electrophoretic transfer cell against defects in materials and workmanship for 1 year. If any defects occur in the instrument during this warranty period, Bio-Rad Laboratories will repair or replace the defective parts free. The following defects, however, are specifically excluded:

1. Defects caused by improper operation.
2. Repair or modification done by anyone other than Bio-Rad Laboratories or an authorized agent.
3. Use of fittings or other spare parts supplied by anyone other than Bio-Rad Laboratories.
4. Damage caused by accident or misuse.
5. Damage caused by disaster.
6. Corrosion due to use of improper solvent or sample.

For any inquiry or request for repair service, contact Bio-Rad Laboratories after confirming the model and serial number of your instrument.

# Table of Contents

<b>Section 1</b>	<b>Introduction</b> .....	<b>1</b>
1.1	Specifications .....	1
1.2	Safety Instructions .....	2
<b>Section 2</b>	<b>Trans-Blot Cell Assembly</b> .....	<b>4</b>
2.1	Assembly of the Unit for Standard Transfers .....	4
2.2	High Intensity Field Option .....	7
2.3	Acidic Transfers .....	8
<b>Section 3</b>	<b>Transfer Conditions</b> .....	<b>8</b>
3.1	General Guide for Transfer Buffers and Running Conditions .....	8
3.2	Notes on Electrophoretic Transfer Conditions .....	9
3.3	Buffer Formulation .....	11
<b>Section 4</b>	<b>Strategies for Optimizing Electro-Elution</b> .....	<b>12</b>
4.1	Optimizing Protein Transfer .....	12
4.2	Optimizing DNA and RNA Transfer .....	13
<b>Section 5</b>	<b>Choice of Blotting Membranes</b> .....	<b>14</b>
5.1	Protein Blotting Membranes .....	14
5.2	DNA and RNA Blotting Membrane .....	14
<b>Section 6</b>	<b>Troubleshooting Guide</b> .....	<b>15</b>
<b>Section 7</b>	<b>Product Information</b> .....	<b>22</b>
<b>Section 8</b>	<b>References</b> .....	<b>22</b>

## Section 1 Introduction

Blotting was first performed by Southern<sup>1</sup> in 1975 with the transfer of DNA from agarose gels to nitrocellulose membranes. Blotting has subsequently been applied to RNA<sup>2-4</sup> and protein<sup>5,6</sup> from both agarose and polyacrylamide gels. To overcome the inefficiencies observed in various capillary transfers, electric current was used to elute protein from polyacrylamide gels, as first described by Towbin<sup>7</sup> *et al.* in 1979. Since that time, electrophoretic transfer has been applied to DNA and RNA.<sup>8-14</sup> Numerous publications have dealt with the topic of protein electrophoretic transfer technique.<sup>15-26</sup> There have also been several reviews summarizing the literature on electrophoretic blotting.<sup>27-33</sup>

The Trans-Blot cell is available with standard electrode cards or with plate electrodes, which consist of a platinum-coated titanium anode and a stainless steel cathode. The Trans-Blot cell allows electrophoretic transfers to be performed in the standard configuration, with either set of electrodes positioned 8 cm apart, using up to three gel holders positioned between the electrodes. Removable electrodes can be positioned 4 cm apart with one gel holder between. This allows generation of a high intensity electrical field, when used in combination with the PowerPac 200 Power Supply. An optional Super Cooling Coil is available for temperature control during routine blots and high power transfers.

### 1.1 Specifications

#### Construction

Trans-Blot tank	Molded polysulfone
Electrode cards	Cathode 18" Anode 25"
Anode plate electrode	Platinum-coated titanium
Cathode plate electrode	Stainless steel
Gel holder	Acrylic

**Overall dimensions** 24 x 18 x 9.5 cm

**Gel holder dimensions** 21 x 17 cm

**Buffer capacity** 3 L

**Transfer capacity** 3 gel holders


**Cleaning** Use mild soap and warm water to clean the electrodes, cassettes, and buffer tank. Use special care when cleaning the electrode cards or plate electrodes. Avoid stretching or breaking the platinum wires. Avoid scratching or marring the platinum plate. Do not use abrasives or strong detergents. The cathode plate (stainless steel) can be cleaned with a mild abrasive to remove salt that may be deposited during normal operation. Rinse the fiber pads under hot water and then in distilled, deionized water.

**Chemical compatibility** The Trans-Blot cell components are not compatible with chlorinated hydrocarbons (e.g., chloroform), aromatic hydrocarbons (e.g., toluene, benzene), or acetone. **Use of organic solvents voids all warranties.**

## 1.2 Safety Instructions



Power to the Trans-Blot cell is supplied by an external DC voltage power supply. This power supply must be ground isolated in such a way that the DC voltage output floats with respect to ground. All of Bio-Rad's power supplies meet this important safety requirement. Regardless of which power supply is used, the maximum specified operating parameters for the cell are:

<b>300 VDC</b> 	Maximum voltage limit
<b>200 Watts</b>	Maximum power limit
<b>50 °C</b>	Maximum ambient temperature limit



Current to the cell, provided from the external power supply, enters the unit through the lid assembly, providing a safety interlock to the user. Current to the the cell is broken when the lid is removed. **Do not attempt to circumvent this safety interlock, and always turn the power supply off before removing the lid, or when working with the cell in any way.**

### Important

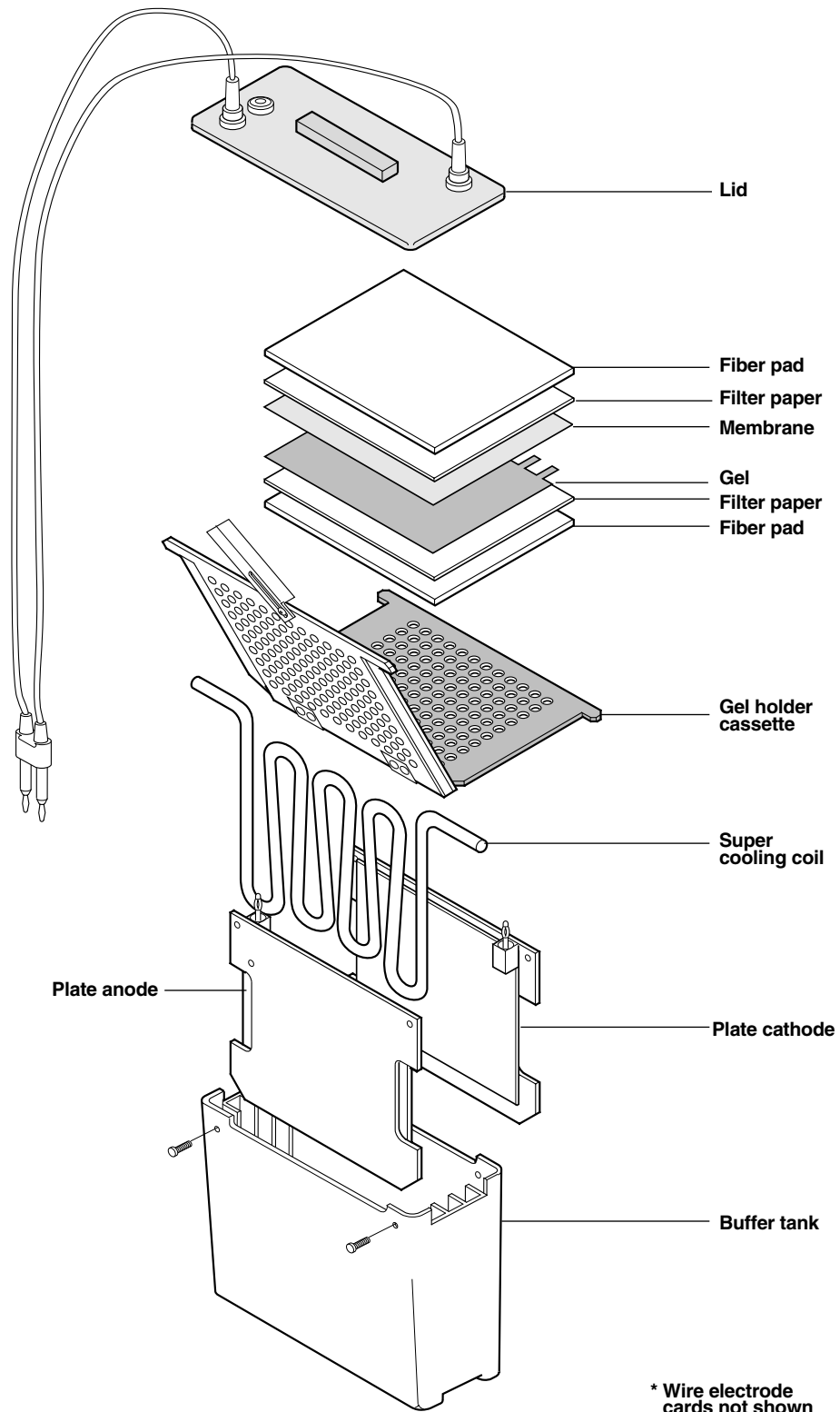
This Bio-Rad instrument is designed and certified to meet IEC 1010-1\* safety standards. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will:

- **Void the manufacturer's warranty**
- **Void the IEC1010-1 safety certification**
- **Create a potential safety hazard**

Bio-Rad is not responsible for any injury or damage caused by the use of this instrument for purposes other than for which it is intended or by modifications of the instrument not performed by Bio-Rad or an authorized agent.

