

# Trichomonas Foetus DNA Test Kit

VetMAX™-Gold Trich Detection Kit

Catalog Number 4483869

Pub. No. 4485468 Rev. B

**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](http://thermofisher.com/support).

## Product information

### Name, intended use, and principle of the procedure

The Applied Biosystems™ Trichomonas Foetus DNA Test Kit (VetMAX™-Gold Trich Detection Kit, Cat. No. 4483869) is a highly sensitive, qualitative, real-time PCR assay for detection of *Trichomonas foetus* DNA isolated from enriched smegma samples.

Bovine trichomonosis is a sexually transmitted infection caused by *T. foetus* that results in significant monetary losses to the cattle industry in the United States and other parts of the world where open range management and natural breeding are practiced. *T. foetus* is a flagellated protozoan found in bovines that colonizes the vaginal, uterine, oviduct, and preputial epithelium, resulting in embryonic death, abortion, and infertility in the female. Although bulls are the main carriers of *T. foetus*, they remain asymptomatic for their entire lives. The VetMAX™-Gold Trich Detection Kit enables detection of venereal trichomonosis in carrier bulls.

The VetMAX™-Gold Trich Detection Kit can be used to test pools of up to 5 samples for the presence of *T. foetus*. Pool size should be determined by the testing laboratory, based on the prevalence of bovine trichomonosis in the area from which the samples were collected. Pooled samples yielding a positive result should then be tested individually to determine the infection status of each animal in the positive pool.

The assay is a single-well/tube, real-time PCR in which *T. foetus* DNA and Xeno™ DNA Control targets are amplified and detected in real time using fluorescent TaqMan® probes (hydrolysis probe chemistry). The kit includes:

- T. foetus-Xeno™ Control DNA Mix, to serve as a positive control for the real-time PCR components, and it is also used to set the cycle threshold (C<sub>t</sub>) for evaluating test results.
- Xeno™ DNA Control, to serve as an internal control for the DNA isolation process, and it is also used to monitor for the presence of PCR inhibitors.
- T. foetus Primer Probe Mix, optimized for multiplex PCR amplification of both Xeno™ DNA Control and *T. foetus* targets.

### Limitations

- Handle samples as recommended in Table 2 to prevent degradation of any *T. foetus* DNA that is present.
- Prepare pooled samples from no more than 5 individual samples.
- Pooling samples may cause loss of sensitivity of detection of infected animals with an individual C<sub>t</sub> of ≥35.
- DNA extraction methods should yield DNA free of PCR inhibitors, which can prevent amplification of target DNA.
- Follow “Good laboratory practices for PCR and RT-PCR” on page 5 to prevent false positive amplifications due to contamination of test samples with PCR products.

### Kit contents and storage conditions

Reagents for 100 25-µL real-time PCR tests are supplied.

Table 1 VetMAX™-Gold Trich Detection Kit

| Component                                         | Amount   | Storage        |
|---------------------------------------------------|----------|----------------|
| 2X qPCR Master Mix                                | 1.375 mL | –30°C to –10°C |
| T. foetus Primer Probe Mix                        | 110 µL   |                |
| Xeno™ DNA Control (10,000 copies/µL)              | 250 µL   |                |
| T. foetus-Xeno™ Control DNA Mix (1,000 copies/µL) | 80 µL    |                |
| Nuclease-free Water                               | 1.75 mL  | –30°C to 25°C  |

