

Monarch[®] RNA Cleanup Kit Protocol Card

NEB #T2030, #T2040, and #T2050

To download the full manual or view a detailed protocol (including guidance on RNA gel extraction or fractionation into small and large RNA pools), please visit:

Monarch RNA Cleanup Kit (10 µg): www.neb.com/T2030

Monarch RNA Cleanup Kit (50 µg): www.neb.com/T2040

Monarch RNA Cleanup Kit (500 µg): www.neb.com/T2050

BEFORE YOU BEGIN:

- Add 4 volumes of ethanol ($\geq 95\%$) to one volume of RNA Cleanup Wash Buffer.
- If a precipitate has formed in the RNA Cleanup Binding Buffer, warm to room temperature to re-dissolve before use.
- All centrifugation steps should be carried out at room temperature at $16,000 \times g$ (~13,000 RPM).
- The standard protocol outlined below will purify RNA ≥ 25 nt. A simple modification in step 2 can allow for the purification of RNA as small as 15 nt.

PROTOCOL STEPS:

1. **Add 100 µl RNA Cleanup Binding Buffer to the 50 µl sample.** A starting sample volume of 50 µl is recommended. For smaller samples, nuclease-free water can be used to adjust the volume. For samples larger than 50 µl, scale buffer volumes accordingly. Samples with a starting volume > 150 µl will require reloading of the column during step 3.
2. **Add 150 µl (1 volume) of ethanol ($\geq 95\%$) to your sample and mix by pipetting or flicking the tube. Do not vortex.** This will enable the binding of RNA ≥ 25 nt. If you wish to bind RNA as small as

15 nt, add 2 volumes (300 µl) of ethanol to your sample instead of 1 volume (150 µl). The addition of 2 volumes of ethanol shifts the cutoff size of RNA binding from 25 nt down to 15 nt.

- 3. Insert column into collection tube, load sample onto column and close the cap. Spin for 1 minute, then discard flow-through.** For diluted samples > 900 µl, load a portion of the sample, spin, and then repeat as necessary.
- 4. Re-insert column into collection tube. Add 500 µl RNA Cleanup Wash Buffer, spin for 1 minute, then discard the flow-through.**
- 5. Repeat step 4.**
- 6. Transfer column to an RNase-free 1.5 ml microfuge tube (not provided).** Use care to ensure that the tip of the column does not come into contact with the flow-through. If in doubt, re-spin for 1 minute to ensure traces of salt and ethanol are not carried over.

- 7. Elute in nuclease-free water according to the table below.** The eluted RNA can be used immediately or stored at -70°C.

KIT	ELUTION VOLUME**	INCUBATION TIME	SPIN TIME
T2030	6-20 µl	n/a	1 minute
T2040	20-100 µl	n/a	1 minute
T2050*	50-100 µl	5 minutes (room temp)	1 minute

* When cleaning up large amounts of RNA (> 100 µg, #T2050), some precipitation may occur following the addition of the Monarch RNA Cleanup Binding Buffer and ethanol to the sample (Steps 1 and 2). A pellet containing the RNA of interest may form on the side of the column following the first binding spin (Step 3). To maximize recovery of this RNA, a second elution is recommended.

** Yield may slightly increase if a larger volume is used, but the RNA will be less concentrated.

Questions?

Our tech support scientists would be happy to help.
Email us at info@neb.com

This product is covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information about commercial rights, please email us at glb@neb.com. The use of this product may require you to obtain additional third party intellectual property rights for certain applications.

New England Biolabs is an ISO 9001, ISO 14001 and ISO 13485 certified facility.

© Copyright 2019, New England Biolabs, Inc.; all rights reserved.

This card is made with FSC certified 100% post-consumer fiber. Please recycle.

V2.0 - 7.19 #102961

