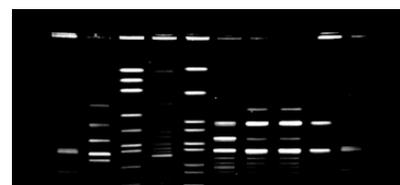


# Pulsed Field Gel Electrophoresis



## Pulsed Field Gel Electrophoresis



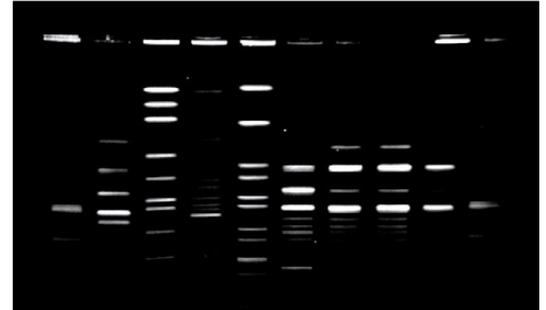
Map It Out



# The Direct Route to Your Large Molecule Applications

## PFGE Is Used in Many Research Areas

Pulsed field gel electrophoresis (PFGE) has enabled progress in cancer research, food safety, public health, quality control, and genome mapping. It is widely used in molecular epidemiology for strain typing and it has been adopted by PulseNet, a global network of health and food regulatory agency laboratories coordinated by the Centers for Disease Control and Prevention (CDC).



**Electrophoretic profile of *Salmonella enteritidis*.**  
Courtesy of Kara L. Cooper, Centers for Disease Control and Prevention, Atlanta, Georgia.

### Molecular Epidemiology

#### Strain Typing in Public Health and Food Safety

PFGE is used for epidemiological studies of pathogenic organisms such as *Escherichia coli* O157:H7, *Salmonella*, *Shigella*, *Listeria*, *Campylobacter*, or *Vibrio cholerae*. When epidemiologists need to precisely identify the strain variants from a sample, genetic fingerprinting is the method of choice. Rare cutting restriction enzymes yield large DNA fragments which are analyzed using the CHEF system; the variant-specific electrophoretic gel pattern is then compared to the PulseNet database. The PulseNet database, which contains thousands of patterns, is used by member laboratories to identify and track foodborne infections worldwide.

#### Food Quality Control

PFGE is in widespread use as a quality control method in the food industry. For example, the beer and wine industries use it to monitor the genetic stability of organisms in fermentation processes.

### Cancer Research

#### DNA Damage and Repair Studies

Research efforts using PFGE are focused on better understanding the factors mediating the damage to DNA caused by ionizing radiation and chemical treatment. It is important to quantitatively measure the dsDNA breaks due to these treatments. DNA from treated cells is subjected to PFGE and the density of the DNA in different molecular weight regions indicates the integrity of DNA and the extent of its repair.

#### Apoptosis Assays

Apoptotic DNA fragmentation is a key characteristic of programmed cell death. Analysis of the fragmentation that

occurs in the apoptosis process demonstrates either a "ladder" pattern at ~200 bp intervals (200–600 bp) or the formation of larger fragments (50–300 kb). Both of these size ranges can be visualized on one gel using field inversion gel electrophoresis (FIGE), available on the CHEF Mapper® XA system, enabling easier assessment of the fragmentation process.

### Genomics Applications

#### Generation of Artificial Chromosome Libraries

Cloning large DNA (100 kb–1 Mb) is the first step in sequencing complex genomes. FIGE is often used to separate and isolate the large digested DNA fragments, which are then cloned into artificial chromosomes to generate yeast, bacterial, human, and mammalian libraries.

