

CUTANA™ CUT&RUN Library Prep Kit

QUICK-START CARD



Scan for
full manual

Section I: End Repair (~75 min)

1. Transfer 5 ng CUT&RUN-enriched DNA to an **8-strip tube**. Adjust volume to 25 μL with **0.1X TE Buffer**.
2. Prepare an **End Repair Master Mix** in a 1.5 mL tube on ice. Per reaction, combine 4.2 μL (●) **End Prep Buffer** and 1.8 μL (●) **End Prep Enzyme**. Gently vortex to mix, quick spin, and return to ice.
3. Add 5 μL **End Repair Master Mix** per reaction. Pipette to clear tips, gently vortex to mix, and quick spin.
4. Place reactions in a thermocycler with heated lid set to $\geq 75^\circ\text{C}$ and run the following program:

STEP #	BLOCK TEMP	TIME	CYCLES	NOTES
1	20°C	20 minutes	1	Reaction temperature
2	65°C	30 minutes	1	Enzyme inactivation
3	4°C	∞		Hold temperature

5. Quick spin 8-strip tubes and place directly on ice or in a pre-chilled aluminum block on ice.

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

6. Prepare a **Ligation Master Mix** in a 1.5 mL tube on ice. Per reaction, combine 16.5 μL (●) **Ligation Mix** and 0.55 μL (●) **Ligation Enhancer**. Gently vortex (setting #7), quick spin, and return to ice.
7. To 8-strip tubes on ice, add 1.25 μL /reaction (●) **Adapter for Illumina®** and 15.5 μL /reaction **Ligation Master Mix**. Vortex tubes thoroughly to mix, quick spin, and return to ice.
8. Incubate 15 min in a thermocycler set to 20°C, without a heated lid.
9. Place tubes in a room temperature (RT) rack and add 1 μL /reaction (●) **U-Excision Enzyme**. Pipette up and down to clear tips, gently vortex to mix, and quick spin.
10. Incubate 15 min in a thermocycler set to 37°C, with heated lid at $\geq 47^\circ\text{C}$.
11. 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 from this point forward.

12. Make 900 μL **FRESH** 85% Ethanol (EtOH) per reaction; if pausing after **Section III**, make 450 μL /reaction.
13. Vortex **SPRIselect** reagent (manufactured by Beckman Coulter, Inc.*) thoroughly to resuspend.
14. Slowly add 47.75 μL /reaction **SPRIselect** reagent in 8-strip tubes.
15. Gently vortex tubes and quick spin. Incubate 5 min at RT.
16. Place 8-strip tubes on a magnet for 2 min at RT. Remove supernatant.
17. Keeping tubes on magnet, add 180 μL /reaction 85% EtOH. Remove supernatant.
18. Repeat the previous step one time.
19. Remove tubes from magnet, quick spin, and return to magnet. Remove residual supernatant.
20. Remove tubes from magnet, leave caps open, and air dry beads for 2 min.
21. Add 12 μL /reaction **0.1X TE Buffer** to elute target DNA.
22. Gently vortex to resuspend beads and quick spin. Incubate 2 min at RT.
23. 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 for future processing.



Section IV: Indexing PCR (~30 min)

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

24. 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).
25. To each 10.5 μL reaction add the following reagents individually and in order. Visually inspect tips to confirm that the correct volume was aspirated and change tips between each addition to avoid cross-contamination.
- 1 μL assigned (•) **i7 primer**
 - 1 μL assigned (o) **i5 primer**
 - 12.5 μL (•) **Hot Start 2X PCR Master Mix** (mix well before using)
26. Vortex to mix, then quick spin. Run PCR using the following parameters, with heated lid set to 105°C:

STEP #	BLOCK 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	∞		Hold temperature

Section V: PCR Cleanup (~15 min)

27. Follow **Section III: DNA Cleanup** using 25 μL /reaction **SPRIselect** reagent. Results in 10.5 μL of each library.

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

28. Use 1 μL of each CUT&RUN library for quantification on the Qubit™ fluorometer.
29. Prepare 5 μL of each library at 10 ng/ μL . Run 1 μL on the Agilent BioAnalyzer® or Tapestation® for quality analysis. Obtain library concentration (200-700 bp range) and confirm fragment size distribution (~300 bp).
30. Store prepared CUT&RUN sequencing libraries at -20°C.

Primer Tracking Table

Mark consumed primer combinations in the table.

Included with both 14-1001 & 14-1002



***IMPORTANT:** Each sequencing run must contain at least one i5 Sequential Primer Set (grey dashed lines) AND one i7 Sequential Primer Set (red dashed lines) to ensure proper index diversity for Illumina® systems. See **Appendix 2** for a Primer Selection Guide. For index sequences, visit epicypher.com/14-1001 and download the Excel spreadsheet under Documents and Resources.

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.