Kit v3 - Manual v3.0

CUTANA[™] CUT&TAG Kit

QUICK-START CARD



DAY 1 (~2 HOURS)

1. Prepare buffers as outlined below. Recipes contain 20% excess - no overage is needed.

BUFFER	COMPONENTS	1X	8X	16X	STORAGE
Nuclei Extraction Buffer	Pre-Nuclei Extraction Buffer	235 µL	1.9 mL	3.8 mL	lce for use on Day 1
	25X Protease Inhibitor	9.8 µL	78.4 μL	157 µL	
	1 M Spermidine	0.13 µL	1.0 µL	2.0 µL	
Wash Buffer 1	Pre-Wash Buffer	1.3 mL	10.4 mL	20.8 mL	4°C for use on Day 2
	25X Protease Inhibitor	56 µL	448 µL	896 µL	
	1 M Spermidine	0.7 µL	5.6 µL	11.2 µL	
	5% Digitonin	2.8 µL	22.4 µL	44.8 µL	
Wash Buffer 2	Wash Buffer 1	600 µL	4.8 mL	9.6 mL	4°C for use on Day 2
	4.5 M NaCl	20.7 µL	166 µL	331 µL	
Antibody Buffer	Wash Buffer 1	60 µL	480 µL	960 µL	lce for use on Day 1
	0.5 M EDTA	0.25 µL	2 µL	4 µL	

- 2. Resuspend ConA Beads and transfer 11 μ L/reaction to a 1.5 mL tube. Place tube on a magnetic rack, allow slurry to clear, and pipette to remove supernatant.
- 3. Resuspend beads in 100 μ L/reaction cold **Bead Activation Buffer**. Return to magnet, allow slurry to clear, and remove supernatant. Repeat one time.
- 4. Resuspend beads in 11 μ L/reaction cold **Bead Activation Buffer**. Aliquot 10 μ L/reaction activated ConA beads into **8-strip Tubes**. Keep on ice.
- 5. Harvest 100,000 cells/reaction. Spin 600 x g, 3 min, room temperature (RT). **High-quality** sample prep is crucial see kit manual or <u>support.epicypher.com</u> for guidance!
- 6. Remove supernatant and resuspend cells in 100 μ L/reaction cold Nuclei Extraction Buffer.
- 7. Incubate 10 min on ice. Spin 600 x g for 5 min at 4°C. Pipette to remove supernatant.
- 8. Resuspend nuclei in 105 μ L/reaction cold **Nuclei Extraction Buffer**. Add 100 μ L nuclei to 10 μ L ConA beads in 8-strip tubes. Gently vortex to mix and quick spin to collect liquid.
- 9. Incubate 10 min at RT. Place tubes on magnet, allow slurry to clear, and remove supernatant.
- 10. Resuspend in 50 μ L/reaction cold Antibody Buffer.
- 11. Quick spin the K-MetStat Panel tube and mix by pipetting (do NOT vortex stock). Add K-MetStat Panel to reactions designated for H3K4me3, H3K27me3, and IgG Control Antibodies. Add 2 μL if using 100,000 nuclei/reaction. For lower nuclei numbers, decrease K-MetStat Panel per manual instructions. Gently vortex tubes to mix and quick spin.
- Add 0.5 μg primary antibody to each reaction. For control reactions, add 1 μL of respective H3K27me3, H3K4me3, or IgG Control Antibody. Gently vortex and quick spin.
- Incubate overnight on a nutator at 4°C, gently rocking tubes with caps elevated. Do NOT rotate tubes, as this will result in sample loss.



DAY 2 (~7 HOURS)

Mixing, steps 17-25: Carefully pipette, avoid bead loss in tip, expel all material back into tubes.

Before and after all incubation steps: Gently vortex tubes ~5 sec and quick spin to collect liquid.

