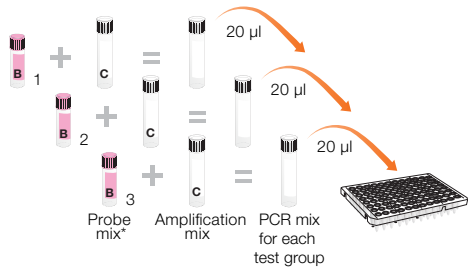
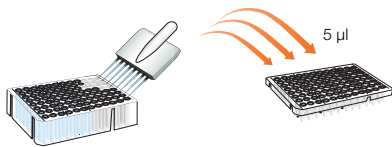


# iQ-Check STEC SerO II, 12013174

## Using CFX Manager IDE Software



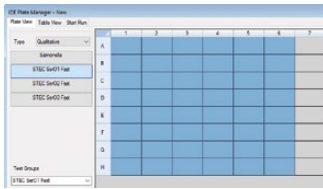
- Prepare the PCR mix for each Test Group (see PCR Mix Calculation Guide on back)
- Use probe vial B1 for Test Group **STEC SerO1 Fast**, probe vial B2 for Test Group **STEC SerO2 Fast**, and probe vial B3 for Test Group **STEC SerO3 Fast**
- Distribute 20 µl/well of each PCR mix in a microplate



- Add 5 µl each of positive and negative controls to designated wells
- Transfer 5 µl/well of extracted DNA (from iQ-Check STEC VirX DNA extraction preparation) to the PCR microplate for each Test Group

**Do not vortex before collecting the sample.**  
Check there are no bubbles.

- Seal the PCR microplate



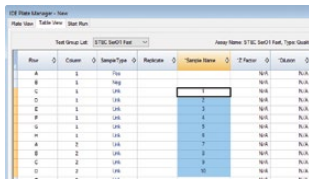
- Launch CFX Manager IDE Software to get the Plate Manager window
- Select **STEC SerO1 Fast** from the list of tests on the left side of the screen
- Select the required wells for the Test Group **STEC SerO1 Fast** assay



- Apply sample types to the wells by selecting Positive Control, Negative Control, or Unknown
- Enter sample name
- Enter Lot Number (for DNA extraction kit and/or PCR kit)



- Repeat for Test Group **STEC SerO2 Fast** and Test Group **STEC SerO3 Fast**



- To duplicate the sample names for all Test Groups, select the Table View tab
- With **STEC SerO1 Fast** selected in the Test Group dropdown, select the Sample Name column and then copy (Ctrl + C)
- In the Test Group List dropdown, move to **STEC SerO2 Fast**, then select the Sample Name column and paste (Ctrl + V)
- Repeat for **STEC SerO3 Fast**



- Start the amplification by clicking Start PCR Run

\* Probe solutions: B1, *E. coli* O157:H7 and O111; B2, *E. coli* O26, O103, and O145; B3, *E. coli* O45 and O121.

## PCR Mix Calculation Guide for the iQ-Check STEC SerO II Kit

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 B1, B2, or B3 and reagent C in the table.

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

Visit [bio-rad.com/iqcheck](http://bio-rad.com/iqcheck) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. IQ-CHECK is a trademark of Bio-Rad Europe GmbH in certain jurisdictions. All trademarks used herein are the property of their respective owner.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

