

DCode™ Universal Mutation Detection Systems, Reagents, and Accessories

| Catalog # | Description |
|-----------------------|---|
| DCode Systems* | |
| 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 |
| Electrophoresis Reagents and DNA Control Reagents | | | |
| 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 |
| DCode Miscellaneous Accessories | | | |
| 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 |
| | | Literature Available | |
| | | 2069 | DCode Universal Mutation Detection System Brochure |
| | | 5479 | DCode System CD (includes WinMelt/MacMelt Software and DCode Literature Library) |
| | | | 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 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