



**Bio-Rad**  
Laboratories, Inc.

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD

Australia 01 877 89 01 Belgium 09 385 55 11  
 Austria 01 877 89 01 Canada 905 364 3435 China 86 21 6426 0808  
 Brazil 55 21 3237 9400 Denmark 44 52 10 00  
 Czech Republic 420 241 430 532 France 01 47 95 69 65 Germany 089 318 84 0  
 Finland 09 804 22 00 Hong Kong 852 2789 3300  
 Greece 30 210 777 4396 India 91 124 4029300 Israel 03 963 6050  
 Hungary 36 1 455 8800 Japan 03 6361 7000 Korea 82 2 3473 4460  
 Italy 39 02 216091 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

08-1033 0109 Sig 0308

Bulletin 5479 Rev B US/EG

DCode System CD — Literature Library | Version 1.1

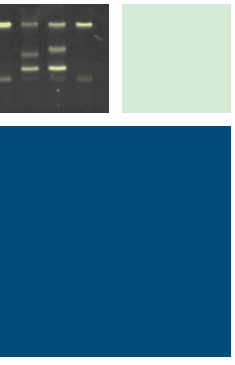


**DCode™ System CD**  
Literature Library

Unravel the Code



Mutation Analysis



Foam  
Hub