\*IEC 1010-1 is an internationally accepted electrical safety standard for laboratory instruments.

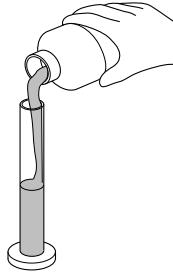
## Trans-Blot Cell Description & Assembly of Parts



## Section 2 Trans-Blot Cell Assembly

### 2.1 Preparation for Blotting

1. Prepare the transfer buffer. (See Section 3.3 for buffer formulation. Using buffer chilled to 4 °C will improve heat dissipation.)



2. Cut the membrane and the filter paper to the dimensions of the gel. Always wear gloves when handling membranes to prevent contamination. Equilibrate the gel and soak the membrane, filter paper, and fiber pads in transfer buffer (15 min—1 hour depending on gel thickness).
3. Prepare the gel sandwich.

Place the cassette, with the gray side down, on a clean surface.

Place one pre-wetted fiber pad on the gray side of the cassette.

Place a sheet of filter paper on the fiber pad.

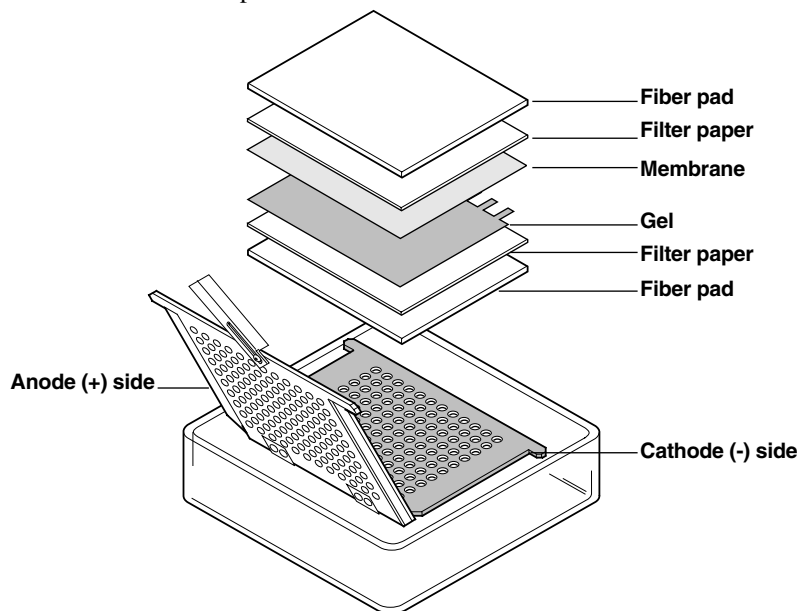
Place the equilibrated gel on the filter paper.\*

Place the pre-wetted membrane on the gel.\*

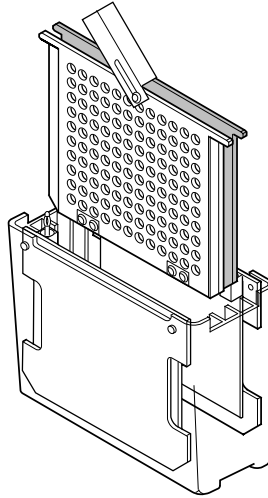
Complete the sandwich by adding a piece of filter paper on the membrane.\*

\* Removing any air bubbles which may have formed is very important for good results. Use a glass tube to gently roll air bubbles out.

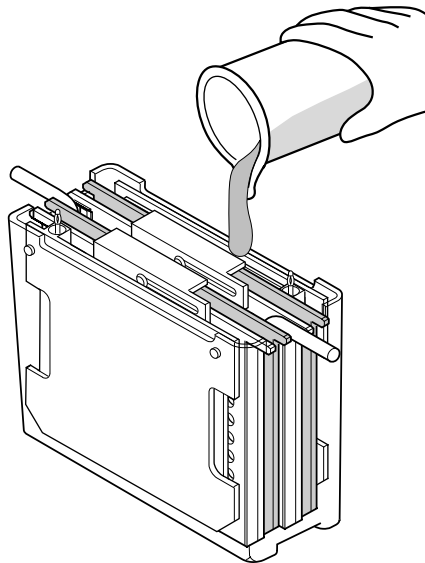
Add the last fiber pad.



4. Close the cassette firmly, being careful not to move the gel and filter paper sandwich. Lock the cassette closed with the white latch.
5. Place the cassette in tank. Repeat for any other cassettes you may want to run. Completely fill the tank with buffer.



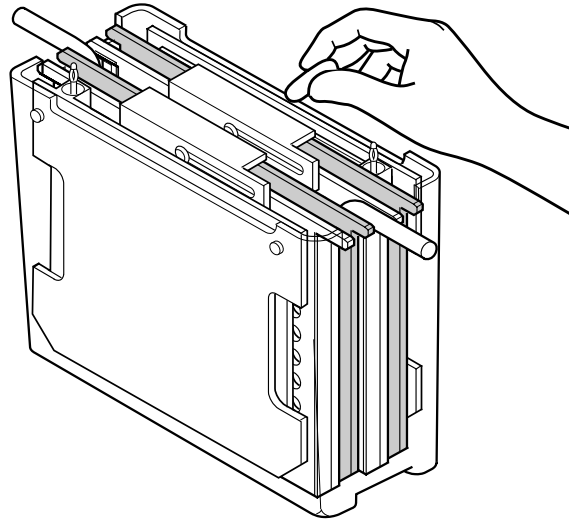
6. If running high intensity or high voltage with plate electrodes, add the super cooling coil in any available slot.



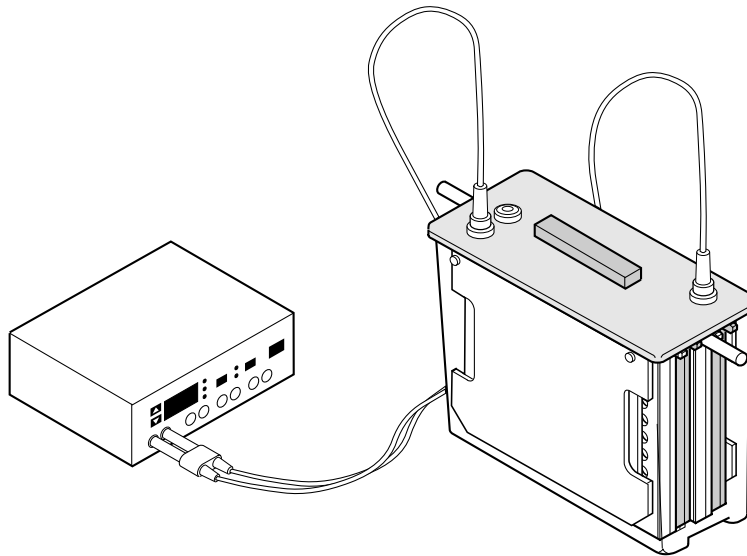
**Note:** For high intensity runs (4 cm inter-electrode distance) only one cassette can be placed in the tank. The Super Cooling Coil must be used. It is placed outside the electrodes. See Section 2.2, high intensity field option. Two cassettes can be run with the Super Cooling Coil in the standard field with the electrodes 8 cm apart, and three cassettes can be run at low voltage without the super cooling coil.



