

Monarch® Spin PCR & DNA Cleanup Kit (5 µg)

NEB #T1130S/L

50/250 preps

Version 1.0 06.24

Table of Contents

| | |
|--|-------------------|
| Kit Contents and Storage | 2 |
| Storage recommendations | 2 |
| Intended Use | 2 |
| Safety Information | 2 |
| Quality Control | 2 |
| Introduction | 3 |
| Features | 3 |
| Sustainability and Recycling Information | 3 |
| Overview | 3 |
| Properties | 5 |
| Applications & Usage | 6 |
| Important Notes | 6 |
| DNA Size | 6 |
| Monarch DNA Cleanup Protocols Overview and Comparison | 6 |
| General Guidelines | 7 |
| Equipment and Reagents Required & Supplied by User | 7 |
| Buffer Preparation | 8 |
| Protocol | 8 |
| Protocol: PCR & DNA Cleanup Kit (5 µg) Using Centrifugation | 8 |
| Protocol: PCR & DNA Cleanup Kit (5 µg) Using Vacuum Manifold | 9 |
| Protocol: Oligonucleotide Cleanup Protocol | 9 |
| Troubleshooting | 9 |
| Ordering Information | 10 |
| Revision History | 10 |

Kit Contents and Storage

| Component | NEB # | Application/Usage | T1130S 50 preps | T1130L 250 preps | Storage Temperature |
|-------------------------------|-------|---|--------------------|---------------------|------------------------|
| Monarch Buffer BZ | T1114 | Binding buffer concentrate (1.42X) | 42 ml | 168 ml | 15-25°C |
| Monarch Buffer WZ | T1115 | Wash buffer concentrate (5X) | 5 ml | 26 ml | 15-25°C |
| Monarch Buffer EY | T1116 | Elution buffer | 3 ml | 7 ml | 15-25°C |
| Monarch Spin Columns S1A | T2037 | Spin column for nucleic acid purification | 50 columns | 250 columns | 15-25°C |
| Monarch Spin Collection Tubes | T2118 | Collection tube | 50 tubes | 250 tubes | 15-25°C |

Storage Recommendation

- All kit components should be stored at room temperature.
- Always keep reagent bottles tightly closed.
- Keep columns sealed in the enclosed bag.
- See individual component labels for specific storage guidance.

Intended Use

The Monarch Spin PCR & DNA Cleanup Kit is developed for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

Safety Information

- Monarch Buffer BZ contains guanidine hydrochloride, which can form highly reactive compounds when combined with bleach. Do not add bleach or acidic solutions directly to the buffers or to the sample preparation waste.
- For more information regarding the composition of buffers, please consult the Safety Data Sheets available on our website www.neb.com/T1130.
- Proper laboratory safety practices should be employed using this kit, including the use of lab coats, gloves, and eye protection.

Quality Control

To help ensure consistent quality and performance, each lot of this kit is tested for predetermined quality control and functional specifications.

Introduction

The Monarch Spin PCR & DNA Cleanup Kit (5 µg) is a rapid and reliable method for the purification and concentration of up to 5 µg of high-quality, double-stranded and single-stranded DNA from enzymatic reactions such as PCR, restriction digestion, ligation, and reverse transcription. Designed with sustainability in mind, these kits use significantly less plastic than other kits on the market.

Features of this kit include:

- **High Performance:** Achieve high yields (up to 5 µg) and high purity in the purification, cleanup, and concentration of DNA. This kit has the capability to remove short primers, detergents, and other low-molecular-weight reaction components (e.g., nucleotides, DMSO, betaine).
- **High Concentration:** Elute in very small volumes, as little as 5 µl, allowing for highly concentrated DNA.
- **Powerful Flexibility** – Purify small DNA fragments, including oligonucleotides, using the modified protocol provided.
- **Time Savings:** The protocol takes only 5 minutes to complete the bind, wash, and elution steps, with minimal incubation and spin times.
- **Unique Column Design:** Spin column features a unique design that enables elution in low volumes and minimizes buffer retention and contaminant carryover.
- **Optimized Buffers:** Buffer system is optimized, without the need to adjust pH.
- **Application Compatibility:** Purified DNA is ready for downstream molecular applications, such as restriction digests, DNA sequencing, ligation, amplification, and other enzymatic reactions.