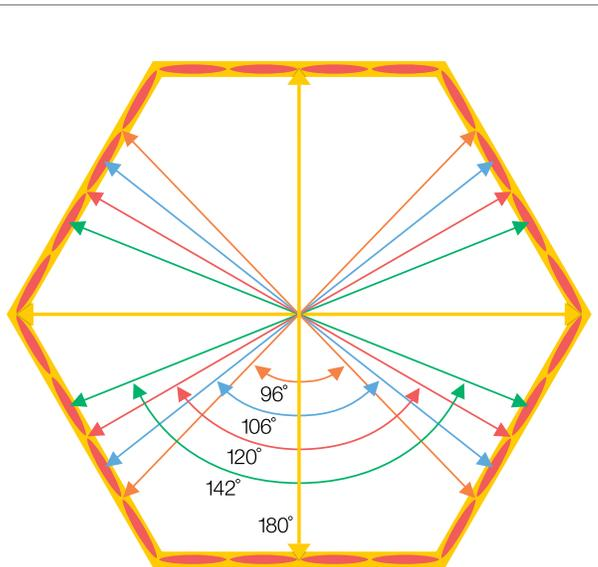
#### Genome Mapping

PFGE is still the benchmark for mapping applications. The libraries constructed using FIGE can also be used for mapping applications and specific assays used in research areas, including mapping specific disease loci, identifying chromosome rearrangements, and RFLP and DNA fingerprinting.

#### DNase I Hypersensitivity Assay

PFGE is used in mapping the genome for DNase I hypersensitive sites, which involves identifying different types of regulatory domains, such as active promoters and enhancers, where DNA-binding proteins are bound within nuclear chromatin. Conformation of chromatin at these sites causes them to be sensitive to DNase I cleavage. These sites are often located near active genes and play a role in eukaryotic gene regulation.

# The Leader in PFGE Technology



**Multistate switching capability.** Use this feature of the CHEF Mapper XA system to select vectors to dramatically speed up your separations and improve resolution.

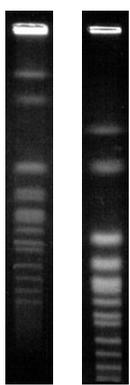
Bio-Rad is the leader in PFGE, offering the exclusive clamped homogenous electrical field (CHEF) technology (Chu et al. 1986) for pulsed field electrophoresis that has been used in mapping genomes since 1988.

Conventional electrophoresis can effectively separate fragments up to 20 kb; DNA fragments larger than 20 kb in a sample comigrate and when imaged appear as a large fuzzy band at the top of the gel. In 1984, Schwartz and Cantor invented PFGE to overcome this problem. PFGE resolves DNA by alternating the electrical field between spatially distinct pairs of electrodes, causing DNA molecules as large as several megabases to reorient and move at different speeds through the pores in an agarose gel.

## Technologies Used in Bio-Rad's PFGE Systems

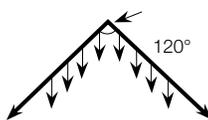
To achieve straight runs and good resolution in PFGE it is necessary to create homogenous electrical fields. There have been multiple approaches to PFGE but the combination of CHEF, PACE, and DR technologies used in Bio-Rad PFGE systems works best in creating the homogenous electrical fields that ensure consistency and run-to-run reproducibility.

- CHEF (clamped homogeneous electrical field) technology resolves DNA over a wide range of molecular weights in a straight lane; it employs the principles of contour-clamped electrophoresis to generate homogenous electrical fields
- PACE (programmable autonomously controlled electrodes) technology allows users to select the angle of electrophoretic pulsing optimal for the desired size range
- DR (dynamic regulation) is the electronics design by which each of the 24 electrodes is regulated; CHEF-DR® systems are capable of compensating for changes in buffer conductivity or gel size, preventing these changes from affecting the reproducibility of results
- FIGE (field inversion gel electrophoresis) is used for rapid sample resolution in the 100 bp–250 kb size range; in FIGE the electrical field is fixed at 1 angle (180°) and is inverted in the forward and reverse directions
- AFIGE (asymmetric field inversion gel electrophoresis) is a further refinement of the FIGE technology; AFIGE applies a different voltage to the forward direction electrical field than to the reverse direction electrical field, which optimizes the sample resolution in the FIGE size range



Multistate mode, 20 hr run,  
120° included angle  
60–120 sec switch-time ramp  
6 V/cm (200 V), 0.5x TBE, 14°C  
1.0% Certified™ molecular  
biology agarose

Secondary pulses  
6 V/cm (200 V), 0° angle  
3 sec switch time  
4 pulses/min



**Increased separation with secondary pulsed field electrophoresis.** *Saccharomyces cerevisiae* chromosomes separated under two-state conditions (left) and under two-state conditions with secondary pulses (right).

