

User Guide

Catalog No. CHP2

CHP2 Experienced User Protocol

sigma-aldrich.com/life-science/epigenetics/chip2.html

SIGMA-ALDRICH®

SIGMA® Where *bio* begins™
Life Science

CHP2 Experienced User Protocol

This protocol has been written for the experienced **CHP2 kit** user who has read & successfully performed ChIPs following the procedure as described in the CHP2 kit user guide. This abbreviated procedure is written for the advanced (repeat) user who has all the required buffers reconstituted & is familiar with the pre-requisites for each section.

Index:

O/N - overnight

H - hour(s)

RT - room temperature

min - minute(s)

PIC – protease inhibitor cocktail

1. Blocking ChIP Next Gen Seq Sepharose®:

- Rehydrate each vial of ChIP Next Gen Seq Sepharose with 200 μ L of water & transfer 100 μ L to 0.5 mL tube, store remaining 100 μ L ChIP Next Gen Seq Sepharose at 4 °C & use within 1 month
- Add 5 μ L BSA (**A7638**, 20 mg/mL), 1 μ L 0.1 M AEBSF & 10 μ L salmon sperm DNA (omit DNA for ChIP-Seq)
- Rotate end-over-end O/N at 4 °C (or 2 H minimum at RT)
- Transfer to 1.5 mL tube
- Spin 14K RCF at 4 °C, 3 min
- Discard supernatant, resuspend pellet with 1 mL Dialysis buffer & spin 14K RCF at 4 °C, 3 min
- Discard supernatant & repeat wash as above
- Resuspend pellet in 100 μ L Dialysis buffer & add 1 μ L 0.1 M AEBSF
- Store at 4 °C & use within 2 weeks

2. Cross-Link Cells:

- Grow healthy cell cultures in 500 cm² dishes
- Remove media, rinse cells with 35 mL HEPES Buffered Salt Solution
- Add 9 mL Trypsin-EDTA, incubate 37 °C, 5 min, dislodge attached cells
- Add 31 mL complete medium / 500 cm² dish & resuspend cells, transfer to 50 mL tube
- Add 1.2 mL 37 % Formaldehyde/ 40 mL cell suspension
- Rotate end-over-end exactly 10 min at RT
- Add 4.1 mL cold 1.25 M Glycine/ 40 mL cell suspension
- Rotate end-over-end exactly 5 min at RT
- Spin 200 RCF at 4 °C, 10 min
- Discard supernatant, resuspend pellet with 50 mL cold 1 \times Phosphate Buffered Saline
- Spin 200 RCF at 4 °C, 10 min
- Prepare chromatin or snap freeze cell pellet in liquid N₂ & store at -80 °C

Complete Cell Lysis Buffer:

Cell Lysis Buffer	0.9 mL
+ 10 % IGEPAL	0.1 mL
+ PIC (P8340)	10 μ L
+ 0.1 M AEBSF	10 μ L
Total	1 mL

Complete Nuclei Lysis Buffer:

Nuclei Lysis Buffer	0.5 mL
+ PIC (P8340)	5 μ L
+ 0.1 M AEBSF	5 μ L
Total	0.5 mL

IP Dilution Buffer:

IP Dilution Buffer	1 mL
+ PIC (P8340)	10 μ L
+ 0.1 M AEBSF	10 μ L
Total	1 mL

3. Chromatin Prep:

- Resuspend cross-linked cells in fresh Complete Cell Lysis Buffer (1 mL/ 50 \times 10⁶ cells)
- Chill on ice, 20 min
- Dounce 1 mL aliquots in 2 mL glass Dounce homogenizer with 15 strokes of B pestle, transfer lysate to 1.5 mL tubes
- Spin 2.5 K RCF at 4 $^{\circ}$ C, 5 min
- Discard supernatant, resuspend nuclei pellet in fresh Complete Nuclei Lysis Buffer (0.5 mL/ 50 \times 10⁶ cells)
- Chill on ice, 10 min
- Sonicate nuclear lysate using pre-determined optimum conditions for each cell line & instrument. (For BioRuptor: Transfer to 15 mL polystyrene tubes, 30s ON + 30s OFF, 3 \times 10 min for ChIP-q-PCR & 4-5 \times 10 min for ChIP-Seq experiments)
- Transfer to 1.5 mL tubes
- Spin 14 K RCF at 4 $^{\circ}$ C, 10 min
- Transfer supernatant to new 1.5 mL tube(s) & aliquot
- Proceed for ChIP or snap freeze in liquid N₂ & store at -80 $^{\circ}$ C

4. Pre-clearing:

- Pre-clear the required amount of chromatin with blocked ChIP Next Gen Seq Sepharose, 10 μ L per 10⁷ cells
- Rotate end-over-end exactly 15 min at 4 $^{\circ}$ C
- Spin 14 K RCF at 4 $^{\circ}$ C, 5 min
- Transfer supernatant (pre-cleared chromatin) to new 1.5 mL tube(s), 125 μ L/ tube

5. Ab Binding:

- Remove 12.5 μ L chromatin for "10 % Input" & keep aside on ice
- Dilute 125 μ L chromatin with 375 μ L IP Dilution buffer (+ protease inhibitors) + Ab (specific/control)
 - For positive control ChIP: Add 0.5 μ L Mu IgG (**I5381**)
 - For negative control ChIP: Add 1 μ L RNAP II Ab (**WH0005430M1**)
- Rotate end-over-end O/N at 4 $^{\circ}$ C (or for abundant target rotate end-over-end for 2 H at 4 $^{\circ}$ C)

