

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° Intracommunautaire FR 62 449 990 712 - Siret 449 990 712 00019



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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