

# MagMAX™ CORE Mechanical Lysis Module

## USER GUIDE

Automated purification of high-quality nucleic acid from *Mycobacterium avium* subspecies *paratuberculosis* (MAP) and other difficult-to-lyse samples

for use with:

MagMAX™ CORE Nucleic Acid Purification Kit

KingFisher™ Flex Purification System

MagMAX™ Express-96 Deep Well Magnetic Particle Processor

KingFisher™ Duo Prime Purification System

KingFisher™ mL Purification System

Catalog Numbers A32836, A37489, A37487, A37488

Publication Number MAN0015945

Revision A.0



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

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. MAN0015945

| Revision | Date             | Description  |
|----------|------------------|--------------|
| A        | 16 November 2017 | 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.

**TRADEMARKS:** All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2017 Thermo Fisher Scientific Inc. All rights reserved.

# Contents

|   |                   |   |    |
|---|-------------------|---|----|
| ■ | <b>CHAPTER 1</b>  | <b>Product information</b>  | 5  |
|   |                   | Product description   | 5  |
|   |                   | Contents and storage  | 5  |
|   |                   | Required materials not supplied   | 6  |
|   |                   | Workflow  | 8  |
| ■ | <b>CHAPTER 2</b>  | <b>Before you begin</b>   | 9  |
|   |                   | Procedural guidelines   | 9  |
|   |                   | Before first use of the kit   | 9  |
|   |                   | <i>(Optional)</i> Determine the optimal bead mill homogenizer settings          | 9  |
|   |                   | Download the script   | 10 |
| ■ | <b>CHAPTER 3</b>  | <b>Methods</b>  | 11 |
|   |                   | Set up the processing plates  | 11 |
|   |                   | Prepare Lysis/Binding Mix   | 12 |
|   |                   | Prepare the sample  | 13 |
|   |                   | High-input workflow   | 13 |
|   |                   | Low-input workflow  | 15 |
|   |                   | Combine the lysate with Proteinase K and Lysis/Binding Mix                      | 16 |
|   |                   | Process samples on the instrument   | 16 |
| ■ | <b>APPENDIX A</b> | <b>Troubleshooting</b>  | 17 |
| ■ | <b>APPENDIX B</b> | <b>Purification with the KingFisher™ Duo Prime or KingFisher™ mL instrument</b> | 18 |
|   |                   | Required materials not supplied   | 18 |
|   |                   | Purification procedure  | 19 |
| ■ | <b>APPENDIX C</b> | <b>Safety</b>   | 20 |
|   |                   | Chemical safety   | 21 |
|   |                   | Biological hazard safety  | 22 |

|  |           |
|--|-----------|
| <b>Documentation and support .....</b> | <b>23</b> |
| Related documentation .....            | 23        |
| Customer and technical support .....   | 23        |
| Limited product warranty .....         | 23        |



# Product information

---

**IMPORTANT!** Before using this product, read and understand the information in the “Safety” appendix in this document.

---

## Product description

The MagMAX™ CORE Mechanical Lysis Module is a supplemental module for use with the MagMAX™ CORE Nucleic Acid Purification Kit (Cat. Nos. A32700 and A32702). The module is designed for processing tough-to-lyse bacteria in complex sample matrices such as feces. For example, DNA from *Mycobacterium avium* subspecies *paratuberculosis* (MAP), known to cause Johne's Disease in cattle, can be isolated from bovine feces using the procedures that are described in this guide.

The MagMAX™ CORE Nucleic Acid Purification Kit uses a simple magnetic separation process in preparation for downstream molecular analysis, and has been optimized for use with KingFisher™ instruments.