Sustainability and Recycling Information

Monarch DNA and RNA Purification Kits are designed for sustainability and developed for performance. Learn more about Monarch sustainability at www.neb.com/monarchsustainability.

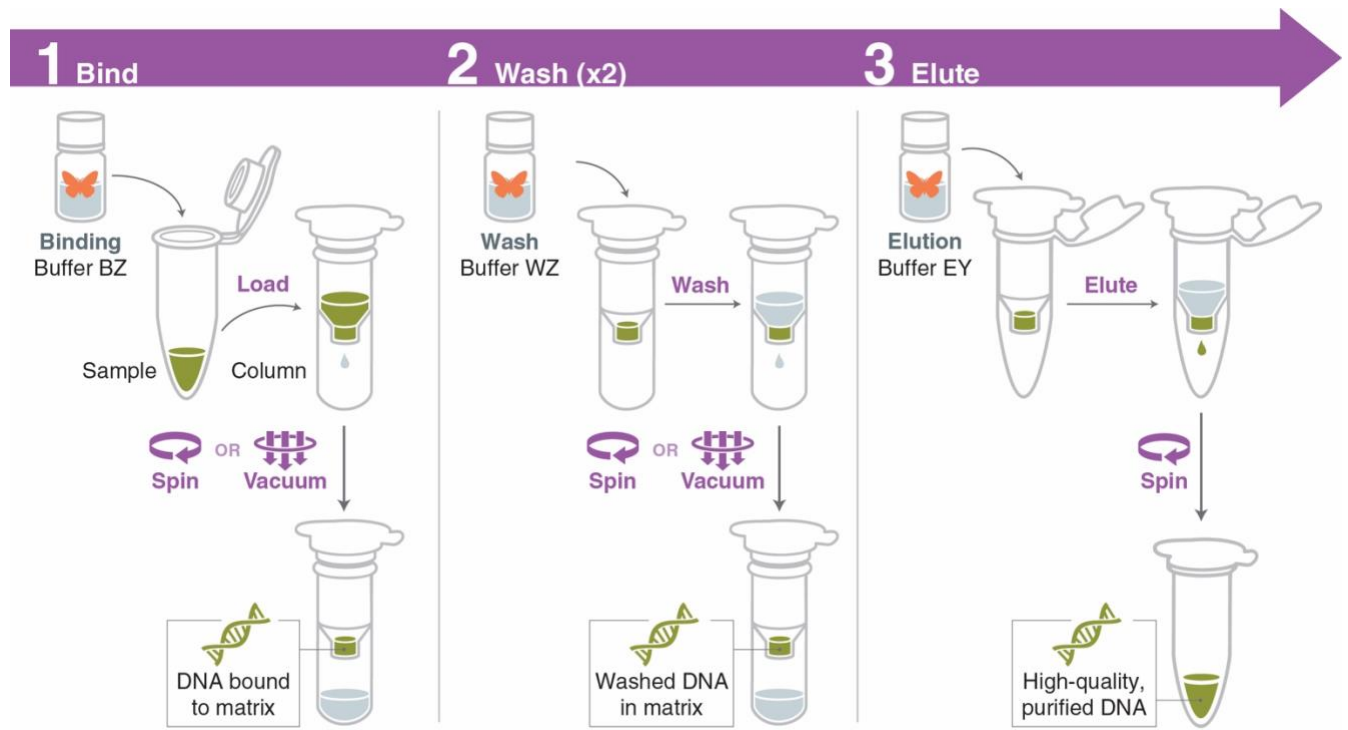
- **Sustainable performance:** Significantly less plastic is used in spin columns, bottles, and other plastic parts, compared to similar kits from other leading suppliers.
- **Thinner-walled columns:** Reduction in total plastic without affecting performance.
- **Flexible purchasing options:** Columns also available separately. Purchase only what you need and avoid wasted materials.
- **Same performance, design, and formulations:** Standalone products are the same components and formulations that are included in complete kits.
- **Streamlined packaging:** Sturdy, reusable boxes at just the right size with concise protocol cards that replace printed manuals.
- **Sustainable and recyclable packaging:** Packaging printed with less ink using eco-friendly practices and powered by sustainable sources such as wind, where possible. Packaging is sourced for recyclability and recycled paper is used where possible to make the kit boxes, inserts, and paper materials.

