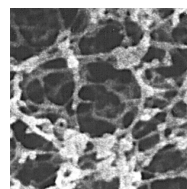
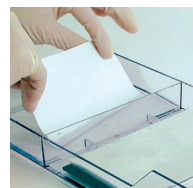


Electrophoresis



Blotting Membranes



Simplify Western Blot Preparation

BIO-RAD

Supports That Mean Results

To meet multiple blotting requirements and experimental designs, Bio-Rad offers a convenient and comprehensive line of blotting membranes, including different grades of nitrocellulose, PVDF, and Zeta-Probe® nylon membranes. All blotting membranes are sold in rolls and precut sheets to fit Bio-Rad's line of blotting equipment, including the Trans-Blot® cell, the Trans-Blot® Plus cell, the Trans-Blot SD semi-dry cell, the Criterion™ blotter, and the Mini Trans-Blot® cell, as well as the Bio-Dot® SF (slot format) apparatus and the Mini-PROTEAN® II multiscreen apparatus.

Bio-Rad also offers a line of precut membrane sandwiches (nitrocellulose or PVDF) assembled for your convenience. Two sheets of thick filter paper and a blotting membrane are preassembled into a blotting membrane/filter paper sandwich. These gel-sized membrane sandwiches save time and simplify western blot preparation.



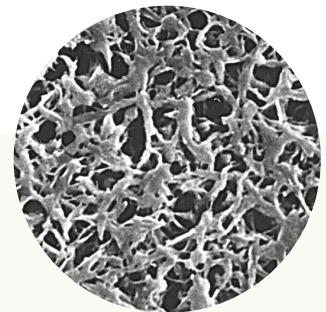
Bio-Rad's blotting membranes are available in rolls, precut sheets, or convenient membrane/filter paper sandwiches.

For more information, contact your local Bio-Rad representative, visit us on the Web at discover.bio-rad.com, or call toll free at 1-800-4BIORAD (1-800-424-6723).

PVDF Membranes

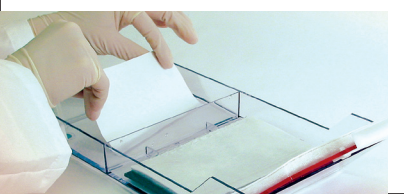
Ideal for:

- Protein blotting (Immun-Blot® PVDF membrane)
- Protein sequencing (Sequi-Blot™ PVDF membrane)



4,000x

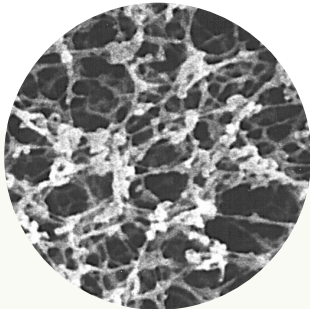
- PVDF blotting membrane is a naturally hydrophobic support ideal for western blotting, protein sequencing, and protein binding assays
- Mechanical strength of PVDF membranes allows multiple reprobing of blots
- Immun-Blot PVDF membrane for chemiluminescent or colorimetric western blots ensures high signal with low background
- Sequi-Blot PVDF membrane offers the highest protein binding capacity and is chemically resistant to solvents for protein sequencing
- PVDF membranes are available in convenient sandwich formats, precut sheets, and rolls



Nitrocellulose Membranes

Ideal for:

Protein blotting
DNA/RNA blotting
Dot/slot blot screening
Colony/plaque screening



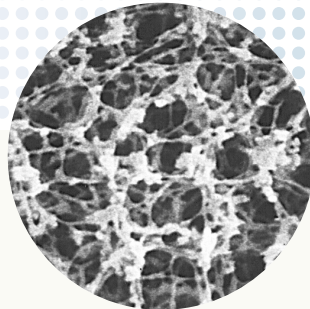
4,000x

- Pure 100% nitrocellulose is the most frequently used membrane support for the transfer of both proteins and nucleic acids
- Nitrocellulose is compatible with all standard detection methods and exhibits superior binding capacity without background interference
- Two pore sizes are available — 0.45 μm for most analytical protein transfers and for DNA and RNA >500 bp, and 0.2 μm for the transfer of low molecular weight proteins and nucleic acids
- Membranes are available in convenient sandwich formats, precut sheets, and rolls

