



Innovating Epigenetics Solutions



WELCOME TO DIAGENODE

ChIP Workshop

Jessica Apulei & Juri Kazakevych

8-9 December 2020

OBJECTIVES

- CHIP overview
 - Chip-qPCR vs. ChIP-Seq
- Chromatin preparation
 - Fixation, Cell lysis and Chromatin shearing
- Setting up IP
 - Antibodies, Replicates, inputs, controls
- ChIP-qPCR
- ChIP-seq: library prep & sequencing
- ChIP-seq: analysis
- Overview of Alternate methods

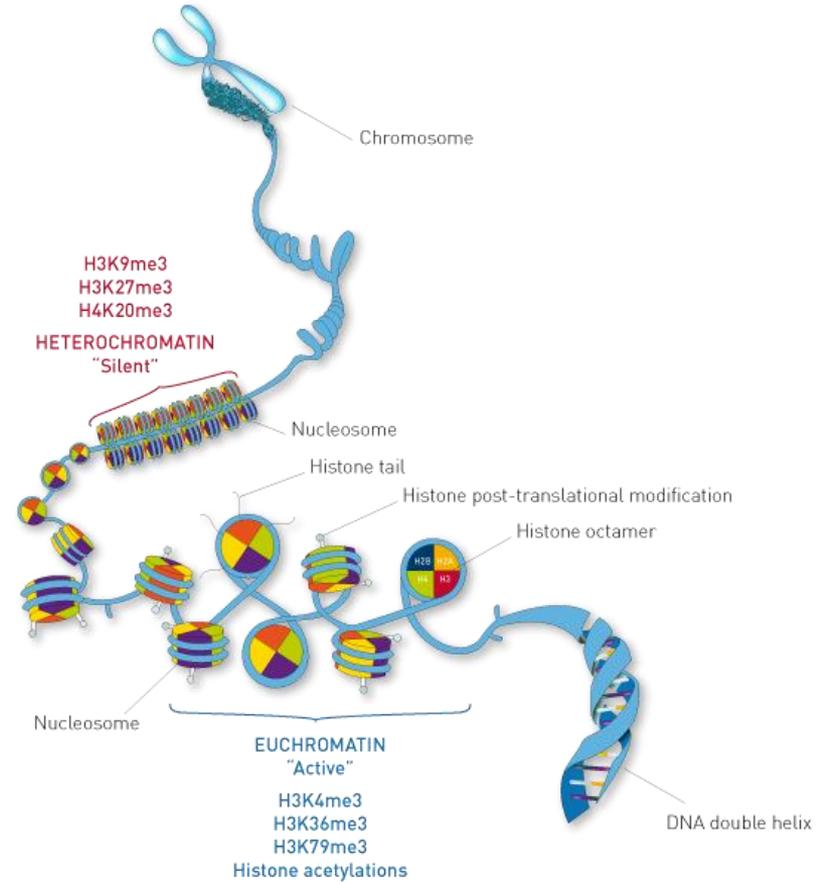




INTRODUCTION

What is Chromatin?