### Required materials not supplied

| Item                                                                                                                                                                                | Source <sup>[1]</sup>                                                          |
|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------------------------------------------------------------------------------|
| <b>Real-time PCR instrument and accessories, one of the following:</b>                                                                                                              |                                                                                |
| 7500 Fast Real-Time PCR System (96-well), running SDS Software v1.4                                                                                                                 | Contact your local sales office.                                               |
| 7500 Fast Precision Plate Holder, for 0.1 mL Tube Strips (Cat. No. A29252), or equivalent                                                                                           |                                                                                |
| QuantStudio™ 5 Real-Time PCR System, 96-well, 0.1-mL                                                                                                                                | Contact your local sales office.                                               |
| <b>Equipment</b>                                                                                                                                                                    |                                                                                |
| Microcentrifuge                                                                                                                                                                     | MLS                                                                            |
| Laboratory mixer (vortex or equivalent)                                                                                                                                             | MLS                                                                            |
| Nuclease-free pipettors                                                                                                                                                             | MLS                                                                            |
| 2 ice buckets: <ul style="list-style-type: none"> <li>• One for the PCR setup area where the master mix is prepared</li> <li>• One for the area where DNA may be present</li> </ul> | MLS                                                                            |
| <b>Plates or tubes and caps</b>                                                                                                                                                     |                                                                                |
| MicroAmp™ Fast Optical 96-Well Reaction Plate with Barcode, 0.1 mL                                                                                                                  | 4366932 (200 plates), 4346906 (20 plates), or equivalent                       |
| MicroAmp™ Fast Optical 96-Well Reaction Plate, 0.1 mL                                                                                                                               | 4346907                                                                        |
| MicroAmp™ Optical Adhesive Film                                                                                                                                                     | 4311971 (100 covers), 4360954 (25 covers), or equivalent                       |
| MicroAmp™ Fast 8-Tube Strip, 0.1 mL                                                                                                                                                 | 4358293, or equivalent                                                         |
| MicroAmp™ Optical 8-Cap Strips                                                                                                                                                      | 4323032, or equivalent                                                         |
| <b>Additional consumables and reagents</b>                                                                                                                                          |                                                                                |
| Filtered pipette tips                                                                                                                                                               | <a href="http://thermofisher.com/pipettetips">thermofisher.com/pipettetips</a> |
| Nuclease-free reagent tubes for preparing master mixes                                                                                                                              | MLS                                                                            |
| 1X Phosphate Buffered Saline (PBS), pH 7.4                                                                                                                                          | MLS                                                                            |
| InPouch™ TF Bovine                                                                                                                                                                  | BioMed Diagnostics 11-1003, 11-1010                                            |

<sup>[1]</sup> Unless otherwise indicated, all materials are available through [thermofisher.com](http://thermofisher.com). MLS: Fisher Scientific ([fisherscientific.com](http://fisherscientific.com)) or other major laboratory supplier.

## Isolate DNA from samples

**Table 2** Sample handling recommendations

| Step or process                                                    | Recommendation                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
|--------------------------------------------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Transport/store samples                                            | <ol style="list-style-type: none"> <li>1. Transport inoculated InPouch™ media to the diagnostic lab, according to the manufacturer's instructions.</li> <li>2. Incubate samples at 37°C for 24 hours.</li> </ol>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
| Prepare cultured smegma samples                                    | <ol style="list-style-type: none"> <li>1. Thoroughly mix the InPouch™ contents.</li> <li>2. Transfer 1 mL from the pouch to a 1.5-mL tube.</li> <li>3. Vigorously vortex the tube.</li> <li>4. Use 300 µL of the cultured sample for DNA isolation.</li> </ol>                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| (Optional) Prepare pooled cultured smegma samples                  | <p>Pool up to 5 individual cultured smegma samples as described below.<sup>[1]</sup></p> <ol style="list-style-type: none"> <li>1. Thoroughly mix the individual cultured smegma sample, and create a 300-µL pool of cultured samples by combining equal volumes of individual samples. For example, transfer 60 µL of 5 individual samples to create one pool.</li> <li>2. Briefly vortex the pooled cultures (properly sealed to prevent cross-contamination) to ensure sufficient mixing, and centrifuge briefly to collect the contents in the lower portion of the tube/plate.</li> <li>3. Use 300 µL of the pooled smegma samples for each DNA purification, and process in the same way as individual samples.</li> </ol> |
| Prepare mock-purified samples (for use in extraction control PCRs) | Prepare duplicate mock-purified samples, using 1X PBS as the starting material. Process mock-purified samples with the same DNA isolation method that is used for test samples.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Proposed DNA isolation method                                      | 5X MagMAX™ Pathogen RNA/DNA Kit (Cat. No. 4462359), or an equivalent DNA isolation method.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                       |
| Required modifications to the DNA isolation method                 | <ul style="list-style-type: none"> <li>• Add 2 µL of undiluted Xeno™ DNA Control (20,000 copies) per purification to the lysis solution used for DNA purification.</li> <li>• Add carrier RNA to the lysis solution according to the manufacturer's recommendation. Add carrier RNA in addition to Xeno™ DNA Control.</li> </ul> <p>Carrier RNA is provided in the 5X MagMAX™ Pathogen RNA/DNA Kit.</p>                                                                                                                                                                                                                                                                                                                          |