7. Addition of a standard stir bar will help to maintain even buffer temperature and ion distribution in the tank. Set the speed as fast as possible to keep ion distribution even.



8. Put on the lid, plug the cables into the power supply, and run the blot. Refer to Section 3 for run times and voltage setting with various buffers.

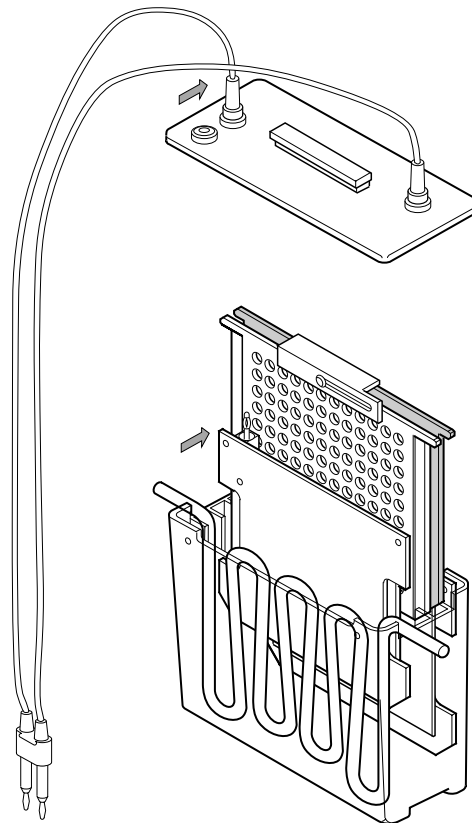


9. Upon completion of the run, disassemble the blotting sandwich and remove the membrane for development. Clean the cell, fiber pads, and cassettes with laboratory detergent and rinse well with deionized water.

## 2.2 High Intensity Field Option

The electrode cards of the Trans-Blot cell can be moved closer together to allow higher field strengths (V/cm), and thus more efficient, more rapid transfer. To decrease the inter-electrode distance, the anode electrode card is moved to the center gel holder position.

1. Remove the screws at the top corners of the anode (red) electrode card.
2. Set up the gel holder with the gel, pads, and membrane following the standard Trans-Blot cell operation instructions outlined in Section 2.1. Insert the sandwich into the gel holder slot closest to the cathode (black) electrode panel.
3. Pull the anode card out of the cell, and move it into the middle gel holder slot. The gel holder will keep the electrode card firmly positioned in the slot.
4. Completely fill the buffer tank to the bottom of the red electrode disc.
5. Place the Super Cooling Coil into the free gel holder slot and place a stir bar in the bottom of the tank. The Super Cooling Coil must be used for adequate heat dissipation of generated in high-intensity transfer. Set stirring speed as high as possible.
6. Remove the white plug from the electrode cable hole in the Trans-Blot cell lid. Unscrew the two red screws anchoring the red power cable to the lid. Loosen the screws 4 turns. Move the cable to its alternative position, aligning shallow holes in the cable with the anchor screws, and retighten all screws.
7. Begin electrophoretic blotting.



**The anode electrode is removed from the standard position and placed 4 cm apart from the cathode. The cable is also moved to line up with electrodes.**

## 2.3 Acidic Transfers

If transferring under acidic conditions, switch the gel and membrane in the set up instructions. This will place the membrane on the cathode side of the gel. Under acidic conditions, proteins will transfer in the opposite direction going toward the negative cathode. **Do not reverse the electrodes themselves.** This will cause damage to the instrument.

## Section 3 Transfer Conditions

### 3.1 General Guidelines to Transfer Buffers and Running Conditions

Tables 3.1 to 3.4 provide guidelines for power conditions using different buffers. Power conditions are provided for various run times. Where multiple conditions are displayed, the higher the voltage, the less time required for the run. Note all “requires cooling” warnings.

**Table 3.1. SDS-PAGE Gels**

	Standard Field Overnight		Standard Field 8 cm electrode distance		High Intensity Field 4 cm electrode distance *COOLING REQUIRED	
	Buffer A or B	Buffer C	Buffer A or B	Buffer C	Buffer A or B	Buffer C
<b>Trans-Blot cell with wire electrodes</b>	30 V, 0.1 A	10 V, 0.1 A	60 V, 0.21 A	20 V, 0.25 A	100–200 V 0.3–0.8 A	30–60 V 0.4–1.0 A
	☞Overnight		☞5 hours approximately		☞30 minutes–4 hours *	
<b>Trans-Blot cell with plate electrodes</b>	10 V 0.1 A	5 V 0.1 A	100–150 V 1–1.5 A	30–50 V 1–1.6 A	50–100 V 0.7–1.6 A	20–40 V 0.7–1.6 A
	☞Overnight		☞1–5 hours *		☞30 minutes–1 hour *	

Buffer A: 25 mM Tris, pH 8.3, 192 mM glycine, with or without 20% MEOH and .025%–0.1% SDS.  
 B: 48 mM Tris, pH 9.2, 39 mM glycine, with or without 20% MEOH and .025%–0.1% SDS.  
 C: 10 mM NaHCO<sub>3</sub>, 3 mM NaCO<sub>3</sub>, pH 9.9, with or without 20% MEOH and .025%–0.1% SDS.

**Table 3.2. DNA and RNA**

	<b>Standard Field Overnight</b>	<b>Standard Field 8 cm electrode distance</b>		<b>High Intensity Field 4 cm electrode distance</b> *COOLING REQUIRED	
<b>Trans-Blot cell with wire electrodes</b>	TAE or TBE 30 V, 0.2 A  ☞Overnight	TAE 60 V, 0.6 A	TBE 80 V, 0.55 A  ☞5 hours approximately	TAE 80–100 V 0.8–1.4 A	TBE 100–130 V 0.7–1.1 A  ☞30 minutes–4 hours *
<b>Trans-Blot cell with plate electrodes</b>	TAE or TBE 15 V, 0.2 A  ☞Overnight	TAE 20–35 V, 0.8–1.5 A	TBE 75–100 V 0.8–1.1 A  ☞1–5 hours *	TAE 15–25 V 0.8–1.4 A	TBE 50–90 V 0.7–1.1 A  ☞30 minutes–1 hour *

1x TAE: 40 mM Tris-Acetate, 1.0 mM EDTA

1x TBE: 90 mM Tris-Borate, 2.0 mM EDTA

**Table 3.3. Native Gels**

	<b>Standard Field Overnight</b>	<b>Standard Field 8 cm electrode distance</b>	<b>High Intensity Field 4 cm electrode distance</b> *COOLING REQUIRED
<b>Trans-Blot cell with wire electrodes</b>	30 V, 0.1 A  ☞Overnight	60 V, 0.21 A  ☞5 hours approximately	100–200 V 0.3–0.8 A  ☞30 minutes–4 hours *
<b>Trans-Blot cell with plate electrodes</b>	10 V 0.1 A  ☞Overnight	100–150 V 1–1.5 A  ☞1–5 hours *	50–100 V 0.7–1.5 A  ☞30 minutes–1 hour *

Buffer: 25 mM Tris, pH 8.3, 192 mM glycine. No methanol.

**Table 3.4. Isoelectric Focusing, Native Gels, Basic Proteins, Acid Urea Gels**

	<b>Standard Field Overnight</b>	<b>Standard Field 8 cm electrode distance</b>	<b>High Intensity Field 4 cm electrode distance</b> *COOLING REQUIRED
<b>Trans-Blot cell with wire electrodes</b>	30 V, 0.2 A  ☞Overnight	70 V, 0.5 A  ☞5 hours approximately	100–150 V, 0.55–0.85 A  ☞30 minutes–4 hours *
<b>Trans-Blot cell with plate electrodes</b>	15 V, 0.2 A  ☞Overnight	40–70 V, 0.6–1.0 A 1–5 hours *	30–60 V, 0.6–1.0 A  ☞30 minutes–1 hour *

Buffer: 0.7% Acetic Acid

## 3.2 Notes on Electrophoretic Transfer Conditions

**These variables will change total resistance and thus the current readings:**

- Alterations in buffer make-up, *i.e.*, addition of SDS, or changes in ion concentration due to addition of acid or base to adjust the pH of the buffers.
- Gel pH, ionic strength, and percentage of acrylamide, especially if the gel has not been properly equilibrated.
- Number of gels; current increases slightly as the number of gels increases.
- Volume of buffer; current increases when volume increases.
- Platinum mass; current increases when mass increases.
- Transfer temperature; current increases when temperature increases.
- Time in transfer at which reading was taken; current normally increases as the buffering capacity diminishes with progress of the run.

