

PRODUCT INFORMATION

Thermo Scientific MagJET Plant RNA Kit

#K2771, #K2772

Read Storage information (p. 4) upon receipt and store kit components appropriately!

www.thermoscientific.com/onebio

#K2771 Lot 00000000 Exp. 00.0000

CERTIFICATE OF ANALYSIS

Thermo Scientific[™] MagJET[™] Plant RNA Kit is qualified by isolation of total RNA from 50 mg of plant tissue following the protocol outlined in the manual. The quality of isolated RNA is evaluated spectrophotometrically and by agarose gel electrophoresis. The purified RNA has an A₂₆₀/A₂₈₀ ratio of 2.0±0.2 and the RNA integrity number (RIN) of ≥7.

Quality authorized by:

Jurgita Žilinskienė

Rev.1

CONTENTS	page
COMPONENTS OF THE KIT	4
STORAGE	4
DESCRIPTION	4
PRINCIPLE	4
IMPORTANT NOTES	5
ADDITIONAL MATERIALS AND EQUIPMENT REQUIRED	6
AVOIDING RIBONUCLEASE CONTAMINATION	6
RNA HANDLING AND STORAGE	6
STARTING MATERIAL HANDLING AND SAMPLE HOMOGENIZATION	7
PROTOCOL SELECTION GUIDE	9
TOTAL RNA PURIFICATION PROTOCOLS AND PIPETTING INSTRUCTIONS	10
Protocol A. Instructions for total RNA purification from up to 50 mg plant sample using KingFisher Flex 96 and Microtiter deep well 96 plates	10
Protocol B. Instructions for total RNA purification from up to 50 mg plant sample using KingFisher Duo with 12-pin magnet head and Microtiter deep well 96 plate	12
Protocol C. Instructions for manual plant total RNA purification from up to 50 mg plant sample	14
TROUBLESHOOTING	15
SAFETY INFORMATION	16

COMPONENTS OF THE KIT

MagJET Plant RNA Kit	#K2771 96 preps	#K2772 384 preps
Lysis Buffer for MagJET Plant RNA Kit	72 mL	2 × 140 mL
MagJET Magnetic Beads	2 × 1.4 mL	10.6 mL
DNase I (lyophilized)	1 vial	4 vials
DNase I Reconstitution Buffer	1 mL	2 × 1 mL
2X DNase I Buffer	12 mL	45 mL
Manganese Chloride Solution	3 × 1 mL	9 × 1 mL
Wash Buffer 1 (conc.) for MagJET Plant RNA Kit	110 mL	3 × 110 mL
Wash Buffer 2 (conc.) for MagJET Plant RNA Kit	50 mL	3 × 50mL
Rebinding Buffer for MagJET Plant RNA Kit	20 mL	70 mL
Water, nuclease free	30 mL	125 mL

STORAGE

When the kit is delivered, remove DNase I (lyophilized), DNase I Reconstitution Buffer and Manganese Chloride Solution and store at -20°C. Reconstituted DNase I should be stored at -20°C. MagJET Magnetic Beads should be stored at 4°C. Other components of the kit should be stored at room temperature (15-25°C).

DESCRIPTION

The MagJET RNA Kit is designed for automated high throughput or manual purification of total RNA from a wide variety of plant species and tissue types.

The kit utilizes paramagnetic bead technology enabling high yields and robust performance. High binding capacity, uniform particle size, and rapid magnetic response of MagJET magnetic beads makes the technology ideal for high throughput automatic nucleic acid purification, as well as for manual nucleic acid purification by low sample throughput users.

The resulting high quality purified RNA is free of proteins, nucleases and other contaminants or inhibitors, and can be used in a wide range of downstream applications, such as RT-PCR, RT-qPCR and other enzymatic reactions. See Table 1 for typical total RNA yields from various sources.

PRINCIPLE

The MagJET Plant RNA Kit uses the highly efficient MagJET magnetic particle-based technology for nucleic acid purification. The nucleic acid isolation process combines the simple steps of sample lysis, RNA binding to the magnetic beads, DNA digestion, washing and elution.

Purification protocols optimized for automated KingFisher instruments utilize a high throughput magnetic bead transfer technique, where magnetic beads are transferred through different reagent plates containing lysis, binding, washing and elution reagents. This enables high throughput nucleic acid purification and eliminates multiple pipetting steps.

