

Monarch[®] Genomic DNA Purification Kit Protocol

NEB #T3010

We recommend that first-time users of this kit review the product manual at neb.com/T3010 before starting; it provides additional information to consider at various steps. This quick protocol is meant for experienced users. The website and manual also contain protocols for gDNA cleanup and extraction of gDNA from additional sample types, including bacteria, yeast, saliva and buccal swabs.

BEFORE YOU BEGIN

- Store RNase A and Proteinase K at -20°C.
- Add ethanol (≥ 95%) to the gDNA Wash Buffer concentrate as indicated on the bottle label.
- Set a thermal mixer (e.g. ThermoMixer[®]) or, if not available, a heating block to 56°C for sample lysis.
- Set a heating block to 60°C. Preheat the appropriate volume of elution buffer to 60°C (35-100 µl per sample).

GENOMIC DNA PURIFICATION CONSISTS OF TWO STAGES:

Part 1: Sample Lysis

Part 2: Genomic DNA Binding and Elution

Part 1: SAMPLE LYSIS

☐ Cultured Cells:

1. **Start with a cell pellet containing 1×10^4 – 5×10^6 cells (typical starting amount is 1×10^6 cells).**

Frozen cells: thaw cell pellet slowly on ice and loosen by flicking the tube several times. Resuspend in 100 μ l cold PBS by pipetting up and down.

Fresh cells: pellet by centrifugation at 1,000 x g for 1 minute and resuspend in 100 μ l cold PBS by pipetting up and down.

Ensure pellet is resuspended completely. If using lower cell inputs, the use of carrier RNA may be beneficial; see product manual for details.

2. **Add 1 μ l Proteinase K and 3 μ l RNase A to the resuspended pellet and mix by vortexing briefly to ensure the enzymes are efficiently dispersed.** Do not add the enzymes and the Cell Lysis Buffer simultaneously, as the high viscosity of the lysate will prevent equal distribution of the enzymes. Addition of RNase A can be omitted if a low percentage of co-purified RNA will not affect downstream applications.
3. **Add 100 μ l Cell Lysis Buffer and vortex immediately and thoroughly.** The solution will rapidly become viscous.
4. **Incubate for 5 minutes at 56°C in a thermal mixer with agitation at full speed (~1400 rpm).** Incubation for longer than 5 minutes is not necessary, but will not negatively affect the quality of the purified gDNA. If a thermal mixer is not available, use a heating block and vortex occasionally.
5. **Proceed to step 1 of Part 2: Genomic DNA Binding and Elution.**

☐ Mammalian Whole Blood (non-nucleated)

1. **Transfer 100 μ l of whole blood to a 1.5 ml microfuge tube.** If processing less than 100 μ l of blood, add cold PBS to bring the total volume to 100 μ l. For pre-aliquoted frozen samples, do not thaw; add Proteinase K, RNase A and Blood Lysis Buffer to the frozen sample in step 2.

2. **Add 10 μ l Proteinase K, 3 μ l RNase A and 100 μ l of Blood Lysis Buffer. Mix immediately by vortexing.** For frozen blood, do not thaw; add enzymes and lysis buffer to frozen sample and proceed immediately to step 3. When working with multiple samples, prepare a master mix of the three reagents to save pipetting steps. Addition of RNase A can be omitted if a low percentage of co-purified RNA will not affect downstream applications.
3. **Incubate for 5 minutes at 56°C in a thermal mixer with agitation at full speed (~1400 rpm).** Incubation for longer than 5 minutes is not necessary, but will not negatively affect the quality of the purified gDNA from most species. If a thermal mixer is not available, use a heating block and vortex occasionally.
4. **Proceed to step 1 of Part 2: Genomic DNA Binding and Elution.**

☐ Nucleated Whole Blood (birds, reptiles)

1. **Transfer 10 μ l whole blood to a 1.5 ml microfuge tube.**
2. **Add 90 μ l cold PBS and mix by vortexing.**
3. **Add 10 μ l Proteinase K and 3 μ l RNase A, and mix again by vortexing.** Do not add the enzymes and the Blood Lysis Buffer simultaneously, as the high viscosity of the lysate will prevent equal distribution of the enzymes. Addition of RNase A can be omitted if a low percentage of co-purified RNA will not affect downstream applications.
4. **Add 100 μ l Blood Lysis Buffer and vortex thoroughly.** The solution will rapidly become viscous.
5. **Incubate for 5 minutes at 56°C in a thermal mixer with agitation at full speed (~1400 rpm).** Incubation for longer than 5 minutes is not necessary, but will not negatively affect the quality of the purified gDNA.
6. **Proceed to step 1 of Part 2: Genomic DNA Binding and Elution**

□ Animal Tissue

1. **Cut tissue into small pieces to ensure rapid lysis and high yields. Weigh the appropriate tissue amount and place in a 1.5 ml microfuge tube (see table below for recommended input amounts).** Using more than the recommended amounts will not lead to better yields and/or purity. If using more than recommended is required, split the sample into 2 or more preps. Ensure frozen material remains frozen until samples are mixed with Tissue Lysis Buffer and Proteinase K. Both stabilized and fresh tissue should be kept cold or on ice during preparation.

