

# Quick Guide

## CheckN'Safe<sup>™</sup> E.coli

355-4700/355-4720 • Escherichia coli analysis in recreational bathing waters (sea water, inland water)

#### Start



 Switch-on respectively the uninterruptible power source (UPS), then computer (printer if necessary) and finally the XplOrer64<sup>™</sup> System





- Start the XplOrer64 Manager software by cliquing on its icon
- Menu Start, Start focused instrument
- Set the temperature of incubator(s) at 44°C and the measuring cycle on 10 minutes

NB: Incubator A (top) will be preferentially used if Enterococci analyses are performed in parallel

Warm-up the incubator(s) defined for 30 minutes

### Routine Sample Preparation



 Prepare the necessary CheckN'Safe<sup>™</sup> tests into a rack and keep them at ambient temperature at least 1 hr before use



- Prepare the necessary sterile caps and open the seals of CheckN'Safe
- Place a membrane filter composed of cellulose esters, diameter
  Ø 55 mm, nominal pore of 0.45 µm, on a membrane filters apparatus
- Filter 100 ml of sample



 Rinsing by filtration of 50-100 ml of sterile deionised or distilled water



- Using sterile tweezers, fold twice the membrane in a cone shape and inoculate it tip downside into the CheckN'Safe test NB: The entire membrane has to be immersed inside the culture medium broth.
   If necessary, rock the vial gently avoiding the formation of foam
- Add a sterile cap on each CheckN'Safe test

#### Sample Insertion

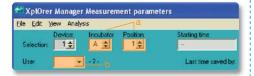


- Place CheckN'Safe E.coli into incubator(s) warm-up at 44°C
- · Verify that each CheckN'Safe test is correctly inserted
- Warming-up of tests inserted start for 1 hr

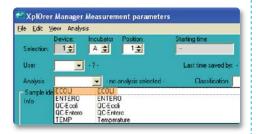
Please read the instruction manual for complete and detailed instructions.



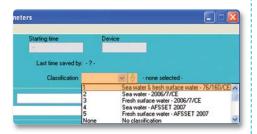
#### Measurement Parameters



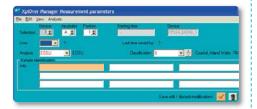
- In the main window XplOrer64 Manager Measurement Parameters, select the option View and Parameters:
  - a. Select an Incubator to parameter and the position 1
  - b. Select an User



 Select the ECOLI analysis (QC ECOLI calibration is dedicated for the internal monitoring of the culture medium performances)



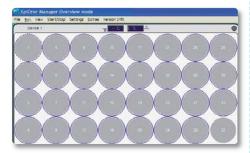
- Select a Classification for the results interpretation, according to the origin of the water sample and the directive to follow
- Confirm rapidly these settings by clicking on the button



- Fill up at minimum the INFO1 identification field. This step could be performed later at once for all samples, within the Edit table function of the Edit menu
- Confirm these settings by clicking on the 
   ✓ button
- Close this window and confirm by Yes



- To set a complete incubator: in the window XplOrer64 Manager Measurement Parameters, click into the Spot 1 with the mouse
- In the menu, select Edit and Copy position parameter



 Press simultaneously shift button on the keyboard and click into the Spot 31

(NB: spot 32 will be dedicated to the CheckN'Safe Temperature Control cell)

- Edit and Insert all parameters
- Edit and Clear copy/paste buffer to clear the buffer. Settings are therefore registered

#### Results Reading







Analysis on going

• From the end of the warming-up, samples are analyzed in continue each 10 minutes, for 8 hrs

#### Approximative Detection Time:

E. coli/100 ml	Detection time		
< 41	≥ 8.00 hr		
100	6.30 hr = <b>6 hr 18 min</b>		
250	5.55 hr = <b>5 hr 33 min</b>		
500	5.09 hr = <b>5hr05 min</b>		
1,000	4.69 hr = <b>4 hr 42 min</b>		
1,800	4.39 hr = <b>4 hr 23 min</b>		
2,000	4.34 hr = <b>4 hr 20 min</b>		
Maximum: 1.59.10°	0.1 hr = <b>6 min</b>		

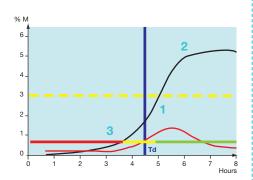
 Real-time analysis moving forward can be followed at any times by simple observation of the spots:

CheckN'Safe <sup>™</sup> <i>E. coli</i> analysis (in germs/100 ml)						
Spot appearance in XplOrer64 Manager	in real time					
	at the end of analysis					
Interpretation	Water quality	Polluted water	Suspicious pollution	Good quality	Excellent quality	
	E. coli/100 ml	≥ Imperative value	]Imperative value; guidance value[	≤ Guidance value	< Detection limit	
76/160/CEE Directive	Coastal and inland water	≥ 2,000	[100-2,000[	< 100	< 41	
2006/7/CE Directive	Coastal (sea water)	< 500	[250-500[	< 250	< 41	
	Inland water	≥ 1,000	[500-1,000[	< 500	< 41	
AFSSET 2007	Coastal (sea water)	≥ 1,000	-	< 1,000	< 41	
	Inland water	≥ 1,800	-	< 1,800	< 41	



 At the end of the cycle of analysis and in all cases before discarding the analyzed tests, finalize the samples identifications with the Edit table function of the Edit menu

#### Positive Signal Confirmation



#### Rules to confirm a positive signal:

- 1 The typical profil of the original impedance curve is a sigmoid curve, reaching the 3% threshold during the analysis
- 2 This original curve has not to decrease before to reach the "stationnary" phase
- 3 If the Td is determined within the 3 first hours of the cycle, the original curve has to reach its 3% threshold (discontinued yellow line) within 5 hrs. In this case, the software will automatically delete the measured points before these 3 hrs, and the result will be considered as negative



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 1 800 268 0213 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 1-800-180-1224 Israel 03 951 4127 Italy 39 02 216 091 Japan 03 5811 6270 Korea 82 2 3473 4460 Latin America 305 894 5960 Moxico 52 555 488 7670 The Netherlands 31 318 540 666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 9P Ortugal 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