# Complete Solutions for PFGE

## Molecular Imager® Gel Doc™ XR+ System

The Molecular Imager Gel Doc XR+ instrument is an easy-to-use gel documentation system that documents and analyzes fluorescent gels, and produces publication-quality output at a fraction of the cost of film. It combines a compact darkroom, UV transilluminator workstation, high-resolution CCD camera, and powerful, user-friendly software for unsurpassed flexibility.

## FPQuest™ and InfoQuest™ FP Software

FPQuest and InfoQuestFP modular software packages offer customizable applications to meet a variety of laboratory informatics requirements.

- FPQuest software offers advanced analysis and statistical tools for analyzing banding patterns and multiple fingerprints in gels
- InfoQuestFP software includes all the functionality of FPQuest software, with the ability to analyze many other data types for more comprehensive studies of biological relationships

## CHEF Genomic DNA Plug Kits

Three DNA plug kits are available for preparing bacterial (lysozyme-sensitive) or mammalian genomic DNA and yeast chromosomes (YACs). Each kit contains sufficient enzymes, reaction buffers, and restriction digest-qualified CleanCut™ agarose for 100 plugs. Disposable molds and screened caps simplify plug preparation.



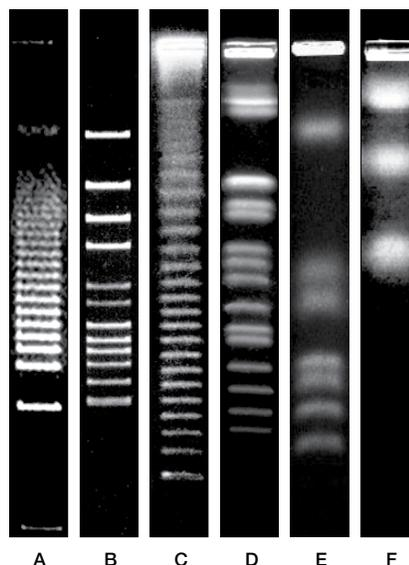
## Pulsed Field Quality Agarose

The type of agarose and the amount (percentage) used in an experiment play a crucial role in achieving optimal fragment resolution. Bio-Rad's Certified agaroses for PFGE are 100% pure and GQT grade (genetic quality tested).

- Pulsed field Certified agarose has an optimal separation range of 1 kb–2 Mb; its running conditions are a preset selectable method of the CHEF Mapper XA system auto-algorithm
- Certified megabase agarose has an optimal separation range of 1 kb–5 Mb and has high gel strength, a high exclusion limit, and high electrophoretic mobility; gels are easy to handle, even at 0.3%, allowing shorter run times

## Standards and Markers

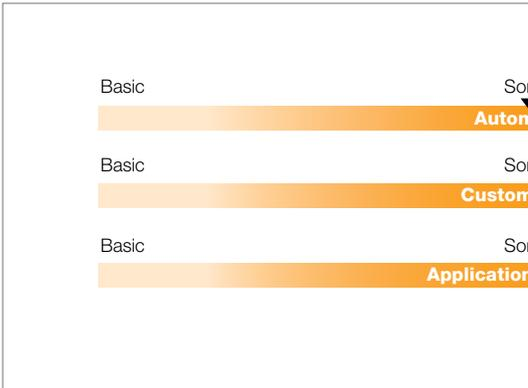
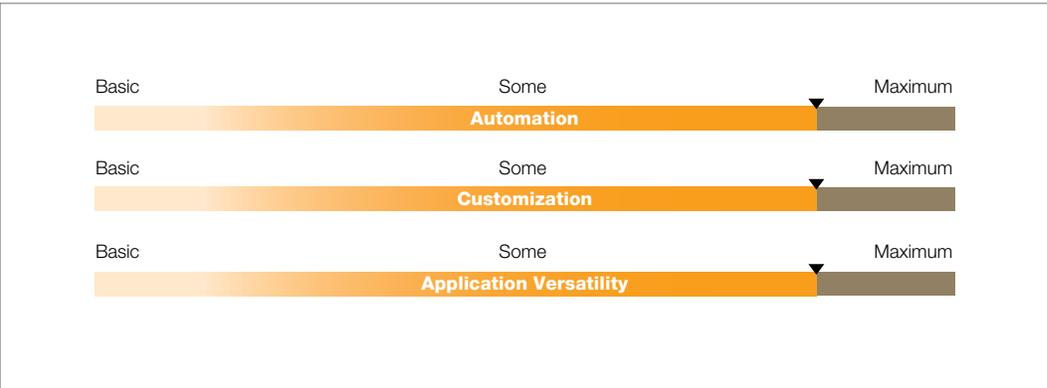
Bio-Rad offers standards for all PFGE applications, from FIGE separation of cosmid inserts to the largest chromosomal separations. Higher molecular weight standards are prepared in low-melt agarose blocks that can be cut to fit any well dimensions.