A complex of DNA and proteins found in eukaryotic cells

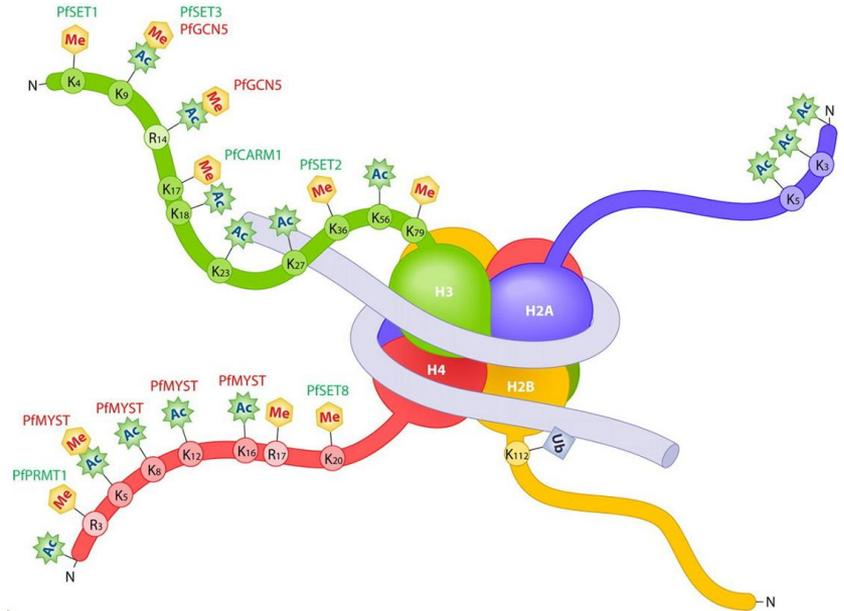




INTRODUCTION

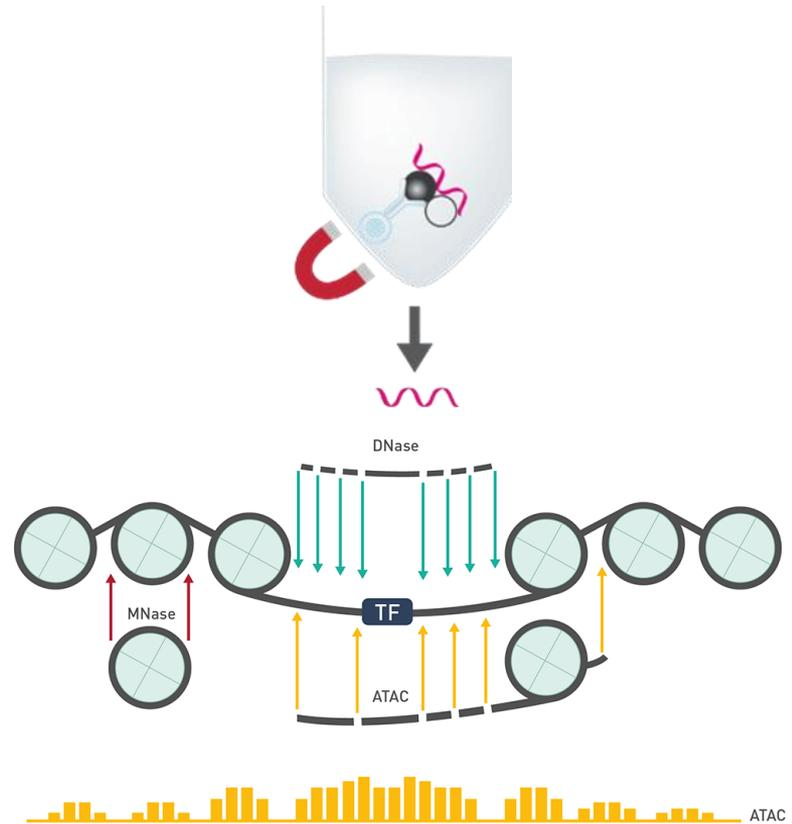
What is Chromatin?

- Identifying genome-wide DNA binding sites for histones, transcription factors and other proteins
- Defines transcription factor binding sites
- Reveals gene regulatory networks in combination with RNA sequencing and methylation analysis



CHROMATIN ANALYSIS

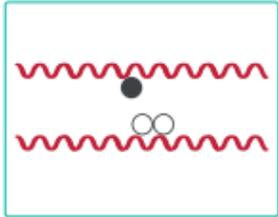
- Interaction between proteins and DNA (immuno- assays)
 - ChIP-qPCR
 - ChIP-Seq
 - Cut&Run and Cut&TAG
- Methods to study chromatin accessibility (non-immuno assays)
 - ATAC-seq





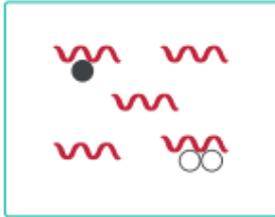
Workflow: Chromatin ImmunoPrecipitation (ChIP):

Step 1



Cross link to fix proteins to DNA

Step 2



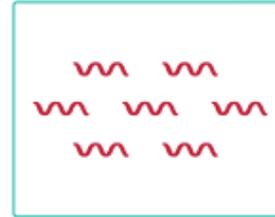
Shear chromatin

Step 3



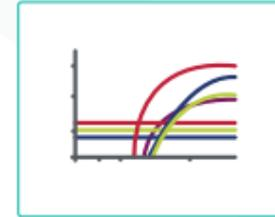
Immunoprecipitate with antibody and magnetic beads

