



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** (800) 480RAD **Australia** (02) 9914 2600 **Austria** (01) 877 89 01 **Belgium** (09) 396 95 11 **Brazil** (51) 507 6 191
Canada (905) 712-2777 **China** (86-21) 63302355 **Czech Republic** (420) 2-4141 0530 **Denmark** 45 44 52 1000 **Finland** 358 9 804 2200
France (33) 1 69 11 20 00 **Germany** (49) 91 25 10 10 **India** (91) 44 43 01 4000 **Italy** (39) 02 21 66 91 **Japan** (81) 3 53 51 1111 **Korea** (82) 2 34 73 4600 **Latin America** 305-894-5950 **Mexico** 52 5 554 2352 6 54
The Netherlands (31) 6-56066 **New Zealand** (64) 9-4152280 **Norway** 47 22 36 41 30 **Poland** (48) 22 8 26 672 **Portugal** 351 21 473 7700
Spain (34) 91 48 41 00 **Sweden** (46) 8 55 12 700 **Switzerland** (41) 05 62 11 555
United Kingdom (44) 1200 811154

BuMkt 2026 US Rev B

00-306 0002 Sig 0402

Multiplex Applications Using the iCycler iQ™ Real-Time PCR Detection System

Marni Brisson, Robert Park, Kent Persson, Luis Uguzzoli, David Chinn, and Keith Hamby

Bio-Rad Laboratories, Inc.
Life Science Group, 2000 Alfred Nobel Drive, Hercules, CA 94547 USA

