



# DNA Extraction from Filamentous Fungi using the GenFind V3 reagent kit

Researchers working on filamentous fungi e.g. *Aspergillus* and *Penicillium* may use this protocol. Please reference the current GenFind V3 protocol for product information. (Part Number C34880, C34881)

## Purpose

Filamentous fungi have tough cell walls that can make lysis and DNA extraction difficult. The most common method for lysing filamentous fungi is bead beating, which is labor intensive and not amenable to automation. This protocol provides two other methods for lysis that are more automation friendly and we show that they produce results that are very similar to the industry standard of bead beating.

Materials Used:

Material	Part Number	Supplier
7 Bar magnet for 96-well plate	771MWZM-1ALT	V&P Scientific
2 mL 96-well plate	609681	Beckman Coulter Life Sciences
Lysozyme from chicken egg white	L6876	Sigma-Aldrich
100% Ethanol (Molecular Grade)	AB00138	AmericanBio
2xYT	AB15063-01000	AmericanBio
PBS, pH 7.4	10010023	ThermoFisher Scientific
Nuclease-free water (Molecular Grade)	AM9932	ThermoFisher Scientific
0.5 mm glass beads	Z250465	Sigma-Aldrich

## Protocol

This GenFind V3 supplemental protocol provides sample DNA isolation with three different protocols. Protocol choice is dependent on workflow needs and the particular fungi strain(s) used.

### Fungal growth should be fresh

- Fungal growth can be done either on a plate or in an overnight culture of a rich media such as 2xYT
- If using fungal growth from a plate, combine fungi with 200  $\mu$ L of 1 x PBS and vortex to mix

### Isolation from up to 200 $\mu$ L of fresh culture using a freeze/thaw lysing method

#### 1. Lysis

- Transfer **200  $\mu$ L** of **fungal growth** to sample wells in a 2 mL 96-well plate
- Add **500  $\mu$ L** of **Lysis (LBB)** to the sample wells
- Add **30  $\mu$ L** of **Proteinase K** to each sample well
- Add **4  $\mu$ L** of **lysozyme (10mg/mL)** to each sample well
- Incubate** the plate for **10 minutes** at **-80°C**
- Mix** by pipetting up and down 10 times, or until thoroughly mixed
- Incubate** at **room temperature** for **5 minutes** or until samples are completely thawed
- Incubate** the samples for **30 minutes** at **37°C**

## 2. Bind

- a. Vortex the bottle of **Bind (BBB)** to fully resuspend the beads
- b. Add **300  $\mu$ L** of **Bind (BBB)** to each sample
- c. Add **100  $\mu$ L** of **80% ethanol** to each sample
- d. **Mix** by pipetting up and down 10 times, or until thoroughly mixed  
**Note: If this step is skipped not all of the beads will settle.**
- e. **Incubate** for **5 minutes** at **room temperature**
- f. Place the plate on a **magnet** for **15 minutes** (or until the supernatant is clear)
- g. Remove and discard the supernatant without disrupting the beads
- h. Remove the plate from the magnet

## 3. Wash

- a. Add **800  $\mu$ L** of **Wash (WBB)** to each sample well
- 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 the 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 **1600  $\mu$ L** of **Wash (WBC)** to each sample well
- 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 the 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  $\mu$ L** of **Nuclease-free water** to each sample well
- b. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- c. **Incubate** for **2 minutes** at **room temperature**
- d. Place the plate on a **magnet** for **2 minutes** (or until the supernatant is clear)
- e. Remove and **Save** the supernatant without disrupting the beads

## Isolation from up to 200 $\mu\text{L}$ of fresh culture by lysing at 65°C for 1 hour

### 1. Lysis

- a. Transfer **200  $\mu\text{L}$**  of **fungal growth** to a 2 mL 96-well plate
- b. Add **500  $\mu\text{L}$**  of **Lysis (LBB)** to each sample well
- c. Add **30  $\mu\text{L}$**  of **Proteinase K** to each sample well
- d. Add **4  $\mu\text{L}$**  of **lysozyme (10mg/mL)** each sample well
- e. **Incubate** the plate for **1 hour** at **65°C**
- f. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- g. **Incubate** the plate for **1 hour** at **65°C**

### 2. Bind

- a. Vortex the bottle of **Bind (BBB)** to fully resuspend the beads
- b. Add **300  $\mu\text{L}$**  of **Bind (BBB)** to each sample well
- c. Add **100  $\mu\text{L}$**  of **80% ethanol** to each sample well
- d. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- e. **Note: If this step is skipped not all of the beads will settle.**
- f. **Incubate** for **5 minutes** at **room temperature**
- g. Place the plate on a **magnet** for **15 minutes** (or until supernatant is clear)
- h. Remove and discard the supernatant without disrupting the beads
- i. Remove the plate from the magnet

