

PrepSEQ™ Sample Preparation Kits for *Mycoplasma*, MMV, and Vesivirus

Manual sample preparation protocols for *Mycoplasma*, MMV, and Vesivirus detection

Catalog Numbers 4452222 and 4443789

Pub. No. 4465875 Rev. C

Note: For safety and biohazard guidelines, see the “Safety” appendix in the *PrepSEQ™ Sample Preparation Kits for Mycoplasma, MMV, and Vesivirus User Guide* (Pub. No. 4465957). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

- Product description 1
- Kit applications 1
- Before you begin 2
- Prepare materials 2
- 1-2-3 manual protocol for *Mycoplasma* and/or MMV detection 2
- 3-in-1 manual protocol for *Mycoplasma*, MMV, and/or Vesivirus detection 4
- Large-scale manual protocol for *Mycoplasma* detection—Option 1: Direct sample testing 5
- Large-scale manual protocol for *Mycoplasma* detection—Option 2: Process pooled cell culture media and mammalian cells 6
- Large-scale protocol for *Mycoplasma* detection—Options 1 and 2: Extract the DNA 8

Product description

The PrepSEQ™ Sample Preparation Kits use Magnetic Particle-based separation technology for manual extraction of DNA and/or RNA from *Mycoplasma* cells or viral particles. A variety of starting material can be used, such as infected cell cultures or *Mycoplasma* liquid cultures. The kits described in this document are:

- PrepSEQ™ 1-2-3 Nucleic Acid Extraction Kit (Cat. No. 4452222)
- PrepSEQ™ *Mycoplasma* Nucleic Acid Extraction Kit (Cat. No. 4443789)

Kit applications

Organisms	Sample volume	For use with kit	Protocol
<i>Mycoplasma</i> and MMV	100 µL (up to 10 ⁶ cells)	PrepSEQ™ 1-2-3 Nucleic Acid Extraction Kit	“1-2-3 manual protocol for Mycoplasma and/or MMV detection” on page 2
<i>Mycoplasma</i> , MMV, and Vesivirus	100 µL (up to 10 ⁶ cells)	PrepSEQ™ 1-2-3 Nucleic Acid Extraction Kit	“3-in-1 manual protocol for Mycoplasma, MMV, and/or Vesivirus detection” on page 4
<i>Mycoplasma</i>	Up to 11 mL (up to 2 × 10 ⁸ cells)	PrepSEQ™ <i>Mycoplasma</i> Nucleic Acid Extraction Kit	“Large-scale manual protocol for Mycoplasma detection—Option 1: Direct sample testing” on page 5 <i>OR</i> “Large-scale manual protocol for Mycoplasma detection—Option 2: Process pooled cell culture media and mammalian cells” on page 6

Before you begin

Before you begin, review the *PrepSEQ™ Sample Preparation Kits for Mycoplasma, MMV, and Vesivirus User Guide* (Pub. No. 4465957) for required materials and procedural guidelines.

Note: This Quick Reference provides brief procedures for sample preparation. For detailed procedures, see the User Guide.

Prepare materials

- (Before first use of the kit)* Prepare the following reagents:
 - **Binding Solution**—Add 30 mL of 100% isopropanol to the empty Binding Solution bottle. Mark the bottle label to indicate that isopropanol has been added.
 - **Wash Buffer**—Add 74 mL of 95% non-denatured ethanol to the Wash Buffer Concentrate bottle, then mix well. Mark the bottle label to indicate that ethanol has been added.
- Power on a heat block to 37°C.
- Incubate the Magnetic Particles suspension at 37°C for a minimum of 10 minutes with intermittent vortexing at setting #7, or until the particles are completely suspended.
- (Large-scale protocols only)* Power on the refrigerated centrifuge to allow it to cool to 2–8°C before use.

1-2-3 manual protocol for *Mycoplasma* and/or MMV detection

Use this protocol to process 100 µL of sample (up to 10⁶ cells) for the detection of *Mycoplasma* and/or MMV.

