



Tissue Sample Preparation

1. For each sample, label a 5 cm petri dish and place on wet ice, along with a labeled 15 mL conical tube containing 5 mL ice-cold PBS.
2. Transfer each sample to its corresponding dish, mince with a razor blade, and transfer to the corresponding 15 mL conical tube containing ice-cold PBS using a 1 mL pipette tip (the tip can be cut to widen the bore in order to avoid clogging the tip).
3. Centrifuge tubes at 500 x *g* for 5 minutes at 4°C. Remove supernatant.
4. Add 1 mL ATAC Lysis Buffer and using a 1 mL wide bore pipette tip, transfer each sample to a 1 mL dounce homogenizer. Slowly dounce for 30 strokes.
5. Filter each sample through a 40 µm mesh strainer and collect in a fresh 1.5 mL microcentrifuge tube.
6. Immediately after filtration, take a 10 µL aliquot for cell counting using Trypan Blue cell-viability staining. Aliquot 50,000-100,000 cells into a new tube.
7. Centrifuge tubes at 500 x *g* at 4°C for 5 minutes. During this time prepare **Tagmentation Master Mix**, see table below. After centrifuge, remove supernatant and proceed to the Tagmentation Reaction and purification steps.

Cell Sample Preparation

1. Count cells and aliquot 50,000 to 100,000 cells into a fresh 1.5 mL centrifuge tube for each sample.
2. Centrifuge cells at 500 x *g* for 5 minutes at 4°C. Remove supernatant. If there is no pellet, do additional centrifuge at 1000 x *g* for 5 minutes at 4°C.
3. Add 100 µL cold PBS. Centrifuge at 500 x *g* at 4°C. Remove supernatant.
4. Resuspend pellet in 100 µL ice-cold ATAC Lysis Buffer. Centrifuge at 500 x *g* at 4°C for 10 minutes.
5. Transfer sample to a PCR tube on ice. Centrifuge at 500 x *g* for 10 minutes at 4°C. During this time prepare **Tagmentation Master Mix**. After centrifuge, remove supernatant and move directly to tagmentation.

Tagmentation Master Mix (Per sample)

Reagents	Volume
2X Tagmentation Buffer	25 µL
10X PBS	2 µL
1.0% Digitonin	0.5 µL
10% Tween 20	0.5 µL
H ₂ O	12 µL
Assembled Transposomes	10 µL

Tagmentation Reaction and Purification

1. Add 50 µL of tagmentation master mix to each sample, pipetter to resuspend. Incubate at 37°C for 30 minutes in a thermomixer at 800 rpm.
2. Immediately following the tagmentation reaction, transfer each sample to a clean 1.5
3. Add 250 µL DNA Purification Binding Buffer and 5 µL 3 M sodium acetate to each sample.
4. If the color of the sample is anything other than bright yellow, add additional 3 M sodium acetate in 5 µL increments until the proper color is achieved.

5. Mix each sample and transfer to a labeled DNA purification column. Centrifuge columns at 17,000 x *g* for 1 minute.
6. Discard the flow-through, return the collection tube to the column, then add 750 μ L of wash buffer. Centrifuge at 17,000 x *g* for 1 minute.

Note: 100% ethanol must be added to the Wash Buffer before the first use, to a final concentration of 80%. (40 mL of 100% Ethanol to the DNA Purification Wash Buffer bottle)

7. Discard the flow-through. With the column cap open, centrifuge at 17,000 x *g* for 2 minutes to remove residual Wash Buffer from the column.
8. Add 35 μ L of DNA purification elution buffer to the center of the column matrix. Incubate at RT for 1 minute. Centrifuge column at 17,000 x *g* for 1 minute.
9. Purified DNA can be stored at -20°C, or you may proceed directly with the PCR Amplification of Tagmented DNA steps that follow.

PCR Amplification of Tagmented DNA

PCR Reaction Master Mix (Per sample)

Reagents	Volume
Tagmented DNA	33.5 μ L
i7 Indexed Primer (25 μ M)	2.5 μ L
i5 Indexed Primer (25 μ M)	2.5 μ L
dNTPs (10 mM)	1 μ L
5X Q5 Reaction Buffer	10 μ L
Q5 Polymerase (2 U/ μ L)	0.5 μ L

1. Set up the PCR reactions by adding the components in the order shown above. If libraries are to be multiplexed for sequencing on the same flow cell, ensure that a unique i5 and/or i7 index combination is used for each.
2. Perform PCR using the following program on a thermal cycler (with a heated lid):
 - 72°C 5 minutes
 - 98°C for 30 seconds
 - 10 cycles of: 98°C for 10 seconds, 63°C for 30 seconds, 72°C for 1 minute
 - Hold at 10°C.
3. Perform SPRI clean-up with 60 μ L SPRI bead solution (1.2X the sample volume), eluting in 20 μ L DNA Purification Elution Buffer. Have 400 μ L of freshly-prepared 80% ethanol ready per sample:
 - A. Add 60 μ L well-mixed, RT SPRI Beads to each sample.
 - B. Vortex briefly to mix and incubate for 5 minutes at RT to allow beads to bind. Apply magnet to collect beads.
 - C. Once the solution is clear, aspirate the supernatant.
 - D. Perform two washes, each with 180 μ L of 80% ethanol, leaving the beads on the magnet for both washes.
 - E. Allow tubes to sit at RT so that residual ethanol can evaporate. Once the beads transition from shiny to matte (2-5 minutes), proceed to the next step.
 - F. Once beads are dry, add 20 μ L of DNA purification Elution Buffer and vortex. Leave at RT for 5 minutes. Apply magnet to collect beads.
 - G. Once the solution is clear, transfer each supernatant containing the eluted DNA to a fresh tube.
4. At this stage, libraries are ready for quantification and sequencing. Use a library quantification kit for next-generation sequencing to quantify the library (e.g. Kapa Biosystems, Catalog No. KR0405). PCR amplified libraries can also be analyzed to assess size distribution with a Bioanalyzer, TapeStation, or similar instrument to assess size distribution.