

# MagMAX™ mirVana™ Total RNA Isolation Kit

High-throughput isolation of RNA (including small RNA) from tissue samples

Catalog Number A27828

Pub. No. MAN0011132 Rev. G



**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 description

The Applied Biosystems™ MagMAX™ mirVana™ Total RNA Isolation Kit (Cat. No. [A27828](#)) is designed for isolation of total RNA, including microRNA, from a wide variety of sample matrices. The kit uses MagMAX™ magnetic-bead technology, ensuring reproducible recovery of high-quality RNA that is suitable for a broad range of applications, including TaqMan™ miRNA Detection Assays.

This protocol describes isolation of RNA from tissue samples. It is optimized for use with the MagMAX™ Express-96 Deep Well Magnetic Particle Processor, the KingFisher™ Flex Purification System with 96 Deep-Well Head (96-well deep well setting), the KingFisher™ Apex Purification System with 96 Deep-Well Head, and the KingFisher™ Duo Prime Purification System (12-well deep well setting).

## Contents and storage

**Table 1** MagMAX™ mirVana™ Total RNA Isolation Kit (Cat. No. [A27828](#), 96 reactions)

Contents	Amount	Storage
<b>Box 1 of 2</b>		
Proteinase K <sup>[1]</sup>	0.48 mL	-25°C to -15°C
Lysis/Binding Enhancer	0.96 mL	
TURBO DNase™, 20 U/μL	0.2 mL	
<b>Box 2 of 2</b>		
Lysis Buffer	115 mL	15–25°C
PK Digestion Buffer <sup>[1]</sup>	4.4 mL	
RNA Binding Beads <sup>[2]</sup>	2 mL	
Wash Solution 1 Concentrate <sup>[3]</sup>	20 mL	
Wash Solution 2 Concentrate <sup>[3]</sup>	60 mL	
Rebinding Buffer	4.8 mL	
MagMAX™ TURBO DNase™ Buffer	4.8 mL	
Elution Buffer	9.6 mL	
Processing Plate <sup>[1]</sup>	1	
Elution Plates	2	
Plate Covers	4	

<sup>[1]</sup> Not used for RNA isolation from tissue samples.

<sup>[2]</sup> Do not freeze the RNA Binding Beads.

<sup>[3]</sup> Final volume; see “Prepare wash solutions” on page 2.

## 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.

Catalog numbers that appear as links open the web pages for those products.

Item	Source
<b>Instrument, one of the following:</b>	
MagMAX™ Express-96 Deep Well Magnetic Particle Processor	— <sup>[1]</sup>
KingFisher™ Flex Purification System with 96 Deep-Well Head <sup>[2]</sup>	<a href="#">5400630</a>
KingFisher™ Apex Purification System with 96 Deep-Well Head <sup>[2]</sup>	<a href="#">5400930</a>
KingFisher™ Duo Prime Purification System <sup>[2]</sup>	<a href="#">5400110</a>
<b>Other equipment</b>	
Thermo Scientific™ Digital Microplate Shaker, or equivalent titer shaker	<a href="#">88882005</a>
Vortex mixer	<a href="#">MLS</a>
PRO250 Homogenizer with 7 × 95 mm Saw Tooth Bottom Generator Probe	PRO Scientific <a href="#">01-01250</a> and <a href="#">02-07095</a>
Adjustable micropipettors	<a href="#">MLS</a>
Multi-channel micropipettors	<a href="#">MLS</a>
(Optional) Magnetic Stand-96	<a href="#">AM10027</a>
<b>Plates and combs<sup>[3]</sup></b>	
Deep-well plates, one of the following:	
KingFisher™ Flex Microtiter Deep-Well 96 plate, sterile	<a href="#">95040460</a>
KingFisher™ 96 Deep-Well Plate, v-bottom, polypropylene	<a href="#">95040450</a>
KingFisher™ 96 Deep-Well Plate, Barcoded	<a href="#">95040450B</a>
Standard well plate:	
KingFisher™ 96 KF microplate	<a href="#">97002540</a>
One of the following tip combs, depending on the instrument used:	
KingFisher™ 96 tip comb for deep-well magnets	<a href="#">97002534</a>
KingFisher™ 12-tip comb, for 96 deep-well plate <sup>[4]</sup>	<a href="#">97003500</a>
<b>Other consumables</b>	
MicroAmp™ Clear Adhesive Film	<a href="#">4306311</a>
Nonstick, RNase-free Microfuge Tubes, 1.5 mL	<a href="#">AM12450</a>
Nonstick, RNase-Free Microfuge Tubes, 2.0 mL	<a href="#">AM12475</a>
5-mL Culture tubes	<a href="#">MLS</a>
Conical Tubes (15 mL)	<a href="#">AM12500</a>
Aerosol-resistant pipette tips	<a href="#">MLS</a>
Reagent reservoirs	<a href="#">MLS</a>
<b>Reagents</b>	
Isopropanol, 100% (molecular grade or higher)	<a href="#">MLS</a>
Ethanol, 200 proof (absolute)	<a href="#">MLS</a>
2-Mercaptoethanol	<a href="#">MLS</a>
(Optional) Chloroform	<a href="#">MLS</a>

