

A Geno Technology, Inc. (USA) brand name

GET™ Viral RNA-Mag

Genomic Efficient Technology for Viral RNA purification with silica magnetic beads (Cat. # 786-1735, 786-1736)



INTRODUCTION	3
ITEMS SUPPLIED	4
STORAGE CONDITIONS	5
ADDITIONAL ITEMS NEEDED	5
IMPORTANT INFORMATION	5
PREPARATION BEFORE USE	5
PROTOCOL	6
SERUM, PLASMA OR CELL CULTURE SUPERNATANT VIRAL SAMPLES	6
SWAB (VIRAL SAMPLE) FROM NOSE, THROAT OR OTHER TEST AREA	7
TROUBLESHOOTING	8
RELATED PRODUCTS	9

INTRODUCTION

Get™ Viral RNA-Mag Kit belongs to our series of kits based on Genomic Efficient Technology (GET) for purification of nucleic acids from diverse sample.

GET™ is based on a highly efficient Genomic lysis buffer that liberates nucleic acid from cellular protein complexes, making nucleic acids free and available for purification in pure form. Free nucleic acids are immobilized, in the presence of high concentration of chaotropic agents, on silica surface. Following the capture of nucleic acid on the silica membrane or beads, a series of washing steps removes interfering impurities. In the final step, pure nucleic acid is eluted in concentrated form with GET GET Elution Buffer.

The Get™ Viral RNA- Mag is based upon the principle of binding of RNA on silica beads with magnetic core in presence of chaotropic salts. RNA is bound to the silica magnetic beads under high concentrations of chaotropic salts and the impurities are removed during the wash steps and nucleic acid eluted with GET GET Elution Buffer (Fig:1). The RNA extraction is rapid with less than 15-30 minutes hands- on time.

Furthermore, the kit is optimized to obtain high quality viral RNA with minimum to no loss of RNA. The purified RNA fraction obtained is concentrated and is suitable for RT-PCR and qRT-PCR studies. These RT-PCR studies are used for virus detection, virus load, and virus genotyping.

Get™ Viral RNA-Mag is available in 50 and 100 prep sizes with the maximum biological sample volume of 200 µl per prep.

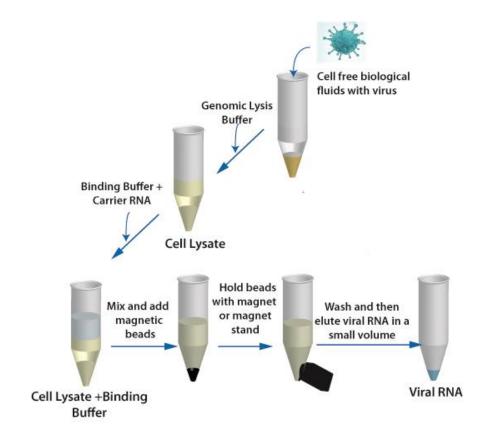


Fig:1

ITEMS SUPPLIED

Description	Cat. # 786-1735 50 preps	Cat. # 786-1736 100 preps
Genomic Lysis Buffer	30 ml	2 x 30 ml
GET Binding Buffer	50	2 x 50 ml
Poly A [Carrier RNA]	1 vial	2 vials
Longlife™ Proteinase K	0.5 ml	2 x 0.5 ml
GET Wash I	30 ml	2 x 30 ml
GET Wash II	20 ml	2x 20 ml
GET Elution Buffer	10 ml	10 ml
Silica Magnetic Beads	2.5 ml	2 x 2.5 ml

STORAGE CONDITIONS

The kit is supplied at ambient temperature. Upon receipt store the kit components as indicated on labels. If stored and used as directed, the kit is stable for 1 year.

ADDITIONAL ITEMS NEEDED

- 1. Magnetic stand (Cat. # 786-888) or a magnet
- Cell-free virus containing sample such as serum, plasma, whole blood, cell culture supernatant, nasal or throat swabs
- 3. Absolute ethanol.
- 4. 1.5 ml nuclease-free (sterile) microfuge tubes
- (Optional) RNAseOUT™ (Cat. #786-70) for cleaning working bench.
- 6. Nuclease free tips for the pipettes.

