

NEBNext® Poly(A) mRNA Magnetic Isolation Module

NEB #E7490S/L

24/96 reactions

Version 11.0_2/24

Table of Contents

Section 1: Isolate mRNA using the NEBNext Poly(A) mRNA Magnetic Isolation Module (Express Protocol).....	2
Section 2: Isolate mRNA using the NEBNext Poly(A) mRNA Magnetic Isolation Module (Standard Protocol)	4
Appendix A.....	7
Kit Components	8
Revision History	9

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

NEBNext Tris Buffer (10 mM Tris HCl, pH 7.5)

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. The new protocol for poly(A) mRNA enrichment (**Section 1, Express Protocol**) enables comparable performance to our current protocol (**Section 2, Standard Protocol**) using a substantially faster and more streamlined workflow.

Please refer to the product page on NEB.com for FAQs about this product.

Where larger volumes, customized or bulk packaging are required, we encourage consultation with the Customized Solutions team at NEB. Please complete the NEB Custom Contact Form at www.neb.com/CustomContactForm to learn more.

Application

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

Section 1

Isolate mRNA Using the NEBNext Poly(A) mRNA Magnetic Isolation Module (Express Protocol)

Note: When using NEBNext Poly(A) mRNA Magnetic Isolation Module E7490 with NEBNext library preparation kits, follow the protocol in the respective library prep kit manual or see Appendix A. Modifications to elution buffer and volumes will need to be made at Step 1.26 to make the enriched mRNA directly compatible with the different library preparation workflows.

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 Integrity

Assess the quality of the Input RNA by running the RNA sample on an Agilent Bioanalyzer RNA 6000 Nano/Pico Chip. For Poly(A) 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 library preparation workflows. Please refer to the respective library prep manual for the appropriate instructions to follow. Modifications to elution buffer and volumes will need to be made here to make the enriched mRNA directly compatible with the different NEBNext library preparation workflows. Please refer to Appendix A and consult the appropriate manual for more information.

- 1.1. 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.
- 1.2. Add 20 µl NEBNext Oligo d(T)₂₅ beads per reaction to a 1.5 ml tube. If preparing multiple libraries, beads for up to 24 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 NEBNext RNA Binding Buffer (2X) does not have to be diluted for this step.
- 1.3. Place the tube on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
- 1.4. Remove and discard all of the supernatant from the tube. Take care not to disturb the beads.
- 1.5. Remove the tube from the magnetic rack.
- 1.6. Add 50 µl NEBNext RNA Binding Buffer (2X) to the beads and wash by pipetting up and down 6 times. If preparing multiple libraries, add 50µl RNA Binding Buffer per sample. The Binding Buffer does not have to be diluted.
- 1.7. Place the tube on the magnet and incubate at room temperature until the solution is clear (~2 minutes).
- 1.8. Remove and discard the supernatant from the tube. Take care not to disturb the beads.
- 1.9. Remove the tube from the magnet and add 50 µl NEBNext RNA Binding Buffer (2X) 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. The Binding Buffer does not have to be diluted.
- 1.10. Add 50 µl beads to each RNA sample from Step 1.1. Mix thoroughly by pipetting up and down 6 times. This first binding step removes most of the non-target RNA.
- 1.11. Heat the sample to denature the RNA and to facilitate binding of the poly(A) RNA to the beads. Place in a thermal cycler, with the heated lid set to ≥ 90°C, and run the following program:
2 minutes at 80°C
5 minutes at 25°C
Hold at 25°C
- 1.12. Remove tubes from the thermal cycler when the temperature reaches hold at 25°C.
- 1.13. 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.