Step 4



Reverse crosslinks and purify

Step 5



Analyze by qPCR

Step 6



Prepare the libraries for NGS

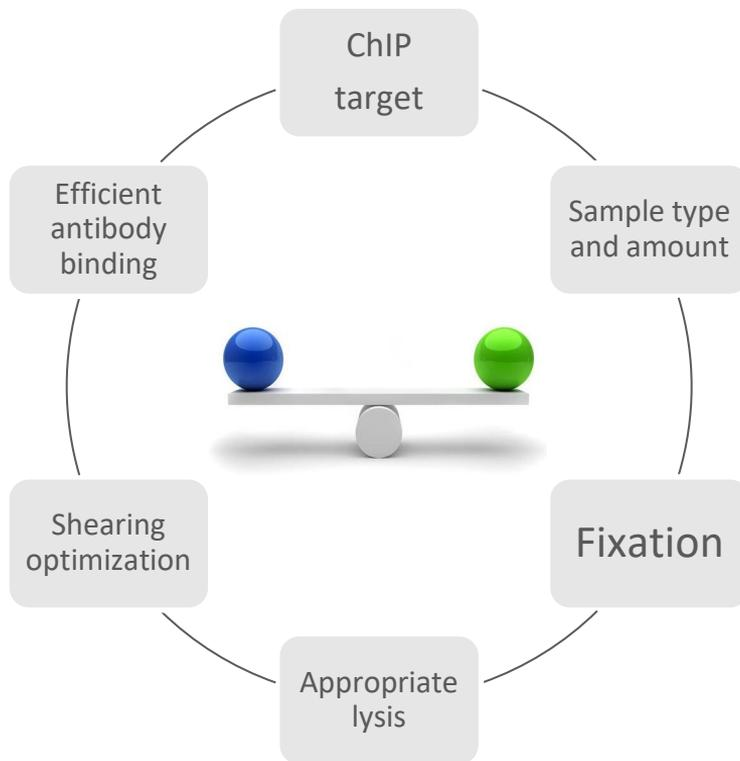


ChIP-qPCR or ChIP-Seq?

ChIP-qPCR	ChIP-Seq
<p>Single-locus data</p> <p>QC step for ChIP-seq</p>	<p>Genome-wide data</p> <p>suitable for exploratory analysis</p>
<p>Low-cost</p> <p>Fast</p>	<p>High sequencing costs</p> <p>Longer protocol</p> <p>High sequencing turnaround time</p>



Summary – Tips to Prepare Good Chromatin



Guide for successful chromatin preparation using the **Bioruptor® Pico**



Starting material: Cells and Tissues

- **Cells**

- ChIP: 1 million/IP for histones, 4 million/IP for TF (less depending on histone/TF)
 - low-input ChIP: 10k/IP for histones

- **Tissues**

- Amount – 20-30 mg/IP

- Dounce homogenization for soft tissues (e.g. liver or brain)

- Bead beater like TissueLyser for hard fibrous frozen tissues (e.g. muscles)

- **FFPE tissue**

- Challenging due to extensive crosslinking

- Heptane instead of xylene for de-paraffinization -> easier, non-toxic workflow



Fixation

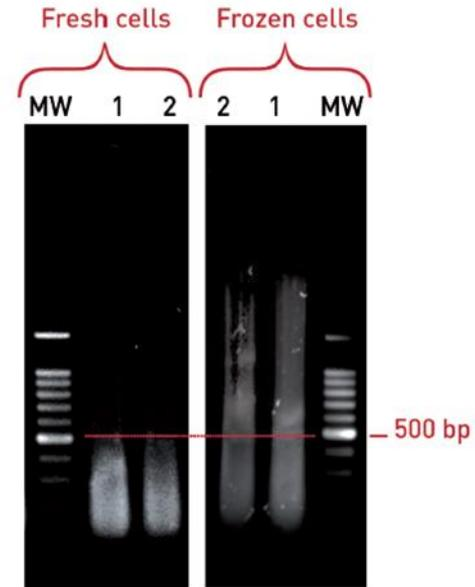
- Covalent stabilization of protein-DNA interactions; Reversible
 - Directly in medium for weak or rare protein-DNA interaction
 - For histone marks, cells can be resuspended by trypsinization before fixation
- Common fixative: Formaldehyde
 - Fresh
 - Methanol-free not mandatory

Target	Fixator	Formaldehyde	ChIP Cross-link Gold C01019021
Histones		Yes (8-10 min)	No need
Transcriptional factors directly bound to DNA		Yes (10-20 min)	No need
Indirect higher order and/or dynamic interactions		Yes (10-15 min)	Yes (30-45 min)



