

MagMAX™ CORE Mastitis & Panbacteria Module

USER GUIDE

Automated and manual purification of high-quality nucleic acid
from all bacteria, including species known to cause mastitis in
cattle

for use with:

KingFisher™ Flex Purification System

KingFisher™ Duo Prime Purification System

Catalog Number A40289

Publication Number MAN0017800

Revision B.0



Life Technologies Ltd | 7 Kingsland Grange | Woolston, Warrington WA1 4SR | United Kingdom

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN0017800

| Revision | Date | Description |
|----------|--------------|---|
| B.0 | 28 July 2023 | <ul style="list-style-type: none">• The MagMAX™ CORE Nucleic Acid Purification Kit was added to the Contents and Storage table (see "Contents and storage" on page 5).• Bulk tank milk samples were added and the procedure was edited accordingly (see Chapter 3, "Methods").• Troubleshooting significant bead coagulate in the eluate was added (see Appendix A, "Troubleshooting"). |
| A.0 | 20 June 2018 | New document for MagMAX™ CORE Mastitis & Panbacteria Module. |

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.

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.

©2023 Thermo Fisher Scientific Inc. All rights reserved.

Contents

| | | | |
|---|-------------------|--|----|
| ■ | CHAPTER 1 | Product information | 5 |
| | | Product description | 5 |
| | | Contents and storage | 5 |
| | | Required materials not supplied | 6 |
| | | Workflow | 7 |
| ■ | CHAPTER 2 | Before you begin | 8 |
| | | Procedural guidelines | 8 |
| | | Before first use of the kit | 8 |
| | | Determine the maximum plate shaker setting | 8 |
| | | Download and install the script | 9 |
| ■ | CHAPTER 3 | Methods | 10 |
| | | Set up the processing plates | 10 |
| | | Prepare the Lysis/Binding/Bead Mix | 10 |
| | | Prepare the sample | 11 |
| | | Combine the sample with MagMAX™ CORE Mastitis Panbacteria solution, then add Proteinase K | 11 |
| | | Process samples on the instrument and add the Lysis/Binding/Bead Mix | 12 |
| ■ | APPENDIX A | Troubleshooting | 13 |
| ■ | APPENDIX B | Manual purification method | 15 |
| | | Required materials not supplied | 15 |
| | | Manual purification procedure | 16 |
| | | Prepare the Lysis/Binding/Bead Mix | 16 |
| | | Prepare the sample | 16 |
| | | Combine the sample with MagMAX™ CORE Mastitis Panbacteria solution, then add Proteinase K | 17 |
| | | Combine the sample mixture with the Lysis/Binding/Bead Mix | 17 |
| | | Wash the sample | 17 |
| | | Dry the beads | 18 |
| | | Elute the nucleic acid | 18 |

- **APPENDIX C Safety 19**
 - Chemical safety 20
 - Biological hazard safety 21
- Documentation and support 22**
 - Related documentation 22
 - Customer and technical support 22
 - Limited product warranty 22



Product information

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

Product description

The MagMAX™ CORE Mastitis & Panbacteria Module is a supplemental module for use with the MagMAX™ CORE Nucleic Acid Purification Kit (Cat. No. [A32700](#)). It is optimized for processing fresh, frozen, or preserved milk samples for the detection of pathogens that are known to cause mastitis in cattle, including gram-positive and gram-negative bacteria. The module uses magnetic bead-based separation, and it is compatible with both automated (using the KingFisher™ Flex Purification System) and manual purification methods.

Contents and storage

Table 1 MagMAX™ CORE Mastitis & Panbacteria Module (Cat. No. [A40289](#))

| Contents | Amount | Storage |
|--|-------------|----------------------------|
| MagMAX™ CORE Nucleic Acid Purification Kit | 1 kit | 15–30°C (room temperature) |
| MagMAX™ CORE Mastitis Panbacteria solution | 4 × 1375 µL | –20°C ^[1] |

^[1] This product is stable for one year when stored as indicated.

