

Urinary Tract Infection (UTI) Multiplex PCR Bacteria Profiling Kit

Code: PCR-UTI-1

Introduction

RayBio® Urinary Tract Infection (UTI) Multiplex PCR Bacteria Profiling Kit is a ready to use real-time PCR assay for the detection of UTI Bacteria DNA in a liquid sample. The kit is based on Taqman™ DNA Polymerase amplification and fluorochrome-labeled specific-probe reporting. This kit detects 21 common bacteria genomic DNA along with an internal control, RNase P across 7 multiplex real-time PCR panels with 4 different fluorescent probes in each panel.

Purpose

This kit is used for the qualitative in vitro detection of Urinary Tract Infection (UTI) Bacteria nucleic acid from purified DNA samples.

Kit Components (10-Sample Kit)

| Component | Catalog Number | Ingredients | Specification | Quantity |
|-----------------------|----------------|---|--------------------------------|----------|
| 2x PCR Master Mix | PCR-UTI-1-MM | Buffer, dNTP's, enzyme | 1350 µl / tube | 1 tube |
| Primers and Probe Mix | PCR-UTI-1-PP-1 | 7 sets Primer and Probe Mix | 40 µl / tube, 1 for each panel | 1 tube |
| | PCR-UTI-1-PP-2 | | | 1 tube |
| | PCR-UTI-1-PP-3 | | | 1 tube |
| | PCR-UTI-1-PP-4 | | | 1 tube |
| | PCR-UTI-1-PP-5 | | | 1 tube |
| | PCR-UTI-1-PP-6 | | | 1 tube |
| | PCR-UTI-1-PP-7 | | | 1 tube |
| Amplification Control | PCR-UTI-1-AC-1 | Target sequences of UTI bacteria for each panel & RNase P | 12 µl / tube, 2 for each panel | 2 tubes |
| | PCR-UTI-1-AC-2 | | | 2 tubes |
| | PCR-UTI-1-AC-3 | | | 2 tubes |
| | PCR-UTI-1-AC-4 | | | 2 tubes |

| | | | | |
|----------------------|----------------|---------------------|---------------|---------|
| | PCR-UTI-1-AC-5 | | | 2 tubes |
| | PCR-UTI-1-AC-6 | | | 2 tubes |
| | PCR-UTI-1-AC-7 | | | 2 tubes |
| Non-Template Control | PCR-UTI-NTC | Nuclease-free water | 500 µl / tube | 2 tubes |

Note: Do not mix reagents from different lots.

Storage

The kit can be stored at -20 °C for a period of 12 months prior to opening. After opening, the reagents are valid for at least 6 months if stored at -20 °C.

Required Materials (NOT INCLUDED)

1. Purified human urine DNA: DNA should be purified according to manufacturer protocols
2. DNA preservation solution
3. Fluorescence PCR instrument capable of reading FAM channel (494 nm maximum absorption, 518 nm maximum emission), VIC channel (520 nm maximum absorption, 558 nm maximum excitation), ROX channel (580 nm maximum absorption, 623 nm maximum excitation) and Cy5 (640 nm maximum absorption, 682 nm maximum excitation).
4. Vortex Mixer
5. Microcentrifuge
6. Pipettes
7. Sterile nuclease-free pipette tips (barrier tips recommended) and microfuge tubes
8. Compatible PCR Plate
9. Contact our technical support team for questions about compatibility: techsupport@raybiotech.com

Sample Requirements

1. All human urine specimens should be regarded as potentially infectious and handled with extreme caution. Measures should be taken to minimize the risk of laboratory transmission based on risk assessment when testing any samples. These precautions, at a minimum, should include proficiency and competency testing, appropriate PPE, avoiding aerosols, and using effective disinfectants (quaternary ammonium compounds and 0.5 % bleach).
2. We recommend centrifuging 1 ml of urine at 5000 g in a microcentrifuge for 10 minutes at room temperature and removing 800 µl of the supernatant. The remaining supernatant and pellet should be used for DNA extraction**. The final extracted DNA should be eluted in 100 µl of DNase, RNase-free water. The real-time PCR procedure requires 10 µl of extracted DNA per reaction.
3. Extracted human urine DNA is the starting material for this kit. The DNA should be purified separately from the PCR assay set-up according to manufacturer protocols with a DNA preservation solution for inactivation of bacteria and preservation of the DNA.

**The data shown in this manual was based on DNA extraction using a MagMAX™ Viral/Pathogen Nucleic Acid Isolation Kit from Thermo Fisher Scientific.