[1] Not available for sale.

[2] See "(If needed) Download, then install the program on the instrument" on page 2.

[3] KingFisher™ Duo Combi Pack (Cat. No. 97003530) includes plates and combs for the KingFisher™ Duo Prime instrument.

[4] For use with the KingFisher™ Duo Prime instrument only.

## Sample collection and storage

- Process tissue immediately, store them in RNA<sup>later</sup>™ Stabilization Solution, or freeze them in liquid nitrogen and store at  $-80^{\circ}\text{C}$ .
- For ease of processing, we recommend pre-weighing and storing the tissue in pieces of incremental sizes, according to the following guidelines:
  - Increments of  $\leq 50$  mg for tissues containing low or normal level of cellular RNase.
  - Increments of  $\leq 30$  mg for tissues containing high level of cellular RNase, such as spleen or pancreas.

## Procedural guidelines

- Perform all steps at room temperature ( $20\text{--}25^{\circ}\text{C}$ ) unless otherwise noted.
- When mixing samples by pipetting up and down, avoid creating bubbles.
- Cover the plate during the shaking steps to prevent spill-over and cross-contamination. The same Plate Cover can be used throughout the procedure, unless it becomes contaminated.
- If you use a titer plate shaker other than the Thermo Scientific™ Digital Microplate Shaker, ensure that:
  - The plate fits securely on your titer plate shaker.
  - The recommended speeds are compatible with your titer plate shaker. Ideal speeds should allow for thorough mixing without splashing.
- Volumes for reagent mixes are given per well. We recommend that you prepare master mixes for larger sample numbers. To calculate volumes for master mixes, refer to the per-well volume and add 5% overage.
- Lysed samples can be stored in Lysis Binding Mix at  $-20^{\circ}\text{C}$  for up to 4 days before adding the Binding Beads Mix. Thaw frozen samples to room temperature before use.

## Before first use of the kit

(If needed) Download, then install the program on the instrument

The appropriate program for the MagMAX™ mirVana™ Total RNA Isolation Kit must be installed on the instrument before first use.

## Lyse the tissue samples (all methods)

- Determine the size of tissue (in mg) to be homogenized.
- Determine the amount of Lysis Binding Mix that is needed to homogenize the tissue, according to the following ratios:
  - For tissues containing low to normal levels of cellular RNase (for example, brain, heart, or liver), use 20  $\mu\text{L}$  of Lysis Binding Mix for 1 mg of tissue (1:20 ratio).  
For example, use 200  $\mu\text{L}$  of Lysis Binding Mix for 10 mg of tissue.
  - For RNA<sup>later</sup>-stored tissues, and tissues containing high levels of cellular RNase (for example, spleen or pancreas), use 40  $\mu\text{L}$  of Lysis Binding Mix for 1 mg of tissue (1:40 ratio).  
For example, use 400  $\mu\text{L}$  of Lysis Binding Mix for 10 mg of tissue.

**Note:** Most mechanical homogenizers require a minimum volume of 200  $\mu\text{L}$ . Therefore, we recommend processing no less than 10 mg when homogenizing.

- Prepare sufficient Lysis Binding Mix, according to the following table.

