

## DCode™ Universal Mutation Detection Systems, Reagents, and Accessories

Catalog #	Description
<b>DCode Systems*</b>	
170-9080	DCode System for DGGE, 120 V, for 16 cm gels with single prep well (1 mm), includes comb gasket, 2 sets of clamps, Model 475 gradient former, all parts required to cast gradient gels
170-9081	DCode System for DGGE, 220/240 V, for 16 cm gels with single prep well (1 mm)
170-9082	DCode System for DGGE, 100 V, for 16 cm gels with single prep well (1 mm)
170-9083	DCode System for DGGE, 120 V, for 10 cm gels with 2 prep wells (1 mm)
170-9084	DCode System for DGGE, 220/240 V, for 10 cm gels with 2 prep wells (1 mm)
170-9085	DCode System for DGGE, 100 V, for 10 cm gels with 2 prep wells (1 mm)
170-9086	DCode System for CDGE, 120 V, for 16 cm gels with 20 wells (1 mm)
170-9087	DCode System for CDGE, 220/240 V, for 16 cm gels with 20 wells (1 mm)
170-9088	DCode System for CDGE, 100 V, for 16 cm gels with 20 wells (1 mm)
170-9089	DCode System for TTGE, 120 V, for 16 cm gels with 20 wells (1 mm)
170-9090	DCode System for TTGE, 220/240 V, for 16 cm gels with 20 wells (1 mm)
170-9091	DCode System for TTGE, 100 V, for 16 cm gels with 20 wells (1 mm)
170-9092	DCode System for SSCP, 120 V, for 20 cm gels with 20 wells (0.75 mm), includes cooling tank adaptor for use with external cooling bath, control reagents for SSCP
170-9093	DCode System for SSCP, 220/240 V, for 20 cm gels with 20 wells (0.75 mm)
170-9094	DCode System for SSCP, 100 V, for 20 cm gels with 20 wells (0.75 mm)
170-9105**	Complete DCode System, 120 V, PC, for all gel sizes and types described above, includes software, standard and cooling tanks, Model 475 gradient former, sandwich clamps, pressure clamp, comb gasket and holder, fittings required for gradient gels
170-9106**	Complete DCode System, 220/240 V, PC
170-9107**	Complete DCode System, 100 V, PC
170-9102**	Complete DCode System, 120 V, Mac
170-9103**	Complete DCode System, 220/240 V, Mac
170-9104**	Complete DCode System, 100 V, Mac

### Training Guide

This self-training guide is an interactive CD-ROM that helps you learn about techniques for screening mutations using the DCode system. It includes the following:

- Principles of DGGE, CDGE, TTGE, and SSCP
- Videos on setting up and using the DCode system
- WinMelt™ software tutorial program, DCode application notes, instruction manual, and other literature

To obtain your self-training guide, request catalog #170-9241. The CD-ROM is PC-compatible only.

### DCode Adaptor Kits\*\*\*

170-9125	DGGE Kit, for 16 cm gels with single prep well (1 mm), includes sandwich clamps, pressure clamp, comb gasket and holder, fittings required for gradient gel casting
170-9126	DGGE Kit, for 10 cm gels with 2 prep wells (1 mm)
170-9127	CDGE/TTGE Kit, for 16 cm gels with 20 wells (1 mm)
170-9128	Complete SSCP Kit, for 20 cm gels with 20 wells (0.75 mm), includes sandwich clamps, cooling finger adaptor for use with external chiller
170-9129	Basic SSCP Kit, for 20 cm gels with 20 wells (0.75 mm), includes sandwich clamps

### DCode Accessories

170-9034	MacMelt™ Software, Mac
170-9240	WinMelt Software, PC/Windows
170-9042	Model 475 Gradient Delivery System, includes cam-operated manual gradient former, 2 each of 10 and 30 ml syringes, all accessories required to cast gradient gels
170-9140	Electrophoresis Cooling Tank, with cooling adaptor for hookup to laboratory recirculating chiller

\* Each system includes electrophoresis/temperature control module, sandwich core, kit to cast gels of indicated size and type (2 sets of plates, 2 sets of clamps and spacers, 2 combs), control reagents for indicated application(s).

\*\* For PC, includes WinMelt software; for Mac, includes MacMelt software.

\*\*\* Each kit includes 2 sets of plates, 2 sets of spacers, 2 combs.

