

PrepSEQ™ Residual DNA Sample Preparation Kit

USER GUIDE

Catalog Number 4413686

Publication Number MAN1001457

Revision A



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: MAN1001457 A (English)

Revision	Date	Description
A	20 March 2025	MAN1001457 Rev A is updated from 4469838 Rev L. Update the wording around Magnetic Particles and update Troubleshooting.
L	28 July 2022	Revision of 4469838, previous part number for <i>PrepSEQ™ Residual DNA Sample Preparation Kit User Guide</i> . Removed the instructions for the PrepSEQ™ Nucleic Acid Sample Preparation Kit (Cat. No. A50485). These instructions will now be found in the <i>PrepSEQ™ Nucleic Acid Sample Preparation Kit User Guide</i> (Pub. No. MAN0026641).
K	16 March 2022	Revision of 4469838, previous part number for <i>PrepSEQ™ Residual DNA Sample Preparation Kit User Guide</i> . Updates for the ViralSEQ™ Lentivirus Physical Titer Kit (Cat. No. A52597).
J	16 September 2021	Revision of 4469838, previous part number for <i>PrepSEQ™ Residual DNA Sample Preparation Kit User Guide</i> . Added the resDNASEQ™ Quantitative E1A DNA Fragment Length Kit (Cat. No. A51969) and the resDNASEQ™ Quantitative Synthetic Vero DNA Kit (A53242).

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.

©2021-2025 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■	CHAPTER 1	Product information	5
		Product description	5
		Contents and storage	5
		Required materials not supplied for manual protocols	6
		Required materials not supplied for automated protocols	7
		Workflow	8
■	CHAPTER 2	Prepare the reagents and samples	10
		Prepare the reagents: before first use of the kit	10
		Magnetic Particles	10
		Binding Solution	10
		Wash Buffer Concentrate	10
		Prepare reagents: before each use of the kit	10
		Proteinase K (PK) mix	10
		Lysis solution	11
		Guidelines for optimal yields	11
		Sample preparation guidelines	12
		Sample dilution (if necessary)	12
		Triplicate extractions	12
		(Kanamycin resistance and synthetic Vero assays only) Add Yeast tRNA	13
		Extraction control guidelines	13
■	CHAPTER 3	Manual protocol for DNA/RNA extraction	15
		Digest the test samples and controls	15
		Bind the DNA/RNA	15
		Wash the DNA/RNA	16
		Elute the DNA/RNA	17

- **CHAPTER 4 Automated protocol for DNA/RNA extraction 18**
 - Before each use of the kit 18
 - Ensure that you have the correct plates 18
 - Prepare the plates 18
 - Prepare the lysis plate 19
 - Process samples on the instrument 19
- **APPENDIX A Troubleshooting 21**
- **APPENDIX B Good laboratory practices 22**
 - Good laboratory practices for PCR and RT-PCR 22
 - Avoiding false positives due to cross-contamination 22
 - Plate layout suggestions 22
- **APPENDIX C Safety 23**
 - Chemical safety 24
 - Biological hazard safety 25
- Documentation and support 26**
 - Related documentation 26
 - Customer and technical support 26
 - Limited product warranty 26



Product information

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

Product description

The PrepSEQ™ Residual DNA Sample Preparation Kit uses chemical lysis and Magnetic Particles to extract DNA from diverse sample types, including samples that contain high protein and low DNA concentration. The kit extracts residual genomic DNA from products that are produced in cell lines such as CHO, *E. coli*, E1A, HEK293, Human, MDCK, NS0, *Pichia*, Sf9 and Baculovirus, and Vero or residual plasmid DNA for Kanamycin resistance. For quantification of residual DNA, we recommend using the resDNASEQ™ Quantitative DNA Kits as described in the *resDNASEQ™ Quantitative DNA Kits User Guide* (Pub. No. 4469836) or *resDNASEQ™ Quantitative E1A DNA Fragment Length Kit User Guide* (Pub. No. MAN0025643).

To help ensure accurate quantitative results, extract each sample in triplicate and perform a single PCR reaction for each extraction.

Contents and storage

The kits contain reagents sufficient for 100 extractions.

