



**AmpLight™ Chemiluminescent  
Western Detection Kit**

**Catalog Number  
170-8234**

**AmpLight Fluorescent  
Western Detection Kit**

**Catalog Number  
170-8232**

**Instruction Manual**

For Technical Service  
Call Your Local Bio-Rad Office or  
in the U.S. Call **1-800-4BIORAD**  
**(1-800-424-6723)**



# Table of Contents

<b>Section 1</b>	<b>Preparation.....</b>	<b>1</b>
1.1	Introduction.....	1
1.2	Method Overview.....	2
1.3	Kit Components.....	2
1.4	Product Storage and Stability.....	3
1.5	Safety Instructions.....	4
<b>Section 2</b>	<b>Protocol.....</b>	<b>4</b>
2.1	Experimental Strategy and General Considerations.....	4
2.2	Reagent Preparation.....	7
2.3	Quick Guide.....	8
2.4	Detailed Protocol.....	10
<b>Section 3</b>	<b>Alternative Protocol .....</b>	<b>15</b>
<b>Section 4</b>	<b>Troubleshooting Guide.....</b>	<b>17</b>
<b>Section 5</b>	<b>Product Information.....</b>	<b>24</b>

# Section 1 Preparation

## 1.1 Introduction

The AmpLight chemiluminescent and fluorescent western detection kits have been designed for instrument-based detection of western blots on nitrocellulose membranes. The kits are optimized for use following application of an antibody conjugated to horseradish peroxidase (HRP), but may also be used following incubation in a biotinylated antibody (see Section 3). The technique employs a proprietary reaction system which amplifies the detection signal. Detection with the ChemiDetect™ chemiluminescent substrate can be documented by CCD-based systems (*e.g.*, Fluor-S™ MultiImager system, Gel Doc™ 1000 video gel documentation system, Insta Doc™ system). In addition, the chemiluminescent detection kit significantly improves detection with the Molecular Imager® light-sensitive phosphor screens. The fluorescent moiety of the FluorDetect™ reagent, fluorescein, may be excited by a broadband ultraviolet signal (centered around 300 nm) or by an argon laser (488 nm); maximum emission is at 520 nm. The kits provide reagents for amplification and detection on 2,500 cm<sup>2</sup> of membrane.

## 1.2 Method Overview

The first step in western blotting is the transfer of antigen onto a solid support membrane by one of several methods. The transfer can be done electrophoretically, following separation of the antigen in a polyacrylamide or agarose gel, passively by directly spotting the antigen onto a membrane, or by vacuum filtration using a microfiltration apparatus. Following antigen binding, the remaining protein binding sites on the membrane surface are blocked with the provided blocking reagent.

The membrane with bound protein is then incubated with a primary antibody specific to the antigen of interest. The blot is washed to remove unbound antibody and incubated with a secondary antibody linked to HRP. The blot is washed to remove unbound secondary antibody and then incubated in the Bio-Rad amplification reagent (BAR). The blot is washed to remove excess BAR, incubated in streptavidin-HRP, and then washed again. At this point the blot is ready for detection with either the ChemiDetect or FluorDetect substrates.

## 1.3 Kit Components

### Reagents provided with the kit

Blocking reagent, 20 grams

10x phosphate buffered saline (PBS), powder to make 1 liter

4x BAR, 53 ml

2x amplification diluent, 105 ml (chemiluminescent kit), 210 ml (fluorescent kit)

Streptavidin-HRP, 0.5 ml

ChemiDetect substrate, 160 ml (chemiluminescent kit)

ChemiDetect oxidizer, 160 ml (chemiluminescent kit)

4x FluorDetect reagent, 53 ml (fluorescent kit)

### Additional required items not provided in the kit

Nitrocellulose membrane

Primary antibody

Secondary antibody conjugated with horseradish peroxidase

Bovine serum albumin, 7.5 g, sufficient for 2,500 cm<sup>2</sup>

Tween 20, 10 ml, sufficient for 2,500 cm<sup>2</sup>

Dimethyl sulfoxide (DMSO), 500 ml, sufficient for 2,500 cm<sup>2</sup>

## 1.4 Product Storage and Stability

The kit is shipped at 4 °C. Store the unopened kit at 4 °C. Powdered blocker, powdered PBS and PBS solutions may be stored at room temperature. After being put into solution, the blocker should be stored at 4 °C. All kit components are guaranteed for 1 year from the time of receipt.