## Contents and storage

Table 1 MagMAX™ CORE Mechanical Lysis Module (Cat. No. A32836)

| Contents                         | Amount <sup>[1]</sup> | Storage |
|----------------------------------|-----------------------|---------|
| MagMAX™ CORE Clarifying Solution | 45 mL                 | 15–30°C |
| MagMAX™ CORE Bead Beating Tubes  | 100 tubes (2 mL)      |         |

<sup>[1]</sup> Also available in combination with the MagMAX™ CORE Nucleic Acid Purification Kit (Cat. No. A37487)

## Required materials not supplied

Unless otherwise indicated, all materials are available through **thermofisher.com**.  
MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

**Table 2** MagMAX™ CORE Nucleic Acid Purification Kit

| Contents                             | Cat. No. A32700<br>(100 reactions) <sup>[1]</sup> | Cat. No. A32702<br>(500 reactions) | Storage                       |
|--------------------------------------|---|------------------------------------|-------------------------------|
| MagMAX™ CORE Lysis Solution          | 50 mL   | 275 mL                             | 15–30°C<br>(room temperature) |
| MagMAX™ CORE Binding Solution        | 45 mL   | 220 mL                             |                               |
| MagMAX™ CORE Wash Solution 1         | 60 mL   | 300 mL                             |                               |
| MagMAX™ CORE Wash Solution 2         | 60 mL   | 300 mL                             |                               |
| MagMAX™ CORE Elution Buffer          | 12 mL   | 55 mL                              |                               |
| MagMAX™ CORE Magnetic Beads          | 2.2 mL  | 11 mL                              |                               |
| MagMAX™ CORE Proteinase K (20 mg/mL) | 1.25 mL   | 5 mL                               |                               |

<sup>[1]</sup> Also available in combination with the MagMAX™ CORE Mechanical Lysis Module (Cat. No. A37487)

**Table 3** Materials required for the MagMAX™ CORE Nucleic Acid Purification Kit

| Item   | Source                           |
|--|----------------------------------|
| <b>Instrument and equipment</b>  |                                  |
| One of the following instruments: <ul style="list-style-type: none"> <li>KingFisher™ Flex Purification System</li> <li>MagMAX™ Express-96 Deep Well Magnetic Particle Processor</li> </ul> See page 18 for other compatible instruments. | Contact your local sales office. |
| Benchtop microcentrifuge capable of 15,000 × <i>g</i>  | MLS                              |
| Laboratory mixer, Vortex or equivalent   | MLS                              |
| <b>Reagents</b>  |                                  |
| PBS, pH 7.4 <sup>[1]</sup>   | 10010023                         |
| <b>(Optional) Internal positive control (IPC), one of the following:</b>   |                                  |
| VetMAX™ Xeno™ Internal Positive Control DNA  | A29764                           |
| VetMAX™ Xeno™ Internal Positive Control RNA  | A29763                           |
| IPC supplied with your VetMAX™ PCR Kit   | <b>thermofisher.com</b>          |
| <b>Tubes, plates, and other consumables</b>  |                                  |
| Adhesive PCR Plate Foils, or equivalent  | AB0626                           |
| KingFisher™ Flex Microtiter Deepwell 96 plates, 50 plates  | 95040460                         |

| Item  | Source   |
|---|----------|
| KingFisher™ 96 KF microplates (200 µL), 48 plates | 97002540 |
| KingFisher™ 96 tip comb for DW magnets, 100 combs | 97002534 |

[1] Can be used for isolation of host nucleic acid, or when processing easy-to-lyse bacteria.

**Table 4** Materials required for the MagMAX™ CORE Mechanical Lysis Module

| Item  | Source   |
|---|--|
| <b>Instrument and equipment</b>   |  |
| Bead mill homogenizer for bead-beating, one of the following, or equivalent: <ul style="list-style-type: none"> <li>Fisher Scientific™ Bead Mill 24 Homogenizer (recommended)</li> <li>Mixer Mill 400 (Verder 207450001)</li> <li>Precellys™ 24 Homogenizer (Bertin)</li> <li>FastPrep-24™ Instrument (MP Biomedical 116004500)</li> <li>Mini-BeadBeater-96 (Glen Mills)</li> </ul> | <ul style="list-style-type: none"> <li>Fisher Scientific™ 15-340-163</li> <li>Fisher Scientific™ 08 418 241</li> <li>Bertin EQ03119.200.RD000.0</li> <li>Fisher Scientific™ MP116004500</li> <li>Fisher Scientific™ NC0141170</li> </ul> |
| <b>Reagents</b>   |  |
| Nuclease-Free Water (or any deionized water)  | AM9938   |

**Table 5** Additional materials for the High-input workflow

| Item  | Source                        |
|---|-------------------------------|
| 50-mL conical tube  | MLS                           |
| {Optional} Fisherbrand™ Disposable Sterile Spoon          | Fisher Scientific™ 14-375-255 |
| {Optional} MagMAX™ CORE Glass Microbeads <sup>[1,2]</sup> | A37489                        |