Table 1 PrepSEQ™ Residual DNA Sample Preparation Kit (Cat. No. 4413686)

Contents	Amount	Storage
Box 1, PrepSEQ™ Nucleic Acid Extraction Kit		
Lysis Buffer	2 × 50 mL	Room temperature
Binding Solution (Isopropanol), empty bottle	1	
Wash Buffer Concentrate	2 × 26 mL	
Elution Buffer	25 mL	
Proteinase K (PK) Buffer Can be used for existing validated manual protocols.	50 mL	
Proteinase K (PK) Buffer II ^[1] Recommended for new manual protocols. Required for automated protocols.	11 mL	

Table 1 PrepSEQ Residual DNA Sample Preparation Kit (Cat. No. 4413686) (continued)

Contents	Amount	Storage
Box 2, PrepSEQ™ Nucleic Acid Extraction Kit		
Magnetic Particles	2 × 1.5 mL	Room temperature
Box 3, PrepSEQ™ Nucleic Acid Extraction Kit		
Proteinase K, 20 mg/mL	1.25 mL	–20°C or below
PrepSEQ™ Residual DNA Sample Preparation Kit		
Proteinase K, 20 mg/mL	1.25 mL	–20°C or below
Yeast tRNA, 10 mg/mL	0.5 mL	
Glycogen, 5 mg/mL	2 × 1.0 mL	

^[1] Also sold separately (Cat. No. 4415320).

Required materials not supplied for manual protocols

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.

Item	Source
Equipment	
Magnetic stand, 16-position	12321D
Block heater for use with 2-mL tubes. Manual DNA/RNA extraction involves two incubations at different settings, so two heaters may be convenient.	MLS
Benchtop microcentrifuge for 1.5-mL and 2-mL tubes	MLS
Vortex	MLS
Vortex Adapter-60, for use with the Vortex-Genie™	AM10014
Consumables	
Disposable gloves	MLS
Aerosol-resistant micropipette tips	MLS
Pipettors, P1000, P200, P20 and P10: <ul style="list-style-type: none"> Positive-displacement Air-displacement Multichannel 	MLS

Item	Source
Nonstick, RNase-free Microfuge Tubes, 1.5-mL (1 box; 250 tubes/box)	AM12450
Safe-Lock Tube, 2.0-mL	VWR™ 62111-754
Reagents	
Ethanol, 95% IMPORTANT! Do not use denatured ethanol. It contains components that are not compatible with the protocol.	MLS
Isopropanol, 100%	MLS
1X PBS (free of Mg and Ca)	MLS

Required materials not supplied for automated protocols

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 Pharma KingFisher™ Flex Purification System with 96 Deep-Well Head (Cat. No. A31508) and Pharma MagMAX™ Express-96 instrument^[1] accessories

Item	Source
Pharma MagMAX™ Express-96 DW plate	A31540
Pharma MagMAX™ Express-96 Deep-Well Tip Combs	A31537
Pharma KingFisher™ Flex Magnetic Head for 96 Deep-Well Plate	A31542
Pharma MagMAX™ 96 PCR Well Magnetic Head	4472991
Pharma MagMAX™ Express-96 Standard Plates	A31541
(Optional) Pharma Magnetic Stand-96	A31543