- 1.14. Remove and discard all of the supernatant. Take care not to disturb the beads. Do not remove the tubes from the magnetic rack.
- 1.15. While still on the magnet rinse the beads by gently adding 200 μ l of NEBNext Wash Buffer to the tubes to remove unbound RNA.
- 1.16. Remove and discard all of the supernatant from each tube. Take care not to disturb the beads.
- 1.17. Remove the tubes from the magnetic rack.
- 1.18. Add the following to each tube containing mRNA bound beads to allow the poly(A) RNA to re-bind to allow the RNA to re-bind to the same beads. Mix thoroughly by gently pipetting up and down 6 times.

COMPONENT	VOLUME PER ONE LIBRARY
NEBNext Tris Buffer	50 μ l
NEBNext RNA Binding Buffer (2X)	50 μ l
Total Volume	100 μl

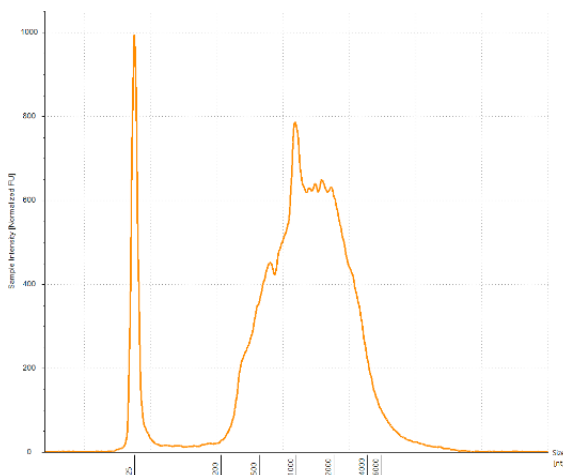
For multiple reactions, a master mix of the reaction components can be prepared on ice before addition to the sample.

- 1.19. Place the tubes in a thermal cycler, with the heated lid set to $\geq 90^{\circ}\text{C}$, and run the following program:
2 minutes at 80°C
5 minutes at 25°C
Hold at 25°C
- 1.20. Remove the tubes from the thermal cycler when the temperature reaches hold at 25°C .
- 1.21. Place the tubes on the magnetic stand at room temperature for 2 minutes or until the solution is clear.
- 1.22. Remove and discard all of the supernatant from each tube. Take care not to disturb the beads. Do not remove the tubes from the magnetic rack.
- 1.23. While still on the magnet, rinse the beads by gently adding 200 μ l of NEBNext Wash Buffer to the tubes to remove unbound RNA.
- 1.24. Remove and discard all of the supernatant from the tubes. Take care to remove all of the NEBNext Wash Buffer and do not disturb the beads that contain the mRNA.

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 NEBNext Wash Buffer. (Caution: Do not disturb beads that contain the mRNA). Avoid letting the beads dry out before adding elution buffer.

- 1.25. Remove the tubes from the magnetic rack.
Note: Modifications to elution buffer and volumes will need to be made here to make the enriched mRNA directly compatible with the different NEBNext library preparation workflows. Please refer to Appendix A and consult the appropriate manual for more information.
- 1.26. Elute the mRNA from the beads by adding 17 μ l of the NEBNext Tris Buffer, mix by pipetting 6 times and incubate the samples at 80°C for 2 minutes, then cool to 25°C . Immediately, place the tubes on the magnetic rack for 2 minutes or until the solution is clear.
- 1.27. 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.
- 1.28. Assess the yield and the size distribution of the purified mRNA. Run 1 μ l on the Bioanalyzer using a RNA Pico Chip or the TapeStation using a High Sensitivity RNA ScreenTape. You may have to dilute your sample before loading.

Figure 1: Example of mRNA distribution on a TapeStation



Section 2

Isolate mRNA Using the NEBNext Poly(A) mRNA Magnetic Isolation Module (Standard Protocol)

Note: When using NEBNext Poly(A) mRNA Magnetic Isolation Module E7490 with NEBNext library preparation kits, follow the protocol in the respective library prep kit manual or see Appendix A. Modifications to elution buffer and volumes will need to be made at Step 2.40 to make the enriched mRNA directly compatible with the different library preparation workflows.

