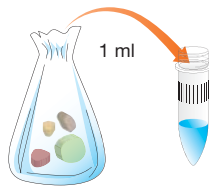
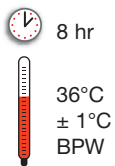


Quick Guide

Simultaneous detection of *E.coli* O157:H7 and *Salmonella* spp. in the same raw ground beef sample

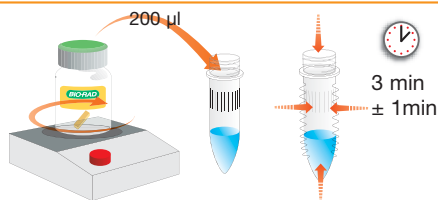
357-8114 • iQ-Check™ *E.coli* O157:H7
357-8123 • iQ-Check™ *Salmonella* II
Standard Extraction II



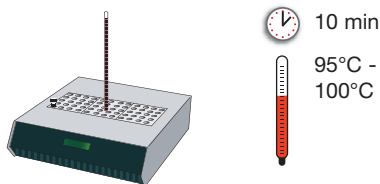
- Enrich the sample in pre-warmed buffered peptone water, 8 hr at 36°C ± 1°C
- Collect 1 ml of enriched sample and place it in a 1.5 ml screwcap tube (*Shake stomacher bag to homogenize, and allow to decant before collecting. Avoid collecting large fragments of food debris*)



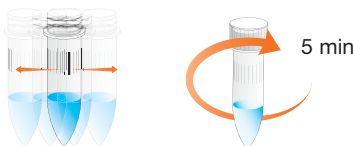
- Centrifuge at 10,000 - 12,000 g for 5 min
- Discard all the supernatant



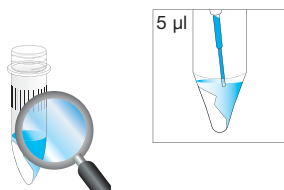
- Add 200 µl of the complete lysis reagent (reagent A + reagent F) to the pellet
- Lysis reagent must be constantly stirring in order to keep it in suspension
- Resuspend the pellet by pipetting the reagent up and down in the tube
- Grind at high speed for 3 min ± 1 min in vortex disruptor



- Place the tube in the heating block
- Incubate at 95 - 100°C for 10 min

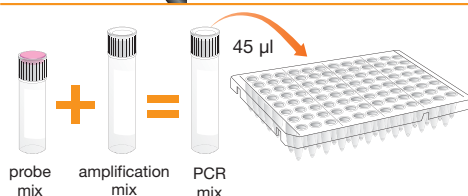


- Vortex at high speed
- Centrifuge at 10,000 - 12,000 g for 5 min



- Use 5 µl of the supernatant obtained for the amplification reaction of iQ-Check *E.coli* O157:H7 and 5 µl of supernatant for the amplification reaction of iQ-Check *Salmonella* II

Do not vortex before collecting 5 µl sample



- Prepare the PCR mix for each iQ-Check test (See PCR mix calculation guide)
- Distribute the PCR mix (45 µl)
- Distribute the samples and the controls (5 µl) for each iQ-Check test
- Check that there are no bubbles
- Seal the microplate



- Set up software for both iQ-Check *E.coli* O157:H7 and iQ-Check *Salmonella* II
- Create the plate setup
- Place the microplate into the thermocycler
- Start the amplification by clicking on "Run"

Please read the kit instruction manual and instrument user guide for complete and detailed instructions.

Quick Guide

iQ-Check™ *E.coli* O157:H7 iQ-Check™ *Salmonella* II PCR Mix Calculation Guide

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 reagent B and reagent C in the table.

