



CFX Manager Software, Industrial Diagnostic Edition (IDE) iQ-Check Prep Sample Setup Protocol Quick Guide



- Turn on the CFX96 Touch Deep Well Real-Time PCR System and open CFX Manager Software, IDE, to launch the IDE Plate Manager window. If necessary, either click the IDE Plate Manager icon on the toolbar, or go to the menu and select View > IDE Plate Manager

Umono Fast
Salmonella

- Select the test you wish to perform from the list of kits on the left side of the screen

	1	2	3	4	5	6	7	8	9	10	11	12
A												
B												
C												
D												
E												
F												
G												
H												

- Select the well(s) for that assay

Positive Control
Negative Control
Unknown (Easy I)
Unknown (Easy II)
Unknown (No extraction)
Clear Selected Wells

- Apply sample types to your wells by selecting Positive Control, Negative Control, or Unknown (Easy I, Easy II, or No extraction)

PCR Kit Lot #	<input type="text"/>
DNA Extraction Lot #	<input type="text"/>

- Type or select lot numbers for PCR Kit and/or DNA Extraction Kit

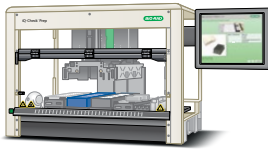
Load	Sample Name
<input type="checkbox"/>	<input type="text"/>

- Type or select sample names: select a well to start, enter this well's name in the Sample Name box, and press the Enter key. The cursor will move to the next well down
- Additional kits can be added by selecting from the list of kits to create new test groups and repeating the steps

continues

Send to iQ-Check® Prep

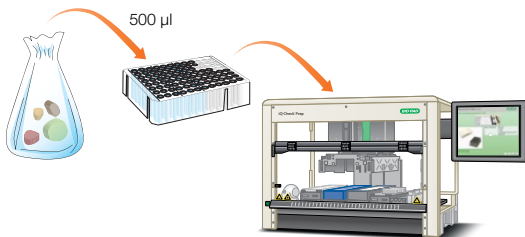
- Click Send to iQ-Check Prep



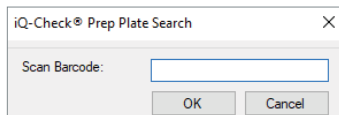
- Turn on the iQ-Check Prep System and open the iQ-Check Prep operating software



- Press or click the Start button
The daily maintenance is required if it has not been performed yet.
- Load the appropriate worklist
- Carefully follow the loading instructions on the screen in the exact order mentioned: PCR plate(s), lysis reagent(s), PCR reagents, sample deep well plate, extraction deep well plate(s), and transfer well plate



- After the enrichment step, transfer 500 µl (minimum 150 µl to maximum 1,000 µl) of sample into a deep well plate
Avoid shaking the sample bag to prevent collecting large fragments of food debris.
- Load the sample deep well plate into the iQ-Check Prep System and start the run



- Remove the PCR plate and seal it properly using optical caps. (If using optical film, a plate sealer is recommended)
- Scan the barcode of the PCR plate into CFX Manager Software, IDE, using the Load Plate File option from the Barcode menu



- Place the PCR plate into the thermal cycler
- Start the amplification by clicking Run

For detailed instructions, review the kit user guide and instrument user manual.

Visit bio-rad.com/foodscience 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 **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