Help keep Monarch sustainable by recycling after using. Learn more on how to recycle Monarch boxes and kit components at www.neb.com/monarchrecycling.

Overview of Monarch Spin PCR & DNA Cleanup Kit (5 µg)

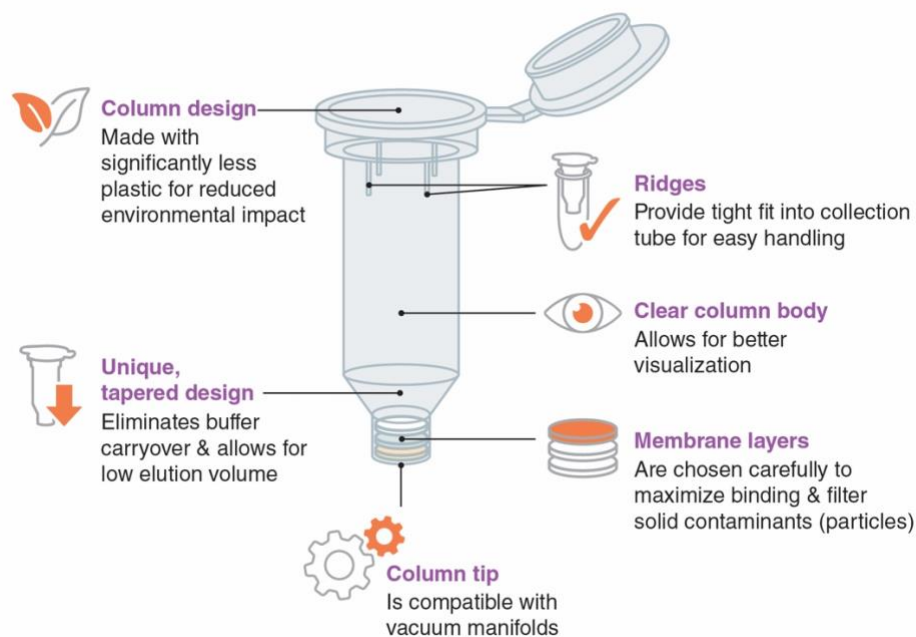
The Monarch Spin PCR & DNA Cleanup kit employs an advanced bind/wash/elute workflow combined with silica-membrane technology. Monarch Buffer BZ (DNA binding buffer) is designed with high salt concentrations, enabling optimized binding of the sample DNA to our unique spin column. Subsequently, the Monarch Buffer WZ (wash buffer) effectively removes enzymes, reaction components, and salts. Elution under low-salt conditions yields highly pure DNA, suitable for diverse downstream applications, including restriction digests, DNA sequencing, ligation, and other enzymatic manipulations. The innovative column design, with precision-engineered silica matrices and layers, enables elution volumes as low as 5 µl with reduced buffer retention and contaminants. Combined with our optimized buffer system, the kit offers rapid and reliable purification and concentration of up to 5 µg of high-quality DNA.

Figure 1: DNA Cleanup Workflow



The Monarch Spin PCR & DNA Cleanup Kit (5 µg) uses the bind/wash/elution method and a unique spin column.

Figure 2. Monarch Column Design



NEB Monarch's unique column design allows high-quality DNA purification with low elution volume. The column is designed and made with significantly less plastic for a reduced environmental impact.