Cell Lysis

- Two step lysis – standard protocol, difficult cells
 - Remove soluble cytosolic proteins first
 - Improves sonication efficiency
 - Reduces background
- One step lysis – for low cell numbers
 - Lyse cells directly with an SDS-containing buffer
- Tips/Tricks/Critical steps:
 - Incubate on ice to start lysis and get narrower fragments size
 - Centrifuge to remove soluble membranes and cytosol
 - Avoid freezing chromatin if possible





Stopping Points

- **Cells**
 - Fix cells, lyse, isolate & shear chromatin -> freeze
 - Fix cells, lyse, isolate chromatin -> freeze
 - Fix cells -> freeze
- **Tissues**
 - Fix tissue, lyse, isolate & shear chromatin -> freeze
 - Freeze prior fixation

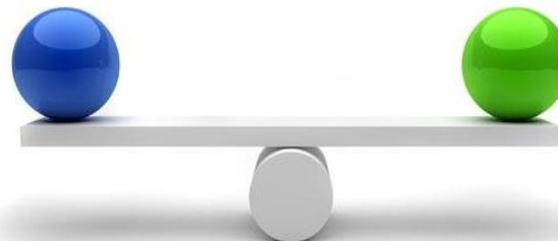


Secrets of ChIP Success

- Prepare “good” chromatin
 - Suitable fragment size and available epitopes
- Use a good antibody at the right concentration
- Optimize for highest specific signal and the lowest background