Component	Volume for 100 $\mu\text{L}$ of Lysis Binding Mix
Lysis Buffer	100 $\mu\text{L}$
2-Mercaptoethanol	0.7 $\mu\text{L}$
<b>Total Lysis Binding Mix</b>	<b>~100 <math>\mu\text{L}</math></b>

- Add tissue to the prepared Lysis Binding Mix.

- On the MagMAX™ mirVana™ Total RNA Isolation Kit web page (at [thermofisher.com](http://thermofisher.com), Cat. No. A27828), go to the **Software & Data Analysis** section.
- Download the appropriate program for your instrument.

**Table 2 Recommended programs**

Instrument	Program name
MagMAX™ Express-96	AM1830DW
KingFisher™ Flex	A27828_FLEX_Tissue_Cells
KingFisher™ Apex	A27828_MagMAX_mirVana_TissueCells
KingFisher™ Duo Prime	A27828_DUO_Tissue_Cells

- See your instrument user guide for instructions for installing the program.

## Prepare wash solutions

- Add 10 mL of isopropanol to Wash Solution 1 Concentrate, then mix.
- Add 48 mL of ethanol to Wash Solution 2 Concentrate, then mix.

Store the prepared wash solutions at room temperature.

## Before each use: Prepare TURBO DNase™ Solution and Binding Beads Mix

- Prepare the TURBO DNase™ Solution as indicated in the following table, mix, then store on ice until use.

Component	Volume per well
MagMAX™ TURBO DNase™ Buffer	48 $\mu\text{L}$
TURBO DNase™	2 $\mu\text{L}$
<b>Total TURBO DNase™ Solution</b>	<b>50 <math>\mu\text{L}</math></b>

- Prepare the Binding Beads Mix as indicated in the following table, mix, then store on ice until use.

Component	Volume per well
RNA Binding Beads	10 $\mu\text{L}$
Lysis/Binding Enhancer	10 $\mu\text{L}$
<b>Total Binding Beads Mix</b>	<b>20 <math>\mu\text{L}</math></b>

- Homogenize the tissue sample using standard homogenization procedures.

**IMPORTANT!** Make sure that the tissue homogenization is complete to ensure maximal RNA recovery. For the tissue input amounts listed previously, we recommend homogenizing for 30 seconds to fully break up the sample and ensure maximal RNA recovery and quality.

Proceed to the appropriate procedures:

- “Isolate RNA using the MagMAX™ Express-96 or KingFisher™ Flex instrument”.
- “Isolate RNA using the KingFisher™ Apex instrument”.
- “Isolate RNA using the KingFisher™ Duo Prime instrument”.

## Isolate RNA from tissues samples

### Isolate RNA using the MagMAX™ Express-96 or KingFisher™ Flex instrument

#### 1 Bind the RNA to the RNA Binding Beads

If samples were frozen at the previous step, thaw them completely to room temperature before proceeding.

- Vortex the lysates, then transfer 100 µL to a separate well in a KingFisher™ 96 Deep-Well Plate.

**Table 3 Recommended tissue mass per 100 µL of Lysis Binding Mix**

Samples	Levels of cellular RNase	
	Low to normal (for example, liver)	High (for example, spleen)
Frozen tissue	up to 5 mg	up to 2.5 mg
Tissue stored in RNAlater™ Stabilization Solution	up to 2.5 mg	up to 2.5 mg

**Note:** Most mechanical homogenizers require at least 200 µL. Therefore, we recommend processing no less than 10 mg of tissue when homogenizing.

- (Optional) Add 10 µL of chloroform to each well.

**IMPORTANT!** The addition of chloroform is required for samples containing high levels of RNase (spleen, pancreas).

- Cover the plate, then shake as indicated.

Time	Speed
5 minutes	1,150 rpm (Speed 10)

During the incubation, set up the processing plates (next section).

- Add 100 µL of isopropanol to each sample, cover the plate, then shake as indicated.

Time	Speed
2 minutes	950 rpm (Speed 7)

- Add 20 µL of the prepared Binding Beads Mix to each sample, then shake as indicated.

Time	Speed
5 minutes	950 rpm (Speed 7)

- Proceed directly to process samples on the instrument (see “Process samples on the instrument” on page 4).

