

Photo Gel Documentation System



The Economical and Compact Film Camera System

The Photo Gel Documentation System provides the conveniences and advantages of the Polaroid® DS-34 fixed, focal-length camera with the compact Mini-Transilluminator. Two different sizes of electrophoresis hoods provide a dark room environment, allowing photography right at the lab bench. The dual intensity setting on the Mini-Transilluminator allows selection for preparative or analytical work, and the UV blocking shield protects the user from exposure to UV rays. The built-in trigger shutter release on the DS-34 camera is easy to use, and the variable shutter speeds and aperture settings allow the user to set the proper exposure for specific documentation. The system comes complete with filter for use with ethidium bromide stained gels, Gel cutter, fluorescent ruler, and a pack of B&W film.

The Photo Gel Documentation System includes

- Mini-Transilluminator
- Darkroom Electrophoresis Hood
- Deep Yellow Filter
- Polaroid DS-34 Camera
- Gel Cutter
- Fluorescent Ruler
- Polaroid 667 B&W Film Pack

Catalog Number	Product Description
170-3742	Standard Documentation System, 120 VAC, includes Mini-Transilluminator, 120 V, Polaroid DS-34 camera, Standard Electrophoresis Hood, Deep Yellow DS-34 camera filter, Polaroid 667 B&W Film Pack, Gel cutter, Fluorescent Ruler
170-3746	Standard Documentation System, 100 VAC
170-3747	Standard Documentation System, 220/240 VAC
170-3743	Wide/Long Documentation System, 120 VAC, includes Mini-Transilluminator, 120 V, Polaroid DS-34 camera, wide/long electrophoresis hood, Deep Yellow DS-34 camera filter, Polaroid 667 B&W Film Pack, Gel cutter, Fluorescent Ruler
170-3748	Wide/Long Documentation System, 100 VAC
170-3749	Wide/Long Documentation System, 220 /240VAC
170-3737	Mini-Transilluminator, 120 VAC
170-3745	Mini-Transilluminator, 100 VAC
170-3738	Mini-Transilluminator, 220/240 VAC
170-3739	Standard Electrophoresis Hood, 5.1" x 7" (13 x 17.8 cm)
170-3740	Wide/Long Electrophoresis Hood, 6" x 9" (15.2 x 22.9 cm)
170-3741	Polaroid DS-34 Camera
170-3744	Deep Yellow DS-34 Camera Filter

Polaroid is a trademark of Polaroid Corp.

Photo Gel Documentation System Specifications

Mini-Transilluminator

Width	13.25"	33.7 cm
Depth	9.50"	24.1 cm
Height	4.75"	12.1 cm
Wavelength	302 nm	
Number of lamps	6 x 8 W	
Intensity	Dual (Analytical and Preparative)	
Filter size	7.9 x 7.9"	20 x 20 cm
Volt/Hz/Amp	120/60/.72	220-240/50/.62
	100/50-60/.85	
Fluorescent Ruler Length	10.25"	25.5 cm

Electrophoresis Hoods

	Standard	Wide/Long
Inside Hood Dimensions	5.1 x 7" 13 x 17.8 cm	5-7/8 x 8-3/4" 14.9 x 22.2 cm
Image Magnification Percentage	0.55 x	0.45 x

Polaroid DS-34 Camera

Lens	105 mm f/4.5
Aperture	4/4.5-4/32
Shutter	Copal Press #0, 1/125-1 sec., plus B
Size	7.25 x 7.25 x 4.75" 18.4 x 18.4 x 12.1 cm
Film size	3.25 x 4.25" 8.3 x 10.8 cm

Deep Yellow Filter

Filter Number	15
Color	Deep yellow
Application	Ethidium Bromide stained gels with UV light



**Bio-Rad
Laboratories**

Life Science Group

2000 Alfred Nobel Drive
Hercules, California 94547
Telephone (510) 741-1000
Fax: (510) 741-1060

Eastern Regional Office, 85A Marcus Dr., Melville, New York 11747 • Phone (516) 756-2575 • Fax (516) 756-2594

European Headquarters, Bio-Rad Laboratories, Dreve du Senechal, 19, B-1180 Brussels • Phone 02 375 59 70 • Fax 02 374 61 62

Australia, Bio-Rad Laboratories Pty Limited, Unit 11, 112-118 Talavera Rd P.O. Box 371, North Ryde, N.S.W. 2113 • Phone 02-805-5000 • Fax 02-805-1920

Austria, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, A-1130 Wien • Phone 0222-877 89 01 • Fax 0222-876 56 29

Belgium, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, B-9810 Nazareth Eke • Phone 091-85 55 11 • Fax 091-85 65 54

Canada, Bio-Rad Laboratories (Canada) Ltd., 5149 Bradco Boulevard, Mississauga, Ontario L4W 2A6 • Phone (416) 624-0713 • Fax (416) 624-3019

China, Bio-Rad Laboratories, Yanshan Hotel Office Tower, #1307, A138 Haidian Road, Beijing • Phone 2563146 • Fax 2564308

France, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94203 Ivry Sur Seine Cedex • Phone 01-49 60 68 34 • Fax 01-46 71 24 67

Germany, Bio-Rad Laboratories GmbH, Heidemannstraße 164, Postfach 45 01 33, D-8000 München 45 • Phone 089-318 84-0 • Fax 089-318 84 100

Italy, Bio-Rad Laboratories S.r.l., Via Cellini, 18A, 20090 Segrate Milano • Phone 02-21609.1 • Fax 02-21609-399

Japan, Nippon Bio-Rad Laboratories, K. K., Sumitomo Seimei Kachidoki Bldg 5-3-6 Kachidoki, Chuo-Ku, Tokyo 104 • Phone 03-3534-7515 • Fax 03-3534-8027

The Netherlands, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 08385-40666 • Fax 08385-42216

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

Scandinavia, Bio-Rad Laboratories, Kanalvägen 10C, 19461 Upplands Väsby, Sweden • Phone 46 (0) 8 590-73489 • Fax 46 (0) 8 590-71781

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

Switzerland, Bio-Rad Laboratories AG, Kanalstrasse, 17, 8152 Glattbrugg • Phone 01-810 16 77 • Fax 01-810 19 33

United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Phone 0800 181134 • Fax 0442 259118

Printed in USA