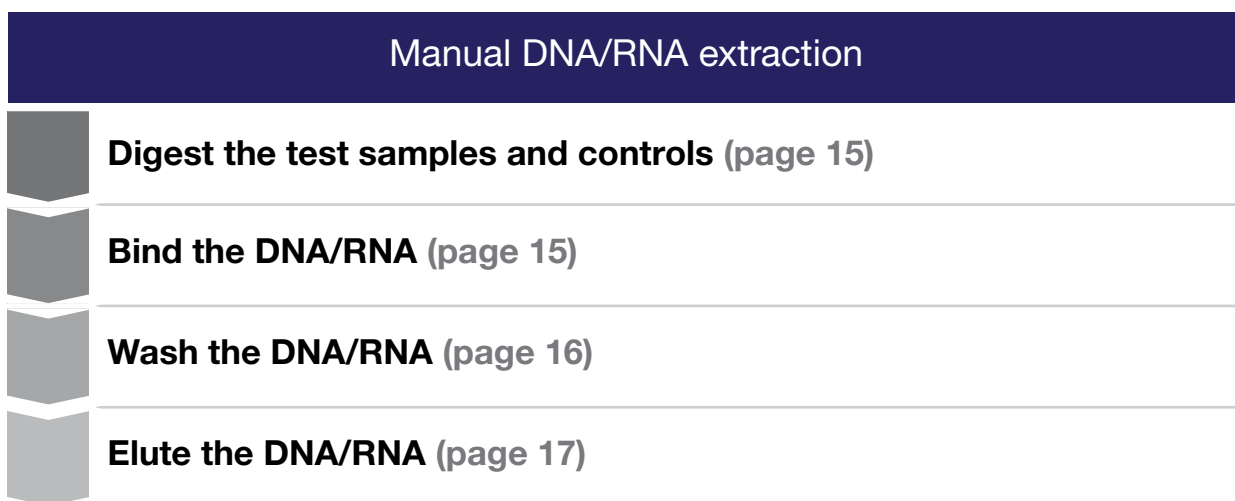
^[1] This instrument is no longer available for purchase

Table 3 Additional materials

Item	Source
Consumables	
Disposable gloves	MLS
Aerosol-resistant micropipette tips	MLS

Item	Source
Pipettors, P1000, P200, P20 and P10: <ul style="list-style-type: none"> • Positive-displacement • Air-displacement • Multichannel 	MLS
Nonstick, RNase-free Microfuge Tubes, 1.5-mL (1 box; 250 tubes/box)	AM12450
Reagents	
Ethanol, 95% IMPORTANT! Do not use denatured ethanol. It contains components that are not compatible with the protocol.	MLS
Isopropanol, 100%	MLS
1X PBS (free of Mg and Ca)	MLS

Workflow



Automated DNA/RNA extraction



Prepare the plates (page 18)



Prepare the lysis plate (page 19)



Process samples on the instrument (page 19)



Prepare the reagents and samples

Prepare the reagents: before first use of the kit

Magnetic Particles

1. Inspect the Magnetic Particle tube before use.
If particle clumping or white precipitates are observed, see Appendix A, “Troubleshooting”.
2. Set the block heater to 37°C.
3. Incubate the Magnetic Particle suspension at 37°C for at least 10 minutes with intermittent vortexing at high speed or 900 rpm until the particles are completely suspended.

Binding Solution

1. Add 45 mL of 100% isopropanol to the Binding Solution bottle.
2. Label the bottle to indicate that it contains isopropanol, then store the bottle at room temperature.

Wash Buffer Concentrate

1. Add 74 mL of 95% ethanol to one bottle of PrepSEQ™ Wash Buffer Concentrate, then mix completely.
2. Label the bottle to indicate that it contains ethanol, then store the bottle at room temperature.

Prepare reagents: before each use of the kit

Proteinase K (PK) mix

- Use Proteinase K (PK) Buffer II for all protocols.

Note: Proteinase K (PK) Buffer is provided in the kit for use by laboratories that had previously validated with this buffer. If required, use exactly as described for Proteinase K (PK) Buffer II.

- Prepare a fresh mix before each use of the kit.
- Include a 10% overage to account for pipetting losses.

Component	Number of extractions				
	1	7	10	13	25
Proteinase K, 20 mg/mL	10 µL	70 µL	100 µL	130 µL	250 µL
Proteinase K (PK) Buffer II	60 µL	420 µL	600 µL	780 µL	1,500 µL

Lysis solution

- Prepare a fresh mixture immediately before use or during Proteinase K incubation.
- Prepare 360 µL (amount required) of lysis solution mix per sample.

Reagent	Volume for 1 extraction	Volume for ~20 extractions
Glycogen, 5 mg/mL	8.32 µL	180 µL
Yeast tRNA, 10 mg/mL	0.18 µL	4 µL
Lysis Buffer	351.5 µL	7,600 µL
Total	360 µL	7,784 µL

Guidelines for optimal yields

- Maintain a homogenous suspension of the Magnetic Particles to maximize the surface area to which the DNA/RNA can bind. The appearance of the mixture should be homogenous after mixing.
- After drying, the DNA/RNA remains bound to the Magnetic Particles. Do not allow the Magnetic Particles to over-dry because this decreases the elution efficiency. Over-dried particles are not easily resuspended.
- During manual elution, vortex every 2 minutes to help elution. This results in better yield during recovery.

