

DCode™ System and Accessories

DCode Universal Mutation Detection System Catalog #170-9080–170-9104

Components

Core and clamps	Tank: molded polycarbonate Core: molded polysulfone Clamps: molded glass-filled polycarbonate
Gradient former	Cast acrylic and aluminum
Lid	Polycarbonate
Electrodes	0.010" diameter platinum
Electrical leads	Flexible, straight
Glass plates	20 x 20 cm (20 cm format inner plate) 22 x 20 cm (20 cm format outer plate) 16 x 20 cm (16 cm format inner plate) 18.3 x 20 cm (16 cm format outer plate) 10.2 x 20 cm (10 cm format inner plate) 12.4 x 20 cm (10 cm format outer plate)
Gel sizes	16 x 20 cm (maximum two per run) 16 x 16 cm (maximum two per run) 16 x 10 cm (maximum two per run) 7.5 x 10 cm (maximum four per run)
Spacers available	0.75, 1.0, 1.5 mm
Combs	16, 20, 25, and 32 (1 mm only) well combs, and one well comb (prep comb for perpendicular gradient gels)
Casting stand	Casts two 16 x 16 cm, two 16 x 10 cm, two 16 x 20 cm, or four 7.5 x 10 cm gels per setup
Heater and control	Temperature control (PID type) $\pm 0.5^{\circ}\text{C}$ variation within gel area, $\pm 0.5^{\circ}\text{C}$ actual in the range of 45–70°C
Maximum set temperature	70.5°C
DC voltage/power limit	500 V DC voltage limit; 50 W power limit
Size and Weight	
Dimensions (W x D x H)	Lid and tank assembly: 39 x 20 x 42 cm
Weight	16 kg (35.3 lb)

Power Requirements

AC power requirement	AC power input: 120 VAC 47–63 Hz, 5 A slow blow fuse (catalog #170-9080/9083/9086/9089/ 9092/9095/9098/9102) AC power input: 100 VAC 47–63 Hz, 5 A slow blow fuse (catalog #170-9082/9085/9088/9091/ 9094/9097/9100/9104) AC power input: 220–240 VAC 47–63 Hz, 2.5 A slow blow fuse (catalog #170-9081/9084/9087/ 9090/9093/9096/9099/9103)
DC power requirement	External DC voltage power supply. Power supply must be ground isolated in such a way that DC voltage output floats with respect to ground
Maximum voltage limit	500 V DC
Maximum power limit	50 W

Environmental Requirements

Storage environment	0–70°C, humidity 0–95% (noncondensing)
Operating environment	0–35°C, humidity 0–95%
Regulatory	Meets requirements of IEC 1010-1 and FCC, Class A

Model 475 Gradient Delivery System Catalog #170-9042

Size and weight	22 x 21 cm (W x H); 2.28 kg (5 lb)
Capacity	50 ml total (25 ml/syringe)

DNA Melting Profile Analysis Software

MacMelt™ Software, Catalog #170-9034

System Requirements	Mac, CD-ROM drive, Mac operating system 7.0 (or later)
---------------------	---

WinMelt™ Software, Catalog #170-9240

System Requirements	PC, 486 MHz or faster CPU, CD-ROM drive, Windows 95 operating system (or later)
---------------------	--

Mac is a trademark of Apple Inc. 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