IMPORTANT INFORMATION

- Get™ Viral RNA is available in 50 prep (Cat. # 786-1735) and 100 prep (Cat. # 786-1736) sizes with sample volume of 200 µl. For more than 200µl sample increase the volume of magnetic beads and buffers added proportionally as per protocol.
- Treat all the viral samples as potentially infectious and use safe lab handling procedures.
- Since the sensitivity and titer of potential pathogen (virus) varies with samples, the end user needs take appropriate safety measures when handling.
- Sterile handling of reagents and sample should be carried out to avoid contamination of reagents and sample with bacteria or nuclease. Frequent change of gloves and cleaning of work area with RNAseOUT™ (Cat. #786-70) is recommended. Avoid touching the mouth of reagent bottles.
- For performing RT-PCR or qRT-PCR experiments using viral RNA, ensure that sample preparation, RT-PCR run, and electrophoresis are carried out in separate work areas to avoid cross-contamination.

PREPARATION BEFORE USE

- Add 18 ml of molecular grade ethanol to 30 ml GET Wash I bottle and check the box on the bottle label to indicate ethanol has been added.
- Add 80ml molecular grade ethanol to the GET Wash II bottle (20 ml) and check the box on the bottle label to indicate ethanol has been added.
- 3. Equilibrate GET Elution Buffer to 60°C.
- 4. Briefly centrifuge the Poly A [Carrier RNA] tube. Add 250 μl of GET Elution Buffer to 1 vial of Poly A [Carrier RNA]. Gently mix with pipette to dissolve. Aliquot 50μl per microfuge vial and store the vials at -20°C as stock solution.
- 5. Store Longlife™ Proteinase K in small aliquots at -20°C for long term use. Before use, Invert the Longlife™ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use

6. Working solution of GET Binding Buffer (*Prepare fresh and do not store after use*): Thaw one vial of Poly A [Carrier RNA] (50 μ l). Transfer the Poly A [Carrier RNA] solution to 10 ml of GET Binding Buffer. Briefly vortex to mix.

PROTOCOL

Serum, plasma or cell culture supernatant viral samples

1. Add 200 μ l of Genomic Lysis Buffer to 200 μ l of viral sample. Vortex to mix. **NOTE**: For sample volume <200 μ l add PBS or 0.9%NaCl to make up the volume to

NOTE: For sample volume <200 μl add PBS or 0.9%NaCl to make up the volume to 200μl.

NOTE: For sample volume >200 μ l increase the volume of reagents added proportionally.

 Add 10 µl Longlife™ Proteinase K suspension into the sample, mix and incubate at 55°C-60°C for 1 hr.

NOTE: Before use, Invert the Longlife[™] Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.

NOTE: Do not exceed 60°C.

- 3. Add 800 μ l of working solution of GET Binding Buffer to the sample and vortex to mix
- 4. Add 50 μ l of Silica Magnetic Beads to sample tube and gently invert 4-5 times to mix. Incubate the sample for 5 minutes with gentle mixing on rotor at room temperature.

NOTE: Do not vortex or shake the sample containing silica magnetic beads vigorously.

- 5. Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end of the tube.
- 6. Gently remove and discard the supernatant without disturbing the beads.
- 7. Remove the tube from magnetic stand or magnet and add 0.5ml GET Wash I to the sample tube and invert the tube gently 4-5 times to mix.
- Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end.
- 9. Gently remove and discard the supernatant without disturbing the beads.
- Remove the tube from magnetic stand or magnet and add 0.5ml GET Wash II to the tube and invert the tube gently 4-5 times to mix.
- 11. Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end.
- 12. Gently remove and discard the supernatant without disturbing the beads.
- 13. Repeat step 10-12.
- 14. Ensure all the liquid is removed from the magnetic beads. Air dry the beads for 5-20 minutes.
- 15. Add 50 μ l of prewarmed GET Elution Buffer to the magnetic beads and resuspend the beads complex by brief vortex or shaking. Incubate for 15 minutes with gentle mixing on rotor at room temperature.

16. Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end of the tube. Transfer the supernatant in a 1.5 ml nuclease free tube.

NOTE: Check the recovery. If recovery is poor, add 25-50 μ l prewarmed (50-60°C) GET Elution Buffer to the magnetic beads and repeat steps 15-16. Combine with previous elution.