Note: Some test samples cause the particles to adhere firmly to the tube wall. Other test samples form loose pellets that detach during the vortex steps. All pellets should dissolve with vortexing during heated elution. If vortexing does not result in full resuspension, then wash the particles off the tube by pipetting.

Sample preparation guidelines

Sample dilution (if necessary)

Test samples from the early purification process often contain levels of DNA/RNA that are above the highest point of the assay standard curve. You must dilute these samples (from 1:100 up to 1:10,000) before PrepSEQ™ sample preparation.

- Dilute test samples before DNA/RNA extraction with a solution of 1X PBS (free of Mg and Ca) or 50 mM Tris, pH 8.0, 0.5 M NaCl.

Note: Diluting samples in water or TE reduces extraction efficiency.

- Use the sample dilution buffer as the negative extraction control instead of water.
- Alternatively, dilute extracted DNA/RNA with elution buffer before running the PCR reaction.

Triplicate extractions

Triplicate extractions are required for post-PCR analysis calculation of mean quantity, standard deviation, and coefficient of variation.

In addition to test samples, we recommend triplicate extractions for the negative control and the extraction/recovery control (ERC).

Perform a single PCR reaction for each extraction.

The table below illustrates the total number of extractions required based on the 1, 2, and 3 samples extracted in a batch.

Table 4 Total number of extractions per batch of test samples

Number of test samples		Total number of extractions for the batch
1	3 extractions required for each:	9
2	<ul style="list-style-type: none"> • Test sample • Test sample extraction/recovery control (ERC) 	15
3	<ul style="list-style-type: none"> • Negative extraction control^[1] 	21

^[1] Optional during routine testing.

(Kanamycin resistance and synthetic Vero assays only) Add Yeast tRNA

(Recommended) Use **Yeast tRNA** with the resDNASEQ™ Quantitative Plasmid DNA - Kanamycin Resistance Gene Kit and the resDNASEQ™ Quantitative Synthetic Vero DNA Kit.

1. Dilute the **Yeast tRNA**.

Table 5 Diluted Yeast tRNA

Component	Volume
Yeast tRNA (10 mg/mL)	5 µL
1X PBS (free of Mg and Ca)	245 µL
Total	250 µL

2. Add 10 µL **Diluted Yeast tRNA** to 740 µL of each test sample before extraction. This is sufficient for triplicate 100 µL extractions.

Extraction control guidelines

We recommend that you use the following extraction controls:

Type of control	Contains	Number to run	Used to
Negative (NEG) ^[1]	1X PBS (free of Mg and Ca)	1 per batch of extractions	Test for cross-contamination of DNA/RNA extraction reagents.
Extraction/recovery (ERC)	Positive control from the kit	3 per sample	<ul style="list-style-type: none"> Evaluate the efficiency of DNA or RNA extraction, recovery, and quantification from test samples. Verify reagent and system performance.

^[1] Optional during routine testing.

For the Extraction/recovery (ERC) :

- Prepare the control standard dilutions as described in the *resDNASEQ™ Quantitative DNA Kits User Guide* (Pub. No. 4469836) or *resDNASEQ™ Quantitative E1A DNA Fragment Length Kit User Guide* (Pub. No. MAN0025643).
- Add a volume of positive control standard dilution to each test sample so that the total DNA/RNA amount is 2–10 times the amount of DNA/RNA measured in the test sample *without* the addition of the DNA/RNA control.

For example (amount):

- The DNA amount measured in a test sample is ≤ 1 pg.
- To prepare a 10 pg ERC for a PCR elution volume of 50 µL, spike samples with 16.7 µL of the 3 pg/µL positive control standard dilution (SD3) = 50 pg spike to yield 10 pg per PCR reaction.