## 1.5 Safety Instructions

Read the entire instruction manual before beginning the protocol.

1. Wear gloves and protective clothing, such as laboratory coats and goggles, when preparing and working with the solutions in the protocol. DMSO is an irritant; it is a colorless liquid which is easily absorbed through the skin and mucous membranes. Avoid skin contact with DMSO and inhalation of DMSO mist. Wash exposed skin thoroughly with soap and water.

**Note:** See Material Data Safety Sheet (MSDS) on DMSO for additional information.

2. Work in well-ventilated areas. Avoid inhalation of vapors when working with solutions containing DMSO.
3. Do not mouth pipet any solution.

## Section 2 Protocol

### 2.1 Experimental Strategy and General Considerations

**Temperature.** All steps are performed at room temperature (22–25 °C).

**Making solutions.** Use only deionized, distilled water to prepare solutions. 1x PBST solutions should be sterile filtered. Do not use azide as a preservative in any solution.

**Membrane selection.** This kit has been optimized for detection on pure nitrocellulose membranes and performs equally well with supported nitrocellulose membranes. However, due to the way in which they are manufactured, the surface of some supported nitrocellulose membranes takes on a ‘wavy’ appearance that can result in less pleasing images than the pure nitrocellulose membranes which remain flat throughout the process. The kit may also be used to detect proteins bound to PVDF.

**Primary Antibody.** Generally when serum or tissue culture supernatants are the source of primary antibody, a 1:100–1:1,000 dilution of the primary antibody in buffer is used for detection of antigens on the membrane surface. For chromatographically purified monospecific antibodies, a 1:500–1:10,000 dilution is used for antigen detection. A 1:1,000–1:100,000 dilution is used when ascites fluid is the source of antibody. Optimal dilution factors must be determined experimentally. The optimal antibody concentration is usually considered the greatest dilution of antibody reagent still resulting in a strong positive signal without membrane background or non-specific reactions.

**Secondary Antibody Conjugates.** The protocols in this manual were worked out using Bio-Rad Blotting Grade secondary antibody conjugates diluted as described below. Using an antibody conjugate at a higher concentration may result in an overall increase in background without any improvement in detection sensitivity.

Secondary antibody conjugates from other sources may be used, but the optimal dilution may be different from that of Bio-Rad antibody conjugates.

**Washes and Incubations.** Continuous gentle agitation should be used during all incubations and washes. For best results, an orbital shaker should be employed to maintain a uniform exposure of the membrane surface to the solution. Use the smallest possible container to hold the membrane and solutions. When a range of washes is specified, as in 'Wash 2–4x', best results are obtained by doing the maximum number of washes. Acceptable results, though potentially with more background, may be obtained with the minimum number of washes. Blocking and washing steps should be done with 0.25 ml per cm<sup>2</sup> of membrane, *e.g.*, for a mini-blot of 60 cm<sup>2</sup>, use 15 ml of solution for those steps. Antibody incubations should be at 0.1 ml per cm<sup>2</sup>, *e.g.*, 6 ml for a 60 cm<sup>2</sup> miniblot. Use dilute BAR and streptavidin-HRP solutions at 85 µl per cm<sup>2</sup> of membrane, *e.g.*, 5 ml for a 60 cm<sup>2</sup> mini-blot.

**Detergents.** Tween-20 is essential in washing to eliminate overall background and non-specific hydrophobic interactions. At 0.1%, Tween-20 will not disrupt binding of primary antibodies to antigens or antigens to the membrane, but will optimize detection sensitivity by eliminating non-specific interactions. Alternative detergents should not be substituted.