Extracted and fragmented chromatin

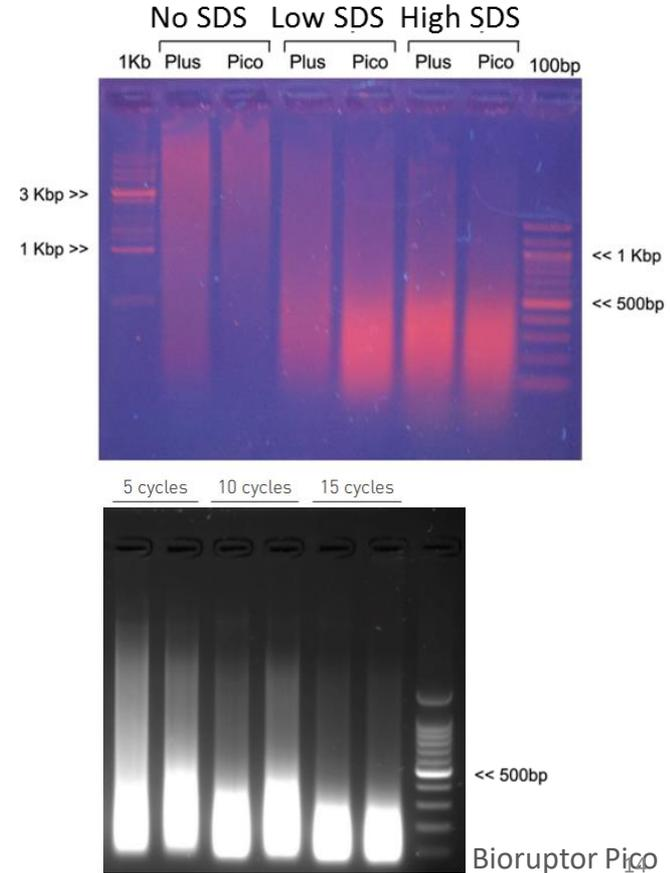
Protein integrity





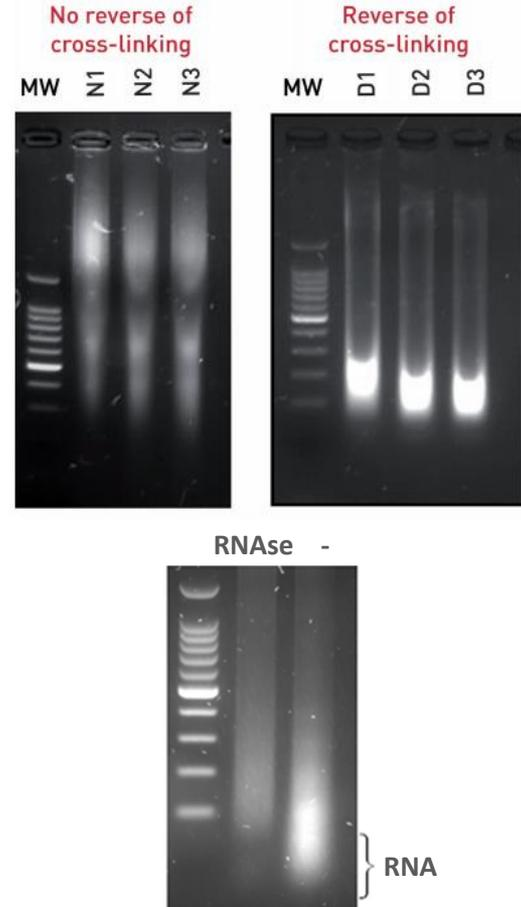
Chromatin Shearing

- 100-800 bp fragments, peak 200-500bp
- Use a good sonicator
 - Gentle - not to dislodge protein
 - Uniform and reproducible energy
 - Temperature control at 4°C
 - Multiplex and easy to use
- Shearing buffer with detergents, preferably SDS
 - Increase sonication efficiency and chromatin yield
 - Improve epitope availability
 - Balance shearing and downstream IP
- Sample concentration
- Select the shortest time resulting in efficient shearing



Analyzing Fragment Size

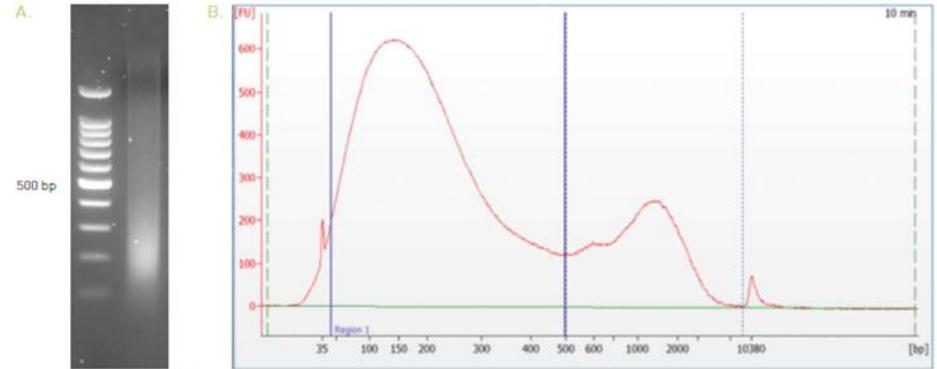
- **De-crosslink**
 - Residual crosslinking retards migration
- **RNase treatment**
 - reduces background
- **DNA purification**
 - IPure beads + DiaMag magnetic rack
 - **Low inputs:** DiaPure columns (eluted in 6 μ l)
- **Electrophoretic analysis**
 - 1.2 - 1.8% agarose gel
 - 300 ng or 60k cells per lane
 - **Low inputs:** FragmentAnalyzer, 2k cells





Analyzing fragment size

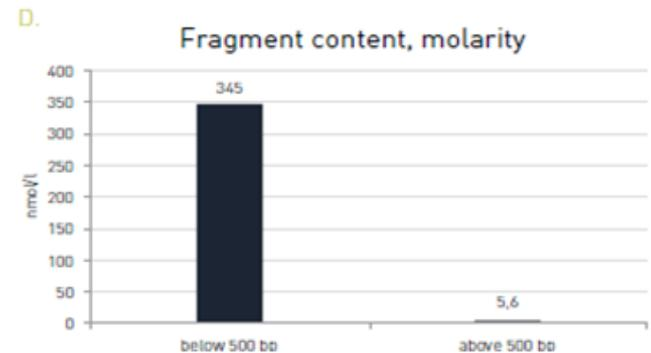
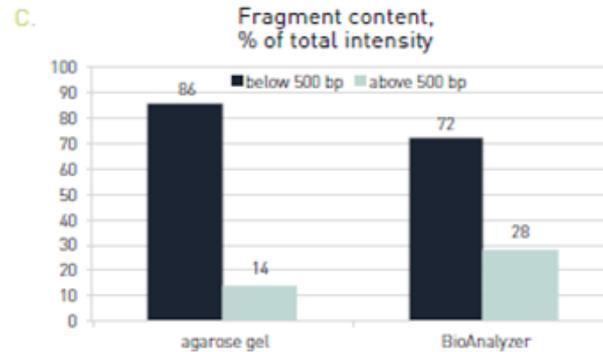
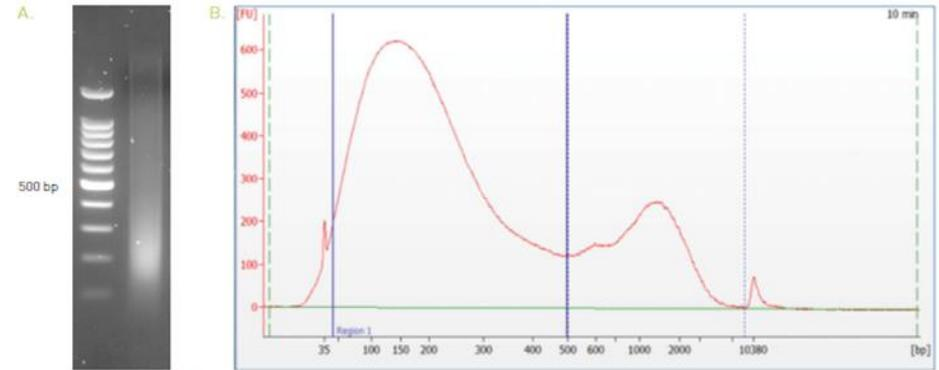
- Use agarose gel or fragment analyzer
- Bioanalyzer or Tapestation:
 - Over-representation of HMW fragments
 - Log-based -> visual misinterpretation of fragment distribution
 - More sensitive to overloading, incomplete reverse crosslinking and residual contaminants