Alternatively, a protocol is provided where buffers and other reagents are transferred in each of the protocol steps, while magnetic beads remain captured on the wall of a tube using a magnetic rack. This allows the kit to be used in various throughput applications using a magnetic rack and manual or automated pipetting equipment.

Plant	Tissue (50 mg)	RNA yield, µg
Arabidopsis thaliana	leaf	18-20
Nicotiana tabacum	leaf	15-20
Tomato	leaf	21-52
Barley seedlings	leaf	25-32
Wheat shoots	leaf	25-36
Spinach	leaf	24-30
Onion	leaf	8-12
Rice	leaf	16-19
Corn	leaf	13-20
	seeds	20-40
	leaf	11-19
Canola (Rape)	stalk	22-28
	seeds	30
Sav	leaf	30-40
Soy	seeds	44
Cucumber	fruit	18
Potato	stalk	10
Sunflower	stalk	4-5
Sulliowei	seeds	40

Table 1. Typical yields of total RNA from various sources.

IMPORTANT NOTES

 Add the indicated volumes of ethanol (96-100%) to Wash Buffer 1 (conc.) and Wash Buffer 2 (conc.) prior to first use:

	#K2771 and #K2772		
	Wash Buffer 1	Wash Buffer 2	
Concentrated buffer	110 mL	50 mL	
Ethanol (96-100%)	110 mL	200 mL	
Total volume:	220 mL	250 mL	

After preparing each solution, mark the bottle to indicate that this step has been completed.

- To prepare the DNase I solution add 0.44 mL of DNase I Reconstitution Buffer to each vial of DNase I (lyophilized) and incubate at room temperature for 5 minutes. Occasional gentle rotation of the vial helps to dissolve the DNase I, but avoid forceful mixing. Do not vortex! Store at -20°C.
- Before each RNA purification, prepare a fresh aliquot of Lysis Buffer supplemented with
 2 M DTT solution. Add 20 μL of 2 M DTT to each 1 mL of Lysis Buffer needed.
- Check all solutions in the kit for salt precipitation before each use. Re-dissolve any
 precipitates by warming the solution at 37 °C, and then equilibrate to room temperature
 (20 ± 5°C).
- Wear gloves when handling the Lysis Buffer, Rebinding Buffer and Wash Buffer 1 as these reagents contain irritants (see page 16 for SAFETY INFORMATION).

ADDITIONAL MATERIALS AND EQUIPMENT REQUIRED

- 2 M DTT (dithiotreitol) solution (store at -20°C).
- polyvinylpyrrolidone (PVP)(for RNA purification from woody, lignified and or polyphenol-rich samples, see p.7-8).
- Sodium chloride (NaCl) (for RNA purification from soy beans, see p.7-8).
- 96-100% ethanol, molecular biology grade.
- Pipettes and RNase-free pipette tips.
- 8 or 12 channel pipettes and pipette tips.
- RNase-free 1.5-2 mL plastic tubes.
- Disposable gloves.
- Thermomixer.
- Centrifuge capable of \geq 16,000 × g for microtubes.
- Centrifuge capable of 3,000-4,000 × g with swinging-bucket rotor for 96-well plates.
- Incubator or water bath capable of 56°C.
- Vortex.
- Mortar and pestle (for manual tissue disruption in liquid nitrogen).
- Liquid nitrogen.
- Commercial homogenizer and stainless steel beads (for mechanical automatic plant tissue disruption).
- Automatic magnetic particle processor and consumables.
- For manual protocol acquire magnetic particle processing rack.

AVOIDING RIBONUCLEASE CONTAMINATION

RNA purity and integrity is essential for downstream applications. RNA can be degraded by RNase A, which is a highly stable contaminant found in any laboratory environment. Care must be taken not to introduce RNases into RNA preparations, especially during wash with Wash Buffer 2 and elution steps. General recommendations to avoid RNase contamination include the following:

- Wear gloves when handling reagents and RNA samples, as skin is a common source of RNases. Change gloves frequently.
- Use sterile, disposable RNase-free pipette tips.
- Use appropriate reagents to remove RNase contamination from nondisposable items (pipettes, centrifuges) and work surfaces.
- Keep all kit components tightly sealed when not in use. After usage close bottles immediately.

RNA HANDLING AND STORAGE

- Keep the RNA on ice after extraction and while working with it.
- Store the extracted RNA at -20°C or -70°C. For long term use store RNA at -70°C.

