

NEBNext® Poly(A) mRNA Magnetic Isolation Module

NEB #E7490S/L

24/96 reactions

Version 10.0_7/22

Table of Contents

Isolate mRNA using the NEBNext Poly(A) mRNA Magnetic Isolation Module (NEB #E7490).....1

Checklist4

Kit Components.....5

Revision History.....6

The NEBNext Poly(A) mRNA Magnetic Isolation Module Includes

The volumes provided are sufficient for preparation of up to 24 reactions (NEB #E7490S) and 96 reactions (NEB #E7490L). All reagents should be stored at 4°C.

- NEBNext Oligo d(T)₂₅ Beads
- NEBNext RNA Binding Buffer (2X)
- NEBNext Wash Buffer
- 10 mM Tris HCl (pH 7.5)
- NEBNext Tris Buffer
- Nuclease-free Water

Required Materials Not Included:

- 96-well 0.2 ml PCR Plates and Microseal® 'B' Adhesive Sealer (Bio-Rad® MSB-1001) or 0.2 ml RNase-free tube, for example Tempassure PCR flex-free 8-tube strips (USA Scientific® #1402-4708)
- Magnetic Rack (NEB#S1515S, Alpaqua®, cat. #A001322 or equivalent)
- 1.5 ml Microcentrifuge tube and NEB #S1506 Magnet stand or equivalent (for washing beads only)
- Thermal cycler or heat block
- Bioanalyzer® (Agilent Technologies, Inc.) or similar instrument and consumables

Description

The NEBNext Poly(A) mRNA Magnetic Isolation Module is designed to isolate intact poly(A)⁺ RNA from previously isolated total RNA. The technology is based on the coupling of Oligo d(T)₂₅ to 1 µm paramagnetic beads which is then used as the solid support for the direct binding of poly(A)⁺ RNA. Thus, the procedure permits the manual processing of multiple samples and can be adapted for automated high-throughput applications. Additionally, magnetic separation technology permits elution of intact mRNA in small volumes eliminating the need for precipitating the poly(A)⁺ transcripts in the eluent. Intact poly(A)⁺ RNA which is fully representative of the mRNA population of the original sample can be obtained in less than one hour.

Application

Isolation of poly(A)⁺ RNA transcript from Total RNA for RNA library preparation and sequencing.

Isolate mRNA Using the NEBNext Poly(A) mRNA Magnetic Isolation Module (NEB #E7490)

Note: If this module is being used with either NEBNext Ultra™ Directional RNA Library Prep Kit for Illumina (NEB #E7420), NEBNext Ultra RNA Library Prep Kit for Illumina (NEB #E7530), NEBNext Ultra II Directional RNA Library Prep Kit for Illumina (NEB #E7760), NEBNext Ultra II Directional RNA Library Prep with Sample Purification Beads (NEB #E7765) or NEBNext Ultra II RNA Library Prep Kit for Illumina (NEB #E7770) or NEBNext Ultra II RNA Library Prep with Sample Purification Beads (NEB #E7775) **do not follow the protocol below**. The correct protocol can be found in NEB #E7420, #E7530, #E7760, #E7765, #E7770 or #E7775 manuals. For use with

NEBNext RNA First Strand Synthesis Module (NEB #E7525 or NEBNext Ultra II RNA First Strand Synthesis Module (NEB #E7771) please follow protocols in the manuals for NEB #E7760 or NEB #E7770.

Note: For best results keep all the reagents used during the Poly(A) isolation except the NEBNext Oligo d(T)₂₅ beads, **on ice** when not in use.

The protocol has been optimized using high quality Universal Human Reference Total RNA.

RNA Sample Requirements RNA Integrity:

Assess the quality of the Input RNA by running the RNA sample on an Agilent Bioanalyzer RNA 6000 Nano/Pico Chip. For PolyA mRNA enrichment, high quality RNA with a RIN score > 7 is required.

RNA Sample Requirements:

The RNA sample should be free of salts (e.g., Mg²⁺, or guanidinium salts, divalent cation chelating agents (e.g., EDTA or EGTA) or organics (e.g., phenol or ethanol). RNA must be free of DNA. gDNA is a common contaminant from RNA preps. It may be carried over from the interphase of organic extractions or when the silica matrix of solid phase RNA purification methods is overloaded. If the total RNA sample may contain gDNA contamination, treat the sample with DNase I to remove all traces of DNA (DNase is not provided in this kit). After treatment with DNase I the enzyme should be removed from the sample. Any residual activity of the DNase I may degrade the oligos necessary for the enrichment. DNase I can be removed from the extraction using phenol/chloroform extraction and ethanol precipitation.