Supported Nitrocellulose Membranes

Ideal for:

Protein blotting
DNA/RNA blotting
Dot/slot blot screening



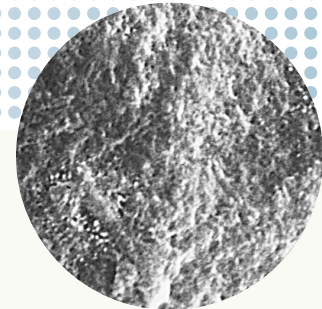
4,000x

- Supported nitrocellulose membranes combine sensitivity with strength
- Pure nitrocellulose is cast on both sides of an inert polyester support material, allowing easier handling during transfers
- Reinforced membrane is ideal for repeated screening and reprobing, and also can be autoclaved
- Two pore sizes are available — 0.45 μm for most analytical protein transfers and for DNA and RNA >500 bp, and 0.2 μm for the transfer of low molecular weight proteins and nucleic acids
- Supported nitrocellulose membranes are available in a variety of precut sheets and rolls

Zeta-Probe Membranes

Ideal for:

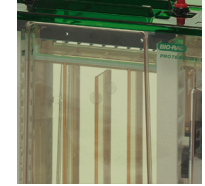
DNA/RNA blotting
Dot/slot blot screening
Colony/plaque screening



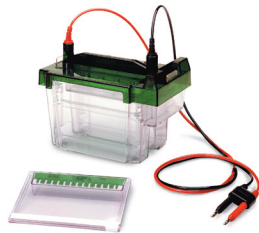
4,000x

- Zeta-Probe and Zeta-Probe GT (genomic DNA-tested) membranes are high-strength nylon membranes that are positively charged with quaternary amine groups
- Ideal for nucleic acid blotting, membranes are resistant to shrinking and tearing during transfer, hybridization, or reprobing
- Membranes can be used in alkaline Southern and northern blotting protocols and in traditional Southern blot transfers
- Only Zeta-Probe GT membranes are pretested for the detection of single-copy sequence in a genomic blot*
- Membranes are available in rolls and precut sheets to fit many agarose gel sizes, including ReadyAgarose™ precast gels

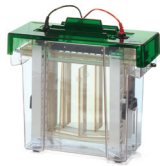
* Each lot is tested to ensure that 3 pg of target single-copy factor VIII human DNA can be detected in 5 μg of total genomic DNA using a radiolabeled probe. A lot-specific result certificate is included.



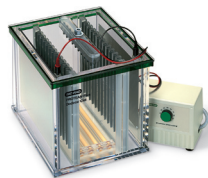
Mini-PROTEAN® Tetra Gel Sizes (W x L)
8.3 x 7.3 cm
8.6 x 6.8 cm



Criterion Gel Size (W x L)
13.3 x 8.7 cm



PROTEAN® II xi/XL Gel Sizes (W x L)
16 x 16 cm
16 x 20 cm
18.5 x 20 cm



PROTEAN® Plus Gel Sizes (W x L)
18.3 x 19.3 cm
19.8 x 17.5 cm
20 x 20.5 cm
25 x 20.5 cm
25.6 x 23 cm

Ordering Information

Catalog #	Description
Immun-Blot PVDF Membranes	
162-0174	8.4 x 7 cm, 10 sheets
162-0175	10 x 15 cm, 10 sheets
162-0176	20 x 20 cm, 10 sheets
162-0177	26 cm x 3.3 m, 1 roll
162-0255	26.5 x 28 cm, 10 sheets

Catalog #	Description
Immun-Blot PVDF Membrane/Filter Paper Sandwiches	
162-0238	13.5 x 8.5 cm, 20 sheets
162-0239	13.5 x 8.5 cm, 50 sheets
162-0218	8.5 x 7 cm, 20 sheets
162-0219	8.5 x 7 cm, 50 sheets

Catalog #	Description
Sequi-Blot PVDF Membranes	
162-0186	8.4 x 7 cm, 10 sheets
162-0180	15 x 10 cm, 10 sheets
162-0181	15 x 15 cm, 10 sheets
162-0182	20 x 20 cm, 10 sheets
162-0185	20 x 20 cm, 3 sheets
162-0184	26 cm x 3.3 m, 1 roll
162-0256	26.5 x 28 cm, 10 sheets

