



DNA extraction from fresh/frozen tissue using GenFind V3

Researchers working on fresh/frozen tissue who want to extract DNA and use the same kit as other sample types may use this protocol. Please reference current GenFind V3 IFU for product information (Part Number C34880, C34881)

Purpose

This provides a solution that allows a user to use one kit to extract DNA from multiple sample types with small modifications in the workflow.

Materials Used

Material	Part Number	Supplier
1.2 mL 96-well plate	AB1127	Thermo Fisher Scientific
100% Ethanol (Molecular Grade)	AB00138	American Bio
Nuclease-free water (Molecular Grade)	AM9932	Thermo Fisher Scientific
2 mL 96-well plate	609681	Beckman Coulter
7 Bar Magnet for 96-Well Plate	771MWZM-1ALT	V&P Scientific
Stainless Steel Beads, 5 mm	69989	Qiagen

Protocol

1. Sample Preparation

- a. Cut off 10 mg of tissue and place in a 1.2 mL 96-well plate

2. Lysis

- a. Add **1 stainless steel** bead to the plate
- b. Add **100 µL** of **Lysis (LBB)** to the plate
- c. Add **100 µL** of **Nuclease-free water** to the plate
- d. Bead beat at **1200 rpm** for **30 seconds**
- e. Rest for **30 seconds**
- f. Bead beat at **1200 rpm** for **30 seconds**
 - i. This step will vary by tissue type and should be optimized for each tissue
- g. Spin sample at **3,000 rpm** for **3 min**
- h. Transfer **200 µL** of the **supernatant** of the **sample** to 2 mL 96-well plate
- i. Add **400 µL** of **Lysis (LBB)** to plate
- j. Add **30 µL** of **Proteinase K (PK)**
- k. **Mix** by pipetting up and down 10 times, or until thoroughly mixed.
- l. **Incubate** the plate for **30 minutes** at **37°C**

3. Bind

- a. Vortex to fully resuspend the **Bind (BBB)**
- b. Add **300 µL** of **Bind (BBB)** to the plate
- c. **Incubate** the plate for **5 minutes** at **room temperature**
- d. Place the plate on a **magnet** for 10 minutes (or until supernatant is clear)
- e. Remove and discard the supernatant without disrupting the beads
- f. Remove the plate from the magnet

4. Wash

- a. Add **800 µL** of Wash (**WBB**) to plate
- b. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- c. Place the plate on a **magnet** for **10 minutes** (or until supernatant is clear)
- d. Remove and discard the supernatant without disrupting the beads
- e. Remove the plate from the magnet
- f. Repeat steps 3.a-3.e for a total of **2 washes**
- g. Add **1.6 mL** of Wash (**WBC**) to plate
- h. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- i. Place the plate on a **magnet** for **6 minutes** (or until supernatant is clear)
- j. Remove and discard the supernatant without disrupting the beads
- k. Remove the plate from the magnet
- l. Repeat steps 3.g-3.k for a total of **2 washes**

5. Elute

- a. Add **40 µL** of **nuclease free water** to plate
- b. **Incubate** the plate for **2 minute** at **room temperature**
- c. Place the plate on a **magnet** for **2 minute** (or until supernatant is clear)
- d. Remove and **Save** the supernatant without disrupting the beads

Example Data

Data shown below resulted from the use of 16.6mg of mouse lung tissue. The three samples represent three technical replicates. DNA yield and purity was assessed on a NanoDrop (Thermo Fisher Scientific) (Table 1). Genomic DNA (gDNA) integrity was assessed on an Agilent Genomic DNA Screen Tape (Agilent) (Figure 1). The DIN scores, which represent the amount of gDNA degradation, were all above 8 indicating low levels of gDNA degradation.

Sample	Conc. (ng/μL)	Yield (μg)	Yield/mg of tissue (μg)	260/280	260/230
A	220.0	8.8	0.53	1.92	2.13
B	186.4	7.5	0.45	1.90	1.94
C	146.7	5.9	0.35	1.92	2.02

Table 1. The concentration, yield, yield per milligram of tissue, and DNA purity of DNA extracted from mouse lung tissue

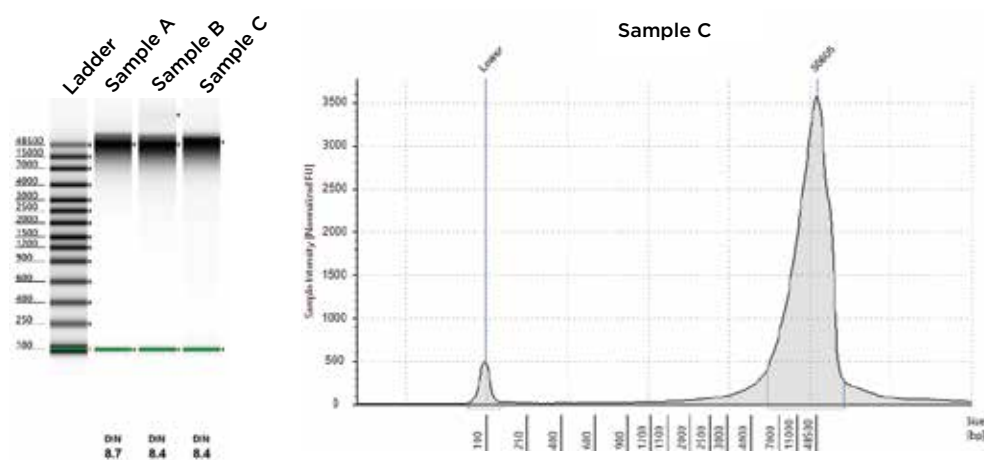


Figure 1. An Agilent Genomic DNA Screen Tape of DNA extracted from 16.6mg of mouse lung tissue. Right is the gel of all three samples and left is electropherogram of Sample A. The DIN scores are at the bottom of the gel.

Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions. This protocol is for demonstration only, and is not validated by Beckman Coulter.

© 2019 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.