Starting Material: 1 -5 µg* of DNA-free total RNA.

* Inputs as low as 10 ng can be used in combination with NEBNext Ultra II RNA Library prep kits (E7770 and E7760). Please refer to the library prep manual for the appropriate instructions to follow.

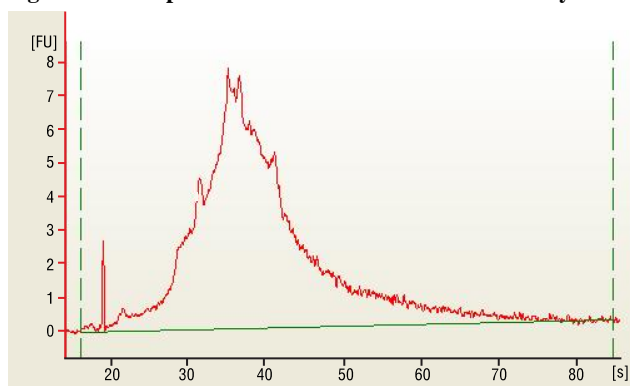
1. Dilute the total RNA with nuclease-free water to a final volume of 50 µl in a nuclease-free 0.2 ml PCR tube and keep on ice.
2. To wash the Oligo dT Beads, add the following to a 1.5 ml nuclease-free tube. If preparing multiple libraries, beads for up to 10 samples can be added to a single 1.5 ml tube for subsequent washes (use magnet NEB #S1506 for 1.5 ml tubes). The purpose of this step is to bring the beads from the storage buffer into the binding buffer. The 2X Binding Buffer does not have to be diluted for this step.

COMPONENT	VOLUME PER ONE LIBRARY
Oligo dT Beads d(T) ₂₅	20 µl
RNA Binding Buffer (2X)	100 µl
Total Volume	120 µl

3. Wash the beads by pipetting up and down six times.
4. Place the tube on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
5. Remove and discard all of the supernatant from the tube. Take care not to disturb the beads.
6. Remove the tube from the magnetic rack.
7. Add 100 µl RNA Binding Buffer to the beads and wash by pipetting up and down six times. If preparing multiple libraries, add 100 µl RNA Binding Buffer per sample. The Binding Buffer does not have to be diluted.
8. Place the tubes on the magnet and incubate at room temperature until the solution is clear (~2 minutes).
9. Remove and discard the supernatant from the tube. Take care not to disturb the beads.
10. Remove tubes from magnet and add 50 µl RNA Binding Buffer to the beads and mix by pipetting up and down until beads are homogenous. If preparing multiple libraries, add 50 µl RNA Binding Buffer per sample.
11. Add 50 µl beads to each RNA sample from Step 1. Mix thoroughly by pipetting up and down six times. This first binding step removes most of the non target RNA.
12. Place the tubes on the thermal cycler and heat the sample at 65°C for 5 minutes and **cool to 4°C with the heated lid set at ≥ 75°C** to denature the RNA and facilitate binding of the mRNA to the beads.
13. Remove tubes from the thermal cycler when the temperature reaches 4°C.
14. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.