For example (copies):

- The number of copies in a test sample is $\leq 2,000$ copies.
- Spike control sample at 1×10^5 copies using appropriate standard control concentration.
- If 1:20 (10 μL out of 200 μL eluate) of the KingFisher™ Flex extraction is used for qPCR, there are 5,000 expected copies of spike control per PCR reaction.
- If 1:5 (10 μL out of 50 μL eluate) of the manual extraction is used for qPCR, there are 20,000 expected copies of spike control per PCR reaction.
- Prepare three separate extractions for each test sample, then add the ERC to each reaction. Do not prepare a large volume of ERC, then aliquot it into three reactions.

Note: To calculate the efficiency of DNA/RNA recovery and quantification from the test samples, subtract the amount of DNA/RNA measured in the sample *without* the addition of DNA/RNA control from the amount of DNA/RNA measured in the ERC sample.



Manual protocol for DNA/RNA extraction

Digest the test samples and controls

1. Set a block heater to 56°C. If available, set a second block heater to 70°C.
2. Label the Safe-Lock Tubes, 2.0-mL:
 - 3 for each sample
 - 3 for each sample + ERC
 - 3 for NEG
3. Add 100 µL of sample, sample + ERC, or 1X PBS (free of Mg and Ca) into each tube.

Note: Ensure that **Diluted Yeast tRNA** was added to each Kanamycin resistance or Vero 2.0 sample. See “(Kanamycin resistance and synthetic Vero assays only) Add Yeast tRNA” on page 13.

4. Add 10 µL of 5 M NaCl and 70 µL Proteinase K/Proteinase K Buffer II mix.
5. Briefly vortex and centrifuge.
6. Incubate at 56°C for 30 minutes.
If only one block heater is available, after this incubation step is complete, reset the block heater to 70°C for the elution step.

Note: For samples with high protein concentration, extending the incubation time to 60 minutes can increase recovery.

7. Cool samples to room temperature.
8. Add 360 µL freshly made Lysis solution mix to each tube.

Bind the DNA/RNA

1. Vortex the Magnetic Particles to resuspend the particles.

Note: The appearance of the mixture should be homogeneous.

2. Add 30 µL of the Magnetic Particles to each tube.

3. Add 400 μL **Binding Solution** to the first sample, then close the cap and invert twice to mix.
4. Repeat step 3 for each additional sample.

Note: Low DNA recovery can be observed if the sample and the **Binding Solution** are not immediately mixed.

5. Vortex all the tubes in the vortex adaptor for 5 minutes at 900 rpm.
6. Briefly centrifuge the tubes for 15 seconds at top speed ($>15,000 \times g$) to collect the Magnetic Particles at the bottom of the tubes.
7. Place the tubes in the magnetic stand with the pellet against the magnet, then let the tubes stand for 5 minutes or until the solution is clear.
8. Without disturbing the Magnetic Particles, remove the supernatant using a pipette or by aspiration.

Wash the DNA/RNA

1. Remove the tube rack (with tubes) from the magnetic stand, then add 300 μL of Wash Solution to the tubes. Vortex the tubes for 5 seconds at room temperature at 900 rpm.
2. Centrifuge the tubes in a microcentrifuge at top speed ($>15,000 \times g$) for a maximum of 20 seconds. Do not centrifuge for >20 seconds.
3. Place the tubes in the magnetic stand, then let the tubes stand for 1 minute.

Note: The Magnetic Particles with the bound DNA/RNA are magnetically captured after about 1 minute.

4. Without disturbing the Magnetic Particles, remove the supernatant by pipette or by aspiration.
5. Remove the tube rack (with tubes) from the magnetic stand, then add 300 μL of Wash Solution to each tube for a second wash. Vortex the tubes for 5 seconds at room temperature at 900 rpm.
6. Centrifuge the tubes in a microcentrifuge at top speed ($>15,000 \times g$) for a maximum of 20 seconds. Do not centrifuge for >20 seconds.
7. Place the tubes in the magnetic stand, then let the tubes stand for 1 minute.

Note: The Magnetic Particles with the bound DNA/RNA are magnetically captured after about 1 minute.

8. Open all tubes, then start the 5-minute timer.
9. Without disturbing the Magnetic Particles, remove the supernatant by pipette or aspiration. Use a P200 to remove the remaining solution from the bottom of the tube.

10. With the tube lid open, air-dry the Magnetic Particles pellet in the magnetic stand for no more than 5 minutes at room temperature.

