

## BLOTTING

# Model 785 Vacuum Blotter

- Fast, accurate transfer of DNA or RNA
- Quick, easy setup
- Controlled vacuum pressure for consistent results
- Molded construction for long-term reliability
- Chemically resistant construction

## Fast Vacuum Transfer System

The Model 785 vacuum blotter is a nucleic acid transfer apparatus that uses low vacuum to transfer DNA or RNA from agarose gels onto nylon membranes in as little as 90 minutes. This vacuum blotter simplifies every step of the blotting process. Its easy assembly reduces setup times for your experiments, while the vacuum regulator accurately controls vacuum pressure to give you consistent results. The Model 785 vacuum blotter has durable molded construction for long-term, heavy use. Detailed protocols for single-copy gene detection are included in the simple, step-by-step instruction manual. The Model 785 vacuum blotter saves valuable experimental time without compromising results.

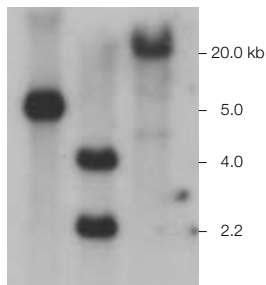
The Model 785 vacuum blotter has several innovative features that help to simplify the nucleic acid blotting process and facilitate complete transfer, including:

- Spring latches make assembly and disassembly quick and easy — no screws or nuts are needed to tighten the sealing frame
- Vacuum stage and porous vacuum plate create a uniform vacuum pressure over the entire transfer area, for even transfers and highly reproducible results
- Blank window gaskets are available for transferring differently sized gels
- Easy-to-see green window gasket, combined with simple base unit and sealing frame, make aligning the gel and membrane quick and easy
- Vacuum regulator accurately controls vacuum pressure, whether you use a Bio-Rad vacuum pump or a house vacuum
- Durable, molded construction withstands long-term, heavy use
- Blotter is chemically resistant to commonly used acids and bases (0.25 N HCl or 0.5 N NaOH)
- Tested protocols for overnight-exposure single-copy gene detection are included in the Model 785 vacuum blotter instruction manual

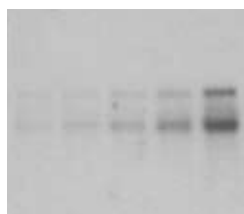


**BIO-RAD**

# Model 785 Vacuum Blotter



**Southern transfer.** Human genomic DNA (5 µg) was cleaved with (left to right) *Bgl*I, *Hind*III, or *Nhe*I, separated on a 1% Certified™ molecular biology agarose gel, and transferred to a Zeta-Probe® GT membrane using a vacuum of 5" Hg for 90 min. The blot was probed with a <sup>32</sup>P-labeled factor VIII fragment and exposed to X-ray film overnight. Probe provided by the Genetics Institute of Boston.



**Northern transfer.** Left to right, 1.0, 2.5, 5, 10, or 20 µg of total RNA from CHO HA-1 cells was separated on a glyoxal gel and transferred to a Zeta-Probe® membrane using a vacuum of 3" Hg for 90 min. The blot was probed with a <sup>32</sup>P-labeled hsp70 cDNA fragment and exposed to X-ray film overnight. Probe and RNA provided by Gloria Lee, University of California, San Francisco, CA USA.

## Specifications

### Dimensions

Overall size (L x W x H)	21 x 12 x 5"
Pump station (L x W x H)	16 x 9 x 8"
Vacuum tubing	1/2" OD x 1/4" ID

### Weight

Basic unit	8.5 lb
Vacuum station	9.5 lb
Vacuum regulator	1.5 lb

### Ratings

Vacuum pump	20" Hg
Vacuum regulator	0–25" Hg
Gel capacity (maximum)	20 x 25 cm

### Required Equipment

Vacuum source	Vacuum pump or house vacuum
Vacuum regulator	Strongly recommended for reproducible results
Liquid trap	Sidearm flask with one-hole stopper (not sold by Bio-Rad Laboratories)
Vacuum tubing	Required for liquid trap assembly

## Ordering Information

Catalog #	Description
165-5000	<b>Model 785 Vacuum Blotter With Regulator</b> includes vacuum regulator, base with vacuum stage, porous vacuum plate, reservoir seal O-ring, sealing frame, assorted window gaskets, lid, instructions
165-5001	<b>Model 785 Vacuum Blotter System, 100/120 V</b> , includes vacuum pump, vacuum regulator, base with vacuum stage, porous vacuum plate, reservoir seal O-ring, sealing frame, assorted window gaskets, lid, instructions
165-5002	<b>Model 785 Vacuum Blotter System, 220/240 V</b>
165-5003	<b>Model 785 Vacuum Blotter Basic Unit</b> , includes base with vacuum stage, porous vacuum plate, reservoir seal O-ring, sealing frame, assorted window gaskets, lid, instructions
165-5004	<b>Vacuum Station 100/120 V</b> , includes vacuum pump, vacuum regulator
165-5005	<b>Vacuum Station, 220/240 V</b>
165-5006	<b>Vacuum Regulator</b>
165-5007	<b>Reservoir Seal O-Ring</b>
165-5008	<b>Blank Window Gaskets, 6</b>
165-5012	<b>Porous Vacuum Plate</b>

### Ultra Pure DNA Agarose

162-0125	<b>High-Strength Analytical Grade Agarose, 100 g</b>
162-0135	<b>Chromosomal Grade Agarose, 25 g</b>
162-0136	<b>Chromosomal Grade Agarose, 100 g</b>

### Blotting Accessories

162-0153	<b>Zeta-Probe Membranes, 9 x 12 cm, 15 sheets</b>
162-0154	<b>Zeta-Probe Membranes, 10 x 15 cm, 15 sheets</b>
162-0155	<b>Zeta-Probe Membranes, 15 x 15 cm, 15 sheets</b>
162-0156	<b>Zeta-Probe Membranes, 15 x 20 cm, 15 sheets</b>
162-0157	<b>Zeta-Probe Membranes, 20 x 20 cm, 15 sheets</b>
162-0158	<b>Zeta-Probe Membranes, 20 x 25 cm, 3 sheets</b>
162-0159	<b>Zeta-Probe Membrane, 30 cm x 3.3 m, 1 roll</b>
162-0165	<b>Zeta-Probe Membrane, 20 cm x 3.3 m, 1 roll</b>
162-0190	<b>Zeta-Probe GT Membranes, 9 x 12 cm, 15 sheets</b>
162-0191	<b>Zeta-Probe GT Membranes, 10 x 15 cm, 15 sheets</b>
162-0192	<b>Zeta-Probe GT Membranes, 15 x 15 cm, 15 sheets</b>
162-0193	<b>Zeta-Probe GT Membranes, 15 x 20 cm, 15 sheets</b>
162-0194	<b>Zeta-Probe GT Membranes, 20 x 20 cm, 15 sheets</b>
162-0195	<b>Zeta-Probe GT Membranes, 20 x 25 cm, 3 sheets</b>
162-0196	<b>Zeta-Probe GT Membrane, 30 cm x 3.3 m, 1 roll</b>
162-0197	<b>Zeta-Probe GT Membrane, 20 cm x 3.3 m, 1 roll</b>
161-0302	<b>SDS (Sodium Dodecyl Sulfate), 1 kg</b>
165-0962	<b>Filter Paper Backing, for stained gels, 35 x 45 cm, 25 sheets</b>
165-0921	<b>Thick Blot Paper, 18 x 34 cm, for Model 224, 443, and 543 slab gel dryers, 25 sheets</b>

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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