STARTING MATERIAL HANDLING AND SAMPLE HOMOGENIZATION

- To minimize RNA degradation, avoid repeated freezing and thawing of the samples and perform extractions from fresh material, or material that has been immediately frozen and stored at **-70°C**.
- Appropriate sample storage is essential for reproducibility and high RNA yields. Yields of RNA may vary depending on sample age, type of sample, and storage conditions.
- Efficient homogenization of the sample material is an essential step before RNA purification in order to gain a good yield of high-quality RNA. The lysis procedure of plant tissue is most effective with well homogenized, powdered samples. Plant tissue can be homogenized using a mortar and pestle in the presence of liquid nitrogen or using commercial homogenizers with steel beads.

Manual plant tissue homogenization

- 1. Place the plant sample in a mortar with liquid nitrogen. Grind the tissue using a pestle.
- Allow the liquid nitrogen to evaporate and transfer the powder sample (10- 50 mg) into a 1.5 mL microcentrifuge tube containing 600 µL Lysis Buffer supplemented with DTT as described on page 5.
 - Transfer the homogenized tissue to the Lysis Buffer as quickly as possible to avoid RNA degradation.
 - All homogenized material must be thoroughly mixed with the Lysis Buffer. RNA degradation can occur in particles that are left to dry on the walls of the tube.
- 3. Vortex for 10-20 s to mix thoroughly.
- 4. Incubate for 5 min at 56°C.
- 5. Centrifuge for 10 min at maximum speed (\geq 16,000 × g).
- 6. Carefully take 400 μL of the cleared supernatant and immediately proceed with the RNA isolation according to protocols **A C**.

Notes for lysis of different plant samples:

- For seed samples, use 10-20 mg starting material into the Lysis Buffer.
- To purify total RNA from woody, lignified and or polyphenol-rich samples such as branches, twigs, needles, wax-coated leaves (such as laurel), supplement the **Lysis Buffer** with polyvinylpyrrolidone (PVP) at a 2% (w/v) final concentration.
- To purify total RNA from soy beans, supplement the Lysis Buffer with sodium chloride (NaCl) at a 2 M final concentration.

Mechanical (automatic) tissue disruption

Plant tissue can be homogenized with commercial homogenizers and grinding mills (with stainless steel beads). High-throughput homogenizers offer a suitable method for handling up to 96 samples simultaneously.

- 1. Place up to 50 mg of plant sample (or up to 20 mg for seeds) into microtube containing a stainless steel bead and 600 µL Lysis Buffer supplemented with DTT as described on page 5.
- 2. Grind the sample according to homogenizer instructions.
- 3. Incubate for 5 min at 56°C.
- 4. Centrifuge for 10 min at maximum speed (≥16,000 × g for microtube or 3,000–4,000 × g for racked 96-well collection microtubes).
- 5. Carefully take 400 µL of the cleared supernatant and immediately proceed with the RNA isolation according to protocols **A C**.

Notes for lysis of different plant samples:

- Use 10-50 mg of fresh plant sample for mechanical automatic disruption (with stainless steel beads) in single microtubes*.
- Use 10-20 mg plant Isample for mechanical automatic disruption (with stainless steel beads) in racked 96-well collection microtubes*.
- Use 10-20 mg of plant seeds as starting material for mechanical automatic disruption (with stainless steel beads) in racked 96-well collection microtubes and/or single mictrotubes*.
- To purify total RNA from woody, lignified and or polyphenol-rich samples such as branches, twigs, needles, wax-coated leaves (e.g. laurel), supplement the Lysis Buffer with polyvinylpyrrolidone (PVP) at a 2% (w/v) final concentration.
- To purify total RNA from soy beans, supplement the Lysis Buffer with sodium chloride (NaCl) at a 2 M final concentration.
- *- Proceed according homogenizer manufacturer recommendations.

PROTOCOL SELECTION GUIDE

The MagJET Plant RNA Kit provides optimized protocols for total RNA purification from different amounts of starting material (10-50 mg). The Kit is compatible with automated and manual sample processing, allowing high and low-throughput nucleic acid purification workflows.

The following selection guide summarizes the available protocols depending on starting sample weight, throughput and sample processing method. Automation protocols are optimized for KingFisher Flex and KingFisher Duo instruments.