A, 5 kb ladder standard;  
B, 8.3–48.5 kb standard;  
C, lambda ladder standard;  
D, *S. cerevisiae* marker;  
E, *H. wingei* marker;  
F, *S. pombe* marker.



# Choose the System You Need for Your Specific



## CHEF Mapper XA System

The CHEF Mapper XA system is the ultimate tool, ideal for all PFGE applications. It offers multistate, secondary pulse, and a combination of CHEF, PACE, DR, FIGE, and AFIGE technologies, making optimal resolution in all size ranges possible.

- Users can achieve optimal resolution of both megabase- and kilobase-sized DNA fragments in 1 lane by selecting any pulse angle (0–360°) and applying asymmetrical angles
- Secondary pulses can be applied to release DNA caught in the gel matrix to further enhance the separation and resolution of very large DNA molecules
- FIGE and AFIGE functions enable enhanced and rapid resolution of small fragments (100 bp–250 kb)

The system is ideal for both the PFGE novice and the expert because it offers two ways to optimize DNA separations:

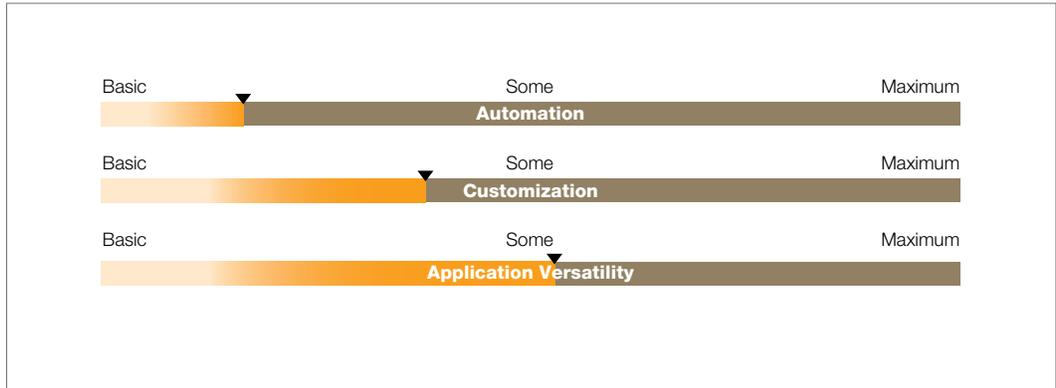
- Built-in auto-algorithm automatically selects and executes optimal separation conditions from only the entered fragment sizes
- Protocols can be refined using the Windows-based interactive algorithm, which allows users to simultaneously specify several run variables to derive optimal separation protocols
- System can store 99 simple programs or 20 complex programs with up to 8 blocks of programming each

## CHEF-DR III Variable Angle System

The CHEF-DR III system is optimized for a 100 bp–10 Mb range using CHEF and PA

- Users enter run conditions, and can optimize fragment resolution by selecting optimal voltage gradient, switch time, and pulse angle (90–120°) for the specific DNA size range
- Users can vary run conditions to obtain accurate size estimations; fine resolution in a complicated digest can be achieved with minimal programming and sample preparation

# Application



rapid separations of DNA in the PFGE technologies.

