

Monarch[®] PCR & DNA Cleanup Kit (5 µg) Protocol Card

NEB #T1030

For a detailed protocol or to download the full manual, visit www.neb.com/T1030.

BEFORE YOU BEGIN:

- Add 4 volumes of ethanol (≥ 95%) to one volume of DNA Wash Buffer.
- All centrifugation steps should be carried out at 16,000 x g (~13,000 RPM).

THERE ARE TWO PROTOCOLS AVAILABLE FOR THIS PRODUCT:

- **DNA Cleanup and Concentration:** for the purification of up to 5 µg of DNA (ssDNA > 200 nt and dsDNA > 50 bp) from PCR and other enzymatic reactions.
- **Oligonucleotide Cleanup:** for the purification of up to 5 µg of DNA fragments ≥ 15 bp (dsDNA) or ≥ 18 nt (ssDNA). Expected recovery is > 70%. When purifying ssDNA of any size, recovery can be increased by using this protocol; however, it is important to note that this protocol shifts the cutoff for smaller fragments to 18 nt (rather than 50 nt for the DNA Cleanup and Concentration Protocol).

DNA CLEANUP AND CONCENTRATION PROTOCOL STEPS:

1. Dilute sample with DNA Cleanup Binding Buffer according to the table below. Mix well by pipetting up and down or flicking the tube. Do not vortex. We recommend a sample volume of 20–100 µl. For smaller samples, adjust the volume with TE. For diluted samples larger than 800 µl, load a portion of the sample, proceed with step 2, and repeat as necessary.

Sample Type	Ratio of Binding Buffer: Sample	Example
dsDNA > 2 kb (plasmids, gDNA)	2:1	200 µl: 100 µl
dsDNA < 2 kb (some amplicons, fragments)	5:1	500 µl: 100 µl
ssDNA* > 200 nt	7:1	700 µl: 100 µl

*Please note that recovery of ssDNA < 200 nts can be increased by using the **Oligonucleotide Cleanup Protocol**, but doing so will shift the cutoff size for DNA binding to 18 nt (versus 50 nt).

2. Insert column into collection tube and load sample onto column. Spin for 1 minute, then discard flow-through.

- 3. Re-insert column into collection tube. Add 200 μ l DNA Wash Buffer and spin for 1 minute.** Discarding flow-through is optional.
- 4. Repeat step 3.**
- 5. Transfer column to a clean 1.5 ml microfuge tube.** 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.
- 6. Add $\geq 6 \mu$ l of DNA Elution Buffer to the center of the matrix. Wait for 1 minute, then spin for 1 minute to elute DNA.** Typical elution volumes are 6–20 μ l. Nuclease-free water (pH 7–8.5) can also be used to elute the DNA. Yield may slightly increase if a larger volume of DNA Elution Buffer is used, but the DNA will be less concentrated. For larger size DNA (≥ 10 kb), heating the elution buffer to 50°C prior to use can improve yield.

Want to use this kit to purify DNA from agarose gels?

Simply purchase the Monarch Gel Dissolving Buffer (NEB #T1021L) and use with this kit. Protocol available at www.neb.com/T1020

Have any questions?

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

One or more of these products are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. For more information, please contact NEB's Global Business Development team at gbd@neb.com. The use of these products 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 2017, New England Biolabs, Inc.; all rights reserved.

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

V2- 12.17

OLIGONUCLEOTIDE CLEANUP PROTOCOL STEPS:

- 1. Add 100 μ l DNA Cleanup Binding Buffer to the 50 μ l sample.** We recommend a sample volume of 50 μ l. For smaller samples, adjust the volume with nuclease-free water.
- 2. Add 300 μ l ethanol ($\geq 95\%$). Mix well by pipetting up and down or flicking the tube. Do not vortex.**
- 3. Insert column into collection tube and load sample onto column. Spin for 1 minute, then discard flow-through.**
- 4. Re-insert column into collection tube. Add 500 μ l DNA Wash Buffer and spin for 1 minute.** Discard flow-through.
- 5. (Optional) Repeat step 4.** This second wash step is optional, but recommended for removal of enzymes that may interfere with downstream applications (e.g., Proteinase K). Please note that if carrying out a second wash step, additional DNA Wash Buffer may be required.
- 6. Transfer column to a clean 1.5 ml microfuge tube.** 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.
- 7. Add $\geq 6 \mu$ l of DNA Elution Buffer to the center of the matrix. Wait for 1 minute, then spin for 1 minute to elute DNA.** Typical elution volumes are 6–20 μ l. Nuclease-free water (pH 7–8.5) can also be used to elute the DNA. Yield may slightly increase if a larger volume of DNA Elution Buffer is used, but the DNA will be less concentrated.