-
- 1 Prepare test samples**
 - Prepare each sample in a new 2-mL microcentrifuge tube:
 - ≤10⁶ total cells—Add 100 µL of sample to the tube.
 - >10⁶ total cells—Spin the sample in a microcentrifuge at 500 × g for 2 minutes, then add 100 µL of supernatant to the tube.
 - (Optional)* For an extraction positive control: Spike the appropriate amount of DPC to 100 µL of the prepared test sample to achieve 200 copies per PCR. Extraction controls can be dual- or triple-spiked with DPCs for each assay to conserve sample. You can also use one of the following in place of test samples: 1X PBS solution, Lysis Buffer, Cell Fractionation Buffer, or plain media.
-
- 2 Prepare sample lysate**

For each sample tube:

 - Add 200 µL of Lysis Buffer, then vortex for ~5 seconds to mix.
 - Add 2 µL of 0.5 M EDTA, 18 µL of RNase Cocktail™ Enzyme Mix, and 2 µL of Proteinase K, then briefly vortex to mix.
 - Incubate at 56°C for a minimum of 15 minutes.
 - Incubate at room temperature for 5 minutes.
 - Add 700 µL of Lysis Buffer.

3 Bind DNA

For each sample lysate tube:

- a. Add 30 μL of Magnetic Particles, add 525 μL of Binding Solution, then immediately invert each tube to mix.
- b. Using a vortex adaptor, vortex the tubes vertically at medium speed for 5 minutes to capture the nucleic acid.
- c. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed (13,000 $\times g$ or greater).
- d. Place in the Magnetic Stand for 5 minutes.
- e. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.

4 Wash DNA

For each tube of Magnetic Particles pellet (bound DNA):

- a. Add 300 μL of Wash Buffer, then vortex for ~5 seconds.
- b. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed (13,000 $\times g$ or greater).
- c. Place in the Magnetic Stand for 1 minute.
- d. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.
- e. Repeat step 4a through step 4d.
Set the timer for 5 minutes just before you remove the wash buffer for the second time. See step 4g.
- f. Use a P200 pipette to aspirate the residual supernatant from the bottom of the tube, then discard the supernatant.
- g. With the cap open, air-dry the Magnetic Particles pellet at room temperature for ≤ 5 minutes to remove any residual ethanol.

IMPORTANT! Set the timer for 5 minutes just before you remove the wash buffer for the second time. Do NOT dry longer than 5 minutes. Overdrying will make the pellets difficult to resuspend in Elution Buffer in the next steps.

IMPORTANT! If you are using an aspirator or processing in a BSC (hood), no additional drying time is required. Aspirate and proceed; do not wait ≤ 5 minutes.

5 Elute DNA

For each sample:

- a. Add 100 μL of Elution Buffer.
- b. Vortex for ~10 seconds.
- c. Incubate at 70°C for 7 minutes.
Vortex 2–3 times during incubation to ensure complete resuspension of the Magnetic Particles. The total time may be 10–12 minutes.
- d. Centrifuge for 5 minutes at top speed (13,000 $\times g$ or greater).
- e. Place in the Magnetic Stand for 3 minutes.
- f. Transfer the eluate to a non-stick 1.5-mL microcentrifuge tube.

3-in-1 manual protocol for *Mycoplasma*, MMV, and/or Vesivirus detection

Use this protocol to process 100 μL of sample (up to 10^6 cells) for the detection of *Mycoplasma*, MMV, and/or Vesivirus.

1 Prepare test samples

- a. Prepare each sample in a new 2-mL microcentrifuge tube:
 - $\leq 1 \times 10^6$ total cells—Add 100 μL of sample to the tube.
 - $> 1 \times 10^6$ total cells—Spin the sample in a microcentrifuge at $500 \times g$ for 2 minutes, then add 100 μL of supernatant to the tube.
 - b. (*Optional*) For an extraction positive control: Spike the appropriate amount of DPC to 100 μL of the prepared test sample to achieve 200 copies per PCR. Extraction controls can be dual- or triple-spiked with DPCs for each assay to conserve sample. You can also use one of the following in place of test samples: 1X PBS solution, Lysis Buffer, Cell Fractionation Buffer, or plain media.
-

2 Prepare sample lysate

For each sample tube:

- a. Add 500 μL of Lysis Buffer, then vortex for ~15 seconds to mix.
 - b. Incubate at 45°C for 10 minutes.
 - c. Vortex for ~10 seconds to mix.
-

3 Bind DNA

For each sample lysate tube:

- a. Add 30 μL of Magnetic Particles.
 - b. Add 330 μL of Binding Solution, then immediately invert each tube to mix.
 - c. Using a vortex adaptor, vortex the tube vertically at medium speed for 5 minutes to capture the nucleic acid.
 - d. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed ($13,000 \times g$ or greater).
 - e. Place in the Magnetic Stand for 5 minutes.
 - f. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.
-