Required materials not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Table 2 Materials required for the MagMAX™ CORE Nucleic Acid Purification Kit

| Item | Source |
|---|----------------------------------|
| Instrument and equipment | |
| KingFisher™ Flex Purification System KingFisher™ Duo Prime Purification System See page 15 for materials required for the manual purification method. | Contact your local sales office. |
| Laboratory mixer, Vortex or equivalent | MLS |
| Tubes, plates, and other consumables | |
| Adhesive PCR Plate Foils, or equivalent | AB0626 |
| KingFisher™ Flex Microtiter Deep-Well 96 plates, 50 plates | 95040460 |
| KingFisher™ 96 KF microplates (200 µL), 48 plates | 97002540 |
| KingFisher™ 96 tip comb for deep-well magnets, 100 combs | 97002534 |
| Reagents | |
| (Optional) Internal positive control (IPC), one of the following: ^[1] | |
| VetMAX™ Xeno™ Internal Positive Control DNA | A29764 |
| VetMAX™ Xeno™ Internal Positive Control RNA | A29763 |
| (Optional) Thermo Scientific Chemicals™ TRIS 1.0 M buffer solution, pH 9.5^[2] | J62084.K2 |

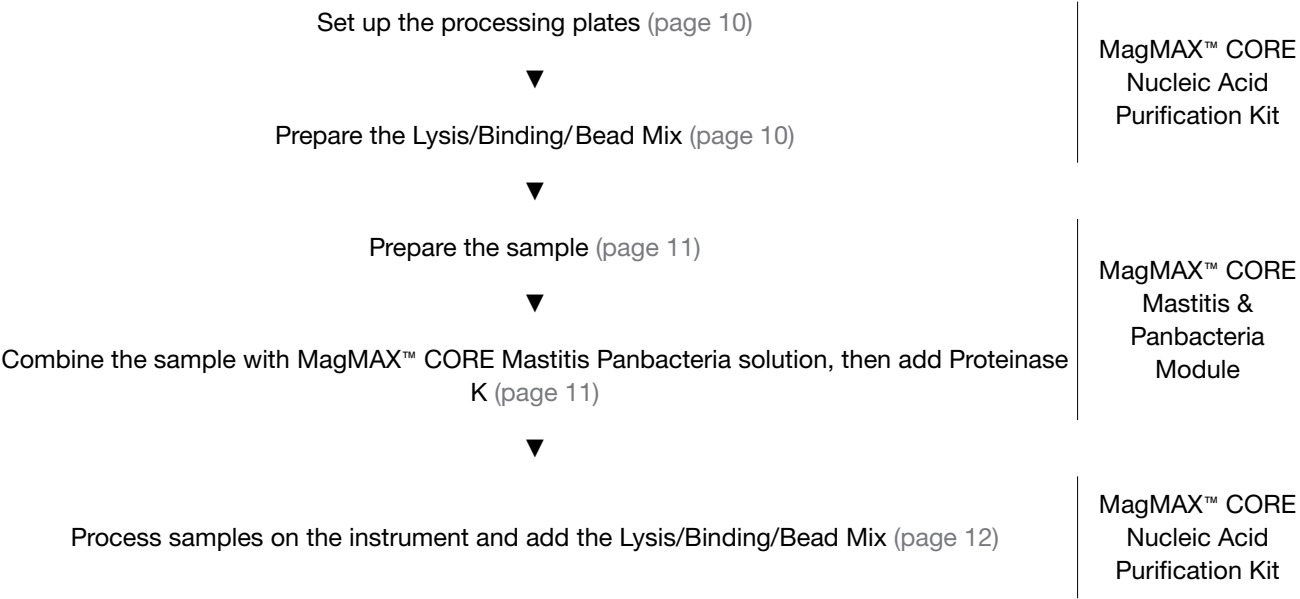
^[1] Not required for VetMAX™ MastiType kits.

^[2] See Appendix A, "Troubleshooting".

Table 3 Materials required for the MagMAX™ CORE Mastitis & Panbacteria Module

| Item | Source |
|--|-------------------------------|
| Boekel Scientific™ 270300 Microplate Shaker, or equivalent | Fisher Scientific™ 15-600-325 |
| (Optional) Magnetic-Ring Stand (96 well) | AM10050 |

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

Determine the maximum plate shaker setting

If a plate shaker is used, use the following steps to determine the maximum setting.

1. Verify that the plate fits securely on your shaker.
2. Add 1 mL of water to each well of the plate, then cover with sealing foil.
3. Determine the maximum setting that you can use on your shaker without any of the water splashing onto the sealing foil.

Download and install the script

The appropriate script for the MagMAX™ CORE Nucleic Acid Purification Kit must be installed on the instrument before first use.

