



DNA extraction from Zeesan@ Saliva DNA sample collection kit using GenFind V3

This method is applicable for scientists who want to extract DNA from Zeesan@ Saliva DNA sample collection kit (SAL-2000L).

Please reference the current GenFind V3 protocol (Part number: C34880 or C34881) for product information.

Purpose

Saliva contains buccal epithelial cells and white blood cells which provide rich genetic data. Due to the easy access and non-invasive character, saliva is an alternative to blood collection. Here we present a high-quality, DNA extraction method from saliva using GenFind V3.

Material Used

Material	Part Number	Supplier
GenFind V3	C34880 or C34881	Beckman Coulter
Proteinase K	C34821 or C34827	Beckman Coulter
Lysis (LBB)	C34822	Beckman Coulter
Bind (BBB)	C34823	Beckman Coulter
96 Square Deep-Well Plate, Polypropylene 2 mL	609681	Beckman Coulter
Magnetic Separation Plate for 96 Deep Well 7 Magnetic Bars	VP-771MWZM-1-ALT	V&P
Nuclease-Free Water	AM9932	Ambion
Ethanol	AB-00138	American Bioanalytical

Protocol

1. Preparation/Lysis

- a. Collect saliva according to manufacturer's instructions
- b. Keep saliva at **room temperature** before DNA extraction
- c. Transfer **300 µL of** saliva to a 2 mL 96-well plate
- d. Add **700 µL LBB** to the plate
- e. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- f. Add **45 µL PK** to the plate
- g. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- h. Incubate the plate for **60 minutes** at **55°C**

2. Bind

- a. Vortex the bottle of BBB to fully resuspend the beads
- b. Add **450 µL** of **BBB** to the plate
- c. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- d. **Incubate** the plate for **5 minutes** at **room temperature**
- e. Place the plate on a **magnet** for **10 minutes** (or until supernatant is clear)
- f. Remove and discard the supernatant without disrupting the beads
- g. Remove the plate from the magnet

3. WBB Wash

- a. Add **1.2 mL** of **WBB** to the 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 the supernatant is clear)
- d. Remove and discard the supernatant without disrupting the beads
- e. Remove the plate from the magnet
- f. Repeat step 3a-e for a total of **2 washes**

4. WBC Wash

- a. Add **1.6 mL** of **Wash buffer WBC** to the 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 the supernatant is clear)
- d. Remove and discard the supernatant without disrupting the beads
- e. Remove the plate from the magnet
- f. Repeat step 4a-e for a total of **2 washes**
- g. Air dry the samples on the magnet for **3 minutes**

5. Elute

- a. Add **150 µL** of **nuclease free water** to the plate
- b. **Mix** by pipetting up and down 10 times, or until thoroughly mixed
- c. **Incubate** for **5 minutes** at **room temperature**
- d. Place the plate on a **magnet** for **4 minutes** (or until supernatant is clear)
- e. Remove and **save** the supernatant without disrupting the beads

Results

Genomic DNA (gDNA) was isolated from three healthy donors using Zeesan@ DNA sample collection kit. DNA yield was measured by Quant-iT™ PicoGreen™ dsDNA Assay (ThermoFisher). An average of 10.3 µg of gDNA was extracted from 300 µL saliva (Fig. 1A). Purity was assessed by NanoDrop (Thermo Fisher Scientific). The average 260/280 is 1.9 and the average 260/230 is greater than 2.0 (Fig. 1B). DNA integrity was assessed using Agilent gDNA ScreenTape Assay. The DIN scores averaged 8.2 and DNA peak size was larger than 55K indicating the gDNA extracted from saliva sample is of high quality (Figure 1C and D).

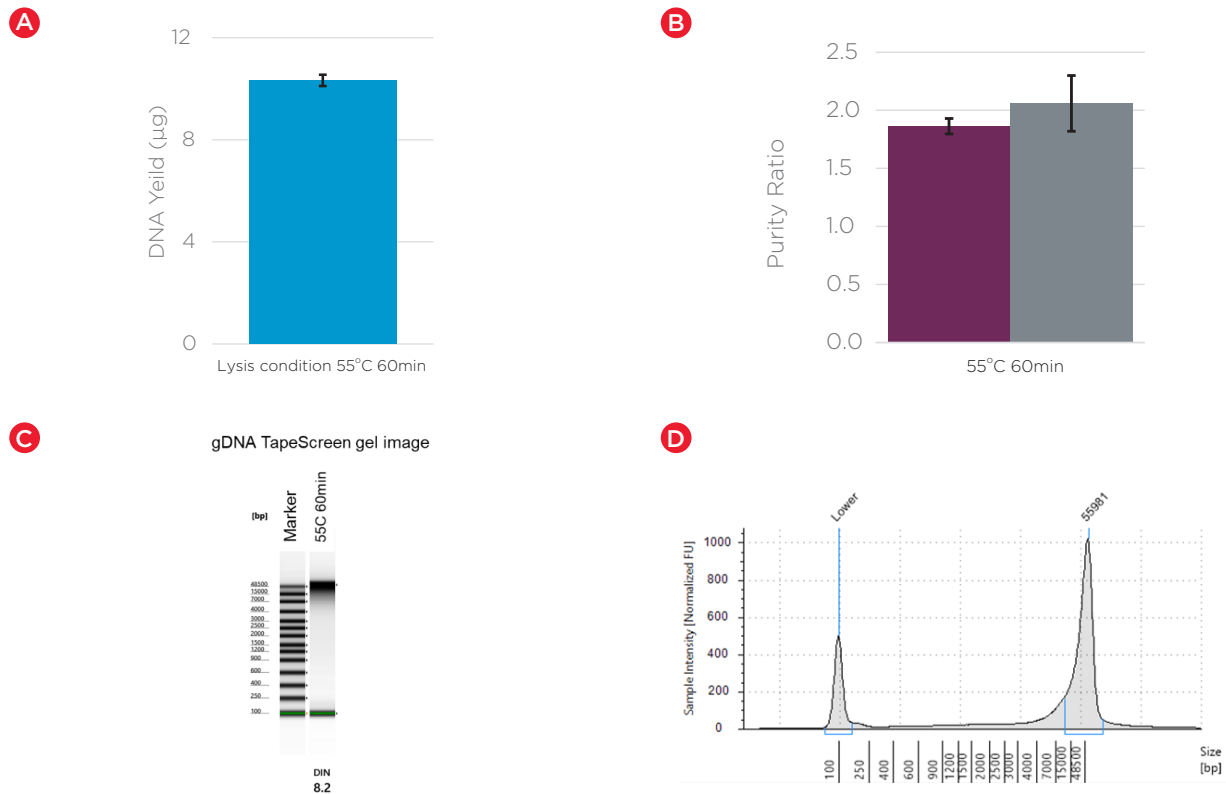


Figure 1. Genomic DNA extraction from 300 µL saliva. DNA was isolated from 3 healthy donors. (A) DNA yield was measured by Quant-iT™ PicoGreen™ dsDNA Assay (ThermoFisher). The error bar represents the standard deviation of three donor replicates. (B) Purity was accessed by NanoDrop (Thermo Fisher Scientific); the average 260/280 is about 1.9 and 260/230 is greater than 2.0. The error bar represents the standard deviation of three donor replicates. (C and D) DNA integrity was analyzed by using Agilent gDNA ScreenTape Assay.

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.