## 2.2 Reagent Preparation

The following reagents should be made upon first receiving the kit.

**10x PBST (Phosphate buffered saline/1% Tween-20).** Pour the contents of the pouch into 950 ml ddH<sub>2</sub>O and stir until dissolved. Add 10 ml Tween-20. Bring final volume to 1 liter with ddH<sub>2</sub>O. Store at room temperature.

**1x PBST.** (Phosphate buffered saline/0.1% Tween-20). Combine 100 ml 10x PBST and 900 ml ddH<sub>2</sub>O. Sterile filter before use. Store at room temperature.

**20% DMSO/PBST Wash.** Combine 100 ml DMSO and 400 ml 1x PBST. This is sufficient for 500 cm<sup>2</sup> of membrane.

**3% Blocker.** While vigorously stirring 665 ml of PBST, very slowly add 20 g powdered blocker. Add the powder a little at a time over a period of 30–45 minutes. Continue stirring another 30–60 minutes after all the powder has been added. Slowly warming the solution to 55 °C will help put the blocker into solution, but do not overheat. This can be accomplished by placing the beaker with the blocker solution inside another beaker on top of the magnetic stirrer. Pour enough room temperature water into the outside container to surround the blocker solution as it stirs. Be careful not to cause the beaker with the blocker to float or tip. Slowly heat the surrounding water to 55 – 60 °C. Do not add azide. Store at 4 °C. Warm to room temperature before use.

Make these solutions on the day of the experiment.

**Antibody dilution buffer (1% BSA in PBST).**

Dissolve 0.75 g of BSA in 75 ml rapidly stirring PBST (sufficient for 4 miniblots). Used for dilution of primary and secondary antibodies and for dilution of streptavidin-HRP.

**Bio-Rad Amplification Reagent.** Prepare 85  $\mu$ l per  $\text{cm}^2$  of membrane. Combine 2 parts 2x amplification diluent, 1 part 4x BAR and 1 part  $\text{ddH}_2\text{O}$ .

**Streptavidin-HRP.** Prepare 85  $\mu$ l per  $\text{cm}^2$  of membrane. Dilute 1:1000 with antibody dilution buffer.

**FluorDetect Substrate.** Prepare 85  $\mu$ l per  $\text{cm}^2$  of membrane. Combine 2 parts 2x amplification diluent, 1 part 4x FluorDetect, and 1 part  $\text{ddH}_2\text{O}$ .

## 2.3 Quick Guide

### Blotting and Amplification

1. Transfer protein to nitrocellulose membrane by electrophoretic blotting, dot blotting, or microfiltration. Allow membrane to air dry.
2. Wet membrane in PBST and then wash 2x for 5 minutes each time in PBST.
3. Block membrane in 3% blocker for 1 hour.
4. Wash 2x with PBST for 3–5 minutes.
5. Incubate in appropriately diluted primary antibody for 1 hour.
6. Wash 2x with PBST 5 minutes each time.

7. Incubate in 1:3,000 – 1:10,000 dilution of GAx -HRP\* secondary antibody for 1 hour.
8. Wash 2x with PBST 5 minutes each time.
9. Incubate membrane in diluted BAR for 10 minutes.
10. Wash 2–4x in 20% DMSO/PBST for 5 minutes each time.
11. Wash 1–2x in PBST for 5 minutes each time.
12. Incubate membrane in diluted streptavidin-HRP for 30 minutes.
13. Wash 2x in PBST for 5 minutes each time.

### Chemiluminescent Detection

1. Mix equal volumes of the ChemiDetect oxidizer and the ChemiDetect substrate. Prepare 0.1 ml per  $\text{cm}^2$  of membrane. Incubate membrane in substrate for 30–60 seconds.
2. Remove from substrate. Place membrane between plastic sheets and image immediately.

### Fluorescent Detection

1. Incubate membrane in diluted FluorDetect substrate for 10 minutes.
2. Wash 2–4x with 20% DMSO/PBST for 5 minutes each time.
3. Wash 1–2x with PBST for 5 minutes each time.
4. Capture image.