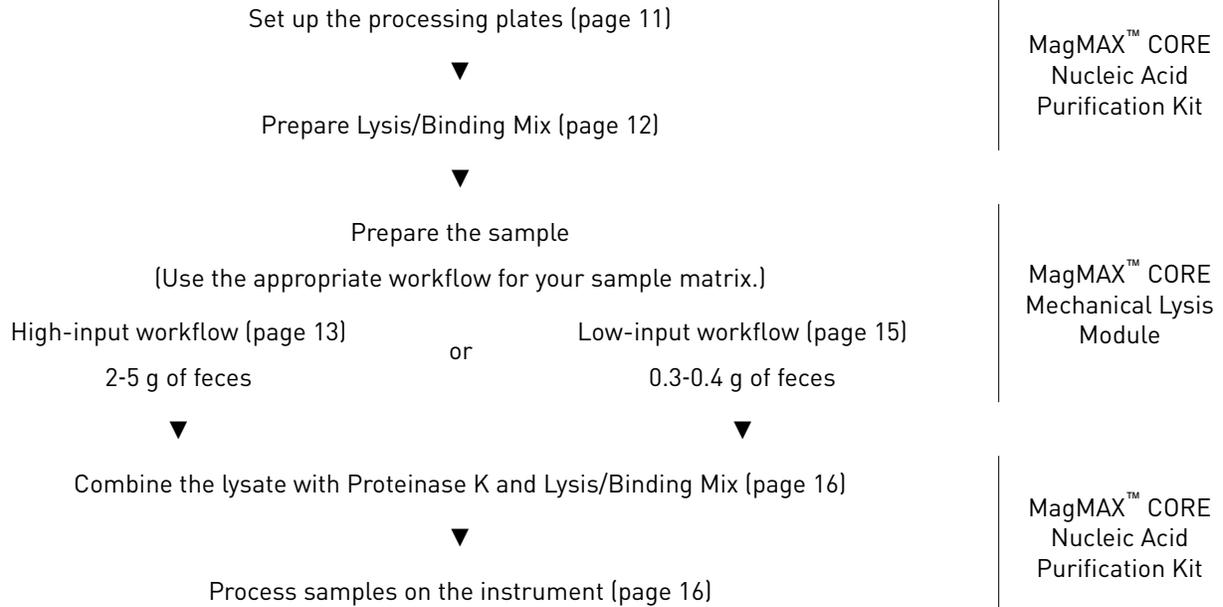
[1] Also available in combination with the MagMAX™ CORE Nucleic Acid Purification Kit and the MagMAX™ CORE Mechanical Lysis Module (Cat. No. A37488)

[2] Required for non-bovine or dried fecal samples.

**Table 6** Additional materials for the Low-input workflow

| Item      | Source |
|-----------|--------|
| 2-mL tube | MLS    |

## Workflow





# Before you begin

## Procedural guidelines

- Before use, invert bottles of solutions and buffers to ensure thorough mixing.
- Mix samples with reagents by pipetting up and down.
- To prevent cross-contamination:
  - Cover the plate or tube strip during the incubation and shaking steps, to prevent spill-over.
  - Carefully pipet reagents and samples, to avoid splashing.
- To prevent nuclease contamination:
  - Wear laboratory gloves during the procedures. Gloves protect you from the reagents, and they protect the nucleic acid from nucleases that are present on skin.
  - Use nucleic acid-free pipette tips to handle the reagents, and avoid putting used tips into the reagent containers.
  - Decontaminate lab benches and pipettes before you begin.

## Before first use of the kit

*(Optional)*  
**Determine the optimal bead mill homogenizer settings**

We recommend using the Fisher Scientific™ Bead Mill 24 Homogenizer for maximum nucleic acid yield. If an alternative instrument is used, follow the manufacturer's guidelines to determine the speed and time settings necessary to achieve sufficient cell lysis.

## Download the script

The scripts for the MagMAX™ CORE Nucleic Acid Purification Kit are not pre-installed on the instruments.

