

DNA extraction from cell culture using GenFind V3

Researchers working with cultured cell lines and primary cells and researchers working on genome editing technologies may use this protocol.

Please reference current GenFind V3 IFU for product information (Part Number C34880, C34881)

Purpose

This protocol provides a method for extracting DNA from cultured cells (≤ 2 million cells). Data shown shows the efficacy of this procedure with 1.4 million Jurkat cells, an immortalized line of human T lymphocyte cells.

Materials Used

Material	Part Number	Supplier
2 ml 96-well plate	609681	Beckman Coulter
PBS, pH 7.4	10010023	Thermo Fisher Scientific
Nuclease-free water (Molecular Grade)	AM9932	Thermo Fisher Scientific
7 Bar Magnet for 96-Well Plate	771MWZM-IALT	V&P Scientific
Microcentrifuge tubes 1.5 mL	357448	Beckman

Protocol

1. Lysis

- Add **200 μ L** of **1 x PBS** to fresh or frozen cell pellet
- Mix by pipetting up and down 10 times, or until thoroughly mixed
- Add **500 μ L** of **Lysis (LBB)**
- Add **30 μ L** of **Proteinase K (PK)**
- Mix** by pipetting up and down 10 times, or until thoroughly mixed
- Incubate** the plate for **15 minutes** at **37°C**

2. Bind

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

3. Wash

- Add **800 μ L** of **Wash (WBB)** to plate
- Mix** by pipetting up and down 10 times, or until thoroughly mixed
- Place the plate on a **magnet** for **10 minutes** (or until supernatant is clear)
- 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**

4. 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 isolated genomic DNA (gDNA) from 1.4 million Jurkat. DNA yield and purity was assessed by a NanoDrop (Thermo Fisher Scientific) (Table 1). The genomic integrity of the three replicates was assessed on an Agilent Genomic DNA Screen Tape (Agilent). The DIN scores, which represent the amount of gDNA degradation, were all above 8.4 indicating highly intact gDNA (Figure 1).

	Cell Number	Conc. (ng/µL)	Yield (µg)	A260/A280	A260/A230
Jurkat cells (Sample A)	1.4 x 10 ⁶	323.8	13.0	1.98	2.16
Jurkat cells (Sample B)	1.4 x 10 ⁶	319.9	12.8	1.97	2.15
Jurkat cells (Sample C)	1.4 x 10 ⁶	281.5	11.3	1.99	2.19
Average	1.4 x 10 ⁶	308.4	12.3	1.98	2.17

Table 1. Concentration and purity of DNA extracted from 1.4 million Jurkat cells. Total average DNA yield from 3 replicates was 12.3 µg. High DNA purity meets requirements for many downstream applications.

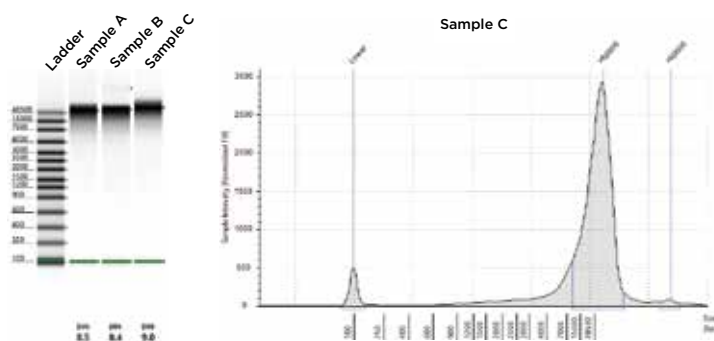


Figure 1. An Agilent Genomic DNA Screen Tape of DNA extracted from 1.4 million Jurkat cells. The gel indicates highly intact DNA predominately >48.5 kb with the DIN scores at the bottom. To the right of the gel is the electropherogram of Sample C.

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.