\* GAx -HRP is Bio-Rad's Goat Anti-Rabbit IgG-HRP (Catalog number 170-6515) or Goat Anti-Mouse IgG-HRP (170-6516) or Goat Anti-Human IgG-HRP (172-1050).

## 2.4 Detailed Protocol

**Note:** Before beginning, read through the entire procedure.

1. **Antigen application.** Apply antigen to the membrane surface using one of the three basic methods described below. A small amount of known antigen or primary antibody dotted on one corner of the membrane prior to blocking will produce a positive reaction if the procedure is successful.
  - a. **Electrophoretic blotting.** The antigens of interest are electrophoretically transferred to the membrane from a gel (*i.e.* SDS-PAGE, IEF, or native gel) using the Trans-Blot®, Mini Trans-Blot®, Trans-Blot Semi-Dry cell or similar device.
  - b. **Microfiltration blotting.** The antigens of interest are transferred by a vacuum device such as the Bio-Dot® or Bio-Dot SF onto the membrane. The membrane should be removed from the apparatus for the blocking and all subsequent steps.
  - c. **Dot blotting.** Cut the membrane sheet to the appropriate size. Draw a grid on the membrane with a pencil. Wet the dry membrane by slowly sliding the membrane at a 45° angle into the PBST. (PVDF membranes must first be wet in 100% methanol; consult membrane instructions for complete information). Remove the thoroughly wetted membrane from the PBST and dry it on filter paper for approximately 5 minutes. Apply

antigen sample to each grid square using a syringe or pipet.

**Note:** Regardless of the method chosen for antigen application, the membrane should be allowed to dry completely before proceeding to the next step.

2. **Wash.** Wet the membrane in PBST and then wash the membrane twice in PBST for 5 minutes each time. These washes help to reduce spotted or blotchy background.
3. **Blocking step.** Immerse the membrane at a 45° angle into the blocking solution. Gently agitate the solution using an orbital shaker platform and incubate for an hour or more.
4. **Wash.** Decant the blocking solution and add PBST to the membrane. Wash for 5 minutes. Repeat the wash with fresh PBST.
5. **Primary antibody incubation.** Decant the PBST and add 0.1 ml of antibody solution per cm<sup>2</sup> of membrane. Dilute the primary antibody in PBST with 1% (w/v) BSA. Incubate 1 to 2 hours with gentle agitation. The optimum conditions of dilution and incubation must be determined experimentally.
6. **Wash.** Decant the antibody solution and add PBST to the membrane. Wash for 5 minutes with gentle agitation and pour off the wash solution. Repeat the wash with fresh PBST.



7. **Secondary antibody incubation.** Decant the PBST and add 0.1 ml of the secondary antibody solution per  $\text{cm}^2$  of membrane. Dilute the secondary antibody 1:3,000 – 1:10,000 with PBST containing 1% (w/v) BSA. Incubate for 30 minutes to 2 hours with gentle agitation. As noted previously, this protocol was developed using Bio-Rad's blotting grade secondary antibodies (Goat Anti-Rabbit IgG-HRP [Catalog number 170-6515] or Goat Anti-Mouse IgG-HRP [170-6516] or Goat Anti-Human IgG-HRP [172-1050]). Secondary antibody HRP conjugates from other sources may be used, but the optimal dilution and incubation time will have to be determined experimentally.
8. **Wash.** Decant the antibody solution and add PBST to the membrane. Wash for 5 minutes with gentle agitation and pour off the wash solution. Repeat the wash with fresh PBST.
9. **Amplification.** Prepare 85  $\mu\text{l}$  of 1x BAR solution per  $\text{cm}^2$  of membrane. Prepare the solution by combining 2 parts 2x amplification diluent, 1 part 4x BAR and 1 part  $\text{ddH}_2\text{O}$ . Pour PBST off the membrane and add the 1x BAR solution. Incubate for 10 minutes with gentle agitation.
10. **Wash.** Pour off the BAR solution and wash the membrane 2–4 times for 5 minutes each time with 20% DMSO/PBST. Follow these washes with 1–2 washes in PBST for 5 minutes each time.