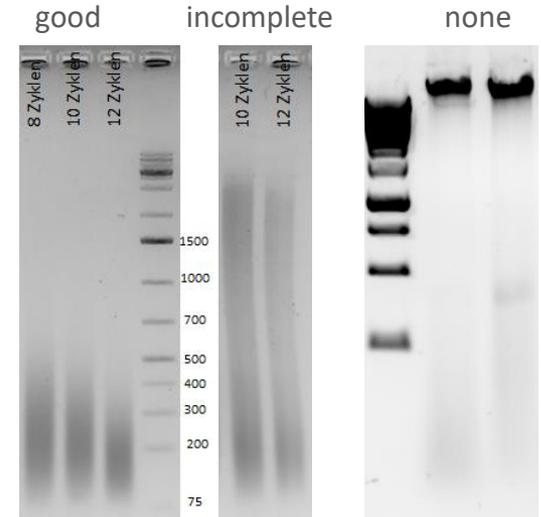
Analyzing fragment size

- Use agarose gel or fragment analyzer
- Bioanalyzer or Tapestation:
 - Over-representation of HMW fragments
 - Log-based -> visual misinterpretation of fragment distribution
 - More sensitive to overloading, incomplete reverse crosslinking and residual contaminants



Troubleshooting Chromatin Shearing

- **No shearing at all**
 - Incomplete lysis – check buffer composition
 - Check instrument efficiency - QC test on sonicator
- **Incomplete shearing**
 - Over-fixation: check fixative and duration
 - Too high cell density
 - Changes in sample require adjustment of shearing protocol
 - Fresh vs. Frozen chromatin
 - Different sample types
 - Wrong consumables (tubes)
 - Sample out of sonication focus
 - droplets on walls/lid of tube
 - Wrong sample volume
 - Wrong temperature (should be 4°C for chromatin)
 - None of the above? -> Check instrument efficiency - QC test on sonicator