15. Place the tubes on the bench and incubate at room temperature for 5 minutes to allow the RNA to bind to the beads.
16. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.
17. Incubate for 5 more minutes on the bench at room temperature to allow the RNA to bind to the beads.
18. Place the tubes on the magnetic rack at room temperature for 2 minutes or until the solution is clear to separate the poly-A RNA bound to the beads from the solution.
19. Remove and discard all of the supernatant. Take care not to disturb the beads.
20. Remove the tubes from the magnetic rack.
21. Wash the beads by adding 200 μ l of Wash Buffer to remove unbound RNA. Pipette the entire volume up and down 6 times to mix thoroughly.
22. Place the tubes on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
23. Remove and discard all the supernatant from each well of the tube. Take care not to disturb the beads.
24. Remove the tubes from the magnetic rack.
25. Repeat Steps 21-24.
26. Add 50 μ l of Tris Buffer (provided in NEB #E7490 kit) to each tube. Gently pipette the entire volume up and down 6 times to mix thoroughly.
27. Place the tubes on the thermal cycler. Close the lid and heat the sample at 80°C for 2 minutes, then cool to 25°C with the heated lid set at $\geq 90^\circ\text{C}$ to do the first elution of the mRNA from the beads.
28. Remove the tubes from the thermal cycler when the temperature reaches 25°C.
29. Add 50 μ l of RNA Binding Buffer to each sample to allow the RNA to re-bind to the same beads. Gently pipette the entire volume up and down 6 times to mix thoroughly.
30. Incubate the tubes on the bench at room temperature for 5 minutes.
31. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.
32. Incubate the tubes on the bench at room temperature for 5 more minutes to allow the RNA to bind to the beads.
33. Place the tubes on the magnetic stand at room temperature for 2 minutes or until the solution is clear.
34. Remove and discard all of the supernatant from each tube. Take care not to disturb the beads.
35. Remove the tubes from the magnetic rack.
36. Wash the beads once with 200 μ l of Wash Buffer. Gently pipette the entire volume up and down 6 times to mix thoroughly. Spin down the tube briefly to collect the liquid from the wall and lid of the tube.
37. Place the tubes on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
38. Remove and discard all of the supernatant from each tube. Take care not to disturb the beads.
Note: It is important to remove all of the supernatant to successfully use the RNA in downstream steps. Spin down the tube. Place the tube on the magnetic rack and with a 10 μ l tip remove all of the wash buffer. (Caution: Do not disturb beads that contain the mRNA). Avoid letting the beads dry out before adding elution buffer.
39. Remove the tubes from the magnetic rack.
40. Elute the mRNA from the beads by adding 17 μ l of the Tris Buffer, mix by pipetting 6 times and incubating the sample at 80°C for 2 minutes, then cool to 25°C to elute the polyA RNA from the beads. Immediately, place the tubes on the magnetic rack for 2 minutes or until the solution is clear.
41. Collect the purified mRNA by transferring 15 μ l of the supernatant to a clean nuclease-free PCR Tube. Place on ice if using immediately. Alternatively, the sample can be placed at -80°C for long term storage.
42. Assess the Yield and the Size Distribution of the purified mRNA. Run 1 μ l on the Bioanalyzer using a RNA Pico Chip. You may have to dilute your sample before loading.

Figure 1: Example of mRNA distribution on a bioanalyzer.



Checklist:

- 1. Dilute RNA to 50 μ l in 0.2 ml tube
- 2. New 0.2 ml tube: aliquot 20 μ l NEBNext Magnetic Oligo d(T)₂₅ Beads
- 3. Add 100 μ l of RNA Binding Buffer and mix 6 times
- 4. Place tube on magnet for 2 minutes
- 5. Remove and discard supernatant
- 6. Remove tube from magnet
- 7. Add 100 μ l of RNA Binding Buffer and mix 6 times
- 8. Place tube on magnet for 2 minutes
- 9. Remove and discard supernatant
- 10. Remove tube from magnet, add 50 μ l RNA Binding Buffer, mix
- 11. Add 50 μ l resuspended beads to 50 μ l total RNA sample mix
- 12. Heat the sample at 65°C for 5 minutes, then hold at 4°C
- 13. Remove tubes when temperature is 4°C
- 14. Resuspend beads by mixing slowly 6 times
- 15. Incubate tubes on bench for 5 minutes
- 16. Resuspend beads by mixing slowly 6 times
- 17. Incubate tubes 5 more minutes on bench
- 18. Place tubes on magnet for 2 minutes
- 19. Remove and discard supernatant
- 20. Remove tube from magnet
- 21. Wash beads with 200 μ l Wash Buffer by mixing 6 times
- 22. Place tubes on magnet for 2 minutes
- 23. Remove and discard supernatant
- 24. Remove tubes from magnet
- 25. Repeat Steps 21 , 22 , 23 and 24
- 26. Add 50 μ l Tris to tubes and mix 6 times
- 27. Place tubes on thermal cycler, heat at 80°C for 2 minutes, then cool to 25°C
- 28. Remove tubes from thermal cycler when temperature is 25°C
- 29. Add 50 μ l RNA Binding Buffer to same beads in tubes and mix 6 times
- 30. Incubate tubes at room temperature for 5 minutes