Properties

| | |
|---|--|
| Purification format | Spin-column |
| DNA Sample* | DNA from PCR and other enzymatic reactions (e.g., restriction digests, kinase reactions, ligations) |
| | ssDNA or dsDNA oligonucleotides from enzymatic reactions can also be purified using the Oligonucleotide Cleanup Protocol |
| Typical Recovery | 70-90% |
| DNA Purity | $A_{260/280} > 1.8$ and $A_{260/230} > 1.8$ |
| Binding Capacity | Up to 5 μg |
| Elution Volume | 5-20 μl |
| DNA Size Range | Standard protocol: 50 bp – 25 kb |
| | Oligonucleotide Cleanup protocol: ssDNA > 16 nt and dsDNA > 12 nt |
| Protocol Time | 5 minutes of spin and incubation time |
| Compatible Downstream Applications | Ligation, restriction digestion, labeling and other enzymatic manipulations, library construction and DNA sequencing |

*See next table for the extended list of DNA samples that can be used with this kit

Applications & Usage

Listed are selected examples of applications and usage. To see the most updated list, refer to the product webpage.

| | |
|--|--|
| PCR Cleanup | DNA from PCR reactions can be purified after amplification to remove polymerases, primers, detergents, dNTPs, etc. |
| Enzymatic reaction cleanup | Restriction enzymes and modifying enzymes such as ligases, kinases, nucleases, phosphatases are efficiently removed, allowing for effective desalting and concentration of the DNA sample. |
| cDNA cleanup | DNA/RNA complexes can be purified post-reverse transcription/ amplification to enable removal of the RT and polymerase as well as nucleotides. |
| Labeling cleanup | Unincorporated radiolabeled or fluorescently labeled nucleotides can be removed from the DNA substrate. |
| Plasmid cleanup | Plasmid preps from unknown sources may contain inhibitors and unwanted contaminants. Purification and concentration can be easily achieved using this kit. |
| Oligonucleotide and ssDNA Purification | ssDNA oligonucleotides (≥ 16 nt) and dsDNA fragments (≥ 12 bp) can be purified using the Oligonucleotide Cleanup Protocol (page 9). |

Important notes before starting

The Monarch Spin PCR & DNA Cleanup Kit (5 μ g) is designed to ensure optimal DNA yield and quality, accommodating for various types of DNA and sizes. Although the kit is optimized for a broad range of conditions, it is crucial to carefully consider these influencing factors to ensure high quality and maximize DNA recovery.

DNA Size

Typically, longer DNA exhibits a stronger affinity to silica in the presence of chaotropic salt, resulting in tight binding. For Monarch Spin PCR & DNA Cleanup Kit (5 μ g), DNA exceeding 15 kb may bind tightly to the silica column and become difficult to elute. If working with a DNA longer than 15 kb, a modified elution method can be employed to increase elution efficiency. For a more detailed procedural guide, we recommend reading the full protocol provided in this manual.

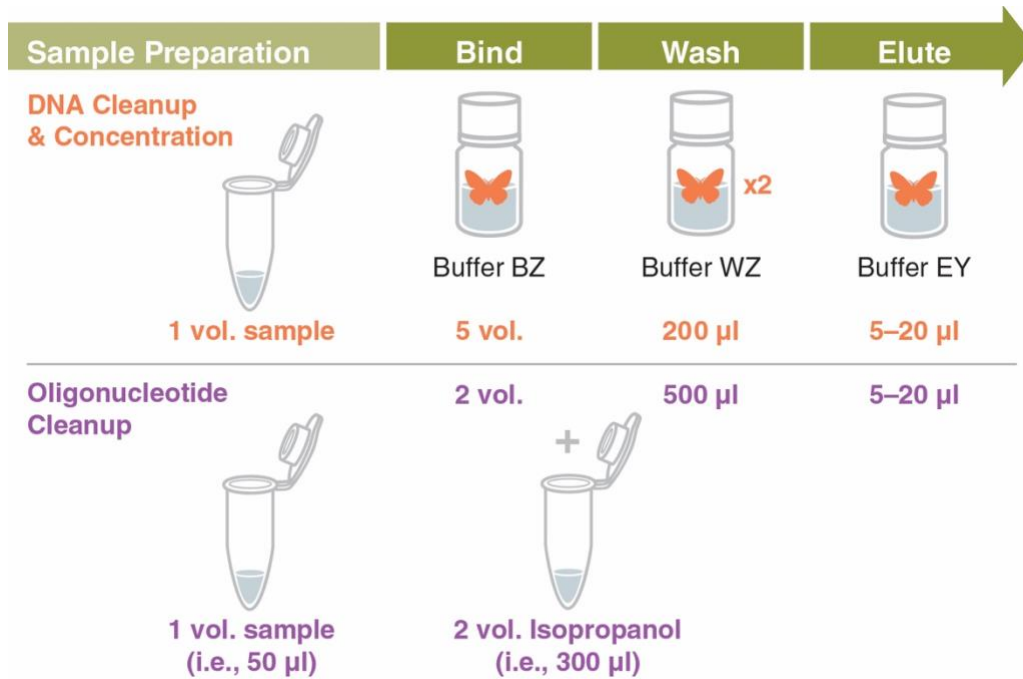
For DNA < 50 bp or oligonucleotide purification, we provide a separate protocol to allow optimal recovery. Please read the next section describing 2 different protocols in detail to choose the best protocol for your samples.