1. On the MagMAX™ CORE Nucleic Acid Purification Kit product web page (at [thermofisher.com](https://www.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 4 Recommended scripts

| Instrument | Script name |
|-----------------------|------------------------------|
| KingFisher™ Flex | MagMAX_CORE_Mastitis.bdz |
| KingFisher™ Duo Prime | MagMAX_CORE_DUO_Mastitis.bdz |

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

- Follow this procedure for automated purification using the KingFisher™ Flex instrument.
- For manual purification, see Appendix B, “Manual purification method”.

Set up the processing plates

1. Set up the processing plates.

Table 5 Plate setup: KingFisher™ Flex instrument

| Plate ID | Plate position ^[1] | 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. | |

^[1] 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 the Lysis/Binding/Bead Mix

1. Combine the following components, in the order indicated, 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 |
| Total Lysis/Binding/Bead Mix (–IPC) | 720 µL |
| (Optional) Internal positive control (IPC), one of the following: ^[1] | |
| VetMAX™ Xeno™ Internal Positive Control DNA | 2 µL ^[2] |
| VetMAX™ Xeno™ Internal Positive Control RNA | 2 µL ^[2] |

(continued)

| Component | Volume per sample |
|--|--|
| Internal positive control (IPC) supplied with your VetMAX™ PCR Kit | As indicated in the instructions for the kit |
| Total Lysis/Binding/Bead Mix (+IPC) | 720 µL + volume of IPC |

^[1] Not required for VetMAX™ MastiType kits.

^[2] Different assays may require different volumes of internal positive control (IPC). See your assay guidelines for IPC recommendations.

2. Vortex the mixture thoroughly.

(Optional) Store the Lysis/Binding/Bead Mix at room temperature for up to 24 hours.

Prepare the sample

Prepare the samples and controls as described.

| Sample type | Action |
|--|---|
| Individual Milk Sample | Proceed with 50 µL of fresh, frozen, or preserved milk sample. |
| Bulk Tank Milk Sample | Proceed with 200 µL of fresh, frozen, or preserved bulk tank milk sample. |
| Negative Extraction Control ^[1] | Proceed with 200 µL of nuclease-free water or 1X PBS. |

^[1] Recommended to detect reagent contamination during sample preparation.

Combine the sample with MagMAX™ CORE Mastitis Panbacteria solution, then add Proteinase K

1. Add 50 µL of the MagMAX™ CORE Mastitis Panbacteria solution to the required wells in the plate.
2. Add the milk sample.
 - 50 µL of the individual milk sample.
 - 200 µL of the bulk tank milk sample.
3. Mix the sample with MagMAX™ CORE Mastitis Panbacteria solution for 5 minutes at room temperature according to your mixing method.
 - **Using a plate shaker**—Shake vigorously for 5 minutes (see “Determine the maximum plate shaker setting” on page 8).
 - **By pipetting**—Pipet up and down several times, then incubate for 5 minutes at room temperature.
4. Add 10 µL of MagMAX™ CORE Proteinase K to the required wells in the plate.
5. Immediately proceed to process samples on the instrument (next section).

Process samples on the instrument and add the Lysis/Binding/Bead Mix

1. Select the appropriate script on the instrument (see “Download and install the script” on page 9).
2. Start the run, then load the prepared plates in the appropriate positions when prompted by the instrument.
The run proceeds for 10 minutes, followed by a pause.
3. While the instrument is paused, remove the sample plate from the instrument.
4. Vortex the tube of Lysis/Binding/Bead Mix to resuspend the beads, then add 720 µL of the Lysis/Binding/Bead Mix to each well.
5. Load the sample plate back on the instrument, then press **Start** to continue the run.
6. When the run is complete, remove the elution plate containing the purified nucleic acid, then cover the plate with adhesive film.

Note: In samples with high DNA content, beads can carry over into the final eluate. If this occurs, place the elution plate on a magnetic ring stand to allow the beads to settle (~1 minute) before using the eluate for PCR amplification.

Store the purified nucleic acid on ice for immediate use, at –20°C for up to 1 month, or at –80°C for long-term storage.