- System will recall the last conditions and use them as the default protocol; it also has a battery-operated backup RAM to recall current run conditions and resume the run without user intervention in the event of a power failure

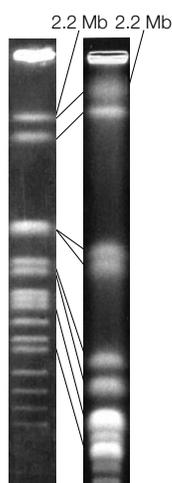
## CHEF-DR II System

The CHEF-DR II system effectively and reliably resolves DNA fragments in the 5 kb–6 Mb range by adjusting the running conditions for low voltage and extended run times.

- System is easy to program and cost effective, allowing enhanced resolution by executing two blocks of running conditions successively; users can input optimal run time, voltage gradient, and switch intervals for each run
- System employs the most common pulse angle for PFGE, 120°; the instrument manual provides examples of run conditions for a variety of size separation ranges for easy startup



Feature	CHEF Mapper XA System	CHEF-DR III Variable Angle System	CHEF-DR II System
<b>Fragment size</b>	100 bp–10 Mb	100 bp–10 Mb	5 kb–6 Mb
<b>Optimal separation size range</b>	100 bp–10 Mb	100 bp–10 Mb	100 kb–2 Mb
<b>Auto-algorithm and interactive algorithm:</b> algorithmic derivation of optimal run conditions	Yes	No	No
<b>Program storage:</b> storage and easy access of run conditions	20 complex programs	Last program run	No
<b>Programming blocks of run conditions:</b> optimized separation of fragments	8 blocks	3 blocks	2 blocks
<b>Battery-operated backup RAM:</b> recalls current run conditions and run progress in the event of power failure	Yes	Yes	No
<b>Pulse angle:</b> selection of different pulse angles optimizes resolution of both chromosomal and plasmid DNA with one system	Any angle from 0 to 360°	Any angle from 90 to 120° (in 1° increments)	Fixed angle of 120°
<b>Asymmetrical angles:</b> further optimizes separation of both chromosomal and plasmid DNA with one system; necessary for difficult samples	Yes	No	No
<b>Nonlinear switch-time ramping:</b> expands linear range of fragment separation to 50–700 kb, thus providing accurate fragment size measurements	Yes	No	No
<b>Multistate separation:</b> optimizes separation of subsets of fragments for enhanced resolution in selected fragment size ranges and faster separation	Yes	No	No
<b>Secondary pulses (voltage interrupts):</b> releases large DNA caught in the gel matrix and enhances separation and resolution of very large DNA molecules	Yes	No	No
<b>FIGE and AFIGE:</b> superior resolution of small fragments in 100 bp–250 kb range	Yes	No	No
<b>Recommended use</b>	<ul style="list-style-type: none"> <li>• Ideal for all PFGE applications</li> <li>• Best resolution in all size ranges</li> <li>• Most accurate results</li> <li>• Most reproducible results</li> <li>• Fastest runs</li> </ul>	<ul style="list-style-type: none"> <li>• Better suited for more advanced separations than CHEF-DR II system</li> <li>• Better separation of DNA fragments &gt;2 Mb</li> </ul>	<ul style="list-style-type: none"> <li>• Suitable for routine separations with the same organism</li> <li>• Separation of DNA fragments &lt;2 Mb</li> </ul>

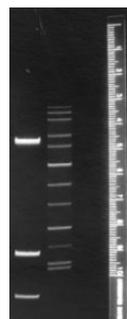


Left, two-state mode,  
24 hr run, 120° included angle  
60–120 sec switch-time ramp  
6 V/cm, 0.5x TBE, 14°C  
1.0% pulsed field Certified agarose

Right, multistate mode, 60 hr run  
State (pulse angle):

1. 90 sec switch time, –60° angle
2. 45 sec switch time, 180° angle
3. 90 sec switch time, 60° angle
4. 90 sec switch time, –60° angle
5. 90 sec switch time, 60° angle
6. 45 sec switch time, 180° angle
7. 90 sec switch time, –60° angle
8. 90 sec switch time, 60° angle

**High-resolution separation with multiple states (pulse angles).**  
*S. cerevisiae* chromosomes separated under two-state conditions (left) and under multistate conditions (right). Notice separation of the comigrating chromosomes under multistate conditions.