<sup>[1]</sup> Pool size is determined by the testing laboratory.

## Perform real-time PCR

1

Determine the quantity of reactions and thaw the reagents

a.

Plan to include the following control reactions on each plate:

- Positive control (prepare duplicate reactions); use 8 μL of T. foetus-Xeno™ Control DNA Mix (1,000 copies/μL) per reaction.
- No-template control (NTC) (prepare duplicate reactions); use 8 μL of Nuclease-free Water in place of sample DNA.

b.

Plan the plate layout so that the wells containing NTCs are located as far as possible from positive controls and test samples to prevent accidental cross-contamination.

c.

Thaw PCR master mix reagents in one ice bucket and controls and samples in a separate ice bucket. Gently vortex each tube to mix the contents thoroughly, then briefly centrifuge to collect the solution at the bottom of the tube. Keep the reagents on ice.

2

Prepare the real-time PCR master mix on ice

Combine the following components for the number of reactions required plus 10% overage.

| Component                                       | Volume per 25 μL reaction |
|-------------------------------------------------|---------------------------|
| 2X qPCR Master Mix                              | 12.5 μL                   |
| T. foetus Primer Probe Mix                      | 1.0 μL                    |
| Nuclease-free Water                             | 3.5 μL                    |
| <b>Total volume of real-time PCR master mix</b> | <b>17.0 μL</b>            |

3

Set up the PCR reactions

a.

Dispense 17 μL of real-time PCR master mix to the appropriate wells of a PCR plate or PCR tubes on ice.

b.

Add the appropriate component for each reaction type, according to the following table:

| Reaction type      | Component                                         | Volume per reaction |
|--------------------|---------------------------------------------------|---------------------|
| Test sample        | Sample DNA                                        | 8.0 μL              |
| NTC                | Nuclease-free Water                               | 8.0 μL              |
| Positive control   | T. foetus-Xeno™ Control DNA Mix (1,000 copies/μL) | 8.0 μL              |
| Extraction control | Mock-purified 1X PBS sample                       | 8.0 μL              |

c.

Seal each reaction vessel, mix, then centrifuge briefly to bring the contents to the bottom.

4

Set up and run the real-time PCR instrument

For detailed information to set up and run the instrument, see the appropriate documentation for your instrument.

a.

Following the manufacturer’s instructions, set up the run using the following parameters:

- Experiment type: Standard curve
- Run mode: Standard
- Reaction volume: 25 μL
- ROX™ passive reference dye: Included in the 2X qPCR Master Mix
- TaqMan® probe reporter dyes and quenchers:

**4** Set up and run the real-time PCR instrument  
(continued)

| Target               | Reporter                | Quencher   |
|----------------------|-------------------------|------------|
| <i>T. foetus</i> DNA | FAM™ dye <sup>[1]</sup> | Eclipse™ Q |
| Xeno™ DNA Control    | VIC™ dye <sup>[2]</sup> | Eclipse™ Q |

<sup>[1]</sup> Absorbance maximum of 495 nm; emission maximum of 520 nm.

<sup>[2]</sup> Absorbance maximum of 540 nm; emission maximum of 552 nm.

- b. Run the thermal cycler program and collect real-time amplification data during stage 2. Use the following thermal cycler settings:

| Stage                |   | Reps. | Temp.          | Time                     |
|----------------------|---|-------|----------------|--------------------------|
| Initial denaturation | 1 | 1     | 95 °C          | 10 minutes               |
| Amplification        | 2 | 40    | 95 °C<br>55 °C | 15 seconds<br>45 seconds |

## Data analysis

See your real-time PCR instrument user guide for instructions on how to analyze your data, using the following method.

