

CUTANA™ CUT&RUN Library Prep Kit

QUICK-START CARD



Scan for
full manual -
read before
first use

Section I: End Repair (~75 min)

1. Thaw kit reagents stored at -20°C and keep on ice during experiment. Mix **End Prep Buffer**; if a white precipitate is observed, bring to room temperature (RT), vortex to dissolve, and place on ice.
2. Transfer 5 ng CUT&RUN-enriched DNA to fresh **8-strip tubes**. Adjust to 25 μL with **0.1X TE Buffer**.
3. Prepare an End Repair Master Mix in a new 1.5 mL tube on ice. Per library prep reaction, combine 4.2 μL (●) **End Prep Buffer** and 1.8 μL (●) **End Prep Enzyme**. Gently vortex to mix, quick spin to collect liquid, and return to ice. This recipe includes excess volume to account for pipetting error.
4. Add 5 μL **End Repair Master Mix** to 25 μL CUT&RUN DNA in 8-strip tubes. Pipette up and down 5X to clear tips, gently vortex to mix, and quick spin.
5. Place reactions in a thermocycler with heated lid set to $\geq 75^{\circ}\text{C}$. Run the program outlined in the table.
6. Quick spin tubes and place on ice. **Note:** It is critical to keep tubes on ice during adapter ligation (below).

| STEP # | TEMP | TIME | CYCLES |
|--------|------------------------|----------|--------|
| 1 | 20°C | 20 min | 1 |
| 2 | 65°C | 30 min | 1 |
| 3 | $4-12^{\circ}\text{C}$ | ∞ | 1 |

Section II: Adapter Ligation and U-Excision (~45 min)

7. Prepare a Ligation Master Mix in a new 1.5 mL tube on ice. Per reaction, combine 16.5 μL (●) **Ligation Mix** and 0.55 μL (●) **Ligation Enhancer**. Gently vortex, quick spin, and return to ice. This recipe includes excess volume to account for pipetting error.
8. Add 1.25 μL (●) **Adapter for Illumina®** and 15.5 μL **Ligation Master Mix** to each reaction, keeping tubes on ice. Vortex tubes thoroughly to mix, quick spin, and return to ice.
9. Place tubes in a thermocycler without a heated lid, with block set to 20°C . Incubate 15 min.
10. Quick spin tubes and place in a RT rack. Add 1 μL /reaction (●) **U-Excision Enzyme**. Pipette up and down 3X to clear tips, gently vortex to mix, and quick spin.
11. Place tubes in a thermocycler with a heated lid set to $\geq 47^{\circ}\text{C}$ and block set to 37°C . Incubate 15 min.
12. Quick spin tubes. Continue to next step or store at -20°C for future processing.

Section III: DNA Cleanup (~15 min)

***NOTE:** Use of multi-channel pipettors is recommended

13. Prepare 1 mL 85% Ethanol (EtOH) per reaction by combining 850 μL 100% EtOH and 150 μL molecular biology grade water. Prepare **fresh**, mix well, and store at RT.
14. Vortex **SPRIselect** reagent (manufactured by Beckman Coulter, Inc.) to thoroughly resuspend beads. Slowly add 47.75 μL /reaction. Ensure pipette tip is free of extra bead droplets.
15. **Mix well** by pipetting and/or vortexing to resuspend. Quick spin to collect liquid. Incubate 5 min at RT.
16. Place tubes on magnet for 2-5 min. Pipette to remove supernatant without disturbing beads.
17. **Keep tubes on magnet.** Add 180 μL /reaction **85% EtOH**. Remove supernatant. Repeat one time.
18. Quick spin with caps facing in to avoid dislodging beads. Return to magnet and remove residual EtOH.
19. Remove tubes from magnet. Air-dry, caps open, 2-3 min at RT. Beads should appear damp matte brown (Figure 1). If beads are crackly/light brown, they are too dry.
20. Add 12 μL /reaction **0.1X TE Buffer** to elute DNA. Pipette/vortex to resuspend. Incubate 2 min, RT.
21. Place tubes on magnet for 2 min at RT. Transfer 10.5 μL eluted DNA to new **8-strip tubes**. Continue or store at -20°C .