- [_] 31. Resuspend beads by mixing slowly 6 times
- [_] 32. Incubate tubes on bench for 5 more minutes
- [_] 33. Place tubes on magnet for 2 minutes
- [_] 34. Remove and discard supernatant
- [_] 35. Remove tubes from magnet
- [_] 36. Wash beads with 200 µl of Wash Buffer by mixing 6 times
- [_] 37. Place tubes on magnet for 2 minutes
- [_] 38. Remove and discard supernatant
- [_] 39. Remove tubes from magnet
- [_] 40. Add 17 µl Tris Buffer to elute mRNA, mix incubate at 80°C for 2 minutes, cool to 25°C, immediately put on magnet for 2 minutes
- [_] 41. Transfer 15 µl of the supernatant (containing the purified mRNA) into a clean nuclease-free PCR Tube. Place on ice if using immediately. Alternatively, the sample can be placed at –80°C for long term storage.
- [_] 42. Run mRNA on a Bioanalyzer RNA Pico Chip.

Kit Components

NEB #E7490S Table of Components

NEB #	PRODUCT	VOLUME
E7499A	NEBNext Oligo d(T) ₂₃ Beads	0.480 ml
E7492A	NEBNext RNA Binding Buffer (2X)	7.2 ml
E7493A	NEBNext Wash Buffer	28.8 ml
E7496A	NEBNext Tris Buffer	6.0 ml
E7495A	Nuclease-free Water	1.2 ml

NEB #E7490L Table of Components

NEB #	PRODUCT	VOLUME
E7499AA	NEBNext Oligo d(T) ₂₃ Beads	1.92 ml
E7492AA	NEBNext RNA Binding Buffer (2X)	28.8 ml
E7493AA	NEBNext Wash Buffer	57.6 ml
E7496AA	NEBNext Tris Buffer	24.0 ml
E7495AA	Nuclease-free Water	4.8 ml

Revision History

REVISION #	DESCRIPTION	DATE
1.2	Added protocol warning note to page 3. Renamed NEBNext Elution Buffer to NEBNext Tris Buffer.	
2.0	Volume of beads increased from 15 µl to 20 µl. Additional mixing and incubation steps were added after each thermocycler incubation.	7/15
3.0	Update the protocol "Isolate mRNA using the NEBNext Oligo d(T)25 Magnetic Beads."	3/16
4.0	Adding clarification to many steps, broke out steps into more individual steps. Added to keep reagents on ice. Adjusted the title of the protocol.	5/16
5.0	Component volume change E7493AA.	11/16
5.1	Added note to protocol.	7/17
6.0	Create "Kit Component – Table of Components" for small and large size kits. Delete individual component information pages.	4/18
7.0	New manual format applied.	1/20
8.0	Update protocol Step 38.	9/20
9.0	Update protocol. Add RNA sample requirements. Add starting materials note.	1/21
10.0	Update protocol and Required materials not included	7/22

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

This product is covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; rather, it indicates where the document was originally developed. For more information about commercial rights, please email us at busdev@neb.com. While NEB develops and validates its products for various applications, the use of this product may require you to obtain additional third party intellectual property rights for certain applications.

ALPAQUA® is a registered trademark of Alpaqua Engineering, LLC.
 B CORPORATION® is a registered trademark of B Lab IP, LLC, Inc.
 BIOANALYZER® is a registered trademark of Agilent Technologies, Inc.
 MICROSEAL® and BIO-RAD® are registered trademarks of Bio-Rad Laboratories, Inc.
 ILLUMINA® and MISEQ® are registered trademarks of Illumina, Inc.
 USA SCIENTIFIC® is a registered trademark of USA Scientific, Inc.

© Copyright 2022, New England Biolabs, Inc.; all rights reserved.



www.neb.com



be INSPIRED
 drive DISCOVERY
 stay GENUINE

New England Biolabs, Inc., 240 County Road, Ipswich, MA 01938-2723 Telephone: (978) 927-5054 Toll Free: (USA Orders) 1-800-632-5227 (USA Tech) 1-800-632-7799 Fax: (978) 921-1350 e-mail: info@neb.com