11. **Streptavidin-HRP.** Prepare 85  $\mu\text{l}$  of a 1:1,000 dilution streptavidin-HRP solution per  $\text{cm}^2$  of membrane. Dilute the streptavidin-HRP in the same 1% BSA/PBST solution used to dilute the antibodies. Pour off the PBST wash and add the diluted streptavidin-HRP solution. Incubate for 30 minutes with gentle agitation.
12. **Wash.** Pour off the streptavidin-HRP solution and add PBST. Wash the membrane for 5 minutes with gentle agitation. Repeat the wash step with fresh PBST.

### Chemiluminescent Detection

1. Prepare 0.125 ml of substrate solution per  $\text{cm}^2$  of membrane by mixing equal volumes of the ChemiDetect substrate and the ChemiDetect oxidizer. Pour off the PBST and add the prepared substrate solution. Incubate for 30–60 seconds. Do not incubate any longer than 60 seconds.
2. Quickly pour off the ChemiDetect solution and place the membrane between two sheets of plastic or wrap it in plastic film.
3. Photography, digitization or exposure to phosphor screens must be initiated immediately after removing the membrane from the substrate solution. For CCD-based or film-based detection, open apertures completely, remove all filters from the lenses and limit exposure time so that the detector is not saturated. A 1- or 2-minute exposure on a CCD-based instrument like the Fluor-S MultiImager is usually sufficient.

Leave the blots on a Molecular Imager CH screen for 5–10 minutes.

### Fluorescent Detection

1. Prepare 0.1 ml of 1x FluorDetect solution per cm<sup>2</sup> of membrane by mixing 2 parts 2x amplification diluent, 1 part 4x FluorDetect reagent and 1 part ddH<sub>2</sub>O. Pour off the PBST and add the diluted FluorDetect solution to the membrane. Incubate for 10 minutes with gentle agitation.
2. Wash. Pour off the FluorDetect solution and wash the membrane 2–4 times for 5 minutes each time with 20% DMSO/PBST. Follow these washes with 1–2 washes in PBST for 5 minutes each time.
3. The fluorescein moiety of the FluorDetect substrate is best detected when the membrane is illuminated from above (epi-illuminated) by either broadband ultraviolet light centered around 300 nm or by a 488 nm laser. Maximum emission is around 520 nm. Strong signals may also be excited by illumination from below the membrane (trans-illumination), but there is a loss of sensitivity when compared to epi-illumination. The signal will persist as long as the excitation source is on. The fluorescein will photobleach after long exposure to the excitation source, *e.g.*, more than 30 minutes cumulative exposure. With the Fluor-S MultiImager system, good sensitivity may be obtained in a 1 to 2 minute exposure using the 530DF60 filter.

## Section 3 Alternative Protocol

An alternative, and potentially more sensitive amplification, may be accomplished by substituting a biotinylated secondary antibody for the HRP-linked secondary antibody. There is an additional incubation in streptavidin-HRP that precedes incubation in the BAR solution. The kit has sufficient streptavidin-HRP to carry out detection over 2,500 cm<sup>2</sup> of membrane, even if you use a biotinylated secondary antibody.

This approach entails two rounds of amplification, so background can be a significant problem and it may not be trivial to work out the best conditions. Background is affected by the dilution of primary antibody, secondary antibody, the BAR and the streptavidin-HRP. It may be necessary to further dilute one or more of these four reagents.

A brief protocol is outlined below.

### Blotting and Amplification

1. Transfer protein to nitrocellulose membrane by electrophoretic blotting, dot blotting or microfiltration. Allow membrane to air dry.
2. Wet membrane in PBST and then wash 2x for 5 minutes each time in PBST.
3. Block membrane in 3% blocker for 1 hour.
4. Wash 2x with PBST for 3–5 minutes.