Monarch Protocol overview and comparison

The Monarch Spin PCR & DNA Cleanup Kit (5 μ g) provides 2 different protocol types for the different applications of 1) PCR and DNA cleanup and 2) oligonucleotide and short DNA fragments cleanup.

| | DNA Cleanup and Concentration | Oligonucleotide Cleanup |
|-----------------------------|---|---|
| Effective size range | dsDNA > 50 bp and ssDNA > 200 nt | dsDNA ≥ 12 bp or ssDNA ≥ 16 nt |
| Typical usage | <ul style="list-style-type: none"> Purification from PCR and other enzymatic reactions Removes primers, nucleotides, enzymes and other enzyme reaction components | <ul style="list-style-type: none"> Purification and recovery of oligonucleotide and ssDNA This protocol will not remove primers from PCR reactions as DNA size cutoff is 12 bp/16 nt. |

Workflow diagram comparing DNA Cleanup and Oligonucleotide Cleanup



See detailed protocol for specific instructions.

General Guidelines for Monarch Spin PCR & DNA Cleanup Kit (5 µg)

- The input amount of DNA to be purified should not exceed the binding capacity of the columns (5 µg).
- A starting sample volume of 20–100 µl is recommended. For smaller samples, nuclease-free water or TE buffer can be used to adjust the volume to the recommended volume range.
- Centrifugation should be carried out at 16,000 x g (~13,000 RPM) in a standard laboratory microcentrifuge at room temperature. This ensures all traces of the buffer are removed at each step.
- The column holds a maximum volume of 800 µl.
- Always keep columns tightly sealed in the provided bag.

Equipment and Reagents Required & Supplied by the User

Equipment

- Benchtop microcentrifuge
- Vacuum manifold (for the vacuum manifold protocol)
- Vacuum pump (for the vacuum manifold protocol)

Reagents/supplies

- Isopropanol (100%)
- Ethanol (≥ 95%)
- 1.5 ml or 2 ml microfuge tubes
- Optional: Nuclease-free water for elution, if provided elution buffer will not be used

Buffer Preparation

Add isopropanol to the Monarch Buffer BZ prior to use (0.43 volume of isopropanol per volume of Buffer BZ).

- For T1130S (50-prep) kit, add 18 ml of isopropanol to Monarch Buffer BZ.
- For T1130L (250-prep) kit, add 72 ml of isopropanol to Monarch Buffer BZ.

Add ethanol to the Monarch Buffer WZ prior to use (4 volumes of $\geq 95\%$ ethanol per volume of Buffer WZ).

- For T1130S (50-prep) kit, add 20 ml of ethanol to Buffer WZ.
- For T1130L (250-prep) kit, add 104 ml of ethanol to Buffer WZ.

Always keep all buffer bottles tightly closed when not actively in use.