4 Wash DNA

For each tube of Magnetic Particles pellet (bound DNA):

- a. Add 300 μL of Wash Buffer, then vortex for ~5 seconds.
- b. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed ($13,000 \times g$ or greater).
- c. Place in the Magnetic Stand for 1 minute.
- d. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.
- e. Repeat step 4a through step 4d.
Set the timer for 5 minutes just before you remove the wash buffer for the second time. See step 4g.
- f. Use a P200 pipette to aspirate the residual supernatant from the bottom of the tube, then discard the supernatant.

4 Wash DNA
(continued)

- g. With the cap open, air-dry the Magnetic Particles pellet at room temperature for ≤ 5 minutes to remove any residual ethanol.

IMPORTANT! Set the timer for 5 minutes just before you remove the wash buffer for the second time. Do NOT dry longer than 5 minutes. Overdrying will make the pellets difficult to resuspend in Elution Buffer in the next steps.

IMPORTANT! If you are using an aspirator or processing in a BSC (hood), no additional drying time is required. Aspirate and proceed; do not wait ≤ 5 minutes.

5 Elute DNA

For each sample:

- a. Add 100 μL of Elution Buffer.
- b. Vortex for ~ 10 seconds.
- c. Incubate at 70°C for 5 minutes.
Vortex 2–3 times during incubation to ensure complete resuspension of the Magnetic Particles. The total time may be 10–12 minutes.
- d. Centrifuge for 5 minutes at top speed ($13,000 \times g$ or greater).
- e. Place in the Magnetic Stand for 3 minutes.
- f. Transfer the eluate to a non-stick 1.5-mL microcentrifuge tube.

Large-scale manual protocol for *Mycoplasma* detection—Option 1: Direct sample testing

Use this protocol to process up to 11 mL of sample (up to 10^8 cells) for the detection of *Mycoplasma*.

1 Prepare test samples

Prepare each sample in a new 50-mL conical tube:

- $\leq 10^8$ cells—Add 11 mL of sample to the tube.
- $> 10^8$ cells—Add 15 mL of sample to the tube, centrifuge at $1,000 \times g$ for 5 minutes to pellet the cells, then transfer 11 mL of supernatant to a new 50-mL conical tube.

2 Separate mammalian cells from cell culture media

- a. Centrifuge each tube at $1,000 \times g$ for 5 minutes at 4°C to pellet the mammalian cells.
- b. Transfer 10 mL of the supernatant to a new 50-mL conical tube, then place on ice.
The supernatant contains free *Mycoplasma*.
- c. Discard the mammalian cell pellet.

3 Treat with RNase and DNase

If the samples have high SYBR™ Green I dye background because of excess cellular nucleic acid in the supernatant, perform RNase treatment.

IMPORTANT! For some high-density samples, both RNase and DNase treatments are needed. In this case, perform the DNase treatment before the RNase treatment.

- a. To treat with DNase:
 1. Add 450 μL of $10\times$ Reaction Buffer and 90 μL of TURBO™ DNase, then gently vortex to mix.
 2. Incubate at 37°C for 30 minutes.

-
- 3** Treat with RNase and DNase (*continued*)
- b. To treat with RNase:
 1. Add 180 μL of 0.5 M EDTA, 225 μL of RNase Cocktail™ Enzyme Mix, and 150 μL of Proteinase K, gently vortex to mix, then briefly spin.
 2. Incubate at 56°C for a minimum of 30 minutes to digest the cellular RNA and proteins.
-
- 4** Process the supernatant to obtain resuspended *Mycoplasma*
- a. Centrifuge the supernatant at 16,000 $\times g$ for 30 minutes at 4°C to pellet the *Mycoplasma*.
 - b. Aspirate and discard the supernatant without disturbing the *Mycoplasma* pellet. Do not decant the liquid and do NOT touch the pellet.
 - c. Add 300 μL of Lysis Buffer, then mix thoroughly by vortexing to resuspend the *Mycoplasma* pellet.
If the pellet is difficult to dislodge, vigorously agitate the tube.
 - d. Transfer the resuspended pellet to a 2-mL microcentrifuge tube.
-
- 5** Treat the resuspended *Mycoplasma*
- Separately process the resuspended *Mycoplasma* in the 2-mL microcentrifuge tube.
- a. Add 2 μL of 0.5 M EDTA, 18 μL of RNase Cocktail™ Enzyme Mix, and 5 μL of Proteinase K, then briefly vortex to mix.
 - b. (*Optional*) Spike with Discriminatory Positive Control (DPC) for optimization or with *Mycoplasma* DNA for lot release validation.
 - c. (*Optional*) For an extraction positive control: Spike the appropriate amount of DPC to 300 μL of the prepared test sample to achieve 200 copies per PCR. Extraction controls can be dual- or triple-spiked with DPCs for each assay to conserve sample. You can also use one of the following in place of test samples: 1X PBS solution, Lysis Buffer, Cell Fractionation Buffer, or plain media.
 - d. Incubate at 56°C for 15 minutes to digest the cellular RNA. Vortex twice during incubation.
 - e. Add 700 μL of Lysis Buffer.
- Proceed directly to “Large-scale protocol for *Mycoplasma* detection—Options 1 and 2: Extract the DNA” on page 8.
-