5. Incubate in appropriately diluted primary antibody for 1 hour.
6. Wash 2x with PBST for 5 minutes each time.
7. Incubate in a dilution of biotinylated goat anti-rabbit IgG secondary antibody (catalog number 170-6401) for 1 hour.
8. Wash 2x with PBST for 5 minutes each time.
9. Incubate in 1:1,000 dilution of streptavidin-HRP for 30 minutes.
10. Wash 2x in PBST for 5 minutes each time.
11. Incubate membrane in diluted BAR for 10 minutes.
12. Wash 2–4x in 20% DMSO/PBST for 5 minutes each time.
13. Wash 1–2x in PBST for 5 minutes each time.
14. Incubate membrane in diluted streptavidin-HRP for 30 minutes.
15. Wash 2x in PBST for 5 minutes each time.

### Chemiluminescent Detection

1. Mix equal volumes of the two ChemiDetect reagents. Prepare 0.125 ml per cm<sup>2</sup> of membrane. Incubate membrane in substrate for 30–60 seconds.
2. Remove from substrate. Place membrane between plastic sheets and image immediately.

### Fluorescent Detection

1. Incubate membrane in diluted FluorDetect for 10 minutes.
2. Wash 2–4x with 20% DMSO/PBST for 5 minutes each time.
3. Wash 1–2x with PBST for 5 minutes each time.
4. Capture image.

## Section 4 Troubleshooting Guide

<b>Problem</b>	<b>Probable Cause</b>	<b>Recommended Solution</b>
1. No reaction or weak signal.	a. Exposure time was too short.	i. Increase the exposure time.
	b. Blot was allowed to dry after incubation with the chemiluminescent substrate.	i. Use heat sealable bags to prevent drying of the membrane.
	c. Chemiluminescent substrate solution is inactive.	i. Store the reagent at the proper temperature, 4 °C.

## Troubleshooting Guide (continued)

<b>Problem</b>	<b>Probable Cause</b>	<b>Recommended Solution</b>
	d. Primary antibody solution is inactive or non-saturating.	<p>i. Store the antibody solution at the proper temperature. Avoid bacterial contamination, heat inactivation, and repeated freeze-thaw cycles.</p> <p>ii. Antibody titer was too low. Increase the concentration of the antibody used in the assay.</p> <p>iii. Tween-20 may affect the reactivity of some antibodies. Eliminate Tween-20 from the assay (except the wash after the blocking step).</p>

## Troubleshooting Guide (continued)

<b>Problem</b>	<b>Probable Cause</b>	<b>Recommended Solution</b>
	e. Conjugate is inactive.	<p>i. Store the conjugate at the proper temperature. Avoid repeated freeze-thaw cycles.</p> <p>ii. The concentration of the conjugate was non-saturating. Increase the concentration of the conjugate used in the assay.</p> <p>iii. Conjugate may be contaminated, causing inactivation of the antibody or enzyme. Tap water may cause inactivation; use only distilled, deionized water to prepare all solutions.</p>

## Troubleshooting Guide (continued)

<b>Problem</b>	<b>Probable Cause</b>	<b>Recommended Solution</b>
	f. Little or no antigen is bound to the membrane.	<p>i. Tween-20 may wash bound antigen from the membrane. Eliminate Tween-20 from the assay (except the wash after the blocking step).</p> <p>ii. Transfer of protein onto the membrane was incomplete. Stain gel to assure transfer of protein. Use Prestained Standards to monitor transfer efficiency. Consult the appropriate instrument manual for proper procedures and recommendations.</p>

## Troubleshooting Guide (continued)

<b>Problem</b>	<b>Probable Cause</b>	<b>Recommended Solution</b>
	g. Primary antibody is not specific or does not recognize denatured antigens (common with monoclonals).	i. Loss of reactivity may have occurred during electrophoretic transfer. Pre-test the reactivity of the antibody against the antigen by a dot blot.
2. High background.	<p>a. Exposure time was too long.</p> <p>b. PBST washes after transfer were omitted or insufficient.</p> <p>c. Blocking was insufficient.</p>	<p>i. Decrease exposure time.</p> <p>i. The washes are critical to reduce spotted or blotchy background development.</p> <p>i. Increase the time of the blocking step and/or the concentration of blocker used.</p>