Troubleshooting

| Observation | Possible cause | Recommended action |
|--|---|--|
| There is significant bead coagulate (dark brown aggregate of magnetic beads) in the eluate. | Milk samples may have excess amounts of fat or proteins that cause the beads to aggregate. | Add 50 μ L of TRIS buffer (pH 9.5, Cat. No. J62084.K2) to 50 μ L of individual milk sample, then vortex thoroughly. Use 100 μ L of the resultant TRIS-milk mixture as the starting sample. |
| 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 PCR reactions. 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. |
| Poor or no DNA signal (that is, the C_t value is higher than expected) In test samples, the C_t value of the Internal Amplification Control (IAC) target is outside of the validated value range (non-compliant IAC C_t value; invalid sample). | Inhibitors are present in the recovered nucleic acid. These workflows yield high-quality nucleic acid for most samples. However, samples that contain exceptionally high amounts of inhibitors can carry over inhibitors at levels sufficient to affect PCR. | <ol style="list-style-type: none"> 1. Dilute the invalid nucleic acid sample 1:10 in 1X TE buffer. 2. Perform a new PCR analysis with the diluted nucleic acid. <ul style="list-style-type: none"> • If the diluted nucleic acid is positive for the target, or if it is negative for the target with a compliant IAC C_t value, the result is validated. • If the diluted nucleic acid is negative for the target with a non-compliant IAC C_t value, the result is not validated. In this case, dilute the original biological sample 1:10 in 1X PBS, then repeat the purification and PCR. If the result is still not validated, then repeat the purification and PCR on a new biological sample. |
| | Samples with high amounts of nucleic acid can saturate the magnetic beads. Bead saturation reduces nucleic acid recovery. | Dilute the samples 1:2, 1:4, 1:8, and 1:16 in 1X PBS, then repeat the purification and PCR. |



Appendix A Troubleshooting

Process samples on the instrument and add the Lysis/Binding/Bead Mix

| Observation | Possible cause | Recommended action |
|--|--|--|
| Well-to-well variation in 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°C. Be sure that you: <ul style="list-style-type: none">• Vortex the magnetic beads thoroughly before preparing a bead mix.• Fully resuspend the bead mix before adding it to the samples. |
| Positive samples are clustered in the PCR plate | High-titer samples (exhibiting a low or early C _t) have contaminated nearby wells. | Repeat the nucleic acid purification of the positive or suspect samples without the high-titer samples. |



Manual purification method

This appendix provides manual purification procedures for the low-throughput processing of nucleic acid from milk samples.

Required materials not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](https://www.fisherscientific.com) or another major laboratory supplier.

Table 6 Materials required for the manual purification method

| Item | Source |
|---|------------------------------|
| Benchtop microcentrifuge capable of 15,000 × <i>g</i> | MLS |
| Laboratory mixer, Vortex or equivalent | MLS |
| Heat block set to 95°C | MLS |
| DynaMag™-2 Magnet, or equivalent 16-tube magnetic stand | 12321D |
| Vortex Adapter-60, or equivalent | AM10014 |
| Eppendorf™ Snap-Cap Microcentrifuge Safe-Lock™ Tubes, or equivalent 1.5 mL tubes | Fisher Scientific™ 05-402-25 |
| (Optional) Thermo Scientific Chemicals™ TRIS 1.0 M buffer solution, pH 9.5 ^[1] | J62084.K2 |

^[1] See Appendix A, "Troubleshooting".

Manual purification procedure

Prepare the Lysis/Binding/Bead Mix

1. Combine the following components, in the order indicated, 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 |
| Total Lysis/Binding/Bead Mix (–IPC) | 720 µL |
| <i>(Optional)</i> Internal positive control (IPC), one of the following: ^[1] | |
| VetMAX™ Xeno™ Internal Positive Control DNA | 2 µL ^[2] |
| VetMAX™ Xeno™ Internal Positive Control RNA | 2 µL ^[2] |
| Internal positive control (IPC) supplied with your VetMAX™ PCR Kit | As indicated in the instructions for the kit |
| Total Lysis/Binding/Bead Mix (+IPC) | 720 µL + volume of IPC |

^[1] Not required for VetMAX™ MastiType kits.

^[2] Different assays may require different volumes of internal positive control (IPC). See your assay guidelines for IPC recommendations.

2. Vortex the mixture thoroughly.

(Optional) Store the Lysis/Binding/Bead Mix at room temperature for up to 24 hours.