Monarch Spin PCR & DNA Cleanup Kit Protocols:

1. Standard Cleanup Protocol using centrifugation
2. Standard Cleanup Protocol using a vacuum manifold
3. Oligonucleotide Cleanup Protocol

Standard Cleanup Protocol using Centrifugation

1. **Add 5 volumes (e.g., 250 μ l) of Monarch Buffer BZ to 1 volume (e.g., 50 μ l) of sample. Mix well by pipetting up and down or flicking the tube. Do not vortex.** Using a sample volume of 20-100 μ l is recommended. For samples less than 20 μ l, adjust the volume with TE or nuclease-free water to 20-100 μ l. For diluted samples larger than 800 μ l, load 800 μ l first, proceed with step 2, and repeat as needed.
2. **Insert the Monarch Spin Column S1A into the Monarch Spin Collection Tube and load the sample onto the column. Spin for 1 minute, then discard the flow-through.**
3. **Re-insert the column into the collection tube. Wash by adding 200 μ l of Monarch Buffer WZ and spin for 1 minute.** Discarding flow-through is optional.
4. **Repeat wash (step 3).**
5. **Transfer the column to a clean 1.5 ml microfuge tube.** Use care to ensure that the tip of the column does not touch the flow-through. If in doubt, re-spin for 1 minute.
6. **Add 5-20 μ l of Monarch Buffer EY to the center of the matrix to elute DNA. Wait for 1 minute, and spin for 1 minute.** Nuclease-free water can also be used to elute the DNA. Yield may slightly increase if a larger volume of Monarch Buffer EY is used, but the DNA will be less concentrated. For larger size DNA (≥ 15 kb), incubate the column with elution buffer at room temperature for 5 minutes to maximize the yield. Alternatively, heating the elution buffer to 50°C prior to use can be used.

Standard Cleanup Protocol using a Vacuum Manifold

1. **Add 5 volumes (e.g., 250 μ l) of Monarch Buffer BZ to 1 volume (e.g., 50 μ l) of sample. Mix well by pipetting up and down or flicking the tube. Do not vortex.** We recommend a sample volume of 20-100 μ l. For smaller samples, adjust the volume with TE.
2. **Insert the column into the vacuum adapter or manifold directly, switch the vacuum on, and load the sample onto the column. Allow the solution to pass through the column, then switch the vacuum source off.** Make sure to follow the manifold manufacturer's instructions to set up the manifold and connect it properly to a vacuum source.
3. **Wash by adding 200 μ l of Monarch Buffer WZ and switch the vacuum on. Allow the solution to pass through the columns, then switch the vacuum source off.**
4. **Repeat wash (step 3).**
5. **(Recommended) Insert the column into the Monarch Spin Collection Tube and centrifuge for 1 minute.** Since vacuum set-ups can vary, centrifugation is recommended before the elution step to ensure no traces of buffer and ethanol are carried over.
6. **Transfer the column to a clean 1.5 ml microfuge tube.** Use care to ensure that the tip of the column does not touch the flow-through. If in doubt, re-spin for 1 minute.
7. **Add 5-20 μ l of Monarch Buffer EY to the center of the matrix to elute DNA. Wait for 1 minute, and spin for 1 minute.** Nuclease-free water can also be used to elute the DNA. Yield may slightly increase if a larger volume of Monarch Buffer EY is used,

but the DNA will be less concentrated. For larger size DNA (≥ 15 kb), incubate the column with elution buffer at room temperature for 5 minutes to maximize the yield. Alternatively, heating the elution buffer to 50°C prior to use can be used.