STARTING MATERIAL	RECOMMENDED INPUT AMOUNT
Rodent tail	up to 25 mg
Brain	up to 12 mg
Fibrous tissue (muscle, heart)	up to 25 mg
Ear clips, skin	up to 10 mg
Liver, lung	up to 15 mg
Spleen, kidney	up to 10 mg

2. **Add Proteinase K (according to the table below) and 200 μ l of Tissue Lysis Buffer to each sample. Mix immediately by vortexing.** Ensure tissue particles are able to move freely in the lysis mix and do not remain stuck on the bottom of the tube. When working with multiple samples, prepare a master mix of Tissue Lysis Buffer and Proteinase K to save pipetting steps.

TISSUE TYPE	PROTEINASE K AMOUNT
Brain, Kidney, Skin, Ear Clips	3 μ l
All other tissues	10 μ l

3. **Incubate at 56°C in a thermal mixer with agitation at full speed (1400 rpm) until tissue pieces have completely dissolved (typically 30-60 minutes).** If time is not limiting, additional incubation up to 3 hours can further improve yields and decrease residual RNA. If an incubator with agitation is not available, use a tube rotator placed within an incubator, shaking water bath or a heating block (vortex samples every 5-15 minutes to speed up lysis).
4. **[Note: The following step can be omitted when working with fresh or frozen tissue amounts <15 mg]. Centrifuge for 3 minutes at maximum speed (>12,000 x g) to pellet debris. Transfer the supernatant to a fresh microfuge tube.** This prevents residual debris from clogging the membrane binding sites and helps to reach maximal yield and purity. It is especially important to perform this step if sample appears turbid, contains residual particles, when working with stabilized tissue, or when working with brain or fibrous tissues.
5. **Add 3 μ l of RNase A to the lysate, vortex thoroughly, and incubate for a minimum of 5 minutes at 56°C with agitation at full speed.** This step can be skipped if a low percentage of co-purified RNA will not affect downstream applications.
6. **Proceed to step 1 of Part 2: Genomic DNA Binding and Elution.**

PART 2: GENOMIC DNA BINDING AND ELUTION

- 1. Add 400 μ l gDNA Binding Buffer to the sample and mix thoroughly by pulse-vortexing for 5-10 seconds.** Thorough mixing is essential for optimal results.
- 2. Transfer the lysate/binding buffer mix (~600 μ l) to a gDNA Purification Column pre-inserted into a collection tube, without touching the upper column area.** Proceed immediately to step 3. Avoid transferring foam that may have formed during lysis.
- 3. Close the cap and centrifuge: first for 3 minutes at 1,000 x g to bind the gDNA (no need to empty the collection tube or remove from centrifuge), and then for 1 minute at maximum speed (>12,000 x g) to clear the membrane. Discard the flow-through and the collection tube.**
- 4. Transfer column to a new collection tube and add 500 μ l gDNA Wash Buffer. Close the cap and invert a few times so that the wash buffer reaches the cap (do not vortex). Centrifuge immediately for 1 minute at maximum speed, then discard the flow through.** The collection tube can be tapped on a paper towel to remove any residual buffer before reusing it in the next step.
- 5. Replace the column into the collection tube. Add 500 μ l gDNA Wash Buffer and close cap. Centrifuge immediately for 1 minute at maximum speed. Discard the collection tube and flow through.**
- 6. Place the gDNA Purification Column in a DNase-free 1.5 ml microfuge tube (not included). Add 35-100 μ l preheated (60°C) gDNA Elution Buffer, close the cap and incubate at room temperature for 1 minute.** Elution in 100 μ l is recommended, but smaller volumes can be used and will result in more concentrated DNA but a reduced yield (20-25% reduction when using 35 μ l). Eluting with preheated elution buffer will increase yields by ~20-40% and eliminates the need for a second elution. The elution buffer (10 mM Tris-Cl, pH 9.0, 0.1 mM EDTA) offers strong protection against enzymatic degradation and is optimal for long term storage of DNA. However, other low-salt buffers or nuclease-free water can be used if preferred.
- 7. Centrifuge for 1 minute at maximum speed (> 12,000 x g) to elute the gDNA.**

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 gd@neb.com. The use of this product may require you to obtain additional third party intellectual property rights for certain applications.

THERMOMIXER® is a registered trademark of Eppendorf AG.

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.

V3.0 - 7.19 #103248

