

# Re-shearing of decrosslinked immunoprecipitated DNA using the Bioruptor<sup>®</sup> Pico

## INTRODUCTION

ChIP-seq is a common tool used to study protein-DNA interactions in cells and tissues on a genome-wide scale. The size of the immunoprecipitated DNA fragments is a key parameter for successful library preparation and reliable ChIP-seq data. Re-shearing of the purified decrosslinked DNA after immunoprecipitation and before library preparation enable enrichment of fragments in the desired optimal size range suitable for next-generation sequencing (Figure 1). This approach eliminates a size-selection step from a conventional library prep workflow and increases the yield of immunoprecipitated DNA available for a library. It improves the sensitivity and/or resolution of standard

ChIP-seq experiments, especially for inactive histone marks (Laczik et al, 2016). The chromatin that represents inactive marks is more resistant to sonication. As a consequence, these regions are relatively more present among the longer fragments after the initial shearing step, and re-shearing shifts these fragments to smaller sizes which are more efficiently incorporated into sequencing libraries. ChIP-seq on active histone marks, transcriptional factors or chromatin modifying proteins benefits from the re-shearing approach by increasing the number of fragments available for library generation for sequencing and permitting reduction of the starting amount of cells.

## REQUIRED MATERIALS AND EQUIPMENT

- Bioruptor<sup>®</sup> Pico (Diagenode, Cat. No. B01060001)
- 0.2ml tube holder (Diagenode, Cat. No. B01200044) and corresponding 0.2 ml microtubes for Bioruptor<sup>®</sup> Pico (Diagenode, Cat. No. C30010020) or 0.1 ml tubes holder (Diagenode, Cat. No. B01200041) and corresponding 0.1 ml microtubes for Bioruptor<sup>®</sup> Pico (Diagenode, Cat. No. C30010015)
- Library preparation Kit (e.g. MicroPlex Kit, Diagenode, Cat. No. C05010012 or iDeal Library Preparation Kit, Diagenode, Cat. No. C05010020).
- (Optional) Fragment Analyzer<sup>™</sup> (Advanced Analytical) or BioAnalyzer (Agilent) with high sensitivity reagents

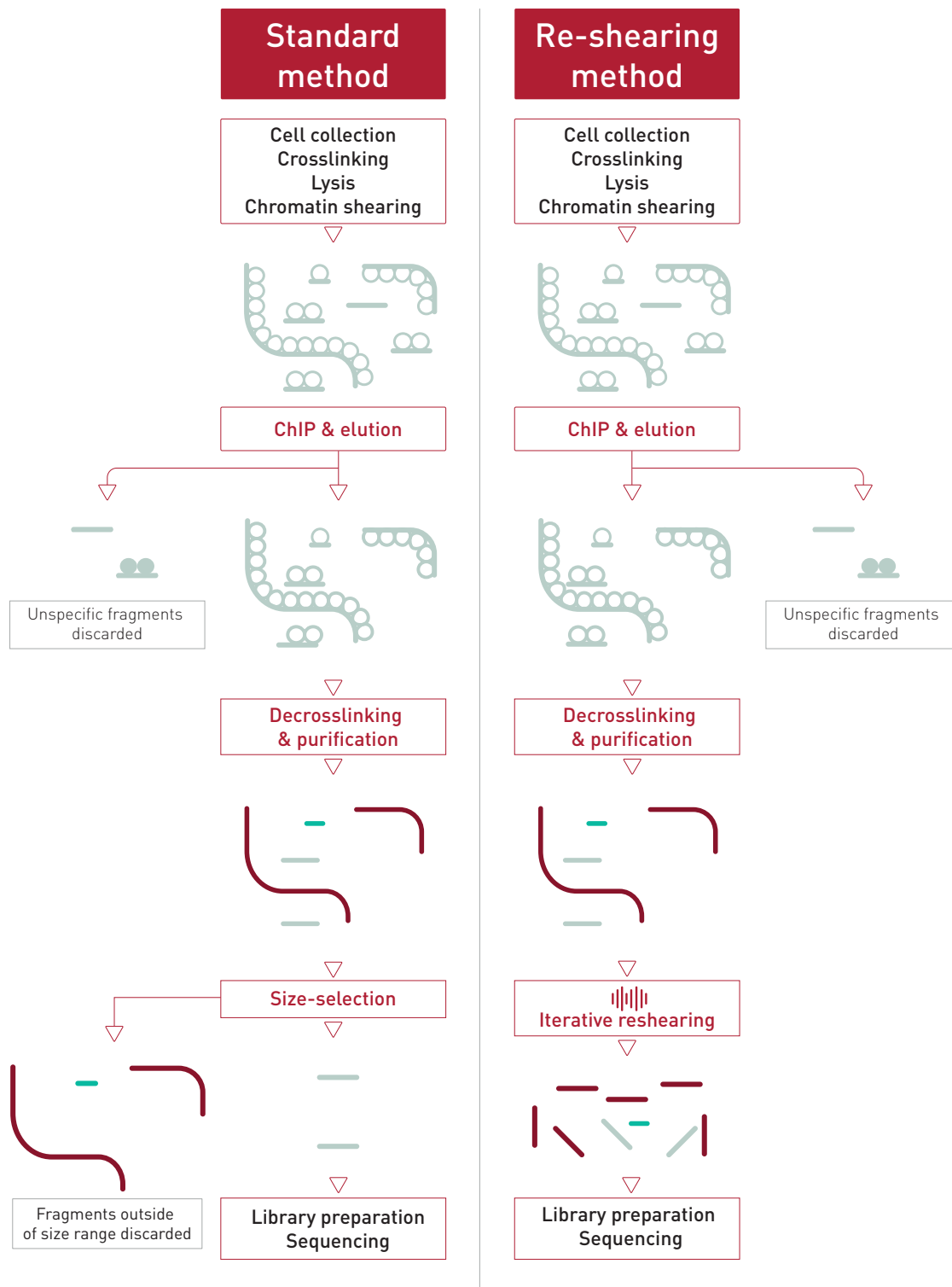


Figure 1. An overview of standard method (left) and re-shearing method (right)