## Troubleshooting Guide (continued)

Problem	Probable Cause	Recommended Solution
	d. Wash stringency was insufficient.	<ul style="list-style-type: none"> <li>i. Tween-20 is necessary in wash steps to reduce background. The concentration can be increased up to 0.3% if background persists.</li> <li>ii. Increase the number and length of washes.</li> </ul>
	e. Second antibody conjugate was used at an excessive concentration.	<ul style="list-style-type: none"> <li>i. Use the recommended dilution, or determine the optimal dilution experimentally.</li> </ul>
	f. Contamination occurred during transfer.	<ul style="list-style-type: none"> <li>i. Refer to the instrument instruction manual for recommendations.</li> </ul>
	g. Insufficient number of DMSO/PBST washes.	<ul style="list-style-type: none"> <li>i. Increase number of work steps following BAR or Fluor Detect.</li> </ul>

## Troubleshooting Guide (continued)

Problem	Probable Cause	Recommended Solution
	h. BAR concentration was too high.	<ul style="list-style-type: none"> <li>i. Dilute BAR further.</li> </ul>
	i. Streptavidin-HRP concentration was too high	<ul style="list-style-type: none"> <li>ii. Dilute Streptavidin-HRP further.</li> </ul>

## Section 5 Product Information

Catalog Number	Product Description
----------------	---------------------

### AmpLight Western Detection Kits

170-8232	AmpLight Fluorescent Western Detection Kit
170-8234	AmpLight Chemiluminescent Western Detection Kit

### Blotting Grade Conjugates

170-6515	Goat Anti-Rabbit IgG-HRP
170-6516	Goat Anti-Mouse IgG-HRP
172-1050	Goat Anti-Human IgG-HRP
170-6401	Biotinylated Goat Anti-Rabbit IgG

### Pure Nitrocellulose Membranes (0.45 micron)

162-0145	Sheets, 7 x 8.4 cm, 10
162-0117	Sheets, 9 x 12 cm, 10
162-0114	Sheets, 15 x 9.2 cm, 10
162-0116	Sheets, 15 x 15, 10
162-0113	Sheets, 20 x 20 cm, 5
162-0115	Roll, 33 cm x 3 m, 1

### Pure Nitrocellulose Membranes (0.2 micron)

162-0146	Sheets, 7 x 8.4 cm, 10
162-0147	Sheets, 13.5 x 16.5 cm, 10
162-0150	Sheets, 20 x 20 cm, 5
162-0112	Roll, 33 cm x 3 m, 1

Catalog Number	Product Description
----------------	---------------------

### Supported Nitrocellulose Membranes (0.45 micron)

162-0090	Sheets, 7 x 8.4 cm, 10
162-0091	Sheets, 10 x 15 cm, 10
162-0092	Sheets, 15 x 15, 10
162-0093	Sheets, 20 x 20 cm, 10
162-0094	Roll, 33 cm x 3 m, 1

### Supported Nitrocellulose Membranes (0.2 micron)

162-0095	Sheets, 7 x 8.4 cm, 10
162-0096	Sheets, 15 x 15 cm, 10
162-0097	Roll, 33 cm x 3 m, 1

### PVDF Membranes (0.2 micron)

162-0186	Sheets, 7 x 8.4 cm, 10
162-0180	Sheets, 10 x 15 cm, 10
162-0181	Sheets, 15 x 15, 10
162-0182	Sheets, 20 x 20 cm, 10
162-0185	Sheets, 20 x 20 cm, 3
162-0184	Roll, 24 cm x 3.3 m, 1

### Miscellaneous

170-6531	Tween-20, 100 ml
----------	------------------