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 Integrity

Assess the quality of the Input RNA by running the RNA sample on an Agilent Bioanalyzer RNA 6000 Nano/Pico Chip. For Poly(A) 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 library preparation workflows. Please refer to the respective library prep manual for the appropriate instructions to follow. Modifications to elution buffer and volumes will need to be made here to make the enriched mRNA directly compatible with the different NEBNext library preparation workflows. Please refer to Appendix A and consult the appropriate manual for more information.

- 2.1. 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.
- 2.2. To wash the NEBNext Oligo d(T)₂₅ 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 NEBNext RNA Binding Buffer (2X) does not have to be diluted for this step.

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

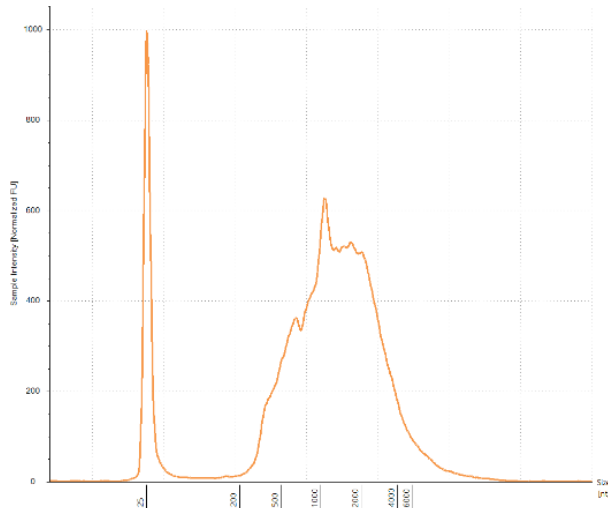
- 2.3. Wash the beads by pipetting up and down 6 times.
- 2.4. Place the tube on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
- 2.5. Remove and discard all of the supernatant from the tube. Take care not to disturb the beads.
- 2.6. Remove the tube from the magnetic rack.
- 2.7. Add 100 µl NEBNext RNA Binding Buffer (2X) to the beads and wash by pipetting up and down 6 times. If preparing multiple libraries, add 100 µl RNA Binding Buffer per sample. The Binding Buffer does not have to be diluted.
- 2.8. Place the tubes on the magnet and incubate at room temperature until the solution is clear (~2 minutes).
- 2.9. Remove and discard the supernatant from the tube. Take care not to disturb the beads.
- 2.10. Remove tubes from magnet and add 50 µl NEBNext RNA Binding Buffer (2X) 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. The Binding Buffer does not have to be diluted.
- 2.11. Add 50 µl beads to each RNA sample from Step 2.1. Mix thoroughly by pipetting up and down 6 times. This first binding step removes most of the non target RNA.