17. Store the eluted RNA at -80°C.

Swab (Viral sample) from nose, throat or other test area

- Take swab (nose, throat or other test area) in a swab tube containing PBS or 0.9
 NaCl. Vortex the tube for 1-2 minutes to release the pathogen (virus) in the solution.
- 2. Remove the swab and centrifuge the tube at 15,000 g for 10 minutes at room temperature.

NOTE: Alternatively, filtration step can be used to remove cells from swabs.

- 3. Transfer the supernatant into a nuclease free microfuge tube.
- 4. Transfer 200 μ l of supernatant in a microfuge tube and add 200 μ l Genomic Lysis Buffer. Briefly vortex to mix.
- 5. Add 10 μl Longlife™ Proteinase K suspension into the sample, mix and incubate at 55°C-60°C for 1 hr.

NOTE: Before use, Invert the Longlife $^{\text{m}}$ Proteinase K tube 3-4 times to mix the enzyme suspension, then remove an aliquot for use.

NOTE: Do not exceed 60°C.

- 6. Add 800 μ l of working solution of GET Binding Buffer to the sample and vortex to mix.
- 7. Add 50 μ l of Silica Magnetic Beads to sample tube and gently invert 4-5 times to mix. Incubate the sample for 5 minutes with gentle mixing on rotor at room temperature.

NOTE: Do not vortex or shake the sample containing silica magnetic beads vigorously.

- Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end of the tube.
- 9. Gently remove and discard the supernatant without disturbing the beads.
- 10. Remove the tube from magnetic stand or magnet and add 0.5ml GET Wash I to the sample tube and invert the tube gently 4-5 times to mix.
- Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end.
- 12. Gently remove and discard the supernatant without disturbing the beads.
- 13. Remove the tube from magnetic stand or magnet and add 0.5ml GET Wash II to the tube and invert the tube gently 4-5 times to mix.
- Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end.
- 15. Gently remove and discard the supernatant without disturbing the beads.
- 16. Repeat step 13-15 once.

- 17. Ensure all the liquid is removed from the magnetic beads. Air dry the beads for 5-20 minutes.
- 18. Add 50 μ l of prewarmed GET Elution Buffer to the magnetic beads and resuspend the beads complex by brief vortex or shaking. Incubate for 15 minutes with gentle mixing on rotor at room temperature.
- 19. Place the sample tube on magnetic stand or use magnet to immobilize the beads at one end of the tube. Transfer the supernatant in a 1.5 ml nuclease free tube.
 NOTE: Check the recovery. If recovery is poor, add 25-50µl prewarmed (50-60°C)
 GET Elution Buffer to the magnetic beads and repeat steps 15-16. Combine with previous elution.
- 20. Store the eluted RNA at -80°C.

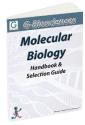
TROUBLESHOOTING

Issue	Suggested reason	Possible solution
Low RNA yield or purity	Kit components are not stored properly	Store kit components as indicated in the label.
	Ethanol not added to the GET Wash I and II Buffers	Add absolute ethanol to GET Wash I and II Buffers before using
	Reagent and samples not properly mixed	Mix the sample tube well after addition of each reagent.
Low RNA yield	High levels of RNAse activity	Create RNase-free work environment. Use RNase OUT to clean the working bench Process starting material immediately or store at - 80°C until it is processed Use eluted RNA directly for
		downstream application or store at -80°C for later use
	Incomplete Proteinase K digestion	Thaw Longlife™ Proteinase K on ice and Resuspend Proteinase K solution by inverting tube 3-4 times for uniform suspension before adding to the sample. Incubate for longer time if necessary.
Poor elution of RNAs	GET Elution Buffer provided in kit is not used	Use the GET Elution Buffer provided in the kit
		p. oaca iii ciic kie

If using own water or GET
Elution Buffer, ensure the pH
of buffer is same as that of
the GET Elution Buffer
provided

RELATED PRODUCTS

Download our Molecular Biology Handbook.



http://info2.gbiosciences.com/complete-molecular-biology-handbook

For other related products, visit our website at www.GBiosciences.com or contact us.

Version1: 5/28/20



www.GBiosciences.com