## **2. Pre-equilibration of gels**

All electrophoresis gels should be pre-equilibrated in transfer buffer prior to electrophoretic transfer. Pre-equilibration will facilitate the removal of contaminating electrophoresis buffer salts and neutralization salts (salts resulting from the denaturation of nucleic acids prior to transfer). If the salts are not removed, they will increase the conductivity of the transfer buffer and the amount of heat generated during the transfer. Also, low percentage gels (<12%) will shrink in methanol buffers. Equilibration allows the gel to adjust to its final size prior to electrophoretic transfer.

## **3. Current limits**

The PowerPac 200 Power Supply is capable of a 200 watt output. Unless a current limit is set, uncontrolled conductivity changes may result in full power being delivered to the Trans-Blot cell. The gel holder and electrode cards may warp, and the transfer buffer may boil and evaporate (further increasing conductivity). This would result in a potential safety hazard. Refer to the PowerPac 200 Power Supply Instruction Manual for setting current limits and run times.

## **4. Polarity of transfer**

Do not reverse polarity with the plate electrodes. This will result in corrosion and rusting of the stainless steel cathode. If this should occur, the stainless steel should be cleaned with a mild abrasive cleanser to remove the rust.

## **5. Heat dissipation with the Super Cooling Coil - Use of the Super Cooling Coil with the Plate Electrodes**

For high power applications, such as high intensity transfers and transfers performed with the plate electrodes, efficient heat removal is only obtained using the Super Cooling Coil connected to a refrigerated recirculating bath. Placing the Trans-Blot cell in the cold room is an inadequate means of controlling transfer buffer temperature. The plastic tank of the Trans-Blot cell is an effective thermal insulator, thus limiting the efficient dissipation of heat. Use of the Super Cooling Coil is necessary with all experiments using the plate electrodes and with high intensity transfers using the standard platinum wire electrodes.

## **6. Use of a stir bar during transfer**

For all blotting applications a stir bar must be placed inside the Trans-Blot cell, so that the transfer buffer is stirred during the course of the experiment. This will help to maintain uniform conductivity and temperature during electrophoretic transfer. Failure to properly control transfer buffer temperature results in poor transfer of macromolecules and poses a potential safety hazard.

## **7. Transfer buffer pH**

Do not adjust the pH of transfer buffers unless specifically indicated. Adjustments of the pH of transfer buffers, when not indicated, will result in increased buffer conductivity. This is manifested by a higher than expected initial current output and a decreased resistance. It is recommended that the buffer conductivity and resistance be checked with the PowerPac 200 Power Supply before starting each transfer.

## 8. Transfer buffer recommendations

Use only high quality, reagent grade methanol. Contaminated methanol can result in increased transfer buffer conductivity, as well as poor transfer of macromolecules. Do not reuse transfer buffers or dilute transfer buffers below recommended levels. Reuse of transfer buffers is not advised, since these buffers have most likely lost their ability to maintain a stable solution pH during transfer. Dilution of transfer buffers below their recommended levels is also not advised, since this will decrease their buffering capacity.

## 9. Voltage limits

Do not increase voltage settings beyond those indicated in Tables 3.1–3.4 for overnight operation. Buffer conductivity must be close to the current listed and a current limit should be set on the power supply. If overnight transfers at low voltages are ineffective for your application, and higher voltages are necessary, transfer times must also be decreased. Failure to do so may result in a potential safety hazard.

## 3.3 Buffer Formulation

All formulas provided below are for a total volume of 1 liter of buffer. Four liters of buffer are required for the Trans-Blot cell.

**DO NOT ADD ACID OR BASE TO ADJUST pH OF THE FOLLOWING BUFFERS.** Methanol should be analytical reagent grade, as metallic contaminants in low grade methanol will plate on the electrodes.

**Note:** Some pH electrodes will not perform a proper measurement for the pH of Tris buffers. If the pH of the buffer is off, check to make sure the electrode is designed to work with Tris buffers. If the pH electrode functions properly for Tris buffers and the pH is below 8.0, remake the buffer.

1. 25 mM Tris, 192 mM glycine, 20% v/v methanol, pH 8.3  
Mix 3.03 g Tris, 14.4 g glycine, and 200 ml of methanol; add distilled deionized water (dd H<sub>2</sub>O) to 1 liter.
2. 25 mM Tris, 192 mM glycine, pH 8.3  
Mix 3.03 g Tris and 14.4 g glycine; add dd H<sub>2</sub>O to 1 liter.
3. 48 mM Tris, 39 mM glycine, 20% v/v methanol, pH 9.2  
Mix 5.82 g Tris and 2.93 g glycine in ddH<sub>2</sub>O, add 200 ml methanol  
Bring to 1 liter with ddH<sub>2</sub>O
4. 48 mM Tris, 39 mM glycine, pH 9.2  
Mix 5.82 g Tris and 2.93 g glycine  
Add ddH<sub>2</sub>O to 1 liter
5. 10 mM NaHCO<sub>3</sub>, 3 mM NaCO<sub>3</sub>, 20% methanol, pH 9.9  
Mix 0.84 g NaHCO<sub>3</sub> and 318 g Na<sub>2</sub>CO<sub>3</sub> in ddH<sub>2</sub>O, add 200 ml methanol  
Bring to 1 liter with ddH<sub>2</sub>O

6. 1.0x TBE (Tris-Borate EDTA) pH 8.3  
90 mM Tris-Borate 1 mM EDTA  
  
5x stock solution  
54 g Tris base  
27.5 boric acid  
20 ml 0.5 m EDTA (pH 8.0)  
  
Add 200 ml 5x stock to 800 ml ddH<sub>2</sub>O to make 1.0x working solution.
7. 1x TAE (Tris-Acetate EDTA)  
40 mM Tris-Acetate 1 mM EDTA  
  
50-x stock solution  
242 g Tris base  
57.1 ml glacial acetic acid  
100 ml 0.5 m EDTA (pH 8.0)  
  
1x working solution, add 20 ml 50x stock to 980 ml ddH<sub>2</sub>O

## Section 4 Strategies for Optimizing Electro-Elution

### 4.1 Optimizing Protein Transfer

Generally, quantitative elution of denatured high molecular weight proteins is difficult. The following tactics, alone or in combination, will increase transfer efficiency.

1. Vary gel composition
  - a. Gradient gels are often more effective than single gel concentrations for elution of a wide range of molecular weight proteins.
  - b. Lower the total monomer to create a more porous gel.
  - c. Increase or decrease the percentage of crosslinker. A 5.26% C gel will contain the smallest pore size of all gels no matter what the concentration of acrylamide. An increase or decrease in %C will make gels more porous with little loss in resolution.

$$\%C = \frac{\text{grams bis}}{\text{grams bis} + \text{grams acrylamide}} \times 100$$

2. Increase transfer time. An initial control should be performed to determine the time required for complete transfer.<sup>17, 24</sup> Times may vary from as little as 30 minutes to as long as overnight. Remember all overnight applications should be performed at 30–50 volts to minimize heating problems. (For long transfers at elevated voltages use the standard Trans-Blot cell with the super cooling coil.)
3. Increase the power. Initial controls should be performed to evaluate the efficiency of increasing the V/cm as well as its effects on the temperature of transfer. The temperature increase may change buffer resistance and subsequent power delivered, as well as the state of protein denaturation, thus affecting transfer efficiency.
4. Reduce buffer strength. Dilution of transfer buffer results in lower current at any given voltage. This will allow the use of higher voltages without excessive heating.