6. Bridging Ab & IP:

- Add 1 μ L Rb anti-Mouse Ab (**M7023**) for ChIPs with monoclonal antibodies (or appropriate bridging antibody)
- Rotate end-over-end 1 H at 4 $^{\circ}$ C
- Add 10 μ L blocked ChIP Next Gen Seq Sepharose (for chromatin from > 2 \times 10⁶ cells) or 5 μ L blocked ChIP Next Gen Seq Sepharose (for chromatin from \leq 2 \times 10⁶ cells)
- Transfer to new 1.5 mL tube
- Spin 14K RCF at 4 $^{\circ}$ C, 4 min

Dialysis Wash Buffer:

Dialysis Wash Buffer	1 mL
+ 0.1 M AEBSF	10 μ L
Total	1 mL

IP Wash Buffer:

IP Wash Buffer	1 mL
+ 0.1 M AEBSF	10 μ L
Total	1 mL

IP Elution buffer:

Water	425 μ L
+ 10 % SDS	50 μ L
+ 1 M NaHCO ₃	25 μ L
Total	500 μL

7. Washings:

- Discard supernatant, resuspend pellet with fresh 1 mL Dialysis buffer (+ AEBSF), spin 14 K RCF at 4 °C, 3 min
- Discard supernatant & repeat wash as above
- Discard supernatant, resuspend pellet with fresh 1 mL IP Wash buffer (+ AEBSF), spin 14 K RCF at 4 °C, 3 min
- Discard supernatant & repeat wash three more times as above
- Remove last traces of IP Wash buffer with a quick spin 14 K RCF at 4 °C, 1 min

8. ChIP Elution:

- Add 50 μ L fresh IP Elution buffer to each ChIP Next Gen Seq Sepharose pellet
- Incubate RT 15 min with vortexing on moderate setting (3 on Vortex Genie)
- Spin 14K RCF at RT, 3 min
- Transfer supernatant to new 1.5 mL tube
- Repeat Elution: Add 50 μ L fresh IP Elution buffer to each ChIP Next Gen Seq Sepharose pellet
- Incubate RT 15 min with vortexing on moderate setting (3 on Vortex Genie)
- Spin 14K RCF at RT, 3 min
- Transfer & pool supernatants in same 1.5 mL tube

9. Cross-Link Reversal:

- Spin 14K RCF at RT, 5 min to remove residual ChIP Next Gen Seq Sepharose
- Transfer supernatant to new 1.5 mL tube
- To "10 % Input" sample add 87.5 μ L IP Elution buffer & 1 μ L Proteinase K (P4850) & incubate 65 °C, 1 H
- Add 12 μ L 5 M NaCl to all samples (ChIPs & Input)
- Rotate end-over-end O/N at 67 °C (or boil for 15 min for ChIP-q-PCR only)

10. DNA Purification:

- Chill on ice 5 min, thaw by hand until clear
- Add 1 µL RNase A
- Incubate 37 °C, 15 min
- Assemble one set of DNA purification column in collection tube for each ChIP & Input sample
- Add 500 µL DNA Binding buffer to each prep, vortex, transfer to column
- Spin 12K RCF at RT, 1 min
- Discard flow through, wash column with 200 µL DNA Wash buffer
- Spin 12K RCF at RT, 1 min
- Discard flow through, wash column again with 200 µL DNA Wash buffer
- Spin 12K RCF at RT, 1 min
- Transfer column to fresh 1.5 mL tube
- Add 30 µL DNA Elution buffer for ChIP-q-PCR (or add 6 µL Elution buffer for ChIP-chip or ChIP-Seq)
- Spin 12K RCF at RT, 1 min
- Use purified DNA for PCR, q-PCR, WGA, labeling, quantitation or sequencing

11. Q-PCR Assay:

- q-PCR setup for Actin & ZNF-333-3' (per reaction):
 - 9.5 µL Water
 - 1.0 µL 5 µM Primers mix
 - 12.5 µL 2X PCR Ready-mix (component of [S4438](#))
 - 2.0 µL ChIP'ed DNA or 1 % Input or water
 - 25 µL Total volume
- Thermocycling conditions:

Denature	95 °C	2 min
40 cycles of	95 °C	30 sec
	60 °C	30 sec
	72 °C	1 min, read
- Use the ChIP Q-PCR Data analysis excel shell (<http://www.sigmaaldrich.com/life-science/epigenetics/chip2.html>) to calculate ChIP yield i.e. “% Input” & “Fold Enrichment” with respect to non-specific Ab

For more information, visit
sigma-aldrich.com/life-science/epigenetics/chip2.html

World Headquarters
3050 Spruce St., St. Louis, MO 63103
(314) 771-5765
sigma-aldrich.com

Order/Customer Service (800) 325-3010 • Fax (800) 325-5052
Technical Service (800) 325-5832 • sigma-aldrich.com/techservice
Development/Bulk Manufacturing Inquiries (800) 244-1173

©2013 Sigma-Aldrich Co. All rights reserved. SIGMA, SAF-C, SIGMA-ALDRICH, ALDRICH, FLUKA, SUPELCO and Imprint are trademarks belonging to Sigma-Aldrich Co. LLC. Bioruptor is a trademark of Diagenode SA. Nanodrop is a trademark of Thermo scientific. Sepharose is a registered trademark of GE Healthcare. Sigma brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or