Protocol selection guide:							
Plant homogenization type	Sample weight	Throughput per run	KingFisher Flex Instrument	KingFisher Duo Instrument	Manual processing	MagJET purification protocol	Page
Plant homogenized		Up to 96	•	-	-	Protocol A	10
with liquid nitrogen (see	10-50 mg (10-20 mg of seeds)	Up to 12	-	•	-	Protocol B	12
the detailed instruction on page 7)		variable	-	-	•	Protocol C	14
Plant homogenized with equipment for disrupting plant tissue* (see the detailed instruction on page 8)		Up to 96	•	-	-	Protocol A	10
	10-50 mg (10-20 mg of seeds)	Up to 12	-	•	-	Protocol B	12
		variable	-	-	•	Protocol C	14

Protocol selection guide:

*Sample quantity may depend on the homogenizer type and throughput. Please proceed according homogenizer manufacturer recommendations.

TOTAL RNA PURIFICATION PROTOCOLS AND PIPETTING INSTRUCTIONS

Protocol A. Instructions for total RNA purification from up to 50 mg plant sample using KingFisher Flex 96 and Microtiter deep well 96 plates

Note:

- When using the MagJET Plant RNA Kit for the first time, prepare working solutions of Wash Buffer 1 and Wash Buffer 2 and reconstitute DNase I as described on page 5.
- Transfer the **Plant_RNA_Flex** protocol file to the KingFisher instrument before first use. The instructions for transferring the protocol can be found in Chapter 4: "Using the software" in the Bindlt Software for KingFisher Instruments version 3.2 User Manual. The protocol files for MagJET Plant RNA Kit can be found on the product web page on <u>www.thermoscientific.com/onebio</u>
- 1. Homogenize the plant samples and prepare lysates as described on page 7-8 (SAMPLE HOMOGENIZATION).
- Prepare DNase I master mix for the number of samples to be processed.
 DNase I master mix for purification of one sample (200 μL): 100 μL 2X DNase I Buffer, 4 μL DNase I (reconstituted), 20 μL Manganese Chloride Solution and 76 μL nuclease free water. Aliquot prepared master mix to the DNase I plate (plate number 3, below).
- 3. Obtain six empty Thermo Scientific Microtiter deep well 96 plates and two empty Thermo Scientific KingFisher Flex 96 KF plates.
- 4. Fill the plates as follows:

Plate number	Plate type	Plate name	Content	Volume per well
2	Microtiter deep well 96 plate	Wash 1_1	Wash Buffer 1 (supplemented with ethanol)	700 µL
3*	Microtiter deep well 96 plate	DNase I	DNase I master mix	200 µL
4	Microtiter deep well 96 plate	Wash 1_2	Wash Buffer 1 (supplemented with ethanol)	700 µL
5	Microtiter deep well 96 plate	Wash 2_3	Wash Buffer 2 (supplemented with ethanol)	700 µL
6	Microtiter deep well 96 plate	Wash 2_4	Wash Buffer 2 (supplemented with ethanol)	700 µL
7	KingFisher Flex 96 KF plate	Elution	Water, nuclease free	100 µL
8	KingFisher Flex 96 KF plate	Tip plate	-	-

*For best results, avoid storage of **DNase I** in **1X Reaction Buffer with MgCI**₂ at room temperature for extended periods of time. It is recommended to prepare the **DNase I** plate just prior to use.

5. Prepare the **Sample** plate (Microtiter deep well 96 plate).

Add the following reagents to the **Sample** plate.

Plate number	Plate type	Plate name	Content	Volume per well
	1 Microtiter deep well 96 plate	Sample	Plant tissue lysate	400 µL
1			Magnetic Beads*	25 µL
1			Ethanol (96-100%)	400 µL

*-Resuspend Magnetic Beads well by vortexing before use.

- 6. Place a Thermo Scientific KingFisher Flex 96 tip comb for deep well magnets on a Tip Plate (an empty KingFisher Flex 96 KF plate).
- 7. Start the **Plant_RNA_Flex** protocol with the KingFisher Flex 96 and load the plates according to the KingFisher display. After all the plates have been loaded into the instrument, the protocol will begin.
- When the KingFisher Flex pauses at the dispense step (after the DNase I digestion, approximately 30 minutes after starting the run), remove the DNase I plate from the instrument and add 150 μL of Rebinding Buffer (concentrated) and 400 μL of ethanol (96-100%) per well to rebind the RNA.