- 2.12. 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 $\geq 75^\circ\text{C}$** to denature the RNA and facilitate binding of the mRNA to the beads.
- 2.13. Remove tubes from the thermal cycler when the temperature reaches 4°C.
- 2.14. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.
- 2.15. Place the tubes on the bench and incubate at room temperature for 5 minutes to allow the RNA to bind to the beads.
- 2.16. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.
- 2.17. Incubate for 5 more minutes on the bench at room temperature to allow the RNA to bind to the beads.
- 2.18. 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.
- 2.19. Remove and discard all of the supernatant. Take care not to disturb the beads.
- 2.20. Remove the tubes from the magnetic rack.
- 2.21. Wash the beads by adding 200 μl of NEBNext Wash Buffer to remove unbound RNA. Pipette the entire volume up and down 6 times to mix thoroughly.
- 2.22. Place the tubes on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
- 2.23. Remove and discard all the supernatant from each well of the tube. Take care not to disturb the beads.
- 2.24. Remove the tubes from the magnetic rack.
- 2.25. Repeat Steps 2.21–2.24.
- 2.26. Add 50 μl of NEBNext Tris Buffer (provided in NEB #E7490 kit) to each tube. Gently pipette the entire volume up and down 6 times to mix thoroughly.
- 2.27. 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.
- 2.28. Remove the tubes from the thermal cycler when the temperature reaches 25°C.
- 2.29. Add 50 μl of NEBNext RNA Binding Buffer (2X) to each sample to allow the RNA to re-bind to the same beads. The Binding Buffer does not have to be diluted. Gently pipette the entire volume up and down 6 times to mix thoroughly.
- 2.30. Incubate the tubes on the bench at room temperature for 5 minutes.
- 2.31. Resuspend the beads. Pipette up and down slowly 6 times to mix thoroughly.
- 2.32. Incubate the tubes on the bench at room temperature for 5 more minutes to allow the RNA to bind to the beads.
- 2.33. Place the tubes on the magnetic stand at room temperature for 2 minutes or until the solution is clear.
- 2.34. Remove and discard all of the supernatant from each tube. Take care not to disturb the beads.
- 2.35. Remove the tubes from the magnetic rack.
- 2.36. Wash the beads once with 200 μl of NEBNext 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.
- 2.37. Place the tubes on the magnetic rack at room temperature for 2 minutes or until the solution is clear.
- 2.38. 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 μl tip remove all of the NEBNext Wash Buffer. (Caution: Do not disturb beads that contain the mRNA). Avoid letting the beads dry out before adding elution buffer.
- 2.39. Remove the tubes from the magnetic rack.

Note: Modifications to elution buffer and volumes will need to be made here to make the enriched mRNA directly compatible with the different NEBNext library preparation workflows. Please refer to Appendix A and consult the appropriate manual for more information.

- 2.40. Elute the mRNA from the beads by adding 17 μ l of the NEBNext Tris Buffer, mix by pipetting 6 times and incubate the samples at 80°C for 2 minutes, then cool to 25°C. Immediately, place the tubes on the magnetic rack for 2 minutes or until the solution is clear.
- 2.41. 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.
- 2.42. Assess the Yield and the Size Distribution of the purified mRNA. Run 1 μ l on the Bioanalyzer using a RNA Pico Chip or the TapeStation using a High Sensitivity RNA ScreenTape. You may have to dilute your sample before loading.

Figure 2: Example of mRNA distribution on a TapeStation.



Appendix A

There are different volumes and reagent requirements to adapt the above protocols directly into a library preparation workflow. Please follow the recommendations below for the library preparation methods listed below and immediately proceed with the steps in the respective manuals:

Product	Reagent(s)	Elution Volume (µl)	Transfer Volume into RT (µl)	Manual Step
NEBNext UltraExpress™ RNA Library Prep Kit (NEB #E3330)	1X Fragmentation Mix*	6.5	5	1.1.1 and 1.2.36
NEBNext Ultra™ II Directional RNA Library Prep Kits (NEB #E7760, #E7765) NEBNext Ultra II RNA Library Prep Kits (NEB #E7770, #E7775)	First Strand Synthesis Reaction Buffer and Random Primer Mix*	11.5	10	1.1.1 and 1.2.36

* These are custom mixes made by the user, and not provided as needed for elution in the kits. Please refer to the respective manuals for more instructions on how to make these mixes before proceeding into bead elution.

Kit Components

NEB #E7490S Table of Components

NEB #	PRODUCT	VOLUME
E7499A	NEBNext Oligo d(T) ₂₅ 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) ₂₅ 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 thermal cycler 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
11.0	Added Express Protocol (Section 1). Also updated header and footer.	2/24

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.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed. See www.neb.com/trademarks. The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email busdev@neb.com.

ALPAQUA® is a registered trademark of Alpaqua Engineering, LLC.

B CORPORATION® is a registered trademark of B Lab IP, LLC, Inc.

AGILENT® and 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 2024, New England Biolabs, Inc.; all rights reserved.