1. On the MagMAX™ CORE Nucleic Acid Purification Kit product web page (at [thermofisher.com](http://thermofisher.com), search by catalogue number), scroll to the **Product Literature** section.
2. Right-click the appropriate file to download the latest version of the MagMAX\_CORE script for your instrument.

**Table 7** Recommended scripts

| Instrument                           | Script name                |
|--------------------------------------|----------------------------|
| KingFisher™ Flex                     | MagMAX_CORE_Flex.bdz       |
| KingFisher™ 96<br>MagMAX™ Express-96 | MagMAX_CORE_KF-96.bdz      |
| KingFisher™ Duo Prime                | MagMAX_CORE_DUO.bdz        |
| KingFisher™ mL                       | MagMAX_CORE_mL_no_heat.bdz |

If required by your laboratory, use one of the following scripts, which do not heat the samples during the elution step.

**Table 8** Alternate scripts without heated elution step

| Instrument                           | Script name                   |
|--------------------------------------|-------------------------------|
| KingFisher™ Flex                     | MagMAX_CORE_Flex_no_heat.bdz  |
| KingFisher™ 96<br>MagMAX™ Express-96 | MagMAX_CORE_KF-96_no_heat.bdz |
| KingFisher™ Duo Prime                | MagMAX_CORE_DUO_no_heat.bdz   |
| KingFisher™ mL                       | MagMAX_CORE_mL_no_heat.bdz    |

See your instrument user guide or contact Technical Support for instructions for installing the script.

Follow this procedure if you are using these instruments:

- KingFisher™ Flex
- MagMAX™ Express-96

Follow Appendix B, “Purification with the KingFisher™ Duo Prime or KingFisher™ mL instrument” if you are using these instruments:

- KingFisher™ Duo Prime
- KingFisher™ mL

## Set up the processing plates

1. Set up the processing plates.

**Table 9** Plate setup: KingFisher™ Flex or MagMAX™ Express-96 instrument

| Plate ID     | Plate position <sup>[1]</sup> | Plate type | Reagent                        | Volume per well |
|--------------|-------------------------------|------------|--------------------------------|-----------------|
| Wash Plate 1 | 2                             | Deep Well  | MagMAX™ CORE Wash Solution 1   | 500 µL          |
| Wash Plate 2 | 3                             | Deep Well  | MagMAX™ CORE Wash Solution 2   | 500 µL          |
| Elution      | 4                             | Standard   | MagMAX™ CORE Elution Buffer    | 90 µL           |
| Tip Comb     | 5                             | Standard   | Place a tip comb in the plate. |                 |

<sup>[1]</sup> Position on the instrument.

2. *(Optional)* To prevent evaporation and contamination, cover the prepared processing plates with sealing foil until they are loaded into the instrument.

## Prepare Lysis/Binding Mix

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

| Component  | Volume per sample                            |
|--|--|
| MagMAX™ CORE Lysis Solution  | 350 µL                                       |
| MagMAX™ CORE Binding Solution  | 350 µL                                       |
| MagMAX™ CORE Magnetic Beads  | 20 µL  |
| <b>Total Lysis/Binding Mix (-IPC)</b>                                    | <b>720 µL</b>                                |
| <i>(Optional)</i> Internal positive control (IPC), one of the following: |  |
| VetMAX™ Xeno™ Internal Positive Control DNA                              | 2 µL <sup>[1]</sup>                          |
| VetMAX™ Xeno™ Internal Positive Control RNA                              | 2 µL <sup>[1]</sup>                          |
| Internal positive control (IPC) supplied with your VetMAX™ PCR Kit       | As indicated in the instructions for the kit |
| <b>Total Lysis/Binding Mix (+IPC)</b>                                    | <b>720 µL + volume of IPC</b>                |

<sup>[1]</sup> Different assays may require different volumes of internal positive control (IPC). See your assay guidelines for IPC recommendations.

2. Mix by inverting the tube or bottle at least 10 times.
3. Store Lysis/Binding Mix at room temperature for up to 24 hours. Do not store on ice.