Plate number	Plate type	Plate name	Content	Volume per well
			Rebinding Buffer (conc.)	150 µL
3	Microtiter deep well 96 plate	DNase I	Ethanol (96-100%)	400 µL

- 9. Place the **DNase I** plate back into the instrument and press **Start**. After the pause, the protocol will continue through to completion.
- 10. After the protocol is complete, remove the plates according to the instructions on the KingFisher Flex display and turn off the instrument. Transfer the purified RNA from the Elution plate to a fresh RNase-free microtiter plate. Keep on ice for immediate use in downstream applications or store at -70°C.

Protocol B. Instructions for total RNA purification from up to 50 mg plant sample using KingFisher Duo with 12-pin magnet head and Microtiter deep well 96 plate

Note:

- When using the MagJET Plant RNA Kit for the first time, prepare working solutions of Wash Buffer 1 and Wash Buffer 2 and reconstitute DNase I as described on page 5.
- Ensure you are using the KingFisher Duo 12-pin magnet head and heating block
- Transfer the Plant_RNA_Duo protocol file to the KingFisher instrument before first use. The instructions for transferring the protocol can be found in Chapter 4: "Using the software" in the BindIt Software for KingFisher Instruments version 3.2 User Manual. The protocol files for MagJET Plant RNA Kit can be found on product web page on www.thermoscientific.com/onebio
- 1. Homogenize the plant samples and prepare lysates as described on page 7-8 (SAMPLE HOMOGENIZATION).
- Prepare DNase I master mix for the number of samples to be processed.
 DNase I master mix for purification of one sample (200 μL): 100 μL 2X DNase I Buffer, 4 μL DNase I (reconstituted), 20 μL Manganese Chloride Solution and 76 μL nuclease free water. Aliquot prepared master mix to the Microtiter deep well 96 plate row A (DNase I row).
- 3. Obtain one empty Thermo Scientific Microtiter deep well 96 plate and one Thermo Scientific KingFisher Duo Elution strip.
- 4. Prepare the **Plant RNA plate** (Microtiter deep well 96 plate).

Plate name and type	Row	Row name	Content	Sample/reagent volume per well
	А	DNase I	DNase I master mix	200 µL
	В	Tip	12-tip comb	N/A
			Plant tissue lysate	400 µL
Plant RNA plate Microtiter deep well	C*	Sample	Magnetic Beads	25 µL
			Ethanol (96-100%)	400 µL
96 plate	D	Wash 1_1	Wash Buffer 1	700 µL
	E	Wash 1_2	Wash Buffer 1	700 µL
	F	Wash 2_3	Wash Buffer 2	700 µL
	G	Wash 2_4	Wash Buffer 2	700 µL
	Н	Empty	Empty	Empty

Add the following reagents to the rows. Note that row B is reserved for the tips and should be left empty. Note that row H is left empty.

Note: For best results avoid storing DNase I master mix at room temperature for long-term. It is recommended to prepare the DNase I row A after filling Wash buffers rows D, E, F, G.

*It is recommended to prepare the **Sample row C** last.

5. Fill the KingFisher Duo **Elution strip** as follows.

Elution strip Content		Volume per well
KingFisher Duo Elution strip	Water, nuclease free	100 µL

- 6. Place a Thermo Scientific KingFisher Duo 12-tip comb into row B on the Plant RNA plate.
- 7. Switch on the KingFisher Duo instrument and ensure you are using the KingFisher Duo 12-pin magnet head and heating block. Start the **Plant_RNA_Duo** protocol. Insert the Plant RNA plate and **Elution strip** into the instrument as indicated on the KingFisher Duo display and press **START.** Make sure that the **Elution strip** is placed in the correct direction into the elution block; ensure that the perforated end is facing toward the user.
- When the KingFisher Duo pauses at the dispense step (after the DNase I digest, approximately 30 minutes after starting the run), remove the **Plant RNA plate** from the instrument and add 150 µL of **Rebinding Buffer** (concentrated) and 400 µL ethanol (96-100%) per well to **row A** (DNase I row) to rebind the RNA.

Plate name and type	Row	Row name	Content	Volume per well
Plant RNA plate			Rebinding Buffer (conc.)	150 µL
Microtiter deep well 96 plate	A	DNase I	Ethanol (96-100%)	400 µL

- 9. Place the **Plant RNA plate** back into the instrument and press **OK.** After the pause, the protocol will continue through to completion.
- 10. After the run is complete, remove the **Plant RNA plate** and **Elution strip** according to the instructions on the KingFisher Duo display and turn off the instrument. Transfer the purified RNA from the **Elution strip** to a fresh RNase-free microtiter plate or microcentrifuge tubes. Keep on ice for immediate use in downstream applications or store at -70°C.