#### 2 Set up the processing plates

While the samples are incubating, set up the Wash, DNase, Elution, and Tip Comb Plates outside the instrument as described in the following table.

**Table 4 Processing plates**

Plate ID	Plate position <sup>[1]</sup>	Plate type	Reagent	Volume per well
Wash Plate 1	2	Standard	Wash Solution 1	150 µL
Wash Plate 2	3	Standard	Wash Solution 2	150 µL
DNase Plate <sup>[2]</sup>	4	Deep Well	TURBO DNase™ Solution	50 µL
Wash Plate 3	5	Standard	Wash Solution 2	150 µL
Wash Plate 4	6	Standard	Wash Solution 2	150 µL
Elution Plate	7	Standard	Elution Buffer	50–100 µL <sup>[3]</sup>
Tip Comb	8	Deep Well or Standard	Place a KingFisher™ 96 tip comb for deep-well magnets in a KingFisher™ 96 Deep-Well Plate or in a KingFisher™ 96 KF microplate.	

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

<sup>[2]</sup> The instrument prompts the user to add 50 µL of Rebinding Buffer and 100 µL of isopropanol to the DNase Plate after the DNase treatment step.

<sup>[3]</sup> Use 50 µL for lower-yielding samples, such as brain or heart, or 100 µL for higher-yielding samples, such as liver or spleen.

### 3 Process samples on the instrument

Ensure that the instrument is set up with the deep well magnetic head.

1. Select the appropriate program on the instrument (see “(If needed) Download, then install the program on the instrument” on page 2).
2. Start the run, then load the prepared processing plates in their positions when prompted by the instrument (see Table 4).
3. Load the sample plate (containing the lysate, isopropanol, and Binding Beads Mix) at position 1 when prompted by the instrument.
4. When prompted by the instrument (30–35 minutes after the initial start):
  - a. Remove the DNase Plate from the instrument.
  - b. Add 50 µL of Rebinding Buffer and 100 µL of isopropanol separately to each sample well. Add Rebinding Buffer and isopropanol immediately after the prompt, to prevent excessive drying of any beads that are still captured on the Tip Comb.

**IMPORTANT!** Do not premix the Rebinding Buffer and isopropanol. Add them separately to the samples.

- c. Load the DNase Plate back onto the instrument, then tap **Start**.
5. At the end of the run (about 45 minutes after the initial start), remove the Elution Plate from the instrument, then seal immediately with a new MicroAmp™ Clear Adhesive Film.
  - (Optional) Transfer the eluates to a storage plate after collection.
  - If excess bead residue is seen in the wells, place the Elution Plate on a Magnetic Stand-96 (Cat. No. AM10027) to capture any residue, then transfer the eluates to a new Elution Plate before using the RNA in downstream applications.

**IMPORTANT!** Do not allow the purified samples to sit uncovered at room temperature for more than 10 minutes to prevent evaporation and contamination.

The purified samples are ready for immediate use. Alternatively, store the covered Elution Plate:

- On ice for up to 8 hours.
- At –20°C or –80°C for long-term storage.

## Isolate RNA using the KingFisher™ Apex instrument

### 1 Bind the RNA to the RNA Binding Beads

If samples were frozen at the previous step, thaw them completely to room temperature before proceeding.

1. Vortex the lysates, then transfer 100 µL to a separate well in a KingFisher™ Flex (non-barcoded) or KingFisher™ Apex (barcoded) plate.

Table 5 Recommended tissue mass per 100 µL of Lysis Binding Mix

Samples	Levels of cellular RNase	
	Low to normal (for example, liver)	High (for example, spleen)
Frozen tissue	up to 5 mg	up to 2.5 mg
Tissue stored in RNAwater™ Stabilization Solution	up to 2.5 mg	up to 2.5 mg

**Note:** Most mechanical homogenizers require at least 200 µL. Therefore, we recommend processing no less than 10 mg of tissue when homogenizing.

2. (Optional) Add 10 µL of chloroform to each well.