- 14. Prepare Tagmentation Buffer in a new 1.5 mL tube. Per reaction, combine 59.4 μL Wash Buffer 2 and 0.6 μL 1 M MgCl₂ (10 mM final). Place on ice. Recipe includes 20% excess.
- 15. For step 25, transfer 60 µL/reaction Pre-Wash Buffer to a new tube and place at RT.
- 16. Quick spin reaction tubes, place on a magnet, and allow slurry to clear. Remove supernatant.
- 17. Resuspend in 50 μL/reaction cold **Wash Buffer 1**. Add 0.5 μg/reaction secondary antibody. Use 0.5 μL **Anti-Rabbit Secondary Antibody** for control and rabbit primary antibodies.
- 18. Incubate 30 min on a nutator at RT. Place on magnet, allow slurry to clear, remove supernatant.
- 19. <u>On magnet</u>, add 200 µL/reaction cold **Wash Buffer 1**. Remove supernatant. Repeat one time.
- 20. Resuspend in 50 µL/reaction cold Wash Buffer 2. Add 2.5 µL/reaction pAG-Tn5.
- 21. Incubate 1 hr on a nutator at RT. Place on magnet, allow slurry to clear, remove supernatant.
- 22. Resuspend in 200 µL/reaction cold **Wash Buffer 2**. Place on a magnet, allow slurry to clear, remove supernatant. Repeat one time.
- 23. Resuspend in 50 µL/reaction cold Tagmentation Buffer.
- 24. Incubate 1 hr in thermocycler set to 37°C (lid at 47°C). Place on magnet, allow slurry to clear, remove supernatant.
- 25. Resuspend in 50 μL/reaction RT **Pre-Wash Buffer**. Do **NOT** vortex. Place tubes on magnet, allow slurry to clear. Remove supernatant.
- 26. Add 5 μ L/reaction RT **SDS Release Buffer**. Do **NOT** pipette. Vortex ~10 sec and quick spin.
- 27. Incubate 1 hr in a thermocycler set to 58°C (lid at 68°C).
- 28. Add 15 µL/reaction RT SDS Quench Buffer, carefully pipetting to rinse beads (vortex if viscous). Vortex ~10 sec max speed to mix, quick spin, and keep at RT.
- Assign a unique pair of i5 & i7 Primers per reaction (Appendix 1). To the entire reaction add: 2 μL i5 Primer, 2 μL i7 Primer, and 25 μL Non-Hot Start 2X PCR Master Mix. <u>Mix well</u>, avoid bubbles, and quick spin.
- 30. Perform PCR per table parameters (lid at 105°C). During PCR, prep 500 µL/reaction 85% EtOH.
- 31. Quick spin tubes. Vortex **SPRIselect reagent** (manufactured by Beckman Coulter, Inc.) to fully resuspend and slowly add 65 μL/reaction. <u>Mix well</u> by vortexing and/or pipetting.
- 32. Incubate 5 min at RT. Place tubes on magnet for 2-5 min. Remove supernatant.
- 33. On magnet, add 180 $\mu\text{L/reaction}$ 85% EtOH. Remove supernatant. Repeat one time.
- 34. Quick spin tubes with caps facing in. Return to magnet and pipette to remove residual EtOH.
- 35. Take tubes off magnet. Air-dry, caps open, 2-3 min at RT. Beads should be damp matte brown.
- 36. Add 17 µL/reaction **0.1X TE Buffer**. Pipette/vortex to resuspend beads. Incubate 2 min at RT.
- 37. Quick spin tubes, place on magnet for 2 min. Transfer 15 μ L libraries to new **8-strip Tubes**.
- 38. Quantify libraries and examine fragment size distribution per manual instructions. Proceed to sequencing or store at -20°C.

Beckman Coulter, the stylized logo, and SPRIselect are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

STEP # TEMP TIME 58°C 5 min 1 1 2 72°C 5 min 1 3 98°C 45 sec 1 4 98°C 15 sec 14-21 10 sec 5 60°C 6 72°C 1 min 1 7 4-12°C ω 1