IMPORTANT! Air-dry to remove ethanol from the Wash Solution. After dry, the DNA/RNA remains bound to the Magnetic Particles. Do not over-dry; over-dried particles are not easily resuspended.

Elute the DNA/RNA

1. Add 50 μ L of Elution Buffer to each tube.
2. Vortex the tubes for 20 seconds at high speed, then incubate the tubes at 70°C for 7 minutes. Vortex the tubes two to three times during the incubation to help resuspend the particles.
3. Centrifuge the tubes in a microcentrifuge at top speed ($>15,000 \times g$) for a maximum of 20 seconds. Do not centrifuge for >20 seconds.
4. Place the tubes in the magnetic stand, then let the tubes stand for 1 minute.
5. Without disturbing the Magnetic Particles, transfer the liquid phase containing the eluted DNA/RNA to a new nonstick 1.5-mL microcentrifuge tube.
6. Centrifuge the tube at top speed ($>15,000 \times g$) for 3 minutes to collect the Magnetic Particles at the bottom of the tube, then place the tubes in the magnetic stand for 1 minute.
7. Without disturbing the Magnetic Particles, transfer the liquid phase containing the eluted DNA/RNA to a new nonstick 1.5-mL microcentrifuge tube.

Note: Magnetic Particles can inhibit PCR.

Store eluted DNA/RNA for up to 6 hrs on ice, or up to 24 hrs at -20°C .

4

Automated protocol for DNA/RNA extraction

You can use the KingFisher™ Flex or MagMAX™ Express 96-deep well automation platforms to automate the extraction of DNA/RNA. For all chemicals, read the Safety Data Sheet (SDS) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Before each use of the kit

Ensure that you have the correct plates

The KingFisher™ Flex or the MagMAX™ Express require 5 plates.

Plate name	Plate type
Lysis	96 deep-well plate
Wash 1	96 deep-well plate
Wash 2	96 deep-well plate
Elution	96 deep-well plate
Comb loading plate	96 deep-well tip comb combined with 96 standard plate

Prepare the plates

Prepare the Wash 1, Wash 2, and Elution plates:

Plate name	Plate type	Volume of buffer to add
Wash 1	96 deep-well plate	300 µL of Wash buffer
Wash 2	96 deep-well plate	300 µL of Wash buffer
Elution	96 deep-well plate	200 µL of Elution buffer

Prepare the lysis plate

In all steps that require pipetting, dispense liquid at bottom center of the wells.

1. Add 100 µL to the appropriate wells of the 96 deep-well Lysis plate:

- 3 wells for each sample
- 3 wells for each sample + ERC
- 3 wells for NEG

Note: Ensure that **Diluted Yeast tRNA** was added to each sample when using the resDNASEQ™ Quantitative Plasmid DNA - Kanamycin Resistance Gene Kit. See “(Kanamycin resistance and synthetic Vero assays only) Add Yeast tRNA” on page 13.

2. Add 10 µL of 5 M NaCl to each sample well.
3. Add 70 µL Proteinase K/Proteinase K (PK) Buffer II mix to each sample well.

Process samples on the instrument

1. Select the script for the instrument and kit that you are using:

Kit	Instrument	Script
PrepSEQ™ Residual DNA Sample Preparation Kit	KingFisher™ Flex	PrepSEQ_resDNA_v1
	MagMAX™ Express-96	PrepSEQ_resDNA_2011 PrepSEQ_1hr_resDNA (if installed)

2. Load the plates into the instrument in the order listed below. After loading each plate, press **START** to move the turntable.
 - a. Comb loading plate
 - b. Elution plate with 200 µL of Elution Buffer
 - c. Wash 2 plate with 300 µL of wash buffer
 - d. Wash 1 plate with 300 µL of wash buffer
 - e. Lysis plate
3. Press **START** to begin the PK digestion process.
The instrument mixes the samples for 10 seconds at fast speed, then incubates the samples at 56°C for 30 minutes, mixing at slow speed. When digestion is complete, the instrument pauses and returns the Lysis plate to the loading position.
4. After the digestion step is complete, add additional components to the Lysis plate:
 - a. Remove the Lysis plate from the instrument.
 - b. Add 360 µL of Lysis Solution to each sample well.