**IMPORTANT!** The addition of chloroform is required for samples containing high levels of RNase (spleen, pancreas).

3. Cover the plate, then shake as indicated.

Time	Speed
5 minutes	1,150 rpm (Speed 10)

During the incubation, set up the processing plates (next section).

4. Add 100 µL of isopropanol to each sample, cover the plate, then shake as indicated.

Time	Speed
2 minutes	950 rpm (Speed 7)

5. Add 20 µL of the prepared Binding Beads Mix to each sample, then shake as indicated.

Time	Speed
5 minutes	950 rpm (Speed 7)

6. Proceed directly to process samples on the instrument (see “Process samples on the instrument” on page 5).

## 2 Set up the processing plates

While the samples are incubating, set up the Wash, DNase, Elution, and Tip Comb Plates outside the instrument as described in the following table.

**Table 6 Processing plates**

Plate ID	Plate position <sup>[1]</sup>	Plate type	Reagent	Volume per well
Wash Plate 1	3	Standard	Wash Solution 1	150 µL
Wash Plate 2	4	Standard	Wash Solution 2	150 µL
DNase Plate <sup>[2]</sup>	5	Deep Well	TURBO DNase™ Solution	50 µL
Wash Plate 3	6	Standard	Wash Solution 2	150 µL
Wash Plate 4	7	Standard	Wash Solution 2	150 µL
Elution Plate	8	Standard	Elution Buffer	50–100 µL <sup>[3]</sup>
Tip Comb	1	Deep Well	Place a KingFisher™ 96 tip comb for deep-well magnets into a KingFisher™ 96 Deep-Well Plate.	

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

<sup>[2]</sup> The instrument prompts the user to add 50 µL of Rebinding Buffer and 100 µL of isopropanol to the DNase Plate after the DNase treatment step.

<sup>[3]</sup> Use 50 µL for lower-yielding samples, such as brain or heart, or 100 µL for higher-yielding samples, such as liver or spleen.

## 3 Process samples on the instrument

Ensure that the instrument is set up with the deep well magnetic head.

1. Select the appropriate program on the instrument (see “(If needed) Download, then install the program on the instrument” on page 2).
2. Start the run, then load the prepared processing plates in their positions when prompted by the instrument (see Table 6).
3. Load the sample plate (containing the lysate, isopropanol, and Binding Beads Mix) when prompted by the instrument.
4. When prompted by the instrument (30–35 minutes after the initial start):
  - a. Remove the DNase Plate from the instrument.
  - b. Add 50 µL of Rebinding Buffer and 100 µL of isopropanol separately to each sample well.  
Add Rebinding Buffer and isopropanol immediately after the prompt to prevent excessive drying of any beads that are still captured on the Tip Comb.

**IMPORTANT!** Do not premix the Rebinding Buffer and isopropanol. Add them separately to the samples.

- c. Load the DNase Plate back onto the instrument, then tap **Start**.
5. At the end of the run (about 45 minutes after the initial start), remove the Elution Plate from the instrument, then seal immediately with a new MicroAmp™ Clear Adhesive Film.
    - (Optional) Transfer the eluates to a storage plate after collection.
    - If excess bead residue is seen in the wells, place the Elution Plate on a Magnetic Stand-96 (Cat. No. [AM10027](#)) to capture any residue, then transfer the eluates to a new Elution Plate before using the RNA in downstream applications.

**IMPORTANT!** Do not allow the purified samples to sit uncovered at room temperature for more than 10 minutes to prevent evaporation and contamination.

The purified samples are ready for immediate use. Alternatively, store the covered Elution Plate:

- On ice for up to 8 hours.
- At –20°C or –80°C for long-term storage.

## Isolate RNA using the KingFisher™ Duo Prime instrument

### 1 Bind the RNA to the RNA Binding Beads

If samples were frozen at the previous step, thaw them completely to room temperature before proceeding.

1. Vortex the lysates, then transfer 100 µL to a separate well of row H of a KingFisher™ 96 Deep-Well Plate.

**Table 7 Recommended tissue mass per 100 µL of Lysis Binding Mix**

Samples	Levels of cellular RNase	
	Low to normal (for example, liver)	High (for example, spleen)
Frozen tissue	up to 5 mg	up to 2.5 mg
Tissue stored in RNAlater™ Stabilization Solution	up to 2.5 mg	up to 2.5 mg

**Note:** Most mechanical homogenizers require at least 200 µL. Therefore, we recommend processing no less than 10 mg of tissue when homogenizing.