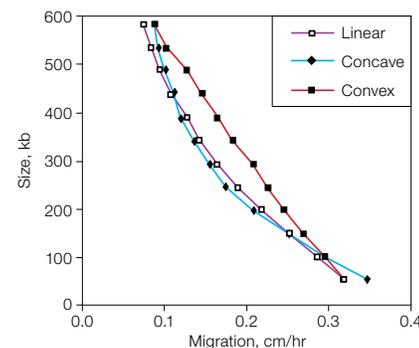


FIGE mode,  
180° angle  
200–800 ms  
switch-time ramp  
9 V/cm forward,  
6 V/cm reverse  
1x TAE, 14°C,  
18 hr run  
Forward  
switch time =  
reverse time

Lane 1: Bio-Rad  $\lambda$  HindIII  
standard (6.6, 9.4,  
23.1 kb)

Lane 2: Bio-Rad 8.3–48.5 kb  
size standard (8.3, 8.6,  
10.0, 12.2, 15.0, 17.1,  
19.4, 22.6, 24.8, 29.9,  
33.5, 38.4, 48.5 kb)

**High resolution of 8.3–48.5 kb  
size standard on the CHEF Mapper  
XA system with AFIGE.**



**Mobility effects of nonlinear switch-time ramps on the CHEF Mapper XA system.**  
Molecular size vs. migration for linear, concave, and convex ramps. The convex ramp results in the widest linear range.