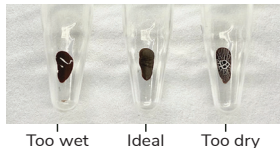


FIGURE 1
Elute DNA when "ideal."

Section IV: Indexing PCR (~30 min)

***NOTE:** Before first use, transfer entire volume of (o) **i5 primers** to a new set of **8-strip tubes**.

- Assign a unique pair of (o) **i5** and (•) **i7 primers** to each reaction, following the primer selection instructions in **Appendix 2**. Mark the combination as consumed in the **Primer Tracking Table** (below).
- Mix the stock tubes containing **Hot Start 2X PCR Master Mix** and (o) **i5** and (•) **i7 primers** and quick spin. To each library prep reaction, add the following reagents individually and in order: 1 μL (•) **i7 primer**, 1 μL (o) **i5 primer**, and 12.5 μL (•) **Hot Start 2X PCR Master Mix**.

Change tips between each addition to avoid contamination. Visually inspect tips to confirm the correct volume was aspirated.

- Mix well by vortexing tubes, avoid bubbles, and quick spin. Place tubes in a thermocycler with a heated lid set to 105°C. Perform PCR using parameters outlined in the table.

| STEP # | TEMP | TIME | CYCLES | NOTES |
|--------|------|----------|--------|----------------------------|
| 1 | 98°C | 45 sec | 1 | Hot start activation |
| 2 | 98°C | 15 sec | 14 | DNA melting |
| 3 | 60°C | 10 sec | | Hybrid annealing/extension |
| 4 | 72°C | 60 sec | 1 | Final extension |
| 5 | 4°C | ∞ | 1 | Hold temperature |

Section V: PCR Cleanup (~15 min)

- Vortex **SPRIselect** reagent to thoroughly resuspend beads. Slowly add 25 μL **SPRIselect** reagent to each indexing PCR reaction. Ensure pipette tip is free of extra bead droplets.
- Proceed with library cleanup following steps 15-21 in **Section III**. The protocol generates 10.5 μL purified CUT&RUN sequencing libraries.

Section VI: Analysis of Library Fragment Size (~1 hr)

- Quantify libraries using the Qubit™ fluorometer. Examine fragment size distribution on the Agilent TapeStation® or Bioanalyzer®. Proceed to sequencing or store at -20°C. Obtain library concentration (200-700 bp range) and confirm fragment size distribution (~300 bp).
- Store prepared CUT&RUN sequencing libraries at -20°C.

Primer Tracking Table

Mark consumed primer combinations in the table.

***IMPORTANT:** Sequencing runs must contain libraries with a sufficient mixture of i5 and i7 indexes to ensure proper diversity for Illumina® systems. At least one i5 Sequential Primer Set (grey dashed lines) AND one i7 Sequential Primer Set (red dashed lines) should be represented per sequencing run. See **Appendix 2** for guidance. For index sequences, visit epicypher.com/14-1001 and download the Excel spreadsheet under Documents and Resources.

Included with both 14-1001 & 14-1002



| | | | | | | | | | |
|--|--|--|--|--|--|--|--|--|--|
| | | | | | | | | | |
| | | | | | | | | | |
| | | | | | | | | | |
| | | | | | | | | | |
| | | | | | | | | | |
| | | | | | | | | | |

Included with 14-1001 Included with 14-1002

i7 Sequential Primer Sets (i7 Sets) **i5 Sequential Primer Sets (i5 Sets)**

*Beckman Coulter, the stylized logo, and SPRIselect are trademarks or registered trademarks of Beckman Coulter, Inc.