Oligonucleotide Cleanup Protocol

1. **Add 100 μ l Monarch Buffer BZ (2 volumes) (ensure that isopropanol has been added, as indicated on the bottle label) to 50 μ l (1 volume) sample.** A sample volume of 50 μ l is recommended. For smaller samples, adjust the volume with nuclease-free water. If using a sample > 90 μ l, reloading the column during step 3 will be required.
2. **Add 300 μ l (2 volumes) isopropanol. Mix well by pipetting up and down or flicking the tube. Do not vortex.**
3. **Insert the column into the collection tube and load the sample onto the column. Spin for 1 minute, then discard the flow-through.**
4. **Re-insert the column into the collection tube. Wash by adding 500 μ l Monarch Buffer WZ and spin for 1 minute.** Discard flow-through. Additionally, a second wash step with Monarch Wash buffer WZ or 80% ethanol can be carried out to ensure the removal of enzymes that may interfere with downstream applications.
5. **Transfer the column to a clean 1.5 ml microfuge tube.** Use care to ensure that the tip of the column does not touch the flow-through. If in doubt, re-spin for 1 minute.
6. **Add 5-20 μ l of Monarch Buffer EY to the center of the matrix to elute DNA. Wait for 1 minute, and spin for 1 minute.** Nuclease-free water can also be used to elute the DNA. Yield may slightly increase if a larger volume of Monarch Buffer EY is used, but the DNA will be less concentrated.

Troubleshooting

| Problem | Common Cause | Suggestions/Solutions |
|--------------------------------|------------------------------------|---|
| No DNA purified | Ethanol not added to wash buffer | Ensure proper amount of ethanol was added to the wash buffer. |
| Low DNA yield | Reagent added incorrectly | Check protocol to ensure correct buffer reconstitution, order of addition of buffers and proper handling of column flow-through and eluates. |
| | Incomplete elution | Ensure the Monarch Buffer EY (elution buffer) is added correctly to the center of the matrix. Larger elution volumes and longer incubation time can increase the yield of DNA, especially when the amount of DNA is close to maximum binding capacity and the DNA size is large (>10 kb). Alternatively, heating the elution buffer to 50°C prior to elution step may also increase the yield. |
| Low DNA purity and performance | Ethanol is carried over | Ensure the final wash spin time is 1 minute for the complete removal of the wash buffer. Carefully transfer the column to a microfuge tube ensuring the tip of the column does not touch the flow-through. |
| | Trace amounts of salt carried over | Carried-over salts will be indicated by a low $A_{260/230}$ ratio. Ensure the column tip does not touch the flow through. |

For more troubleshooting and FAQs, please visit product webpage or reach out to our technical support team at info@neb.com

Ordering Information

Monarch Spin PCR & DNA Cleanup Kit (5 µg)

| PRODUCT | NEB# |
|---|-------|
| Monarch Spin PCR & DNA Cleanup Kit (5 µg) | T1130 |
| Kit components sold separately | |
| Monarch Spin Columns S1A and Tubes | T2037 |
| Monarch Spin Collection Tubes | T2118 |

NEB Companion Products

| PRODUCT | NEB# |
|---------------------------------------|-------|
| Exo-CIP Rapid PCR Cleanup Kit | E1050 |
| Gel Loading Dye, Purple (6x) | B7024 |
| Gel Loading Dye, Purple (6x), no SDS | B7025 |
| Quick-Load® Purple 1 kb DNA Ladder | N0552 |
| Quick-Load Purple 100 bp DNA Ladder | N0551 |
| Quick-Load Purple 1kb Plus DNA Ladder | N0550 |
| T4 DNA Ligase | M0202 |
| Blunt/TA Ligase Master Mix | M0367 |
| Instant Sticky-end Ligase Master Mix | M0379 |

Revision History

| REVISION # | DESCRIPTION | DATE |
|------------|-------------|-------|
| 1.0 | – | 06/24 |

This product is intended for research purposes only. This product is not intended 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. For more information about commercial rights, please email us at busdev@neb.com. While NEB develops and validates its products for various applications, using this product may require the buyer to obtain additional third party intellectual property rights for certain applications.

B CORPORATION™ is a trademark of B Lab, Inc.

© Copyright 2024, New England Biolabs, Inc.; all rights reserved



New England Biolabs, Inc., 240 County Road, Ipswich, MA 01938-2723 | T: (978) 927-5054 F: (978) 921-1350 US Orders (Toll-free): 1-800-632-5227 US Tech Support: 1-800-632-7799 | info@neb.com