## Ordering Information

Catalog #	Description	Catalog #	Description
170-3670	<b>CHEF Mapper XA Chiller System</b> , 120 V, includes CHEF Mapper XA power module, embedded auto-algorithm for protocol optimization, interactive algorithm program disk, electrophoresis cell, cooling module, variable-speed pump, Tygon tubing (12'), 14 x 13 cm (W x L) casting stand, 15-well 1.5 mm comb and comb holder, screened cap, disposable plug molds, leveling bubble, cables, <i>S. cerevisiae</i> DNA size standards, two 0.5 A FB fuses, 5 g pulsed field Certified agarose, 5 g Certified megabase agarose, instructions, for North America	170-3592	<b>CHEF Bacterial Genomic DNA Plug Kit</b> , contains 12 ml cell suspension buffer, 1.3 ml proteinase K, 30 ml proteinase K reaction buffer, 12 ml 2% CleanCut agarose, 60 ml 10x wash buffer, 1.6 ml lysozyme (25 mg/ml), 30 ml lysozyme buffer, screened cap, 2 disposable plug molds, instructions; makes 100 plugs
170-3671	<b>CHEF Mapper XA Chiller System</b> , 100 V, for Japan	170-3593	<b>CHEF Yeast Genomic DNA Plug Kit</b> , contains 12 ml cell suspension buffer, 1.3 ml proteinase K, 30 ml proteinase K reaction buffer, 12 ml 2% CleanCut agarose, 60 ml 10x wash buffer, 1.6 ml lyticase, 25 ml lyticase buffer, screened cap, 2 disposable plug molds, instructions; makes 100 plugs
170-3672	<b>CHEF Mapper XA Chiller System</b> , 220 V, for Asia Pacific/Europe	<b>Agaroses and Size Standards for PFGE</b>	
170-3673	<b>CHEF Mapper XA Chiller System</b> , 240 V, for Asia Pacific/Europe	161-3108	<b>Certified Megabase Agarose</b> , 25 g
170-3700	<b>CHEF-DR III Variable Angle Chiller System</b> , 120 V, includes power module, electrophoresis cell, cooling module, variable-speed pump, 14 x 13 cm casting stand with frame and platform, comb holder, 15-well 1.5 mm thick comb, screened cap, disposable plug molds, 12' Tygon tubing, 2 plugs <i>S. cerevisiae</i> DNA size standards, two 0.5 A FB fuses, 5 g pulsed field Certified agarose, 5 g Certified megabase agarose, instructions, for North America	161-3109	<b>Certified Megabase Agarose</b> , 125 g
170-3702	<b>CHEF-DR III Variable Angle Chiller System</b> , 220/240 V, for Asia Pacific/Europe	161-3110	<b>Certified Megabase Agarose</b> , 500 g
170-3703	<b>CHEF-DR III Variable Angle Chiller System</b> , 100 V, for Japan	161-3100	<b>Certified Molecular Biology Agarose</b> , 25 g
170-3725	<b>CHEF-DR II Chiller System</b> , 120 V, includes electrophoresis cell, drive module, cooling module, control module, variable-speed pump, 14 x 13 cm casting stand with frame and platform, comb holder, 15-well 1.5 mm thick comb, screened cap, disposable plug molds, 12' Tygon tubing, 2 plugs <i>S. cerevisiae</i> DNA size standards, two 0.5 A FB fuses, 5 g pulsed field Certified agarose, 5 g Certified megabase agarose, instructions, for North America	161-3101	<b>Certified Molecular Biology Agarose</b> , 125 g
170-3727	<b>CHEF-DR II Chiller System</b> , 220/240 V, for Asia Pacific/Europe	161-3102	<b>Certified Molecular Biology Agarose</b> , 500 g
170-3728	<b>CHEF-DR II Chiller System</b> , 100 V, for Japan	162-0137	<b>Pulsed Field Certified Agarose</b> , 100 g
170-8190	<b>Molecular Imager Gel Doc XR+ System</b> , PC and Mac	162-0138	<b>Pulsed Field Certified Agarose</b> , 500 g
170-9300	<b>FPQuest Basic Software</b>	170-3594	<b>CleanCut Agarose</b> , 2%, 12 ml; makes 24 ml of sample mixture or 100 plugs
170-9301	<b>FPQuest Cluster Analysis</b>	170-3605	<b>CHEF DNA Size Marker</b> , <i>S. cerevisiae</i> , 0.2–2.2 Mb, 5 agarose blocks, sufficient for 25–40 plugs
170-9302	<b>FPQuest Identification and Library Manager</b>	170-3667	<b>CHEF DNA Size Marker</b> , <i>H. wingei</i> , 1–3.1 Mb, 5 agarose blocks, sufficient for 25–40 plugs
170-9310	<b>InfoQuestFP Basic Fingerprint Types</b>	170-3633	<b>CHEF DNA Size Marker</b> , <i>S. pombe</i> , 3.5–5.7 Mb, 5 agarose blocks, sufficient for 25–40 plugs
170-9314	<b>InfoQuestFP Cluster Analysis</b>	170-3624	<b>CHEF DNA Size Standard</b> , 5 kb ladder, 4.9–120 kb, 20–25 lanes
170-9315	<b>InfoQuestFP Identification and Library Manager</b>	170-3707	<b>CHEF DNA Size Standard</b> , 8.3–48.5 kb, 125 lanes
170-3591	<b>CHEF Mammalian Genomic DNA Plug Kit</b> , contains 12 ml cell suspension buffer, 1.3 ml proteinase K, 30 ml proteinase K reaction buffer, 12 ml 2% CleanCut agarose, 60 ml 10x wash buffer, screened cap, 2 disposable plug molds, instructions; makes 100 plugs	170-3635	<b>CHEF DNA Size Standard</b> , lambda ladder, 0.05–1 Mb, 5 agarose blocks, sufficient for 25–40 plugs
			<b>Premixed Nucleic Acid Electrophoresis Buffers and Stains</b>
		161-0733	<b>10x Tris/Boric Acid/EDTA (TBE)</b> , 1 L bottle
		161-0770	<b>10x Tris/Boric Acid/EDTA (TBE)</b> , 5 L cube
		161-0743	<b>50x Tris/Acetic Acid/EDTA (TAE)</b> , 1 L bottle
		161-0773	<b>50x Tris/Acetic Acid/EDTA (TAE)</b> , 5 L cube
		161-0433	<b>Ethidium Bromide Solution</b> , 10 mg/ml, 10 ml
			<b>Reference</b>
			Chu G et al. (1986). Science 234, 1582–1585. CHEF (U.S. patent 5,549,796 issued to Stanford University) is exclusively licensed to Bio-Rad Laboratories, Inc.
			Tygon is a trademark of Saint-Gobain Performance Plastics Corporation. Windows is a trademark of Microsoft Corporation.



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600 **Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 64 9 415 2280 **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 3170 **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 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000