General Considerations

1. To prevent contamination of PCR reactions, clean and decontaminate all working surfaces, centrifuges, pipettes, and other equipment with 10 % bleach or DNA Away®, followed by 70 % Ethanol before every assay.
2. Conduct sample processing and DNA extraction in a separate area from the PCR assay setup (below termed the “PCR Assay Setup Area”).

- Care should be taken to avoid contamination of samples and reactions with DNA amplicons from used PCR plates. It is thus recommended to place used PCR plates in a sealed bag and discard them appropriately. Never open a used PCR plate or place it in the same area as the PCR instrument.
- To minimize cross-contamination between experimental samples, disposable pipettes and filtered pipette tips are recommended.

Testing Method

1. Assay Assembly (PCR Assay Setup Area)

1.1 Thaw reagents: Remove all components from the kit and fully thaw to room temperature. After thawing, mix gently by pipetting. Briefly centrifuge to collect contents at the bottom of each vial.

1.2 Prepare PCR Reaction Mix: the reaction mix for 1 test for each panel should contain 12.5 µl 2x PCR Master Mix and 2.5 µl Primers and Probe Mix, totally 15 µl. To calculate the total volume necessary for each panel, first calculate the total reaction for each panel (of samples + 1 Amplification Control + 1 Non-Template Control + 2 extra). Then, multiply the volumes of the 2x PCR Master Mix and the Primers and Probe Mix by the number of reactions and mix together to prepare a **PCR Reaction Mix** for each panel.

For example: for 10 samples, the total reactions per panel = 10 + 1 + 1 + 2 = 14 total reactions.

2. Sample Loading (PCR Assay Setup Area)

2.1. Add 15 µl of prepared **PCR Reaction Mix** from 1.2 above to each reaction well of a PCR plate.

2.2. Add 10 µl of extracted DNA (no more than 100 ng recommended), Amplification Control, or Non-Template Control for each panel to each reaction well of the PCR plate and pipette up and down at least 5 times to mix. Refer to Tables 1 and 2 for a summary of reaction components included in each well and an example plate layout.

Table 1: Reaction Components

| Component | Amplification Control Reaction (1 needed for each panel) | Non-Template Control Reaction (1 needed for each panel) | Sample Reaction (1 needed for each panel) |
|-----------------------|---|--|--|
| PCR Reaction Mix | 15 µl | 15 µl | 15 µl |
| Amplification Control | 10 µl | -- | -- |
| Non-Template Control | -- | 10 µl | -- |
| Extracted DNA | -- | -- | 10 µl |
| Total Volume | 25 µl | 25 µl | 25 µl |

Table 2: Example layout on a 96-well plate

| Sample 1 | Sample 2 | Sample 3 | Sample 4 | Sample 5 | Sample 6 | Sample 7 | Sample 8 | Sample 9 | Sample 10 | | |
|----------|----------|----------|----------|----------|----------|----------|----------|----------|-----------|-----|-----|
| Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | Panel 1 | NTC | AC1 |
| Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | Panel 2 | NTC | AC2 |
| Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | Panel 3 | NTC | AC3 |
| Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | Panel 4 | NTC | AC4 |
| Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | Panel 5 | NTC | AC5 |
| Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | Panel 6 | NTC | AC6 |
| Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | Panel 7 | NTC | AC7 |

2.3. Seal the plate tightly.

2.4. Centrifuge the plate to collect the contents at the bottom of the wells and run plate immediately.

Note: The sealed, prepared PCR plate can be stored at 2-8 °C for up to 4 hours before the “PCR Amplification” step 3 below.

3. PCR Amplification (PCR Assay Setup Area)

3.1. Instrument setup: Set the sample number, Non-Template Control, and Amplification Control accordingly to your 96-well plate setup. See Table 2 as reference. Each panel contains up to 4 targets including an Internal Amplification Control (IAC), which amplifies the RNase P gene from human samples.