**Table 3** Data analysis

| Method                                                            | Details                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                      |
|-------------------------------------------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Use the <b>Control-Based Threshold</b> setting for data analysis. | <ol style="list-style-type: none"> <li>1. Select <b>Manual CT</b>.</li> <li>2. Export <math>\Delta R_n</math> values for the positive control samples (<i>T. foetus</i>-Xeno™ Control DNA Mix).</li> <li>3. Average the FAM™ and VIC™ dye values (separately) for the <math>\Delta R_n</math> at cycle 40 for all replicates of the positive control reaction.</li> <li>4. Set the threshold for the <i>T. foetus</i> (Trich) DNA target reactions at 10% of the average maximum fluorescence value of the <i>T. foetus</i> amplification signal in the positive control reactions (containing approximately 8,000 copies of <i>T. foetus</i>-Xeno™ Control DNA Mix per reaction).<br/>Example: If the average maximum fluorescence value for the <i>T. foetus</i> target in the positive control reactions is 0.15, set the <i>T. foetus</i> threshold at 0.015.</li> <li>5. Repeat step 4 for the Xeno™ DNA Control target using a 10% threshold.<br/>Example: If the average maximum fluorescence value for the Xeno™ DNA Control target in the positive control reactions is 2.0, set the Xeno™ DNA Control threshold at 0.2.</li> </ol> |
| Check the raw fluorescence data.                                  | Verify that increased fluorescence seen in the normalized data is also evident without mathematical data processing.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |

## Interpretation of test results

Verify that your real-time PCR run is valid before analyzing test sample results.

**Table 4** Criteria for a valid real-time PCR run

| Reaction type      | C <sub>t</sub> value for <i>T. foetus</i> DNA | C <sub>t</sub> value for Xeno™ DNA Control |
|--------------------|-----------------------------------------------|--------------------------------------------|
| Positive control   | 25–32                                         | 25–31                                      |
| NTC                | 40 (no signal detected) <sup>[1]</sup>        | 40 (no signal detected) <sup>[1]</sup>     |
| Extraction control | 40 (no signal detected) <sup>[1]</sup>        | 27–34                                      |

<sup>[1]</sup> The run is invalid if the C<sub>t</sub> values for either *T. foetus* or Xeno™ DNA Control targets in the NTC are <40. See “Troubleshooting” on page 5.

**Table 5** Interpretation of test results from individual samples

| C <sub>t</sub> value for <i>T. foetus</i> DNA | C <sub>t</sub> value for Xeno™ DNA Control                 | Interpretation                    |
|-----------------------------------------------|------------------------------------------------------------|-----------------------------------|
| <38                                           | ≤40 <sup>[1]</sup>                                         | <i>T. foetus</i> -positive sample |
| 40 (no signal detected)                       | Xeno™ DNA Control C <sub>t</sub> shift ≤2.5 <sup>[2]</sup> | <i>T. foetus</i> -negative sample |
| ≥38                                           | ≤40 <sup>[2]</sup>                                         | Suspect result <sup>[2]</sup>     |

<sup>[1]</sup> High levels of *T. foetus* DNA in the sample can reduce the signal from the Xeno™ DNA Control, resulting in a higher C<sub>t</sub> value for Xeno™ DNA Control in *T. foetus*-positive samples. See “Troubleshooting” on page 5.

<sup>[2]</sup> See Table 6.