## PROTOCOL

1. Perform chromatin immunoprecipitation using an appropriate protocol (e.g. Diagenode iDeal ChIP-seq Kit for Histones (Cat. No. C01010051) or iDeal ChIP-seq Kit for Transcription Factors (Cat. No. C01010055) and ChIP-seq grade antibodies ([www.diagenode.com/categories/antibodies](http://www.diagenode.com/categories/antibodies)). Ensure that an expected enrichment is achieved.
2. Transfer 20  $\mu$ l of immunoprecipitated DNA to Bioruptor<sup>®</sup> microtubes (0.2 ml or 0.1 ml).
3. Vortex and spin down samples and proceed to sonication in the Bioruptor<sup>®</sup> Pico with the following settings:

Sonication cycle: 30 sec ON/30 sec OFF

Total sonication time: 5 -10 cycles

Temperature: 4°C

*Note: To guarantee homogeneity of sonication, the tube holder should be always completely filled with tubes.*

4. (Optional) if DNA concentration is sufficient (at least 2-3 ng/ $\mu$ l), DNA fragment size distribution before preparing libraries can be assessed on the Fragment Analyzer<sup>™</sup> (Advanced Analytical) using the DNF-474 High Sensitivity NGS Fragment Analysis Kit or BioAnalyzer (Agilent) using High Sensitivity DNA Analysis kit.
5. Proceed to the library preparation using re-sheared DNA.

*Note: A size selection step can be skipped.*

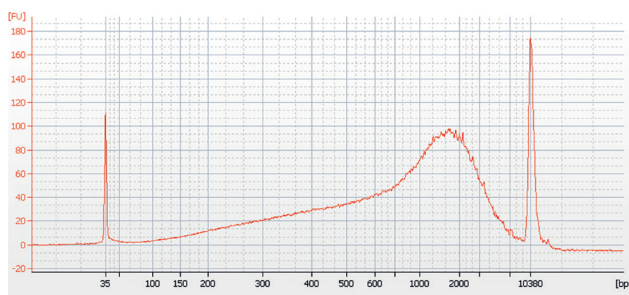


Figure A

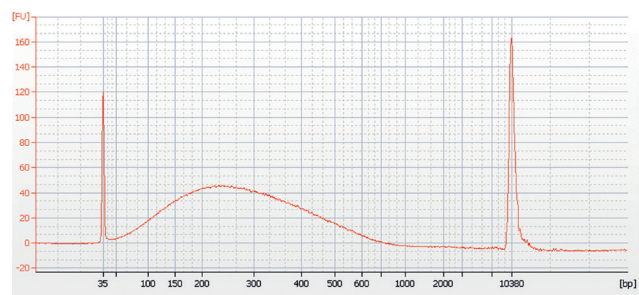


Figure B

**Figure 2. The re-shearing method enriches fragments in an optimal size range suitable for next-generation sequencing.** Shearing profile of immunoprecipitated chromatin before (Figure A) and after (Figure B) re-shearing. Chromatin from HeLa cells was prepared according to the iDeal ChIP-seq Kit for Histones and sheared using the Bioruptor<sup>®</sup> Pico for 8 cycles. ChIP was performed with an H3K27me3 antibody (Diagenode, Cat. No. C15410195). De-crosslinked and purified DNA was analysed on a BioAnalyzer. The observed fragment range is compatible with ChIP-seq library preparation.

## LIST OF DIAGENODE ANTIBODIES SUCCESSFULLY TESTED USING THE RE-SHEARING METHOD:

| Product      | Cat. No.           |
|--------------|--------------------|
| H3K27me3     | Cat. No. C15410195 |
| H3K4me3      | Cat. No. C15410195 |
| H3K4me1      | Cat. No. C15410194 |
| p53          | Cat. No. C15410083 |
| H3K27me3S28p | Cat. No. C15410091 |
| H4K20me3     | Cat. No. C15410057 |
| HDAC2        | Cat. No. C15200201 |
| EZH2         | Cat. No. C15200180 |
| LSD1         | Cat. No. C15410028 |
| CTCF         | Cat. No. C15410210 |

| Product  | Cat. No.           |
|----------|--------------------|
| NFkB     | Cat. No. C15310255 |
| E2F6     | Cat. No. C15410314 |
| RNF2     | Cat. No. C15410313 |
| OCT4     | Cat. No. C15410305 |
| H3K36me3 | Cat. No. C15410058 |
| NFYB     | Cat. No. C15410241 |
| KAP1     | Cat. No. C15410236 |
| H3K9me3  | Cat. No. C15410193 |
| CBX8     | Cat. No. C15410333 |

## REFERENCES

Laczik et al. Iterative Fragmentation Improves the Detection of ChIP-seq Peaks for Inactive Histone Marks. *Bioinformatics and Biology Insights* 2016:10 209–224 doi: 10.4137/BBI.S40628

Mokry M, Hatzis P, de Bruijn E, Koster J, Versteeg R, Schuijers J, et al. (2010) Efficient Double Fragmentation ChIP-seq Provides Nucleotide Resolution Protein-DNA Binding Profiles. *PLoS ONE* 5(11): e15092. doi:10.1371/journal.pone.0015092