5. Vary buffer type and pH
  - a. Maximize charge-to-mass ratio. It appears that alcohols present in SDS transfer buffer strip SDS from proteins. Basic proteins in Tris, glycine, methanol buffer at pH 8.3 may assume a state near isoelectric neutrality and thus transfer poorly. For example, lysozyme exhibits this behavior. Buffers with pH of 9.5 to 10.0 have shown much better elution and binding characteristics for basic proteins such as lysozyme and histones.<sup>41</sup>
  - b. Different buffer types at similar V/cm may yield different efficiencies. Generally Tris buffers allow more efficient transfer than acetate or phosphate buffers.
6. Addition of 0.1% SDS detergent to Tris, glycine, methanol buffer has been reported to increase transfer efficiency.<sup>24</sup> SDS, however, increases relative current, power, and heating. Also, temperatures below 10 °C may precipitate the SDS so the starting buffer temperature will be higher. SDS may also affect the antigenicity of some proteins. SDS will aid in eluting the proteins from the gel, but it may reduce the binding efficiency of those proteins to the nitrocellulose membrane.<sup>42</sup>
7. Eliminate alcohol from the transfer buffer. Alcohol in the transfer buffer improves binding of SDS proteins to nitrocellulose only. Elimination of alcohol results in increased transfer efficiency but diminishes binding to nitrocellulose. Transfer efficiency is increased because alcohol causes gel pores to contract resulting in fixation of large molecular weight proteins within the gel matrix. Use of PVDF membrane for SDS protein transfers eliminates the alcohol requirement, and constitutes a logical strategy for analysis of high molecular weight or difficult-to-transfer proteins.<sup>26, 27</sup>
8. Limited protease treatment. A protocol for protease digestion of protein during transfer has been published.<sup>22</sup> Efficient transfer without loss of immunological reactivity was reported.
9. Alter membrane type. As mentioned in 7, PVDF membrane allows transfer in the absence of alcohol.
10. Alter gel system. If possible, use non-denaturing gradient pore gels for separation of proteins by molecular weight. Isoelectric focusing gels, or native gels, may be considered if separation by molecular weight is not mandatory.
11. Failure of molecules to bind efficiently to the membrane, caused by poor gel-membrane contact, is often confused with inefficient elution. Poor contact is usually due to excess moisture in the gel-membrane interface. Proper technique and the use of a test tube or glass pipet as a “rolling pin” should assure good contact. Proper selection of filter paper spacers will help assure good compression. Gel and membrane equilibration in transfer buffer for 30 minutes to 1 hour prior to transfer will help prevent shrinking of either component during transfer, and will eliminate reactants such as urea or SDS from the gel.

## 4.2 Optimizing DNA and RNA Transfer

Problems with elution of nucleic acids can be solved by altering the gel percentage. It may be somewhat more difficult to quantitatively transfer large amounts of DNA used in genomic blots. The following tactics should be considered for optimizing elution in such transfers.

1. Alter gel composition.
  - a. Lower % total monomer or % crosslinker for polyacrylamide gels.
  - b. Lower % agarose. This allows better elution of high molecular weight DNA.
2. Alter DNA denaturants. It has been found that glyoxal denaturation allows more efficient elution of DNA than NaOH. Boiling polyacrylamide gels to denature DNA has also been found to give excellent results.<sup>11</sup> Base denaturation often causes polyacrylamide gels to weaken and stick to blotting membranes.



## **Section 5**

### **Choice of Blotting Membranes**

#### **5.1 Protein Blotting Membranes**

##### **Nitrocellulose Membrane**

Nitrocellulose membranes have been used extensively for protein binding and detection.<sup>7,20,23,24,27</sup> They can be easily stained for total protein by a dye stain (Amido Black, Coomassie® Blue, Ponceau S, Fast Green FCF, etc.),<sup>27</sup> or the more sensitive Colloidal Gold Total Protein Stain, and also allow either RIA, FIA or EIA.<sup>7</sup> Nitrocellulose has a high binding capacity of 80-100 µg/cm<sup>2</sup>. Nonspecific protein binding sites are easily and rapidly blocked, avoiding subsequent background problems. No pre-activation is required. Low molecular weight proteins (especially <20,000 daltons) may be lost during post transfer washes, thus limiting detection sensitivity.<sup>19</sup> Smaller pore size nitrocellulose membrane (0.2 µm), has been shown to be effective in eliminating this loss.<sup>37</sup> Large proteins ( 100,000 daltons) denatured by SDS may transfer poorly due to the addition of alcohol to the transfer buffer. Alcohol increases binding of SDS-proteins to nitrocellulose, but decreases pore sizes in the gel. Elimination of alcohol from SDS-protein transfers results in considerably diminished binding. Adding SDS (up to 0.1%) to the transfer buffer increases the transfer efficiency of proteins, but reduces the amount of binding to the membrane.<sup>17</sup> Also, SDS increases the conductivity of the buffer and the heat generated during transfer.

##### **PVDF Membrane**

PVDF (Polyvinylidene difluoride) membrane is an ideal support for amino-terminal sequencing, amino acid analysis and immunoassays of blotted proteins. PVDF retains proteins under extreme conditions of exposure to acidic or basic conditions, and in the presence of organic solvents. Greater retention during sequencing manipulations enhances the likelihood of obtaining information from rare, low abundance proteins, by increased initial coupling and higher repetitive yields. In addition, PVDF membrane exhibits better binding efficiency of blotted material in the presence of SDS in the transfer buffer. PVDF must first be wetted in 100% MeOH but can then be used in buffer which does not contain MeOH.

#### **5.2 DNA and RNA Blotting Membrane**

##### **Zeta-Probe® Nylon Membrane**

Nitrocellulose is not a suitable medium for electrophoretic transfer of nucleic acids, as high concentrations of salt ( 10 x SSC) are required for efficient binding.<sup>13</sup> Molecules > 500 bp are not bound at all, even at high salt. Low resistance results when an electric current is passed through a solution of high salt. This causes potentially damaging high currents (and power) at very low voltages. Since V/cm is the eluting force, inefficient transfer occurs under conditions required for proper binding. Zeta-Probe membrane allows efficient binding of all sizes of single stranded DNA and RNA in the presence of low ionic strength buffers.<sup>13</sup> Zeta-Probe membrane is an ideal alternative to nitrocellulose for the analysis of nucleic acids. Binding is more stable through post transfer washes, and reprobing may be performed as many as 10 times.

**Table 5.1 Guide to Protein Blotting Membranes**

A variety of blotting membranes is available for immunoblotting, each with particular advantages depending on the needs of the experiment. The physical properties and performance characteristics of a membrane should be evaluated in selecting the appropriate transfer conditions.

Membrane	Pore Size	Binding Capacity ( $\mu\text{g}/\text{cm}^2$ )	Notes
Nitrocellulose	0.45 $\mu\text{m}$ 0.2 $\mu\text{m}$	80–100	General purpose protein blotting membrane
Supported Nitrocellulose	0.45 $\mu\text{m}$ 0.2 $\mu\text{m}$	80–100	Pure nitrocellulose cast on an inert synthetic support; increased strength for easier handling and for reprobing.
PVDF	0.2 $\mu\text{m}$	170–200	High mechanical strength and chemical stability, used for protein sequencing, and western blotting; enhanced binding in the presence of SDS. Must be wet in alcohol before equilibration in buffer.
Nylon	0.2 $\mu\text{m}$	170	Recommended for nucleic acids.

**Note:** Nucleic acids cannot be transferred to nitrocellulose by electrophoretic blotting. Use Zeta-Probe membrane.

## Section 6 Troubleshooting Guide

### Electrophoretic Transfer

#### *Poor electrophoretic transfer (as detected by staining the gel)—Proteins*

1. Transfer time is too short.
  - Increase the transfer time.
2. Power is too low.
  - Always check the current at the beginning of the run. The current may be too low for a particular voltage setting. If the buffer is prepared improperly, the conductivity may be too low, and not enough power will be delivered to the cell. See the power guidelines for specific applications in Section 3.
  - Remake the buffer or increase the voltage.
  - Try the high intensity blotting option.
3. Transfer apparatus is assembled incorrectly, and the proteins are moving in the wrong direction.
  - The gel/membrane sandwich may be assembled in the wrong order or the cassette is inserted in the tank facing the opposite orientation. Check the polarity of the connections to the power supply.
4. Charge-to-mass ratio is incorrect.
  - Try a more basic or acidic transfer buffer to increase protein mobility. Proteins near their isoelectric point at the pH of the buffer will transfer poorly. (It has been suggested that buffer pH should be 2 pH units higher or lower than the pI of the protein of interest for optimal transfer efficiency.)
5. Protein is precipitating in the gel.
  - Try using SDS in the transfer buffer. SDS can increase transfer efficiency, but can also reduce binding efficiency to nitrocellulose and affect reactivity of some proteins with antibodies.