3.2. Fluorescence Channel Selection:

3.2.1. Select FAM, VIC, and Cy5 (or equivalent channels) and set the targets based on Table 3.

3.2.2. Select ROX (or equivalent channel) and set the target name to “IAC”, this channel will detect the RNase P gene.

3.2.3. Ensure the “reference fluorescence dye” (passive reference) is set to “None”.

Table 3: Panels and fluorescence channel setup

| | FAM | VIC | Cy5 | ROX |
|---------|-------------------------|------------------------------|------------------------|-----|
| Panel 1 | Acinetobacter baumannii | Streptococcus agalactia | Citrobacter freundii | IAC |
| Panel 2 | Morganella morganii | Candida albicans | Klebsiella pneumoniae | IAC |
| Panel 3 | Proteus vulgaris | Klebsiella oxytoca | Enterobacter aerogenes | IAC |
| Panel 4 | Enterobacter cloacae | Neisseria gonorrhoeae | Pseudomonas aeruginosa | IAC |
| Panel 5 | Candida glabrata | Providencia stuartii | Chlamydia trachomatis | IAC |
| Panel 6 | Staphylococcus aureus | Staphylococcus saprophyticus | Proteus mirabilis | IAC |
| Panel 7 | Enterococcus faecium | Escherichia coli | Enterococcus faecalis | IAC |

3.3. Set reaction conditions according to Table 4. Set the reaction volume to 25 µl.

Table 4: PCR Program

| | Step | Temperature (°C) | Time | Temperature Ramp Rate | Number of Cycles |
|---------|-------------------------------------|------------------|-------|-----------------------|------------------|
| UNG | Remove Carry-over Contamination | 50 | 5 min | 1.6 °C/sec | 1 |
| Stage 1 | Initial Denaturation | 95 | 5 min | 1.6 °C/sec | 1 |
| Stage 2 | Denature | 95 | 10 s | 1.6 °C/sec | 40 |
| | Anneal, extend, detect fluorescence | 58 | 40 s | 1.6 °C/sec | |

3.4. Save the file, load the prepared plate, and run the program.

4. Results and Analysis

4.1. Open the data file and click 'save as' to save the file with a new name.

4.2. In the analysis settings, select automatic baseline, or select according to your lab validation.

4.3. Set the appropriate threshold values for each target as validated by your laboratory.

4.4. Determine Ct cutoff values for each target and controls.

Note: an example of established Ct target cutoffs is listed in Table 5 but each laboratory should establish their own cutoffs for the equipment and software in use.

Table 5. Example established Cutoff Ct values using a QuantStudio™ 5 Real-Time PCR machine (Thermo Fisher Scientific, Waltham, MA) and analyzed with Desing & Analysis Software 2.6.0.

| Panel | Target | Cutoff Ct |
|---------|------------------------------|-----------|
| Panel 1 | Acinetobacter baumannii | 33.6 |
| | Streptococcus agalactia | 35.0 |
| | Citrobacter freundii | 35.2 |
| Panel 2 | Morganella morganii | 34.5 |
| | Candida albicans | 35.1 |
| | Klebsiella pneumoniae | 34.6 |
| Panel 3 | Proteus vulgaris | 34.5 |
| | Klebsiella oxytoca | 35.5 |
| | Enterobacter aerogenes | 35.6 |
| Panel 4 | Enterobacter cloacae | 35.8 |
| | Neisseria gonorrhoeae | 35.9 |
| | Pseudomonas aeruginosa | 34.7 |
| Panel 5 | Candida glabrata | 34.5 |
| | Providencia stuartii | 32.8 |
| | Chlamydia trachomatis | 35.5 |
| Panel 6 | Staphylococcus aureus | 35.3 |
| | Staphylococcus saprophyticus | 32.2 |
| | Proteus mirabilis | 35.7 |
| Panel 7 | Enterococcus faecium | 33.9 |
| | Escherichia coli | 35.4 |
| | Enterococcus faecalis | 35.9 |

4.5. Interpretation of PCR reaction results.

Table 6. Interpretation of PCR reaction results for kit QC parameters (AC and NTC)

| Amplification Control (AC) | | Non-Template Control (NTC) | Interpretation |
|------------------------------|-----|------------------------------|----------------|
| Ct < 40 | and | No amplification, or Ct ≥ 40 | Valid result |
| No amplification, or Ct ≥ 40 | or | Ct < 40 | Invalid result |

Note: The Amplification Control and Non-Template Control are considered general controls for the sample preparation process. The PCR reaction results are considered valid if the Amplification Control and Non-Template Controls meet the established criteria listed in Table 6. The PCR reaction result is invalid if 1) the Ct of Amplification Control has no amplification or ≥ 40 , or 2) the Ct of Non-Template Control < 40 . If the reaction is invalid for a panel, the measurement of all DNA samples in this panel should be discarded, and the assay should be repeated. Result of IAC channel is excluded for QC interpretation.