Protocol C. Instructions for manual plant total RNA purification from up to 50 mg plant sample

Following this protocol is based on the transfer of liquids by pipetting through different purification steps rather than magnetic bead transfer as in the KingFisher automatic protocols. This allows the kit to be used in various throughput applications using a magnetic rack and manual or automated pipetting equipment. Protocols for automated pipetting platforms should be optimized for each platform and sample type. To enable protocol optimization, all buffers are available for purchase separately.

Note: When using the MagJET Plant RNA Kit for the first time, prepare working solutions of Wash Buffer 1 and Wash Buffer 2 and reconstitute DNase I as described on page 5.

- 1. Homogenize the plant samples and prepare lysates as described on page 7-8 (SAMPLE HOMOGENIZATION).
- 2. Carefully transfer 400 µL of lysate supernatant to a clean microcentrifuge tube.
- 3. Binding to magnetic beads: add 25 µL MagJET Magnetic beads resuspended well by vortexing and 400 µL 96-100% ethanol. Vortex the tube and briefly spin to remove droplets from the inside of the lid. Place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes (or until the solution clears). Discard the supernatant using a pipette.
- 4. Wash 1: remove the tube from the magnetic stand and add 700 μL Wash Buffer 1 (supplemented with ethanol, see page 5). Resuspend the magnetic beads by vortexing, briefly spin, then place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes. Discard the supernatant using a pipette.
- 5. **Dry:** leave the tube on the magnetic rack for 5 minutes to dry the magnetic beads at room temperature. Remove the tube from the magnetic rack.
- DNase I treatment: add 200 μL of DNase I solution for each purification (mix 100 μL 2X DNase I Buffer, 4 μL DNase I (reconstituted), 20 μL Manganese Chloride Solution and 76 μL nuclease free water. Prepare common mix for all samples to be processed). Mix and incubate in the thermomixer at 37°C and 350 rpm for 15 minutes.
- 7. Rebinding to magnetic beads: after DNase I treatment, add 150 µL of Rebinding Buffer (concentrated) and 400 µL 96-100% ethanol. Resuspend the magnetic beads by vortexing and briefly spin to remove droplets from the inside of the lid. Place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes. Discard the supernatant using a pipette.
- 8. Wash 2: remove the tube from the magnetic stand and add 700 µL Wash Buffer 1 (supplemented with ethanol, see page 5). Resuspend the magnetic beads by vortexing, place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes. Discard the supernatant using a pipette.
- 9. Wash 3: remove the tube from the magnetic stand and add 700 µL Wash Buffer 2 (supplemented with ethanol, see page 5). Resuspend the magnetic beads by vortexing, place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes. Discard the supernatant using a pipette.
- 10. Wash 4: repeat step 9 using 700 μL of Wash Buffer 2. Make sure that all the supernatant is completely removed in the last washing step.
- 11. **Dry:** leave the tube on magnetic rack for 5 min to dry the magnetic beads at room temperature. Remove the tube from the magnetic rack.
- 12. Elution: add 100 μL Nuclease free water. Resuspend the magnetic beads by vortexing, briefly spin, then place the tube on the magnetic stand and let the magnetic beads collect at the magnet for 2-3 minutes.
- 13. While on the magnetic stand, transfer the eluate (which contains the purified RNA) to a new RNase free tube, then close immediately. Keep on ice for immediate use in downstream applications or store at -70°C.