Catalog #	Description
Sequi-Blot PVDF Membrane/Filter Paper Sandwiches	
162-0236	13.5 x 8.5 cm, 20 sheets
162-0237	13.5 x 8.5 cm, 50 sheets
162-0216	8.5 x 7 cm, 20 sheets
162-0217	8.5 x 7 cm, 50 sheets

Catalog #	Description
Nitrocellulose Membranes (0.45 µm)	
162-0113	20 x 20 cm, 5 sheets
162-0114	15 x 9.2 cm, 10 sheets
162-0115	33 cm x 3 m, 1 roll
162-0116	15 x 15 cm, 10 sheets
162-0117	9 x 12 cm, 10 sheets
162-0145	8.4 x 7 cm, 10 sheets
162-0148	11.5 x 16 cm, 10 sheets
162-0251	26.5 x 28 cm, 10 sheets

Catalog #	Description
Nitrocellulose/Filter Paper Sandwiches (0.45 µm)	
162-0234	13.5 x 8.5 cm, 20 sheets
162-0235	13.5 x 8.5 cm, 50 sheets
162-0214	8.5 x 7 cm, 20 sheets
162-0215	8.5 x 7 cm, 50 sheets

Catalog #	Description
Nitrocellulose Membranes (0.2 µm)	
162-0112	33 cm x 3 m, 1 roll
162-0146	8.4 x 7 cm, 10 sheets
162-0147	16.5 x 13.5 cm, 10 sheets
162-0150	20 x 20 cm, 5 sheets
162-0252	26.5 x 28 cm, 10 sheets

Catalog #	Description
Nitrocellulose/Filter Paper Sandwiches (0.2 µm)	
162-0232	13.5 x 8.5 cm, 20 sheets
162-0233	13.5 x 8.5 cm, 50 sheets
162-0212	8.5 x 7 cm, 20 sheets
162-0213	8.5 x 7 cm, 50 sheets

Catalog #	Description
Supported Nitrocellulose Membranes (0.45 µm)	
162-0090	8.4 x 7 cm, 10 sheets
162-0091	15 x 10 cm, 10 sheets
162-0092	15 x 15 cm, 10 sheets
162-0093	20 x 20 cm, 10 sheets
162-0094	30 cm x 3 m, 1 roll
162-0254	26.5 x 28 cm, 10 sheets

Catalog #	Description
Supported Nitrocellulose Membranes (0.2 µm)	
162-0095	8.4 x 7 cm, 10 sheets
162-0096	15 x 15 cm, 10 sheets
162-0097	30 cm x 3 m, 1 roll
162-0253	26.5 x 28 cm, 10 sheets

Catalog #	Description
Zeta-Probe Membranes	
162-0153	9 x 12 cm, 15 sheets
162-0154	15 x 10 cm, 15 sheets
162-0155	15 x 15 cm, 15 sheets
162-0156	15 x 20 cm, 15 sheets
162-0157	20 x 20 cm, 15 sheets
162-0158	25 x 20 cm, 3 sheets
162-0159	30 cm x 3.3 m, 1 roll
162-0165	20 cm x 3.3 m, 1 roll
162-0206	7 x 10 cm, 15 sheets

Catalog #	Description
Zeta-Probe GT Membranes	
162-0190	9 x 12 cm, 15 sheets
162-0191	15 x 10 cm, 15 sheets
162-0192	15 x 15 cm, 15 sheets
162-0193	20 x 15 cm, 15 sheets
162-0194	20 x 20 cm, 15 sheets
162-0195	25 x 20 cm, 3 sheets
162-0196	30 cm x 3.3 m, 1 roll
162-0197	20 cm x 3.3 m, 1 roll
162-0208	7 x 10 cm, 15 sheets

Catalog #	Description
C/P Lift® Membranes	
162-0162	C/P Lift Membrane Disks, 85 mm, 50
162-0163	C/P Lift Membrane Disks, 137 mm, 15
170-3202	Nitrocellulose Membrane Disks, 82.5 mm, 50



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000