6. Power supply circuit is inoperative, or an inappropriate power supply was used.
  - Check the fuse. Be sure the voltage and current output of the power supply match the needs of the blotting instrument.
7. Methanol in the transfer buffer is restricting elution.
  - Reduction of methanol results in increased transfer efficiency of proteins from the gel, but it also diminishes binding to nitrocellulose and PVDF.
8. Gel percentage too high.
  - Reduce %T (total monomer) or %C (crosslinker). A 5% C (with bis as the crosslinker) will produce the smallest pore size gel. Decreasing from this concentration will increase the pore size and increase transfer efficiency.

**Poor transfer—nucleic acid**

1. Gel percentage is too high.
  - Reduce the %T or %C in the acrylamide gel or reduce % agarose in an agarose gel.
  - Prior to transfer, cleave DNA in dilute 0.25 M HCl or RNA in dilute NaOH.
2. Transfer time is too short or power conditions are too low.
  - Increase the transfer time, or try high intensity transfer.
3. DNA or RNA cannot be transferred electrophoretically to nitrocellulose, since high salt concentrations are required for efficient binding.
  - Use Zeta-Probe membrane instead of nitrocellulose.

**Swirls or missing bands; diffuse transfers**

1. Poor contact between the membrane and the gel. Air bubbles or excess buffer remain between the blot and gel.
  - Use a test tube or pipet as a rolling pin, and roll over the membrane carefully in both directions until air bubbles or excess buffer is removed from between gel and membrane, and complete contact is established.
  - Use thicker filter paper in the gel/membrane sandwich.
  - Replace the fiber pads. Pads will compress with time, and will not hold the membrane to the gel.
2. Power conditions are too high.
  - Always check the current at the beginning of the run. The current may be too high for a particular voltage setting. If the buffer is prepared improperly, the conductivity may be too high, resulting in excessive power delivered to the cell. See the power guidelines for specific applications in Section 3.
3. The membrane is not properly wet or has dried out.
  - White spots on the nitrocellulose membrane indicate dry areas where protein will not bind. If wetting does not occur immediately by immersion of the sheet in transfer buffer, heat distilled water until just under the boiling point, and soak the membrane until completely wet. Equilibrate in transfer buffer until ready for use.
  - Because of the hydrophobic nature of PVDF, the membrane must be prewet in methanol prior to equilibration in aqueous transfer buffer. Follow the directions in the product insert.

4. The gel electrophoresis may be at fault.
  - Artifacts of electrophoresis may be produced by poor polymerization, inappropriate running conditions, contaminated buffers, sample overload, etc. Consult your electrophoresis manual for more details.

#### ***Gel cassette pattern transferred to blot***

1. Contaminated or thin fiber pads are used.
  - Replace the fiber pads, or thoroughly clean the contaminated pads.
2. Excessive amounts of protein were loaded on the gel, or too much SDS was used in the transfer buffer. Proteins can pass through the membrane without binding, and recirculate through the tank blotting system.
  - Reduce the amount of protein on the gel, and SDS in the transfer buffer. Add a second sheet of membrane to bind excess protein.
3. The transfer buffer is contaminated.
  - Make fresh solutions.

#### ***Poor binding to the membrane — Nitrocellulose***

1. Nitrocellulose requires 20% methanol in the transfer buffer for optimal protein binding.
  - Make sure the buffer contains the proper amount of methanol.
2. Proteins may be transferring through the nitrocellulose.
  - Use PVDF or nylon (higher binding capacities) or 0.2  $\mu\text{m}$  nitrocellulose (smaller pore size). Decrease the voltage or move the electrodes to the standard position if using the high intensity option.
3. Mixed ester celluloses bind proteins poorly.
  - Use pure nitrocellulose.
4. Proteins <15,000 daltons may show diminished binding to 0.45  $\mu\text{m}$  nitrocellulose, or may be washed from the membrane during assays.
  - To increase stability of binding, proteins can be crosslinked to nitrocellulose with glutaraldehyde.
  - Use PVDF or nylon membrane, which have higher binding capacities.
  - Use Tween-20 detergent in the wash and antibody incubation steps. Reduce or eliminate the more stringent washing conditions.
5. SDS in the transfer buffer will reduce binding efficiency of proteins.
  - Reduce or eliminate the SDS from the transfer buffer.
6. The membrane may not be completely wet.
  - White spots on the membrane indicate dry areas where protein will not bind. If wetting does not occur immediately by immersion of the sheet in transfer buffer, heat distilled water until just under the boiling point, and soak the membrane until completely wet. Equilibrate in transfer buffer until ready for use.

#### ***Poor Binding to the Membrane — PVDF***

1. The membrane may not be completely wet.
  - Because of the hydrophobic nature of PVDF, the membrane must be prewet in alcohol prior to equilibration in aqueous transfer buffer. Follow the directions in the product insert.
2. The membrane may have been allowed to dry during handling.
  - A completely wet membrane has a gray, translucent appearance. White spots will form on the surface of the membrane, indicating that it has been allowed to dry. Since proteins will not bind to the dry spots, rewet the membrane with methanol and re-equilibrate in transfer buffer.

## Immune-Specific Detection

### **Overall High Background**

1. Blocking conditions are inappropriate.
  - Match the blocker to the membrane. For example, nylon and PVDF membranes require more extensive blocking, usually with non-fat dry milk.
  - Increase the concentration or blocking time as necessary.
  - The blocker must be a pure protein. The blocker may be contaminated with material that bind probes non-specifically.
2. Insufficient wash protocols are used.
  - Increase the number, duration, or stringency of the washes. Include progressively stronger detergents in the washes, *e.g.* SDS is stronger than NP-40 which is stronger than Tween-20. Also, include Tween-20 in the antibody dilution buffers to reduce nonspecific binding.
3. The blot is left in the substrate too long.
  - Remove the blot from the substrate solution when the signal-to-noise level is acceptable. Do not overdevelop. Stop the reaction immediately by immersing the blot in dd H<sub>2</sub>O.
4. Contamination occurred during a previous step, *e.g.* electrophoresis or transfer.
  - Discard and remake the gel and transfer solutions.
  - Replace or thoroughly clean contaminated fiber pads, if a tank blotter was used. Excessive amounts of protein were loaded on the gel, or too much SDS was used in the transfer buffer. Proteins can pass through the membrane without binding and recirculate through a tank blotting system. Reduce the amount of protein on the gel or SDS in the transfer buffer. Add a second sheet of membrane to bind excess protein.
5. Primary or secondary antibody is too concentrated.
  - Increase the dilution of the antibodies. Perform a dot-blot experiment to optimize the working concentrations.
6. Incubation trays are contaminated.
  - Clean the trays or use disposable trays.

### **Nonspecific Reactions between Bound Proteins and Probes**

1. Primary or secondary antibody is contaminated with nonspecific or species cross-reactive IgG.
  - Use purified IgG first antibody fractions and affinity-purified blotting grade secondary antibody.
2. Monoclonal antibodies may react non-specifically with SDS denatured proteins.
  - Compare the binding of other monoclonal or polyclonal antibodies.
  - Blot native proteins as a comparison.
3. Nonsense interactions are occurring due to ionic associations. For example, avidin, a glycosylated protein, may bind to more acidic proteins on blots.
  - Increase the ionic strength of the incubation buffers. Increase the number, duration or stringency of the washes. Include progressively stronger detergents in the washes, *e.g.* SDS is stronger than NP-40 which is stronger than Tween-20. Include Tween-20 in the antibody dilution buffers to reduce nonspecific binding.