## Prepare the sample

Select a workflow for your sample matrix.

| Option             | Description         |
|--------------------|---------------------|
| 2-5 g of feces     | High-input workflow |
| 0.3-0.4 g of feces | Low-input workflow  |

### High-input workflow

1. Add the following components to a 50-mL conical tube:

| Component   | Amount               |
|---|----------------------|
| Feces   | 2-5 g <sup>[1]</sup> |
| Deionized water <sup>[2]</sup>                          | 30 mL                |
| [Optional] MagMAX™ CORE Glass Microbeads <sup>[3]</sup> | 3 beads              |

<sup>[1]</sup> Equivalent to one scoop, using a Disposable Sterile Spoon (Fisher Scientific 14-375-255).

<sup>[2]</sup> 1X PBS can be used for isolation of host nucleic acid, or when processing easy-to-lyse bacteria.

<sup>[3]</sup> Recommended when working with non-bovine or dried fecal samples.

2. Vortex vigorously for 30 seconds, or until the sample is suspended.
3. Incubate at room temperature for 10±2 minutes.  
**Note:** Incubate samples on a surface that is free of vibration, to allow the contents to settle before transferring the supernatant.
4. Transfer 1.8 mL of supernatant to a 2-mL MagMAX™ CORE Bead Beating Tube.  
**Note:** If samples are difficult to pipette, the incubation time in step 3 may be increased in 5-minute increments.
5. Centrifuge at 15,000 × g for 10 minutes.
6. Remove the supernatant without disturbing the pellet.  
**Note:** Keep the tube vertical while pipetting to help prevent dislodging the pellet. It is acceptable to aspirate some debris while removing the supernatant.
7. Add 400 µL of MagMAX™ CORE Clarifying Solution to each sample.
8. Disrupt (bead-beat) the samples.

| Option                                      | Settings                  |
|---|---------------------------|
| Fisher Scientific™ Bead Mill 24 Homogenizer | 6 m/s; 3 minutes          |
| Mixer Mill 400                              | 30 Hz; 10 minutes         |
| Precellys™ 24 Homogenizer                   | 6,800 rpm; 2 × 90 seconds |
| FastPrep-24™ Instrument                     | 6.5 M/s; 4 × 45 seconds   |
| Mini-BeadBeater-96                          | 5 minutes                 |

**Note:** If an alternative bead mill homogenizer is used, see “(Optional) Determine the optimal bead mill homogenizer settings” on page 9.

9. Centrifuge at  $15,000 \times g$  for 3 minutes.
10. Proceed with 300  $\mu\text{L}$  of supernatant (clarified lysate) to “Combine the lysate with Proteinase K and Lysis/Binding Mix” on page 16.

Store the clarified lysate at room temperature. Recentrifuge the tubes if not used for  $\geq 1$  hour.

Low-input  
workflow

1. Add the following to a new 2-mL tube:

| Component                      | Amount    |
|--------------------------------|-----------|
| Feces                          | 0.3-0.4 g |
| Deionized water <sup>[1]</sup> | 1 mL      |

<sup>[1]</sup> 1X PBS can be used for isolation of host nucleic acid, or when processing easy-to-lyse bacteria.

2. Vortex vigorously for 3 minutes, or until the sample is suspended.
3. Centrifuge at  $100 \times g$  for 30 seconds.

---

**IMPORTANT!** Do not centrifuge at higher speeds or for longer than 1 minute. Excessive centrifugation can result in bacteria in the pellet instead of in the fecal supernatant.

---

4. Add 400  $\mu\text{L}$  of MagMAX™ CORE Clarifying Solution to the required number of MagMAX™ CORE Bead Beating Tubes.
5. Add 175  $\mu\text{L}$  of the fecal supernatant to the MagMAX™ CORE Bead Beating Tube containing 400  $\mu\text{L}$  of MagMAX™ CORE Clarifying Solution.
6. Disrupt (bead-beat) the samples.

| Option                                      | Settings                         |
|---|----------------------------------|
| Fisher Scientific™ Bead Mill 24 Homogenizer | 6 m/s; 3 minutes                 |
| Mixer Mill 400                              | 30 Hz; 10 minutes                |
| Precellys™ 24 Homogenizer                   | 6,800 rpm; 2 $\times$ 90 seconds |
| FastPrep-24™ Instrument                     | 6.5 M/s; 4 $\times$ 45 seconds   |
| Mini-BeadBeater-96                          | 5 minutes                        |

**Note:** If an alternative bead mill homogenizer is used, see “(Optional) Determine the optimal bead mill homogenizer settings” on page 9.

