



From Start to Finish

## Multiplex Nucleic Acid Solutions

**From flexible assay design to straightforward data analysis, Bio-Rad offers the complete solution for bead-based multiplex nucleic acid analysis.**

- Easy multiplexing of up to 100 targets in one well using the Bio-Plex suspension array system
- Simple single nucleotide polymorphism (SNP) genotyping using the Bio-Plex® SNP Manager™ macro, specifically designed for integrated allelic discrimination using Bio-Plex Manager™ software
- Cost-effective applications in SNP analysis, pathogen detection, strain typing, halotyping, and gene expression
- Easy-to-follow protocols for homebrew assay development
- Dedicated field application specialists and worldwide technical support



For more information, visit us on the Web at [www.bio-rad.com/bio-plex/na-apps/](http://www.bio-rad.com/bio-plex/na-apps/)

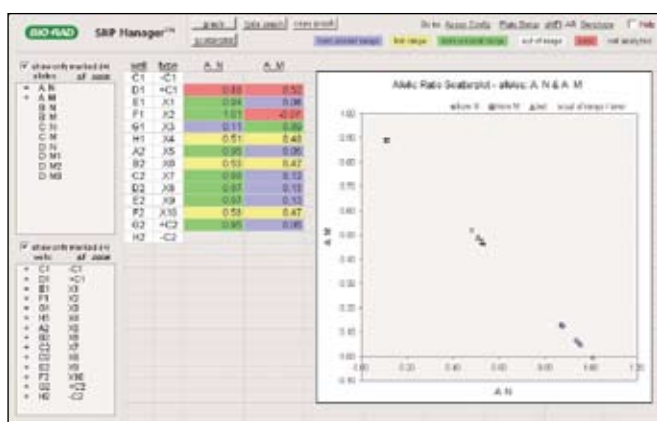
**BIO-RAD**

## Simple SNP Analysis

The Bio-Plex SNP Manager macro facilitates SNP analysis and genotyping for customers performing nucleic acid applications using the Bio-Plex suspension array system. Obtain Bio-Plex SNP Manager at no cost from the Bio-Rad web site at

[www.bio-rad.com/snpmanager/](http://www.bio-rad.com/snpmanager/)

- Automatic allelic ratio calculation and gene calls are based on user-configured tolerance ranges
- User-configured error thresholds are used to flag abnormal results
- All data can be sorted, filtered, and graphed for easy analysis
- Data can be automatically imported directly from Bio-Plex Manager software into SNP Manager
- Graphs can be exported for use in reports and presentations



## Ordering Information

Catalog # Description

### Bio-Plex 200 Systems

- 171-000201 Bio-Plex 200 System, 100–240 V, includes array reader, microplate platform, Bio-Plex Manager 4.1 Standard software with workstation license, PC, monitor, calibration kit, validation kit 4.0, maintenance, calibration, and validation (MCV) plate III, 20 L sheath fluid, spare sample needle
- 171-000205 Bio-Plex 200 System With HTF, same as 171-000201 with high-throughput fluidics (HTF)

### Bio-Plex COOH Beads

- 171-606xxx\* Bio-Plex COOH Beads, (xxx = 000–099)  $1.25 \times 10^7$  beads/ml, 16 ml
- 171-506xxx\* Bio-Plex COOH Beads, (xxx = 000–099)  $1.25 \times 10^7$  beads/ml, 1 ml

### Kits for RNA Isolation

- 732-6800 Aurum™ Total RNA 96 Kit
- 732-6820 Aurum Total RNA Mini Kit

\* A complete selection of 100 spectrally distinct COOH beads is available in 1 ml tubes and 16 ml bottles. For detailed ordering information, request bulletin 2935, or visit us on the Web at [www.bio-rad.com/bio-plex/](http://www.bio-rad.com/bio-plex/)

Catalog # Description

### Kits for Genomic DNA Isolation

- 732-6340 AquaPure™ Genomic DNA Isolation Kit
- 732-6343 AquaPure Genomic DNA Tissue Kit
- 732-6345 AquaPure Genomic DNA Blood Kit

### PCR Reagents

- 172-5302 iProof™ High-Fidelity DNA Polymerase
- 172-5311 iProof HF Master Mix
- 172-5321 iProof GC Master Mix
- 172-5331 iProof High-Fidelity PCR Kit
- 170-8870 iTaq™ DNA Polymerase
- 170-8874 dNTP Mix
- 170-8891 iScript™ cDNA Synthesis Kit, 100 x 20 µl reactions
- 170-8897 iScript™ Select cDNA Synthesis Kit, 100 x 20 µl reactions

### Thermal Cyclers

- PTC-0200G DNA Engine® Thermal Cycler Chassis, does not include Alpha™ unit (requires 1)
- PTC-0220G DNA Engine Dyad® Dual-Bay Thermal Cycler Chassis, does not include Alpha units (requires 2)
- PTC-0240G DNA Engine Tetrad® 2 Thermal Cycler Chassis, does not include Alpha units (requires 4)

### Single-Block Alpha Units With Heated Lid

- ALS-1200G Flat Block™ Alpha Unit With Flat, Heated Inner Lid, includes customizable flat-surface block with 4 screw-down points
- ALS-1238G 384-Well High-Capacity Alpha Unit With Hot Bonnet® Heated Lid, holds 384-well microplates
- ALS-1260G 60-Well Alpha Unit With Hot Bonnet Heated Lid, holds 60 x 0.5 ml tubes
- ALS-1296G 96-Well Alpha Unit With Hot Bonnet Heated Lid, holds one 96-well V-bottom plate or 96 x 0.2 ml tubes

### Accessories

- ALS-0225G Alpha Unit, dummy, fills a bay so the cycler can function with fewer Alpha units than required

### Consumables

- 223-9014 BR-14 Tips, natural, 1,000
- 223-9039 BR-39 Tips, natural, graduated, beveled, 1,000
- 223-9041 BR-41 Tips, natural, 1,000
- MLL-9601 Multiplate™ Low-Profile 96-Well Unskirted PCR Plates, natural, 25 plates
- MLL-9661 Multiplate Low-Profile 96-Well Unskirted PCR Plates, black, 25 plates
- 223-9472XTU 0.2 ml Thin-Wall PCR 8-Cap Strips, natural, 120
- 223-9477XTU 0.2 ml Thin-Wall PCR 12-Cap Strips, natural, 80
- MSF-1001 Microseal® 'F' Foil, package of 100 seals, for long-term storage
- MSA-5001 Microseal 'A' Film, package of 50 seals, for short-term storage
- MSB-1001 Microseal 'B' Adhesive Seals, 100, for PCR
- 224-0096 Costar 96-Well Flat-Bottom EIA Plates, polystyrene, 5 per package, box of 100

### Notice regarding Bio-Rad thermal cyclers.

Purchase of this instrument conveys a limited non-transferable immunity from suit for the purchaser's own internal research and development and for use in applied fields other than Human In Vitro Diagnostics under one or more of U.S. Patents Nos. 5,656,493, 5,333,675, 5,475,610 (claims 1, 44, 158, 160–163 and 167 only), and 6,703,236 (claims 1–7 only), or corresponding claims in their non-U.S. counterparts, owned by Applied Biosystems. No right is conveyed expressly, by implication or by estoppel under any other patent claim, such as claims to apparatus, reagents, kits, or methods such as 5' nuclease methods. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

Costar is a trademark of Corning Costar Corporation.

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.

**BIO-RAD**

**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 905 712 2771 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 5811 6270 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 34 911 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000