**Table 6** Assessment of suspect results (for individual samples only)

| Result                                                                                                                                                                                                                                                                                                                                                                                                                                                              | Action                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                           |                                                           |                                                           |                                                                                                                                                                                                                                                                                                                                                                                                                                                                     |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------------------------------------------------------|-----------------------------------------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Suspect result 1:</b> The sample <i>T. foetus</i> C <sub>t</sub> value is <40 and the C <sub>t</sub> value for one or more NTCs or extraction controls is <40.                                                                                                                                                                                                                                                                                                   | Repeat the extraction or real-time PCR. See “No template control reaction:” on page 5.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                           |                                                           |                                                           |                                                                                                                                                                                                                                                                                                                                                                                                                                                                     |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          |
| <b>Suspect result 2:</b> A <i>T. foetus</i> (TF) sample is considered suspect if the TF C <sub>t</sub> value is 38–40.                                                                                                                                                                                                                                                                                                                                              | <p>Analyze suspect DNA samples for the presence/absence of PCR inhibitors by calculating the Xeno™ DNA Control C<sub>t</sub> Shift:</p> <p><b>Xeno™ DNA Control C<sub>t</sub> Shift = SS – XEC</b>, where:</p> <p><b>SS</b> = C<sub>t</sub> of Xeno™ DNA Control in the suspect sample</p> <p><b>XEC</b> = Average C<sub>t</sub> of Xeno™ DNA Control in the extraction controls</p> <table border="1"> <thead> <tr> <th>Workflow A<br/>Xeno™ DNA Control C<sub>t</sub> shift &lt;2.5</th><th>Workflow B<br/>Xeno™ DNA Control C<sub>t</sub> shift ≥2.5</th></tr> </thead> <tbody> <tr> <td> <ol style="list-style-type: none"> <li>Repeat the DNA purification on triplicate aliquots of the original diagnostic sample.</li> <li>Repeat the real-time PCR with 8 µL of purified DNA from step 1.</li> <li>Determine the number of samples with a C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol> </td><td> <ol style="list-style-type: none"> <li>Repeat the real-time PCR with 2 µL of the suspect DNA sample. (PCR inhibitors may be present in the DNA.) If the <i>T. foetus</i> C<sub>t</sub> value is: <ul style="list-style-type: none"> <li>&lt;38: The sample is <i>T. foetus</i> positive. No further testing is required.<sup>[1]</sup></li> <li>≥38: Continue with steps 2–5 of this workflow.</li> </ul> </li> <li>Dilute the original individual diagnostic sample 1:4 with 1X PBS.</li> <li>Repeat the DNA purification on triplicate aliquots of the 1:4 diluted sample.</li> <li>Repeat the real-time PCR with 8 µL of the purified DNA from step 3.</li> <li>Determine the number of samples with a <i>T. foetus</i> C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol> </td></tr> </tbody> </table> | Workflow A<br>Xeno™ DNA Control C <sub>t</sub> shift <2.5 | Workflow B<br>Xeno™ DNA Control C <sub>t</sub> shift ≥2.5 | <ol style="list-style-type: none"> <li>Repeat the DNA purification on triplicate aliquots of the original diagnostic sample.</li> <li>Repeat the real-time PCR with 8 µL of purified DNA from step 1.</li> <li>Determine the number of samples with a C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol> | <ol style="list-style-type: none"> <li>Repeat the real-time PCR with 2 µL of the suspect DNA sample. (PCR inhibitors may be present in the DNA.) If the <i>T. foetus</i> C<sub>t</sub> value is: <ul style="list-style-type: none"> <li>&lt;38: The sample is <i>T. foetus</i> positive. No further testing is required.<sup>[1]</sup></li> <li>≥38: Continue with steps 2–5 of this workflow.</li> </ul> </li> <li>Dilute the original individual diagnostic sample 1:4 with 1X PBS.</li> <li>Repeat the DNA purification on triplicate aliquots of the 1:4 diluted sample.</li> <li>Repeat the real-time PCR with 8 µL of the purified DNA from step 3.</li> <li>Determine the number of samples with a <i>T. foetus</i> C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol> |
| Workflow A<br>Xeno™ DNA Control C <sub>t</sub> shift <2.5                                                                                                                                                                                                                                                                                                                                                                                                           | Workflow B<br>Xeno™ DNA Control C <sub>t</sub> shift ≥2.5                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                        |                                                           |                                                           |                                                                                                                                                                                                                                                                                                                                                                                                                                                                     |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          |
| <ol style="list-style-type: none"> <li>Repeat the DNA purification on triplicate aliquots of the original diagnostic sample.</li> <li>Repeat the real-time PCR with 8 µL of purified DNA from step 1.</li> <li>Determine the number of samples with a C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol> | <ol style="list-style-type: none"> <li>Repeat the real-time PCR with 2 µL of the suspect DNA sample. (PCR inhibitors may be present in the DNA.) If the <i>T. foetus</i> C<sub>t</sub> value is: <ul style="list-style-type: none"> <li>&lt;38: The sample is <i>T. foetus</i> positive. No further testing is required.<sup>[1]</sup></li> <li>≥38: Continue with steps 2–5 of this workflow.</li> </ul> </li> <li>Dilute the original individual diagnostic sample 1:4 with 1X PBS.</li> <li>Repeat the DNA purification on triplicate aliquots of the 1:4 diluted sample.</li> <li>Repeat the real-time PCR with 8 µL of the purified DNA from step 3.</li> <li>Determine the number of samples with a <i>T. foetus</i> C<sub>t</sub> value &lt;40: <ul style="list-style-type: none"> <li>0 of 3: <i>T. foetus</i> negative</li> <li>1 of 3: Presumptive positive</li> <li>≥2 of 3: <i>T. foetus</i> positive</li> </ul> </li> </ol>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |                                                           |                                                           |                                                                                                                                                                                                                                                                                                                                                                                                                                                                     |                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                          |