- c. Add 30 µL of Magnetic Particle suspension to each sample well.
- d. Add 400 µL of Binding Solution to the first sample well, then immediately pipet up-and-down three times to mix. Repeat for each additional sample well.

Note: Low DNA recovery can be observed if the sample and the **Binding Solution** are not immediately mixed.

- e. Place the plate back into the instrument loading position, then press **START** to begin binding.
5. When DNA/RNA extraction is finished, the instrument returns the Elution plate to the loading position.

Store eluted DNA/RNA for up to 6 hrs on ice, or up to 24 hrs at –20°C.

Upon completion of the sample extraction procedure, see the relevant assay user guide for specific assay instructions.



Troubleshooting

Observation	Possible cause	Recommended action
Poor extraction efficiency (low yields)	Over-drying the sample.	Start the 5-minute timer before removing ~300 μ L from the first 6–8 samples. Then continue removing wash buffer from the remaining samples.
	Magnetic Particles are difficult to resuspend during the elution.	<ul style="list-style-type: none"> Incubate the pellets at 70°C for >7 minutes. Vigorously vortex the tubes three times during incubation to help resuspension. Do not over-dry. If needed, repeat the incubation and vortexing steps.
	PK Buffer was used instead of PKII Buffer.	Use PKII Buffer.
Particles no longer produce consistent results Fine brown sandy particles and brown color are observed in the supernatant.	Samples have low pH.	Determine the pH of the sample and adjust the pH to between 6 and 8.
	Magnetic Particles were stored at –20°C.	Order new materials. Store the new materials at room temperature.
Clumps or white precipitate in the Magnetic Particles tube	The Magnetic Particles were shipped or stored at temperatures below room temperature.	<ol style="list-style-type: none"> Incubate the Magnetic Particles tube at 37°C for 30 minutes. During incubation, vortex intermittently at high speed or 900 rpm. Continue to incubate and vortex until all clumps or white precipitate is dissolved.
		For hard to resuspend particles, Magnetic Particles incubation can be performed at 50°C for 30 minutes. In addition to vortexing, pipette tip displacement is recommended to help ensure particles are in solution before use.



Good laboratory practices

Good laboratory practices for PCR and RT-PCR

- Wear clean gloves and a clean lab coat.
 - Do not wear the same gloves and lab coat that you have previously used when handling amplified products or preparing samples.
- Change gloves if you suspect that they are contaminated.
- Maintain separate areas and dedicated equipment and supplies for:
 - Sample preparation and reaction setup.
 - Amplification and analysis of products.
- Do not bring amplified products into the reaction setup area.
- Open and close all sample tubes carefully. Avoid splashing or spraying samples.
- Keep reactions and components capped as much as possible.
- Use a positive-displacement pipettor or aerosol-resistant barrier pipette tips.
- Clean lab benches and equipment periodically with a 10% bleach solution or decontamination solutions (for example, DNAZap™ PCR DNA Degradation Solutions (Cat. No. [AM9890](#)) and RNaseZap™ RNase Decontamination Solution (Cat. No. [AM9780](#)).

Avoiding false positives due to cross-contamination

To avoid false positives due to cross-contamination:

- Prepare and close all negative control and unknown sample tubes before pipetting the positive control.
- Do not open tubes after amplification.
- Use different sets of pipettors when pipetting negative control, unknown, and positive control samples.

Plate layout suggestions

- For each plate row, dispense in sequence from left to right the: negative controls, unknown samples and ERCs, and positive controls (at the end of the row or column).
- Place positive controls in one of the outer columns (10–12).
- If possible, separate all samples from each other by at least one well; if space is limiting, place at least one well between unknown samples and controls.



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
<https://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>PrepSEQ™ Residual DNA Sample Preparation Kit Quick Reference</i>	4469839
<i>resDNASEQ™ Quantitative DNA Kits User Guide</i>	4469836
<i>resDNASEQ™ Quantitative E1A DNA Fragment Length Kit User Guide</i>	MAN0025643
<i>Thermo Scientific™ KingFisher™ Flex User Manual</i>	MAN0019870
<i>Applied Biosystems™ MagMAX™ Express 96 User Manual</i>	N07849

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.