7. Centrifuge at  $15,000 \times g$  for 3 minutes.
8. Proceed with 300  $\mu\text{L}$  of supernatant (clarified lysate) to “Combine the lysate with Proteinase K and Lysis/Binding Mix” on page 16.

Store the clarified lysate at room temperature. Recentrifuge the tubes if not used for  $\geq 1$  hour.

## Combine the lysate with Proteinase K and Lysis/Binding Mix

1. Add 10  $\mu\text{L}$  of MagMAX™ CORE Proteinase K to the required wells in the plate or tube strip.
2. Add 300  $\mu\text{L}$  of clarified lysate.
3. Mix the clarified lysate with Proteinase K by pipetting up and down several times, then incubate for 2 minutes at room temperature.
4. Add 720  $\mu\text{L}$  of Lysis/Binding Mix to each sample.
5. Immediately proceed to process samples on the instrument (next section).

## Process samples on the instrument

1. Select the appropriate script on the instrument (see “Download the script” on page 10).
2. Start the run, then load the prepared plates or tube strips in their positions when prompted by the instrument.

Store purified nucleic acid on ice for immediate use, at  $-20^{\circ}\text{C}$  for up to 1 month, or at  $-80^{\circ}\text{C}$  for long-term storage.



# Troubleshooting

| Observation   | Possible cause   | Recommended action   |
|---|--|--|
| Poor or no RNA or DNA signal (that is, the $C_t$ value is higher than expected) | There are inhibitors in the recovered nucleic acid.<br><br>These workflows yield high-quality nucleic acid from most samples. However, samples that contain exceptionally high amounts of inhibitors can carry over inhibitors at levels sufficient to affect RT-PCR or PCR. | Include a negative extraction control, to determine the $C_t$ value of the IPC without inhibitors.<br><br>Reduce the amount of RNA used in the RT-PCR or the amount of DNA used in the PCR: Dilute the eluted nucleic acid 10-fold and repeat the RT-PCR or PCR. If a signal is observed using the diluted sample, inhibitors might be present in the eluted nucleic acid. |
|   | Samples with high amounts of nucleic acid, such as tissue, avian blood, and bacterial cultures, can overwhelm the magnetic beads. Overwhelming the beads reduces nucleic acid extraction efficiency.   | Dilute the sample before use, or use less sample.  |
|   | Insufficient lysis resulting in low yields.  | Increase the speed and time of the bead mill homogenizer during the bead-beating step.   |
| Well-to-well variation in RNA/DNA yield from replicate samples                  | The magnetic beads were not fully resuspended/dispersed.   | In general, the magnetic beads disperse more easily when the temperature of the mixture is $> 20^{\circ}\text{C}$ . Be sure that you: <ul style="list-style-type: none"> <li>• Vortex the magnetic beads thoroughly before preparing a bead mix.</li> <li>• Fully resuspend the bead mix before adding it to the samples.</li> </ul>                                       |
| The eluate is light brown in color  | Magnetic beads were carried over into the eluate.  | A small quantity of beads in the sample does not inhibit RT-PCR or PCR reactions.<br><br>Remove the beads from the eluted nucleic acid by placing the plate or tube strip on a magnetic stand (~1 minute), then transfer the nucleic acid solution to a new nuclease-free plate or tube strip.   |



# Purification with the KingFisher™ Duo Prime or KingFisher™ mL instrument

## Required materials not supplied

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.

**Table 10** Materials required for processing on the KingFisher™ Duo Prime instrument

| Item  | Source   |
|---|----------|
| KingFisher™ Duo Prime Purification System   | 5400110  |
| KingFisher™ Duo Combi pack for Microtiter 96 Deepwell plate (tip combs, plates and elution strips for 96 samples) | 97003530 |
| KingFisher™ Duo Elution Strip, 40 pieces <sup>[1]</sup>   | 97003520 |
| KingFisher™ Duo 12-tip comb, for Microtiter 96 Deepwell plate, 50 pieces <sup>[1]</sup>                           | 97003500 |
| KingFisher™ Flex Microtiter Deepwell 96 plates <sup>[1]</sup>   | 95040460 |

<sup>[1]</sup> Included in the KingFisher™ Duo Combi pack (Cat. No. 97003530).