2. (Optional) Add 10 µL of chloroform to each well.

**IMPORTANT!** The addition of chloroform is required for samples containing high levels of RNase (spleen, pancreas).

1 (continued)

3. Cover the plate, then shake as indicated.

Time	Speed
5 minutes	1,150 rpm (Speed 10)

4. Add 100 µL of isopropanol to each sample, cover the plate, then shake as indicated.

Time	Speed
2 minutes	950 rpm (Speed 7)

5. Add 20 µL of the prepared Binding Beads Mix to each sample, then shake as indicated.

Time	Speed
5 minutes	950 rpm (Speed 7)

2 Set up the processing plate

Add processing reagents as indicated in the following table.

**Table 8 Volume of processing reagents and plate location**

Row ID	Plate row <sup>[1]</sup>	Reagent	Volume per well
Elution	A	Elution Buffer	50–100 µL <sup>[2]</sup>
Tip Comb	B	Place a KingFisher™ Duo 12-Tip Comb in Row B.	
Wash 4	C	Wash Solution 2	150 µL
Wash 3	D	Wash Solution 2	150 µL
DNase <sup>[3]</sup>	E	TURBO DNase™ Solution	50 µL
Wash 2	F	Wash Solution 2	150 µL
Wash 1	G	Wash Solution 1	150 µL

<sup>[1]</sup> Row on the MagMAX™ Express-96 Deep Well Plate.

<sup>[2]</sup> Use 50 µL for lower-yielding samples, such as brain or heart, or 100 µL for higher-yielding samples, such as liver or spleen.

<sup>[3]</sup> The instrument prompts the user to add 50 µL of Rebinding Buffer and 100 µL of isopropanol to the DNase Plate after the DNase treatment step.

3 Process samples on the instrument

Ensure that the instrument is set up with the deep well 96–well plates.

1. Select the appropriate program on the instrument (see “(If needed) Download, then install the program on the instrument” on page 2).
2. Start the run, then load the prepared processing plate when prompted by the instrument (see Table 8).
3. When prompted by the instrument (30–35 minutes after the initial start):

- a. Remove the plate from the instrument.
- b. Add 50 µL of Rebinding Buffer and 100 µL of isopropanol separately to each sample well in Row E. Add Rebinding Buffer and isopropanol immediately after the prompt to prevent excessive drying of any beads that are still captured on the Tip Comb.

**IMPORTANT!** Do not premix the Rebinding Buffer and isopropanol. Add them separately to the samples.

- c. Load the plate back onto the instrument, then tap **Start**.

4. At the end of the run (about 45 minutes after the initial start), remove the Elution Plate from the instrument, then transfer the eluted RNA (Row A) to an Elution Plate.
5. Seal immediately with a new MicroAmp™ Clear Adhesive Film.

**IMPORTANT!** Do not allow the purified samples to sit uncovered at room temperature for more than 10 minutes to prevent evaporation and contamination.

The purified samples are ready for immediate use. Alternatively, store the covered Elution Plate:

- On ice for up to 8 hours.
- At –20°C or –80°C for long-term storage.

**Limited product warranty**

Life Technologies Corporation and its affiliates 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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



**Revision history:** Pub. No. MAN0011132 G

Revision	Date	Description
G	13 September 2024	<ul style="list-style-type: none"><li>The Proteinase K concentration was removed from the contents and storage table.</li><li>Copy edits and formatting changes were made to align with current documentation style (throughout the document).</li></ul>
F.0	2 October 2023	Component errors were corrected and another plate type was added to required materials.
E.0	13 September 2022	The Box 2 storage condition was changed from 15–30°C to 15–25°C.
D.0	19 April 2021	Support was added for the KingFisher™ Apex Purification System.
C.0	14 December 2018	Centrifugation speeds were updated.
B.0	14 December 2015	Shaking conditions were added for RNA binding.
A.0	14 May 2015	Initial release.

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:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2015-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a trademark of Roche Molecular Systems, Inc., used under permission and license.