TROUBLESHOOTING

	TROUBLESHOUTING
Problem	Possible cause and solution
Low yield of purified RNA	 RNA degraded during sample storage. Make sure the plant samples were properly stored and make sure the samples were processed immediately after collection or removal from storage. Insufficient homogenization of plant material. To disrupt the cell wall, it is important to homogenize the sample thoroughly until it is ground to a fine powder. Ethanol was not added to the lysate. Ensure that ethanol was added to the lysate during bindind to magnetic beads step. Incomplete resuspension of magnetic particles. Resuspend the magnetic particles by vortexing before use. Ethanol was not added to Wash Buffer 1: add ethanol to Wash buffer 1 as described on page 5. Ethanol was not added to Wash Buffer 2: ensure that ethanol was added to Wash Buffer 2 as described on page 5. Rebinding step was missed: ensure that rebinding buffer and ethanol was added to was added after DNase I treatment. Loss of magnetic beads during purification: use sufficient magnetic beads antivative beads during purification and buffer and ethanol was added to be accepted back adding purification and buffer and ethanol was added and the provide the store adding back adding purification and buffer and ethanol was added after DNase I treatment.
Purified RNA is degraded	 capture time, be careful not remove magnetic beads during procedures. RNase contamination. To avoid RNase contamination wear gloves during the procedure and change gloves frequently. Use sterile, disposable RNase-free pipette tips. Use reagents designed to remove RNase contamination from nondisposable items (pipettes, centrifuges) and work surfaces. RNA degraded during sample storage. Make sure the plant samples were properly stored and make sure the samples were processed immediately after collection or removal from storage. RNA degraded during sample homogenization. To minimize RNA degradation, avoid repeated freezing and thawing of the samples and perform extractions from fresh material or material that has been immediately frozen and stored at -70°C. Transfer the ground tissue to the Lysis Buffer for MagJET Plant RNA Kit as quickly as possible to avoid RNA degradation. All ground material must be thoroughly mixed with the Lysis Buffer for MagJET Plant RNA. RNA degradation can occur in particles that are left to dry on the walls of the tube. Lysis Buffer was not supplemented with DTT. Ensure that 2 M DTT solution has been added to the Lysis Buffer before RNA purification procedure. Purified RNA was not stored properly. Purified RNA should be used immediately in downstream applications or stored at -70°C is recommended.
Inhibition of downstream enzymatic reactions	 Purified RNA contains residual salt. Use the correct order of Wash Buffers steps. Always wash the magnetic beads with Wash Buffer 1 first and then proceed with Wash Buffer 2. Purified RNA contains residual ethanol. It is important to dry magnetic beads before elution step.
Genomic DNA contamination	Inefficient DNase I digestion. Ensure that the temperature of DNase I digestion is 37°C. Prepare DNase I mix according to protocol (see p. 7 and p. 10). Too much sample input: if starting amount of plant sample is too high (> 50 mg), the DNA digestion step may not be effective.
Carryover of the magnetic beads in the elution	Carryover of magnetic beads in the eluted RNA will not affect downstream applications. To remove the carryover from eluted RNA, simply magnetize the magnetic beads and carefully transfer eluate to a new tube or plate.



Lysis Buffer for MagJET Plant RNA Kit

Xn Harmful

Hazard-determining components of labelling: guanidinium thiocyanate

Risk phrases:

R20/21/22: Harmful by inhalation, in contact with skin and if swallowed.

R32-52/53: Contact with acids liberates very toxic gas. Harmful to aquatic organisms, may cause long-term adverse effects in the aquatic environment.

Precautionary statements

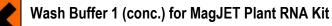
P261 Avoid breathing dust/fume/gas/mist/vapours/spray.

P280 Wear protective gloves/protective clothing/eye protection/face protection.

P305+P351+P338 IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.

P301+P312 IF SWALLOWED: Call a POISON CENTER or doctor/physician if you feel unwell.

P304+P340 IF INHALED: Remove victim to fresh air and keep at rest in a position comfortable for breathing. P501 Dispose of contents/container in accordance with local/regional/national/international regulations.



Xn Harmful

Hazard-determining components of labelling: guanidinium chloride

Risk phrases:

22 Harmful if swallowed.

36/38 Irritating to eyes and skin.

Safety phrases:

23 Do not breathe gas/fumes/vapour/spray.

26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

36/37 Wear suitable protective clothing and gloves.

60 This material and its container must be disposed of as hazardous waste



Rebinding Buffer for MagJET Plant RNA Kit

Xn Harmful

Hazard-determining components of labelling: guanidinium chloride

Risk phrases:

22 Harmful if swallowed.

36/38 Irritating to eyes and skin.

Safety phrases:

23 Do not breathe gas/fumes/vapour/spray.

26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice.

36/37 Wear suitable protective clothing and gloves.

60 This material and its container must be disposed of as hazardous waste

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively for research purposes and *in vitro* use only. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to <u>www.thermoscientific.com/onebio</u> for Material Safety Data Sheet of the product. © 2013 Thermo Fisher Scientific Inc. All rights reserved. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.