<sup>[1]</sup> The Xeno™ DNA Control C<sub>t</sub> value on diluted samples is not considered utilizing this workflow.

**Table 7** Interpretation of test results from pooled samples

| C <sub>t</sub> value for <i>T. foetus</i> DNA | C <sub>t</sub> value for Xeno™ DNA Control  | Interpretation                  | Action                                                                                                                                                                          |
|-----------------------------------------------|---------------------------------------------|---------------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <40 (signal detected)                         | <40 (signal detected)                       | <i>T. foetus</i> -positive pool | Retest individual samples from the positive pool to determine which individual sample or samples caused the positive result. Interpret the retest results according to Table 5. |
| 40 (no signal detected)                       | Xeno™ DNA Control C <sub>t</sub> shift ≤2.5 | <i>T. foetus</i> -negative pool | None required                                                                                                                                                                   |
|                                               | Xeno™ DNA Control C <sub>t</sub> shift ≥2.5 | Suspect result                  | Repeat the DNA purification with the individual cultured smegma samples, then repeat the real-time PCR with each DNA sample. Interpret the retest results according to Table 5. |

## Troubleshooting

| Observation                                                                                                   | Possible cause                                                                                                                                                                                  | Recommended action                                                                                                                                                                                                                                                                                                                                                                |
|---------------------------------------------------------------------------------------------------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Positive control reaction:</b><br><i>T. foetus</i> DNA—no signal<br>Xeno™ DNA Control—no signal            | The <i>T. foetus</i> -Xeno™ Control DNA Mix was improperly handled, resulting in DNA degradation.                                                                                               | Use appropriate precautions against DNase contamination when handling the control DNAs. For example, wear clean gloves and use nuclease-free barrier pipette tips.                                                                                                                                                                                                                |
|                                                                                                               | The 2X qPCR Master Mix was stored or handled improperly and it lost activity.                                                                                                                   | Repeat the real-time PCR with fresh reagents.                                                                                                                                                                                                                                                                                                                                     |
|                                                                                                               | The thermal cycler was not properly set up.                                                                                                                                                     | Check the thermal cycler settings. See “Set up and run the real-time PCR instrument” on page 2.                                                                                                                                                                                                                                                                                   |
|                                                                                                               | The real-time PCR master mix was prepared incorrectly.                                                                                                                                          | Repeat the test with correctly prepared real-time PCR master mix.                                                                                                                                                                                                                                                                                                                 |
| <b>No template control reaction:</b><br>Signal detected [ $C_t$ value is <40]                                 | There was contamination during the PCR.                                                                                                                                                         | <ul style="list-style-type: none"> <li>Repeat the real-time PCR with fresh reagents and freshly decontaminated pipettes.</li> <li>Set up the real-time PCR in an area separate from areas used for DNA isolation and PCR product analysis.</li> </ul>                                                                                                                             |