**Table 11** Materials required for processing on the KingFisher™ mL instrument

| Item   | Source   |
|--|----------|
| KingFisher™ mL Purification System                 | 5400050  |
| KingFisher™ mL Tubes and tip combs for 240 samples | 97002141 |
| KingFisher™ mL Tip comb, 800 pieces                | 97002111 |
| KingFisher™ mL Tube, 20 x 45 pieces                | 97002121 |

## Purification procedure

**Note:** When performing this procedure for processing on the KingFisher™ mL instrument, mix samples by pipetting up and down. Do not use a plate shaker with the large tube strips required by this instrument.

1. Follow the workflow for your sample type, starting with sample lysate preparation through combining the samples with beads and lysis solution.

**Note:** Do not set up processing plates or tubes before preparing samples.

2. Add MagMAX™ CORE Wash Solutions and MagMAX™ CORE Elution Buffer to the indicated positions, according to your instrument.

Load the Tip Comb and all of the plates or tube strips at the same time. The instrument does not prompt you to load items individually.

**Table 12** Plate setup: KingFisher™ Duo Prime instrument

| Row ID                 | Row in the plate                   | Plate type    | Reagent                        | Volume per well  |
|------------------------|------------------------------------|---------------|--------------------------------|------------------|
| Sample                 | A                                  | Deep Well     | Sample lysate/bead mix         | Varies by sample |
| Wash 1                 | B                                  |               | MagMAX™ CORE Wash Solution 1   | 500 µL           |
| Wash 2                 | C                                  |               | MagMAX™ CORE Wash Solution 2   | 500 µL           |
| Elution <sup>[1]</sup> | Separate tube strip <sup>[2]</sup> | Elution strip | MagMAX™ CORE Elution Buffer    | 90 µL            |
| Tip Comb               | H                                  | Deep Well     | Place a tip comb in the plate. |                  |

<sup>[1]</sup> Ensure that the elution strip is placed in the correct direction in the elution block.

<sup>[2]</sup> Placed on the heating element.

**Table 13** Tube strip setup: KingFisher™ mL instrument

| Position ID | Tube strip position | Tube     | Reagent                                      | Volume per well  |
|-------------|---------------------|----------|--|------------------|
| Sample      | 1                   | Standard | Sample lysate/bead mix                       | Varies by sample |
| Wash 1      | 2                   |          | MagMAX™ CORE Wash Solution 1                 | 500 µL           |
| Wash 2      | 3                   |          | MagMAX™ CORE Wash Solution 2                 | 500 µL           |
| Elution     | 4                   |          | MagMAX™ CORE Elution Buffer                  | 90 µL            |
| Tip Comb    | N/A                 | N/A      | Slide the tip comb into the tip comb holder. |                  |

3. Follow “Process samples on the instrument” on page 16.



# Safety



**WARNING! GENERAL SAFETY.** Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
  - Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, etc). To obtain SDSs, see the “Documentation and Support” section in this document.
-



## Chemical safety



---

**WARNING! GENERAL CHEMICAL HANDLING.** To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
  - Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
  - Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
  - Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
  - Handle chemical wastes in a fume hood.
  - Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
  - After emptying a waste container, seal it with the cap provided.
  - Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
  - Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
  - **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.
-



## Biological hazard safety



---

**WARNING! BIOHAZARD.** Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:  
[www.cdc.gov/biosafety/publications/bmb15/BMBL.pdf](http://www.cdc.gov/biosafety/publications/bmb15/BMBL.pdf)
  - World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:  
[www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf](http://www.who.int/csr/resources/publications/biosafety/Biosafety7.pdf)
-

# Documentation and support

## Related documentation

| Document   | Publication number |
|--|--------------------|
| <i>MagMAX™ CORE Nucleic Acid Purification Kit User Guide</i>     | MAN0015944         |
| <i>Thermo Scientific™ KingFisher™ Flex User Manual</i>           | N07669             |
| <i>Thermo Scientific™ KingFisher™ Duo Prime Technical Manual</i> | N16621             |
| <i>Thermo Scientific™ KingFisher™ mL User Manual</i>             | 1508260            |
| <i>Applied Biosystems™ MagMAX™ Express 96 User Manual</i>        | N07849             |

## Customer and technical support

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

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

**Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

## 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).