Large-scale manual protocol for *Mycoplasma* detection—Option 2: Process pooled cell culture media and mammalian cells

Use this protocol to process up to 11 mL of sample (up to 10^8 cells) for the detection of *Mycoplasma*.

- 1** Prepare test samples
- Prepare each sample in a new 50-mL conical tube:
- $\leq 10^8$ cells—Add 11 mL of sample to the tube.
 - $> 10^8$ cells—Add 15 mL of sample to the tube, centrifuge at 1,000 $\times g$ for 5 minutes to pellet the cells, then transfer 11 mL of supernatant to a new 50-mL conical tube.
-
- 2** Separate mammalian cells from cell culture media
- a. Centrifuge each tube at 1,000 $\times g$ for 5 minutes at 4°C to pellet the mammalian cells.
 - b. Transfer 10 mL of the supernatant to a new 50-mL conical tube, then place on ice.
The supernatant contains free *Mycoplasma*.
 - c. Remove residual supernatant from the mammalian cell pellet, then place the cell pellet on ice.

3 Treat with RNase and DNase

If the samples have high SYBR™ Green I dye background because of excess cellular nucleic acid in the supernatant, perform RNase treatment.

IMPORTANT! For some high-density samples, both RNase and DNase treatments are needed. In this case, perform the DNase treatment before the RNase treatment.

- a. To treat with DNase:
 1. Add 450 µL of 10× Reaction Buffer and 90 µL of TURBO™ DNase, then gently vortex to mix.
 2. Incubate at 37°C for 30 minutes.
- b. To treat with RNase:
 1. Add 180 µL of 0.5 M EDTA, 225 µL of RNase Cocktail™ Enzyme Mix, and 150 µL of Proteinase K, gently vortex to mix, then briefly spin.
 2. Incubate at 56°C for a minimum of 30 minutes to digest the cellular RNA and proteins.

4 Process the supernatant to obtain the *Mycoplasma* pellet

- a. Centrifuge the supernatant at 16,000 × g for 30 minutes at 4°C to pellet the *Mycoplasma*.
- b. Carefully remove and discard the supernatant; retain the *Mycoplasma* pellet for use in the next section.

IMPORTANT! Do not decant the liquid and do NOT touch the pellet. Use a P200 pipette to remove the last of the supernatant.

- c. Place the 50-mL tube containing the *Mycoplasma* pellet on ice.

5 Process the mammalian cell pellet to obtain free *Mycoplasma* and combine with the *Mycoplasma* pellet

Perform this procedure during the 30-minute centrifugation step in the previous section.

- a. Add 550 µL of ice-cold Cell Fractionation Buffer to the mammalian cell pellet. Gently vortex or pipet up and down several times with a P1000 pipette to completely resuspend the mammalian cells.
If the pellet is difficult to dislodge, vigorously agitate the tube.
- b. Transfer the mammalian cell suspension to a 2-mL microcentrifuge tube, then place on ice for 5 minutes.
- c. Centrifuge the 2-mL tube at 1,500 × g for 10 minutes at 4°C to pellet the cellular membranes and nuclei.
- d. Carefully transfer 300 µL (two 150-µL aliquots) of the Cell Fractionation Buffer supernatant (mammalian cell lysate) to the *Mycoplasma* pellet obtained in the previous section. Avoid the pellet and viscous material.
- e. Resuspend the *Mycoplasma* pellet in the supernatant by pipetting up and down or by vortexing on medium speed.
- f. Transfer the resuspended *Mycoplasma* pellet to a new 2-mL microcentrifuge tube.