### **No Reaction or Weak Signal**

1. The sample load was insufficient.
  - Increase the amount of protein applied. Concentration of the sample prior to loading may be necessary. Use a more sensitive assay system (see Table 2.10).
2. Insufficient antigen binding to the membrane is occurring.
  - Stain the gel after transfer or use prestained or Kaleidoscope standards to assess transfer efficiency. See the previous section for suggestions on improving transfer related problems.
3. Antigen denaturation is occurring during electrophoresis or transfer. Antibodies, especially monoclonals, may not recognize denatured antigens.
  - Electrophorese and transfer proteins under native conditions. Use the Super Cooling Coil and a refrigerated recirculating bath to transfer heat-sensitive proteins.
4. Primary or secondary antibodies may be inactive or non-saturating.
  - Store the reagents at recommended conditions. Avoid repeated freeze-thaw cycles, bacterial contamination, or heat inactivation.
  - Detergents may affect the activity of some antibodies. Eliminate them from the assay, except for the wash after blocking.
  - If the antibody titer is too low, optimize the concentration using a dot-blot experiment.
  - Increase the antibody incubation times.
5. The enzyme conjugate is inactive or non-saturating.
  - Test the reagent for activity (see below).
  - Store the reagents at recommended conditions. Avoid repeated freeze-thaw cycles, bacterial contamination, or heat inactivation.
  - Sodium azide is a potent inhibitor of horseradish peroxidase. Use Thimerosal as a bacteriostat.
  - Impure water may cause inactivation of the enzyme. Use only distilled, deionized water.
  - If the conjugate concentration is too low, optimize using a dot-blot experiment.
6. Color development reagent is inactive.
  - Test the reagent for activity (see below) and remake if necessary.

### **Tests for Monitoring Reagent Activity**

1. Activity test for the color development solution.
  - Combine 1.0 ml of the color development solution with 10  $\mu$ l of full strength second antibody conjugate. The color reaction should develop immediately. If color fails to develop within a few minutes, the color development solution is inactive. Make up a fresh working solution and repeat the color development assay.
2. Activity test for the conjugate solution.
  - Combine 1.0 ml of the color development solution tested above and 1.0 ml of the 1:3,000 dilution conjugate solution. A light blue tinge should develop within 15 minutes. If color fails to develop within 25 minutes, the conjugate solution is suspect. Repeat the procedure with a freshly prepared dilution of conjugate.

3. Activity test for the first antibody solution.
  - Use an ELISA, RID, Ouchterlony immunodiffusion, or precipitation test to determine reactivity of the antibody with the antigen. If possible, repeat the assay procedure with several dilutions of first antibody solution.

## **Total Protein Detection**

### ***Colloidal Gold Total Protein Stain — High Background***

1. The blocking step is insufficient or omitted.
  - Block with 0.3% Tween-20 in TBS, using three washes of 20 minutes each.
2. The membrane used is not compatible with this stain.
  - Positively charged nylons cannot be used with Colloidal Gold stain. Use the Biotin-Blot Total Protein Detection Kit instead.
3. Contamination of the membrane occurred at a previous step, *i.e.* electrophoresis or transfer.
  - Discard and remake the gel and transfer solutions.
  - Replace or thoroughly clean contaminated fiber pads.
4. Excessive amounts of protein are loaded on the gel, or too much SDS is used in the transfer buffer. Proteins can pass through the membrane without binding and recirculate through a tank blotting system.
  - Reduce the amount of protein on the gel or SDS in the transfer buffer. Add a second sheet of membrane to bind excess protein.
5. Colloidal gold stain solution is contaminated.
  - The stain is a reusable reagent. Be sure to use a separate, clean plastic container to store previously used reagent in the refrigerator. Discard any reagent that has viscous sediment at the bottom of the bottle. If the solution does not have a dark burgundy color, but is a light blue, the stain was contaminated with buffer salts. Buffer salts will react with the gold sol causing non-specific precipitation of the reagent onto the membrane. Discard this solution.

### ***Colloidal Gold Total Protein Stain — Low Sensitivity***

1. Increase the incubation time for detection of low level signals.
  - Overnight incubations are possible, although background staining can increase.
2. Transfer is incomplete.
  - See poor transfer for suggestions on how to enhance transfer efficiency.
3. Stain is exhausted, as evidenced by the loss of the dark burgundy color and longer staining times.
  - Discard the reagent.
4. Buffer salt contamination has occurred. The solution will be light blue instead of dark burgundy.
  - Discard the reagent.
5. The sample load may be too low for the reagent to detect.
  - Use the Gold Enhancement Kit for detection levels as low as 10 pg of protein per band.

### ***Biotin-Blot Total Protein Detection — High Background***

1. Blocking conditions are insufficient.
  - Match the blocker to the membrane. Nylon membranes require the addition of 1-methyl-2-pyrrolidinone (MPO) to several solutions. Consult the Biotin-Blot manual for specific details.
2. Membrane is left in color development solution too long.
  - Remove the membrane from the color development solution when the signal is apparent and the background has not developed. Transfer the blot to distilled water immediately to stop the development.
3. Excessive amounts of protein are loaded on the gel, or too much SDS is used in the transfer buffer. Proteins can pass through the membrane without binding and recirculate through a tank blotting system.
  - Reduce the amount of protein on the gel or SDS in the transfer buffer. Add a second sheet of membrane to bind excess protein.

### ***Biotin-Blot Total Protein Detection — No Reaction or Weak Color Development***

1. Transfer is incomplete.
  - See Poor Transfer for suggestions on how to enhance transfer efficiency.
2. The sample load may be too low for the reagents to detect.
  - Increase the amount of protein loaded on the gel.
3. NHS-biotin solution is inactivated.
  - NHS-biotin hydrolyzes in aqueous solutions. Equilibrate the reagent vial to room temperature before opening to prevent condensation of water inside the container. Use a sterile syringe to remove reagent to prevent contamination.
  - Add the NHS-biotin reagent to the Borate-Tween solution just prior to use.
4. Amine containing buffer salts compete for the biotinylation reagents.
  - Wash the membrane thoroughly in Borate-Tween to remove any residual buffer salts from electrophoresis and transfer.
5. Avidin-HRP conjugate is inactive.
  - Follow the activity test procedures to determine if the reagent is inactive.
6. Color development solution is inactive.
  - Follow the activity test procedures to determine if the reagent is inactive.

### ***Anionic Dyes — High Background***

1. Destaining is insufficient.
  - Increase the number and duration of washes with the destaining solution.
2. Dye solution is too concentrated.
  - Remake the solution.
3. Nylon membranes are not compatible with anionic dyes.
  - Use the Biotin-Blot Protein Detection Kit.

### ***Anionic Dyes — Low Sensitivity***

1. Anionic dye stains do not detect protein bands below ~100 ng.
  - Use a more sensitive stain such as the Colloidal Gold stain or the Biotin-Blot Protein Detection Kit.
  - Increase the sample load to achieve the detection level of the anionic dye stains.



## Section 7 Product Information

Catalog Number	Product Description
170-3939	<b>Trans-Blot Cell with Plate Electrodes, complete</b> , includes 2 Gel Holder cassettes, Super Cooling Coil, cell with lid and power cables, and fiber pads
170-3946	<b>Trans-Blot Cell with Plate Electrodes</b> , includes Gel Holder cassettes, cell with lid and power cables, and fiber pads
170-3910	<b>Trans-Blot Cell with Wire Electrodes</b> , includes Gel Holder, cell with lid and power cables, and fiber pads

### PowerPac 200 Power Supply

165-5052	<b>PowerPac 200 Power Supply</b> , 110/120 V
165-5053	<b>PowerPac 200 Power Supply</b> , 220/240 V

### Trans-Blot Cell Accessories

170-3943	<b>Trans-Blot Platinum Anode Plate Electrode</b>
170-3944	<b>Trans-Blot Stainless Steel Cathode Plate Electrode</b>
170-3945	<b>Trans-Blot Plate Electrode Pair</b> , platinum anode and stainless steel cathode
170-3920	<b>Trans-Blot Standard Wire Electrode Card</b> , cathode
170-3921	<b>Trans-Blot Standard Wire Electrode Card</b> , anode
170-3912	<b>Super Cooling Coil</b> (required for all high intensity transfers)
170-3914	<b>Fiber Pads</b> , 15.5 cm x 20.5 cm x 6 mm, 6
170-3913	<b>Gel Holder</b>
170-3922	<b>Trans-Blot Cell Buffer Tank</b>
170-3923	<b>Trans-Blot Cell Lid with Cables</b>
170-3955	<b>Filter Paper</b> , 14 x 16 cm, 25
170-3956	<b>Filter Paper</b> , 15 x 20 cm, 25

## Section 8 References

1. Southern, E.M, *J Mol. Biol.*, **98**, 503 (1975).
2. Alwine, J. C., Kemp, D. J., Parker, B. A., Reiser, J., Renart j., Stark, G. R. and Wahl, G. W., *Methods Enzymol.*, **68**, 220 (1979).
3. Thomas, P. S., *Proc. Nat. Acad. Sci.*, **77**, 5201 (1980).
4. Seed, B., *Nuc. Acids Res.*, **10**, 1799 (1982).
5. Renart, J., Peiser, J. and Stark, G. R., *Proc. Nat. Acad. Sci.*, **76**, 3116 (1979).
6. Bowen, P., Steinberg, J., Laemmli, U. K. and Weintraub, H., *Nuc. Acids Res.*, **8**, 1 (1980).
7. Towbin, H., Staehelin, T. and Gordon, J., *Proc. Nat. Acad. Sci.*, **76**, 4350 (1970).
8. Bittner, M., Kupferer, P. and Morris, C. R., *Anal. Biochem.*, **102**, 459 (1980).