Catalog #	Description	170-9205	Spacers, 16 cm, 1 mm, 1 pair
<b>Electrophoresis Reagents and DNA Control Reagents</b>			
170-9150	DCode Control Reagent Kit for DGGE/CDGE/TTGE, includes primers (one GC-clamped) and DNA templates for production of wild-type and mutant DNA	170-9206	Spacers, 16 cm, 1.5 mm, 1 pair
170-9151	DCode Control Reagent Kit for SSCP, includes primers and DNA templates for production of wild-type and mutant DNA	170-9207	Spacers, 20 cm, 0.75 mm, 1 pair
170-9170	DCode Electrophoresis Reagent Kit for DGGE, includes 500 ml 40% acrylamide/bis (37.5:1), 2 x 1 L 50x TAE buffer, 225 ml 100% deionized formamide, 10 ml 10 mg/ml ethidium bromide, 10 ml DCode dye solution, 5 ml TEMED, 1 ml 2x gel loading dye, 10 g ammonium persulfate	170-9208	Spacers, 20 cm, 1 mm, 1 pair
170-9171	DCode Electrophoresis Reagent Kit for TTGE, includes 500 ml 40% acrylamide/bis (37.5:1), 1 kg urea, 2 x 1 L 50x TAE buffer, 10 ml 10 mg/ml ethidium bromide, 1 ml 2x gel loading dye, 5 ml TEMED, 10 g ammonium persulfate	170-9209	Spacers, 20 cm, 1.5 mm, 1 pair
170-9172	DCode Electrophoresis Reagent Kit for SSCP, includes 500 ml 40% acrylamide solution, 500 ml 2% bis-acrylamide solution, 100 ml glycerol, 6 x 1 L 10x TBE buffer, 2x SSCP gel loading dye, 5 ml TEMED, 10 g ammonium persulfate	170-9210	Sandwich Clamps, DCode, 10 cm, one pair
<b>DCode Miscellaneous Accessories</b>			
170-9181	Comb, one-well prep, 0.75 mm	170-9211	Sandwich Clamps, DCode, 16 cm, one pair
170-9182	Comb, one-well prep, 1 mm	170-9212	Sandwich Clamps, DCode, 20 cm, one pair
170-9183	Comb, one-well prep, 1.5 mm	170-9213	Glass Plate, inner, 10 cm, 2
170-9184	Comb, two-well prep for 10 cm DGGE, 0.75 mm	170-9214	Glass Plate, outer, 10 cm, 2
170-9185	Comb, two-well prep for 10 cm DGGE, 1 mm	170-9215	Glass Plate, inner, 16 cm, 2
170-9186	Comb, two-well prep for 10 cm DGGE, 1.5 mm	170-9216	Glass Plate, outer, 16 cm, 2
170-9187	Comb, 16-well, 0.75 mm	170-9217	Glass Plate, inner, 20 cm, 2
170-9188	Comb, 16-well, 1 mm	170-9218	Glass Plate, outer, 20 cm, 2
170-9189	Comb, 16-well, 1.5 mm	170-9219	Electrophoresis Core, DCode
170-9190	Comb, 20-well, 0.75 mm	170-9220	Replacement Gasket for Core, DCode
170-9191	Comb, 20-well, 1 mm	170-9221	DCode Lid Stand
170-9192	Comb, 20-well, 1.5 mm	170-9222	DCode Electrophoresis Tank
170-9193	Comb, 25-well, 0.75 mm	170-9223	DCode Casting Stand with 2 Gaskets
170-9194	Comb, 25-well, 1 mm	170-9224	Replacement Gasket for DCode Casting Stand, 2
170-9195	Comb, 25-well, 1.5 mm	170-9225	Replacement Belt for DCode Stirrer
170-9197	Comb, 32-well, 1 mm	170-9226	Pressure Clamp, 10 cm, DCode System for DGGE
170-9198	Comb, 10 cm, grooved for gradient casting, 0.75 mm, 1 pair plus middle spacer	170-9227	Pressure Clamp, 16 cm, DCode System for DGGE
170-9199	Spacers, 10 cm, grooved for gradient casting, 1 mm, 1 pair plus middle spacer	170-9228	Inlet Fitting, 2, with O-rings, for gradient casting
170-9200	Spacers, 10 cm, grooved for gradient casting, 1.5 mm, 1 pair plus middle spacer	170-9229	DCode Gradient Delivery System Tubing and Fitting Kit, includes 3' Tygon tubing, 5 Y-fittings, 4 vent port plugs, 4 female luer fittings, 4 male luer fittings, 2 stopcocks, two 15 ml syringes, two 30 ml syringes
170-9201	Spacers, 16 cm, grooved for gradient casting, 0.75 mm	170-9196	Y-Fittings for Gradient Delivery System
170-9202	Spacers, 16 cm, grooved for gradient casting, 1 mm	170-9230	Comb Gasket Holder, DCode system for DGGE
170-9203	Spacers, 16 cm, grooved for gradient casting, 1.5 mm	170-9231	Comb Gasket for 0.75 and 1.0 mm Combs, DCode system for DGGE
170-9204	Spacers, 16 cm, 0.75 mm, 1 pair	170-9232	Comb Gasket for 1.5 mm Combs, DCode system for DGGE
		170-9233	DCode Alignment Card
		170-9235	Temperature Control Module, to upgrade D GENE™ to do TTGE
		Bulletin #	Title
		<b>Literature Available</b>	
		2069	<b>DCode Universal Mutation Detection System Brochure</b>
		5479	<b>DCode System CD (includes WinMelt/MacMelt Software and DCode Literature Library)</b>
			Mac is a trademark of Apple Inc. Tygon is a trademark of Norton Co. Windows is a trademark of Microsoft Corporation.



**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 905 364 3435 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 6361 7000 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