| Total number of samples & controls | Probes - Reagent B (µl) | Amplification Mix Reagent C (µl) |
|------------------------------------|-------------------------|----------------------------------|
| 1 | 5 | 40 |
| 2 | 11 | 86 |
| 3 | 16 | 130 |
| 4 | 22 | 173 |
| 5 | 27 | 216 |
| 6 | 32 | 259 |
| 7 | 38 | 302 |
| 8 | 43 | 346 |
| 9 | 49 | 389 |
| 10 | 54 | 432 |
| 11 | 59 | 475 |
| 12 | 65 | 518 |
| 13 | 70 | 562 |
| 14 | 76 | 605 |
| 15 | 81 | 648 |
| 16 | 86 | 691 |
| 17 | 92 | 734 |
| 18 | 97 | 778 |
| 19 | 103 | 821 |
| 20 | 108 | 864 |
| 21 | 113 | 907 |
| 22 | 119 | 950 |
| 23 | 124 | 994 |
| 24 | 130 | 1000 |
| 25 | 135 | 1100 |
| 26 | 140 | 1100 |
| 27 | 146 | 1200 |
| 28 | 151 | 1200 |
| 29 | 157 | 1300 |
| 30 | 162 | 1300 |
| 31 | 167 | 1300 |
| 32 | 173 | 1400 |
| 33 | 178 | 1400 |
| 34 | 184 | 1500 |
| 35 | 189 | 1500 |
| 36 | 194 | 1600 |
| 37 | 200 | 1600 |
| 38 | 205 | 1600 |
| 39 | 211 | 1700 |
| 40 | 216 | 1700 |
| 41 | 221 | 1800 |
| 42 | 227 | 1800 |
| 43 | 232 | 1900 |
| 44 | 238 | 1900 |
| 45 | 243 | 1900 |
| 46 | 248 | 2000 |
| 47 | 254 | 2000 |
| 48 | 259 | 2100 |

| Total number of samples & controls | Probes - Reagent B (µl) | Amplification Mix Reagent C (µl) |
|------------------------------------|-------------------------|----------------------------------|
| 49 | 265 | 2100 |
| 50 | 270 | 2200 |
| 51 | 275 | 2200 |
| 52 | 281 | 2200 |
| 53 | 286 | 2300 |
| 54 | 292 | 2300 |
| 55 | 297 | 2400 |
| 56 | 302 | 2400 |
| 57 | 308 | 2500 |
| 58 | 313 | 2500 |
| 59 | 319 | 2500 |
| 60 | 324 | 2600 |
| 61 | 329 | 2600 |
| 62 | 335 | 2700 |
| 63 | 340 | 2700 |
| 64 | 346 | 2800 |
| 65 | 351 | 2800 |
| 66 | 356 | 2900 |
| 67 | 362 | 2900 |
| 68 | 367 | 2900 |
| 69 | 373 | 3000 |
| 70 | 378 | 3000 |
| 71 | 383 | 3100 |
| 72 | 389 | 3100 |
| 73 | 394 | 3200 |
| 74 | 400 | 3200 |
| 75 | 405 | 3200 |
| 76 | 410 | 3300 |
| 77 | 416 | 3300 |
| 78 | 421 | 3400 |
| 79 | 427 | 3400 |
| 80 | 432 | 3500 |
| 81 | 437 | 3500 |
| 82 | 443 | 3500 |
| 83 | 448 | 3600 |
| 84 | 454 | 3600 |
| 85 | 459 | 3700 |
| 86 | 464 | 3700 |
| 87 | 470 | 3800 |
| 88 | 475 | 3800 |
| 89 | 481 | 3800 |
| 90 | 486 | 3900 |
| 91 | 491 | 3900 |
| 92 | 497 | 4000 |
| 93 | 502 | 4000 |
| 94 | 508 | 4100 |
| 95 | 513 | 4100 |
| 96 | 518 | 4100 |



Bio-Rad
Laboratories, Inc.

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 43 (0) 1 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 712 2771 China 86 21 6305 2255 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 33 1 47 95 62 59
Germany 49 (0) 89 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 2398 112/113/114
Israel 03 951 4127 Italy 39 02 216 091 Japan 03 5811 6270 Korea 82 2 3473 4460 Latin America 305 894 5960 Mexico 52 555 200 0520
The Netherlands 31 318 540 666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700
Romania 4021 210 1703 Russia 7 095 721 1404 Singapore 65 6415 3188 South Africa 27 11 442 8508 Spain 34 91 590 5200 Sweden 46 8 555 12700
Switzerland 41 (0) 61 717 95 55 Taiwan 886 2 2578 7189 Thailand 662 651 8311 United Kingdom 44 20 8328 2000 Vietnam 848 823 6757