



To find the correct volumes to use when preparing the PCR mix, add the total number of samples and controls to be analyzed, and find the corresponding volumes of reagents B1, B2, or B3 and reagent C in the table.

| Total number of samples & controls | Probes Reagents B1, B2 or B3 (µl) | Amplification Mix Reagent C (µl) |
|------------------------------------|-----------------------------------|----------------------------------|
| 1                                  | 5                                 | 15                               |
| 2                                  | 11                                | 33                               |
| 3                                  | 16                                | 48                               |
| 4                                  | 22                                | 66                               |
| 5                                  | 27                                | 81                               |
| 6                                  | 32                                | 96                               |
| 7                                  | 38                                | 115                              |
| 8                                  | 43                                | 130                              |
| 9                                  | 49                                | 147                              |
| 10                                 | 54                                | 160                              |
| 11                                 | 59                                | 177                              |
| 12                                 | 65                                | 195                              |
| 13                                 | 70                                | 210                              |
| 14                                 | 76                                | 230                              |
| 15                                 | 81                                | 245                              |
| 16                                 | 86                                | 260                              |
| 17                                 | 92                                | 275                              |
| 18                                 | 97                                | 290                              |
| 19                                 | 103                               | 310                              |
| 20                                 | 108                               | 325                              |
| 21                                 | 113                               | 340                              |
| 22                                 | 119                               | 357                              |
| 23                                 | 124                               | 370                              |
| 24                                 | 130                               | 390                              |
| 25                                 | 135                               | 405                              |
| 26                                 | 140                               | 420                              |
| 27                                 | 146                               | 440                              |
| 28                                 | 151                               | 450                              |
| 29                                 | 157                               | 470                              |
| 30                                 | 162                               | 485                              |
| 31                                 | 167                               | 500                              |
| 32                                 | 173                               | 520                              |

Bio-Rad, S.N.C. au capital de 50 000 000 Euros, Locataire-Gérant, 449 990 712 RCS Nanterre - N° Intracomunautaire FR 62 449 990 712 - Siret 449 990 712 00019



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
 Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00  
 France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300  
 Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Malaysia 60 3 2117 5260 Mexico 52 555 488 7670  
 The Netherlands 0318 540666 New Zealand 64 9 415 2280 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 3170 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 Thailand 66 2 6518311 United Kingdom 020 8328 2000