### 3. Wash

- a. Add **800  $\mu\text{L}$**  of **Wash (WBB)** to each sample well
- 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 **1600  $\mu\text{L}$**  of **Wash (WBC)** to the sample
- 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 the 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  $\mu\text{L}$**  of **Nuclease-free water** to each sample well
- b. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- c. **Incubate** for **2 minutes** at **room temperature**
- d. Place the plate on a **magnet** for **2 minutes** (or until the supernatant is clear)
- e. Remove and **Save** the supernatant without disrupting the beads

## Isolation from up to 200 $\mu\text{L}$ of fresh culture using Bead Beating

### 1. Lysis

- a. Add 0.5 g of 0.5 mm glass beads to each sample well of a 2 mL 96-well plate
- b. Transfer **200  $\mu\text{L}$**  of **fungal growth** to each sample well in the plate
- c. Add **100  $\mu\text{L}$**  of **Lysis (LBB)** to each sample well
- d. Add **100  $\mu\text{L}$**  of **Nuclease-free water** to each sample well
- e. Bead beat at **max speed** ( $\geq 1200\text{rpm}$ ) for **3 minutes**
- f. Let the beads settle and transfer the supernatant to new wells of a 2 mL 96-well plate
- g. Add **30  $\mu\text{L}$**  of **Proteinase K** to each sample well
- h. Add **4  $\mu\text{L}$**  of **lysozyme (10mg/mL)** to each sample well
- i. Add **400  $\mu\text{L}$**  of **Lysis (LBB)** to each sample well
- j. Remove the plate from the magnet
- k. **Incubate** the samples for **30 minutes** at **37°C**

### 2. Bind

- a. Vortex the bottle of **Bind (BBB)** to fully resuspend the beads
- b. Add **300  $\mu\text{L}$**  of **Bind (BBB)** to each sample well
- c. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- d. **Incubate** for **5 minutes** at **room temperature**
- e. Place the plate on a **magnet** for **15 minutes** (or until the supernatant is clear)
- f. Remove and discard the supernatant without disrupting the beads

### 3. Wash

- a. Add **800  $\mu\text{L}$**  of **Wash (WBB)** to each sample well
- 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 the 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 **1600  $\mu\text{L}$**  of **Wash (WBC)** to each sample well
- 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 the supernatant is clear)
- j. Remove and discard the supernatant without disrupting the beads
- k. Remove the plate from magnet
- l. Repeat steps 3.g-3.k for a total of **2 washes**

### 4. Elute

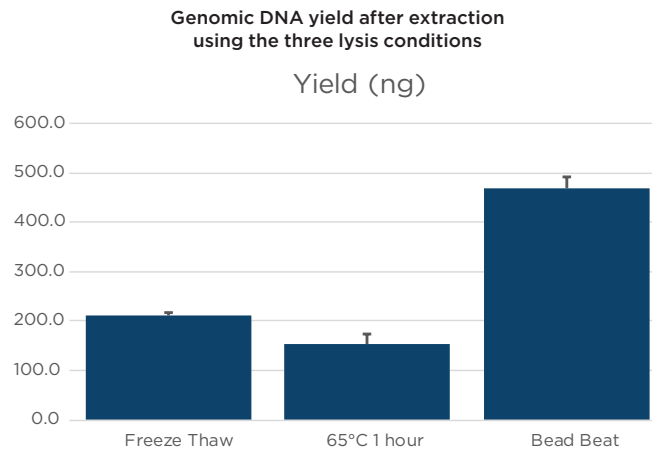
- a. Add **40  $\mu\text{L}$**  of **Nuclease-free water** to each sample well
- b. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- c. **Incubate** for **2 minutes** at **room temperature**
- d. Place the plate on a **magnet** for **2 minutes** (or until the supernatant is clear)
- e. Remove and **Save** the supernatant without disrupting the beads

## Example Data

### Genomic DNA extraction from *A. niger*

Data shown below was generated using *Aspergillus niger* that was streaked onto a YT plate and grown for 2 weeks and kept at 4°C for 2 months. The fungal growth and spores were scraped off the plate into 1.5 mL of 1xPBS and thoroughly vortexed. The fungal PBS mixture was distributed evenly among the three treatment groups.

Bead beating resulted in the highest yield of DNA; this is most likely due to the large amounts of spores that were used in this study. The other two treatment groups, while resulting in about half the total yield, are more permissible to high-throughput applications.



**Figure 1.** The average yield of DNA from *A. niger* using one of the three lysis conditions. Error bars represent the standard deviation of three technical replicates.

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.

© 2024 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.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://beckman.com)  
2023-GBL-EN-102497-v2