| <b>Extraction control reaction:</b><br>Signal detected [ $C_t$ value is <40]                                  | There was contamination during the DNA isolation or PCR.                                                                                                                                        | <ul style="list-style-type: none"> <li>Repeat the DNA isolation or real-time PCR with fresh reagents and freshly decontaminated pipettes.</li> <li>Set up the real-time PCR in an area separate from areas used for DNA isolation and PCR product analysis.</li> </ul>                                                                                                            |
| <b>Test samples</b><br>Xeno™ DNA Control—no or low signal<br><i>T. foetus</i> DNA—high signal                 | The Xeno™ DNA Control primers and probe are at limiting concentrations in the real-time PCR. High levels of <i>T. foetus</i> DNA in a sample can reduce amplification of the Xeno™ DNA Control. | No or low signal from Xeno™ DNA Control is expected in a reaction that has a strong signal for <i>T. foetus</i> DNA.                                                                                                                                                                                                                                                              |
| <b>Test samples:</b><br>Xeno™ DNA Control—no signal<br><i>T. foetus</i> DNA—no signal or suspect-range signal | Poor DNA recovery.                                                                                                                                                                              | <ul style="list-style-type: none"> <li>Check the <math>C_t</math> values of Xeno™ DNA Control in the mock-purified samples.</li> <li>A <math>C_t</math> value <math>\geq 38</math> in individual or pooled samples indicates that Xeno™ DNA Control was omitted or that DNA recovery was poor.</li> <li>Repeat the DNA purification of the original diagnostic sample.</li> </ul> |
|                                                                                                               | The DNA samples contain PCR inhibitors.                                                                                                                                                         | See Table 6.                                                                                                                                                                                                                                                                                                                                                                      |

## Explanation of symbols

The symbols present on the product label are explained in the following table.

|                                                                                    |                |                                                                                     |                              |
|------------------------------------------------------------------------------------|----------------|-------------------------------------------------------------------------------------|------------------------------|
|  | MANUFACTURER   |  | USE BY                       |
|  | CATALOG NUMBER |  | CONSULT INSTRUCTIONS FOR USE |
|  | BATCH CODE     |  | CAUTION                      |
|  | SERIAL NUMBER  |  | TEMPERATURE LIMIT            |

## Good laboratory practices for PCR and RT-PCR

When preparing samples for PCR or RT-PCR amplification:

- Wear clean gloves and a clean lab coat.
  - Do not wear the same gloves and lab coat that you have previously used when handling amplified products or preparing samples.
- Change gloves if you suspect that they are contaminated.
- Maintain separate areas and dedicated equipment and supplies for:
  - Sample preparation and reaction setup.
  - Amplification and analysis of products.
- Do not bring amplified products into the reaction setup area.
- Open and close all sample tubes carefully. Avoid splashing or spraying samples.
- Keep reactions and components capped as much as possible.
- Use a positive-displacement pipettor or aerosol-resistant barrier pipette tips.
- Clean lab benches and equipment periodically with 10% bleach solution or DNA decontamination solution.

## Documentation and support

### Customer and technical support

In the United States, call 1-800-955-6288.

Visit [thermofisher.com/support](http://thermofisher.com/support) for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support
- Order and web support

- Product documentation

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



**Manufacturer:** Life Technologies Corporation | 2130 Woodward Street | Austin, TX 78744

VLN 432 / PCN 5292.80

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** Pub. No. 4485468

| Revision | Date              | Description                                                                                                                                                                                                                 |
|----------|-------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| B        | 2 August 2018     | <ul style="list-style-type: none"><li>• Updated the list of compatible real-time PCR systems.</li><li>• Updated to the current document template, with associated updates to the warranty, trademarks, and logos.</li></ul> |
| A        | 18 September 2013 | Baseline for this revision history.                                                                                                                                                                                         |

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Eclipse is a trademark of Elitech Holding BV. InPouch is a trademark of VWR International, Inc. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license.