Prepare the sample

Prepare the samples and controls as described.

| Sample type | Action |
|--|---|
| Individual Milk Sample | Proceed with 50 µL of fresh, frozen, or preserved milk sample. |
| Bulk Tank Milk Sample | Proceed with 200 µL of fresh, frozen, or preserved bulk tank milk sample. |
| Negative Extraction Control ^[1] | Proceed with 200 µL of nuclease-free water or 1X PBS. |

^[1] Recommended to detect reagent contamination during sample preparation.

Combine the sample with MagMAX™ CORE Mastitis Panbacteria solution, then add Proteinase K

1. Add 50 µL of MagMAX™ CORE Mastitis Panbacteria solution to each 1.5-mL tube.
2. Add the milk sample.
 - 50 µL of the individual milk sample.
 - 200 µL of the bulk tank milk sample.
3. Using a vortex with a microtube adapter, shake at moderate speed for 5 minutes.
4. Centrifuge at $500 \times g$ for 10 seconds to bring the contents to the bottom of the tube.
5. Add 10 µL of MagMAX™ CORE Proteinase K to each tube, then shake at moderate speed for 5 minutes.
6. Incubate at 95°C for 5 minutes.
7. Centrifuge at $500 \times g$ for 10 seconds to collect condensation that can accumulate in the cap of the tube.

Combine the sample mixture with the Lysis/Binding/Bead Mix

Note: A 16-tube magnetic stand is required for the remainder of the manual purification procedure.

1. Add 720 µL of the Lysis/Binding/Bead Mix to each tube, then vortex at moderate-to-high speed for 3 minutes.
2. Centrifuge at $500 \times g$ for 10 seconds.
3. Place the tubes and tube rack on the magnetic stand, then incubate at room temperature for 3 minutes.
4. With the tubes on the magnetic stand, carefully remove, then discard the supernatant.
5. Remove the tube rack from the magnetic stand.

Wash the sample

1. Add 500 µL of MagMAX™ CORE Wash Solution 1 to each tube.
 - a. Vortex at moderate high speed for 1 minute.
 - b. Centrifuge at $500 \times g$ for 10 seconds.
 - c. Place the tubes and tube rack on the magnetic stand, then incubate at room temperature for 1 minute.
 - d. With the tubes on the magnetic stand, carefully remove, then discard the supernatant.
 - e. Remove the tube rack from the magnetic stand.

2. Add 500 μL of MagMAX™ CORE Wash Solution 2 to each tube.
 - a. Vortex at moderate high speed for 1 minute.
 - b. Centrifuge at $500 \times g$ for 10 seconds.
 - c. Place the tubes and tube rack on the magnetic stand, then incubate at room temperature for 1 minute.
 - d. With the tubes on the magnetic stand, carefully remove, then discard the supernatant.

Dry the beads

1. Keep the tubes on the magnetic stand and with the tubes open, dry the beads at room temperature for 5 minutes.
2. Use a 20 μL pipette to remove any remaining wash solution.

Elute the nucleic acid

1. Remove the tube rack from the magnetic stand.
2. Add 90 μL of MagMAX™ CORE Elution Buffer to each tube.
3. Vortex at moderate high speed for 3 minutes.
4. Centrifuge at $500 \times g$ for 10 seconds.
5. Place the tubes and tube rack on the magnetic stand, then incubate at room temperature for 3 minutes.
6. With the tubes on the magnetic stand, transfer the purified nucleic acid to a clean, labeled tube or plate.

Store the purified nucleic acid on ice for immediate use, at -20°C for up to 1 month, or at -80°C for long-term storage.



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, and so on). To obtain SDSs, visit [thermofisher.com/support](https://www.thermofisher.com/support).



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 sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer 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 needed) 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.



WARNING! HAZARDOUS WASTE (from instruments). Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



WARNING! 4L Reagent and Waste Bottle Safety. Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.



Biological hazard safety



WARNING! Potential Biohazard. Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



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)*, 6th Edition, HHS Publication No. (CDC) 300859, Revised June 2020
www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2020-P.pdf
- Laboratory biosafety manual, fourth edition. Geneva: World Health Organization; 2020 (Laboratory biosafety manual, fourth edition and associated monographs)
www.who.int/publications/i/item/9789240011311

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> | MAN0019870 |
| <i>Thermo Scientific™ KingFisher™ Duo Prime Technical Manual</i> | N16621 |
| <i>Thermo Scientific™ KingFisher™ Duo Prime Brief User Manual</i> | N16622 |

Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - 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 at 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.