6 Treat the resuspended *Mycoplasma*

Separately process the resuspended *Mycoplasma* in the 2-mL microcentrifuge tube.

- a. Add 2 μL of 0.5 M EDTA, 18 μL of RNase Cocktail™ Enzyme Mix, and 5 μL of Proteinase K, then briefly vortex to mix.
- b. (Optional) Spike with Discriminatory Positive Control (DPC) for optimization or with *Mycoplasma* DNA for lot release validation.
- c. (Optional) For an extraction positive control: Spike the appropriate amount of DPC to 300 μL of the prepared test sample to achieve 200 copies per PCR. Extraction controls can be dual- or triple-spiked with DPCs for each assay to conserve sample. You can also use one of the following in place of test samples: 1X PBS solution, Lysis Buffer, Cell Fractionation Buffer, or plain media.
- d. Incubate at 56°C for 15 minutes to digest the cellular RNA. Vortex twice during incubation.
- e. Add 700 μL of Lysis Buffer.

Proceed directly to “Large-scale protocol for *Mycoplasma* detection—Options 1 and 2: Extract the DNA” on page 8.

Large-scale protocol for *Mycoplasma* detection—Options 1 and 2: Extract the DNA

1 Bind DNA

For each sample lysate tube:

- a. Add 30 μL of Magnetic Particles, add 525 μL of Binding Solution, then immediately invert each tube to mix.
 - b. Using a vortex adaptor, vortex the tubes vertically at medium speed for 5 minutes to capture the nucleic acid.
 - c. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed (13,000 $\times g$ or greater).
 - d. Place in the Magnetic Stand for 5 minutes.
 - e. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.
-

2 Wash DNA

For each tube of Magnetic Particles pellet (bound DNA):

- a. Add 300 μL of Wash Buffer, then vortex for ~5 seconds.
- b. Spin in a microcentrifuge: Press the short spin button for 15 seconds, or until the microcentrifuge reaches top speed (13,000 $\times g$ or greater).
- c. Place in the Magnetic Stand for 1 minute.
- d. Aspirate the supernatant without disturbing the Magnetic Particles, then discard the supernatant.
- e. Repeat step 2a through step 2d.
Set the timer for 5 minutes just before you remove the wash buffer for the second time. See step 2g.
- f. Use a P200 pipette to aspirate the residual supernatant from the bottom of the tube, then discard the supernatant.

2 Wash DNA (continued)

- g. With the cap open, air-dry the Magnetic Particles pellet at room temperature for ≤5 minutes to remove any residual ethanol.

IMPORTANT! Set the timer for 5 minutes just before you remove the wash buffer for the second time. Do NOT dry longer than 5 minutes. Overdrying will make the pellets difficult to resuspend in Elution Buffer in the next steps.

IMPORTANT! If you are using an aspirator or processing in a BSC (hood), no additional drying time is required. Aspirate and proceed; do not wait ≤5 minutes.

3 Elute DNA

For each sample:

- a. Add 100 µL of Elution Buffer.
- b. Vortex for ~10 seconds.
- c. Incubate at 70°C for 7 minutes.
Vortex 2–3 times during incubation to ensure complete resuspension of the Magnetic Particles. The total time may be 10–12 minutes.
- d. Centrifuge for 5 minutes at top speed (13,000 × g or greater).
- e. Place in the Magnetic Stand for 3 minutes.
- f. Transfer the eluate to a non-stick 1.5-mL microcentrifuge tube.



Manufacturer: Thermo Fisher Scientific | 7 Kingsland Grange | Warrington, Cheshire WA1 4SR | United Kingdom

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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. 4465875

Revision	Date	Description
C	30 January 2018	Update the following protocols: <ul style="list-style-type: none">• 1-2-3 manual protocol for <i>Mycoplasma</i> and/or MMV detection• 3-in-1 manual protocol for <i>Mycoplasma</i>, MMV, and/or Vesivirus detection• Large-scale manual protocol for <i>Mycoplasma</i> detection
B	01 April 2015	Update storage temperature for the Magnetic Particles. Remove Cat. No. 4460627.
A	August 2011	New document.

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.