Table 7. Interpretation of PCR reaction results

| Target Gene Ct (FAM, VIC, CY5) | Interpretation |
|---|----------------|
| $< \text{Cutoff Ct} - 0.5 \text{ Ct}$ | + (Positive) |
| $= \text{Cutoff Ct} \pm 0.5 \text{ Ct}$ | Inconclusive |
| No amplification, or $> \text{Cutoff Ct} + 0.5$ | - (Negative) |

Note: If the PCR reaction result is inconclusive, it is suggested to retest the sample. After retest, if the Ct \leq Cutoff Ct, the PCR reaction result is considered as Positive. If the Ct $>$ Cutoff Ct, PCR reaction result is considered as Negative.

Table 8. Interpretation of individual PCR reaction channels within panels

Note: each channel should be interpreted independently of the other channels in the panel

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 1 | + | + | + | + | Sample contains DNA of Acinetobacter baumannii (FAM), Streptococcus agalactia (VIC), Citrobacter freundii (Cy5) |
| | + | + | + | -** | Sample contains DNA of Acinetobacter baumannii (FAM), Streptococcus agalactia (VIC), Citrobacter freundii (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 2 | + | + | + | + | Sample contains DNA of Morganella morganii (FAM), Candida albicans (VIC), Klebsiella pneumoniae (Cy5) |
| | + | + | + | -** | Sample contains DNA of Morganella morganii (FAM), Candida albicans (VIC), Klebsiella pneumoniae (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |

| | | | | | |
|--|---|---|---|---|----------------|
| | - | - | - | - | Invalid Result |
|--|---|---|---|---|----------------|

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 3 | + | + | + | + | Sample contain DNA of Proteus mirabilis (FAM), Klebsiella oxytoca (VIC), Enterobacter aerogenes (Cy5) |
| | + | + | + | -** | Sample contain DNA of Proteus mirabilis (FAM), Klebsiella oxytoca (VIC), Enterobacter aerogenes (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 4 | + | + | + | + | Sample contain DNA of Enterobacter cloacae (FAM), Neisseria gonorrhoeae (VIC), Pseudomonas aeruginosa (Cy5) |
| | + | + | + | -** | Sample contain DNA of Enterobacter cloacae (FAM), Neisseria gonorrhoeae (VIC), Pseudomonas aeruginosa (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 5 | + | + | + | + | Sample contain DNA of Candida glabrata (FAM), Providencia stuartii (VIC), Chlamydia trachomatis (Cy5) |
| | + | + | + | -** | Sample contain DNA of Candida glabrata (FAM), Providencia stuartii (VIC), Chlamydia trachomatis (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 6 | + | + | + | + | Sample contain DNA of Staphylococcus aureus (FAM), Staphylococcus saprophyticus (VIC), Proteus vulgaris (Cy5) |
| | + | + | + | -** | Sample contain DNA of Staphylococcus aureus (FAM), Staphylococcus saprophyticus (VIC), Proteus vulgaris (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

| | FAM | VIC | CY5 | ROX (IAC) | Interpretation |
|---------|-----|-----|-----|-----------|---|
| Panel 7 | + | + | + | + | Sample contains DNA of Enterococcus faecium (FAM), Escherichia coli (VIC), Enterococcus faecalis (Cy5) |
| | + | + | + | -** | Sample contains DNA of Enterococcus faecium (FAM), Escherichia coli (VIC), Enterococcus faecalis (Cy5) |
| | - | - | - | + | Sample does not contain bacteria DNA in this panel |
| | - | - | - | - | Invalid Result |

Positive Result: the sample contains the target DNA.

Negative Result: the sample does not contain the target DNA.

Invalid Result: the sample should be rerun with fresh samples and controls.

**if the IAC (ROX) control signal is not detected while one or more of the target specific signals (FAM, Cy5, VIC) are detected in the sample, the reaction result should be considered positive and the sample is considered to contain the target DNA. The IAC is designed to be a weak signal to ensure it does not compete with the target and lower the sensitivity of the assay.

Product performance index

- Limit of Detection:** The LOD of the assay is 1~247 copies/μl, or 10~2470 copies per reaction (10 μl of sample volume).
- Repeatability:** Precision testing showed that the coefficient of variation (CV) of the precision Ct values within a kit lot are ≤ 1.2 %. Repeatability between lots of product are ≤ 10 %.

Specificity

For each panel target, no cross reactivity was identified when the kit was evaluated against the other included targets.