9. Stellwag, E. J. and Dahlberg, A. E., *Nuc. Acids Res.*, **8**, 299 (1980).
10. Kutateladze, T. V., Axelrod, B. D., Gorbulev, V. G., Belzhelarshaya, S. N. and Vartikyan, R. M., *Anal. Biochem.*, **100**, 129 (1979).
11. Peudelhuber, T. L., Ball, D. J., Davis, A. H. and Garrard, W. J., *Nuc. Acids Res.*, **10**, 1311 (1982).
12. Danner, D. B., *Anal. Biochem.*, **125**, 139 (1982).
13. Bio-Rad Technical Bulletin 1110 "Zeta-Probe Blotting Membranes" (1982).
14. Holland, L. J. and Wangh, L. H., *Nuc Acids Res.*, **10**, 3283 (1983).
15. Syminton, J., Green, M. and Brackmann, K., *Proc. Nat. Acad. Sci.*, **78**, 177 (1981).
16. Reiser, J. and Wardale, J., *Eur. J. Biochem.*, **114**, 569 (1981).
17. Burnette, W. N., *Anal. Biochem.*, **112**, 195 (1981).
18. Legocki, R. P. and Verma, D. P. S., *Anal. Biochem.*, **111**, 385 (1981).
19. Lin, W. and Kasamatsu, H., *Anal. Biochem.*, **128**, 302 (1983).
20. Anderson, N. L., Nance, S. L., Pearson, T. W. and Anderson, N.G., *Electrophoresis*, **3**, 135( 1982).
21. McLellan, T. and Pamshaw, J. A. M., *Biochem. Genetics*, **19**, 647 (1981).
22. Gibson, W., *Anal. Biochem.*, **118**, 1 (1981).
23. Howe, J. G. and Hershey, J. W. B., *J. Biol. Chem.*, **2566**, 12836 (1981).
24. Erickson, P. G., Minier, L. N. and Lasher, P. S., *J. Immun. Meth.*, **51**, 241 (1982).
25. Tsang, V. C. W., Peralta, J. M. and Simons, A. R., *Meth. Enzymol.*, **92**, 377 (1983).
26. Gershoni, J. M. and Palade, G. E., *Anal. Biochem.*, **124**, 396 (1982).
27. Gershoni, J. M. and Palade, G. E., *Anal. Biochem.*, **131**, 1 (1983).
28. Symington, J., "Two Dimensional Gel Electrophoresis of Proteins: Methods and Applications." Celis, J. E. and Bravo, R., eds. Academic Press, N.Y., (1983).
29. Andrews, A. T., "Electrophoresis: Theory, techniques, and biochemical and clinical application," 2nd ed., Clarendon Press, Oxford, (1986).
30. Beisiegel, V., *Electrophoresis*, **7**, 1 (1986).
31. Bio-Rad Laboratories, unpublished.
32. Gershoni, J. M., in *Advances in Electrophoresis*, Vol. 1. Chrambach, A., Dunn, M. J. and Radola, B. J., eds., VCH, Weinheim, in press.
33. Gershoni, J. M. , in *Methods of Biochemical Analysis*, Vol. 33, Glick, D., ed., Wiley, New York, in press.
34. Bjerrum, O. J. and Schafer-Nielsen, C., *Analytical Electrophoresis*, M. J. Dunn, ed. (VCH, Weinheim), p. 315.
35. Dunn, S. D., *Anal. Biochem.*, **157**, 144 (1986).
36. Zeta-Probe Instruction Manual, Bio-Rad Laboratories, (1988).
37. Polvino, W. J., Saravis, C. A., Sampson, C. E. and Cook, R. B., *Electrophoresis*, **4**, 368 (1983).
39. Bio-Rad Laboratories, Biotin-Blot Total Protein Stain Instruction Manual (1985).
40. LaRochelle, W. J. and Froehner, S. C., *J. Immunol. Meth.*, **92**, 65 (1986).
41. Szewczyk, B. and Kozloff, L. M., *Anal. Biochem.*, **150**, 403 (1985).
42. Perides, G., Plagens, U. and Traub, P., *Anal. Biochem.*, **152**, 94 (1986).

Scotch-Brite® is a registered trademark of 3M Company.

Gel-Bond™ is a trademark of FMC.

Mylar® is a registered trademark of E.I. DuPont de Nemours Co.

Coomassie is a trademark of ICI.



**Bio-Rad  
Laboratories**

**Life Science  
Group**

2000 Alfred Nobel Drive  
Hercules, California 94547  
Telephone (510) 741-1000  
Fax: (510) 741-5800

**Australia**, Bio-Rad Laboratories Pty Limited, Block Y Unit 1, Regents Park Industrial Estate, 391 Park Road, Regents Park, NSW 2143 • Phone 02-9414-2800 • Fax 02-9914-2888  
**Austria**, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, 1130 Wien • Phone (1) 877 89 01 • Fax (1) 876 56 29  
**Belgium**, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, 9810 Nazareth Eke • Phone 09-385 55 11 • Fax 09-385 65 54  
**Canada**, Bio-Rad Laboratories (Canada) Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone (905) 712-2771 • Fax (905) 712-2990  
**China**, Bio-Rad Laboratories, 14, Zhi Chun Road, Hai Dian District, Beijing 100088 • Phone (01) 2046622 • Fax (01) 2051876  
**Denmark**, Bio-Rad Laboratories, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 39 17 9947 • Fax 39 27 1698  
**Finland**, Bio-Rad Laboratories, Business Center Länsikeskus, Pihatörmä 1A SF-02240, Espoo, • Phone 90 804 2200 • Fax 90 804 1100  
**France**, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94 203 Ivry Sur Seine Cedex • Phone (1) 49 60 68 34 • Fax (1) 46 71 24 67  
**Germany**, Bio-Rad Laboratories GmbH, Heidemannstraße 164, D-80939 München/Postfach 450133, D-80901 München • Phone 089 31884-0 • Fax 089 31884-100  
**India**, Bio-Rad Laboratories, C-248 Defence Colony, New Delhi 110 024 • Phone 91-11-461-0103 • Fax 91-11-461-0765  
**Italy**, Bio-Rad Laboratories S.r.l., Via Cellini, 18/A, 20090 Segrate Milano • Phone 02-21609 1 • Fax 02-21609-399  
**Japan**, Nippon Bio-Rad Laboratories, 7-18, Higashi-Nippori 5-Chome, Arakawa-ku, Tokyo 116 • Phone 03-5811-6270 • Fax 03-5811-6272  
**The Netherlands**, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 0318-540666 • Fax 0318-542216  
**New Zealand**, Bio-Rad Laboratories Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097  
**Pacific**, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257  
**Singapore**, Bio-Rad Laboratories (Singapore) Ltd., 221 Henderson Rd #05-19, Henderson Building, Singapore 0315 • Phone (65) 272-9877 • Fax (65) 273-4835  
**Spain**, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98  
**Sweden**, Bio-Rad Laboratories AB, Gärdsvägen 7D, Box 1276, S-171 24 Solna • Phone 46-(0)8-735 83 00 • Fax 46-(0)8-735 54 60  
**Switzerland**, Bio-Rad Laboratories AG, Kanalstrasse 17, Postfach, CH-8152 Glattbrugg • Phone 01-809 55 55 • Fax 01-809 55 00  
**United Kingdom**, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 01442 259118