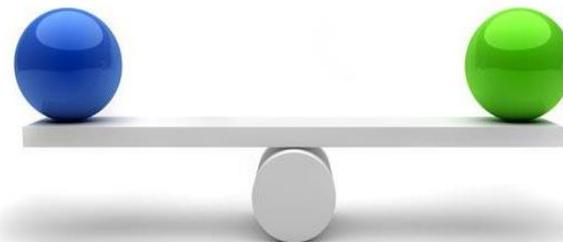


Secrets of ChIP Success

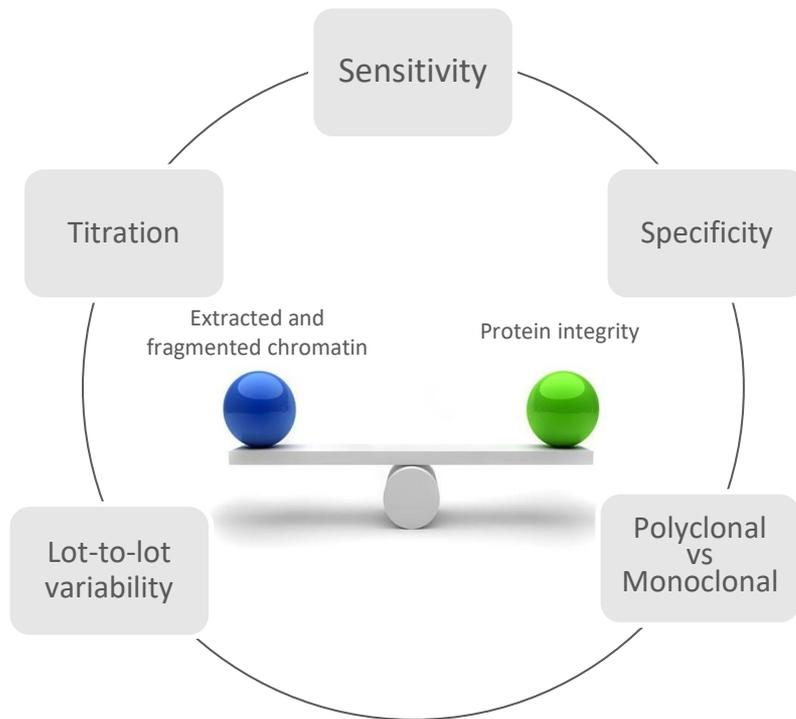
- Prepare “good” chromatin
 - Suitable fragment size and available epitopes
- Use a good antibody at the right concentration
- Optimize for highest specific signal and the lowest background

Extracted and
fragmented chromatin

Protein integrity



Antibodies for ChIP



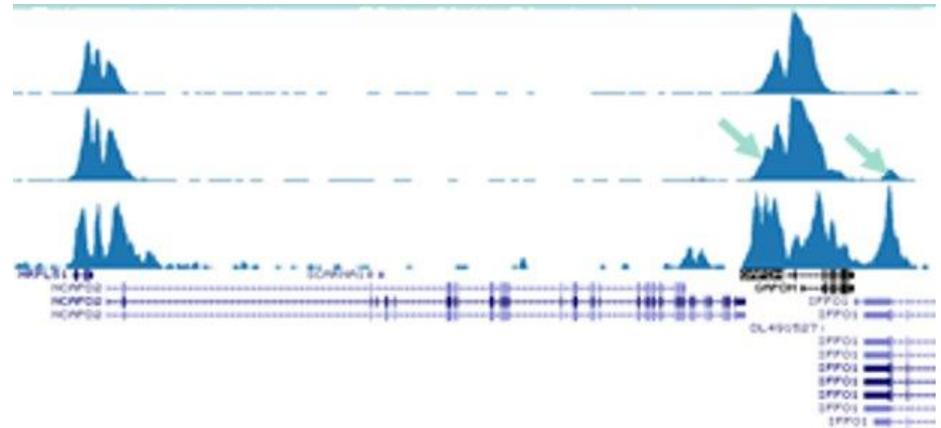
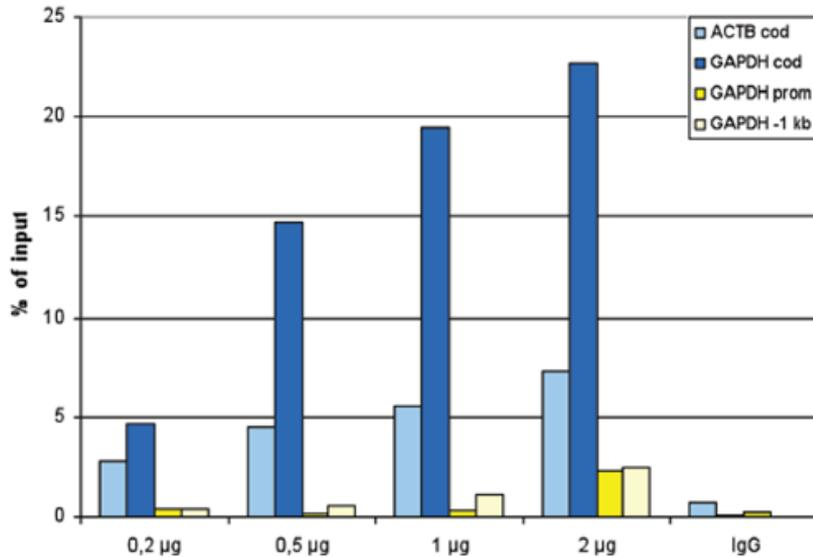


Antibodies for ChIP

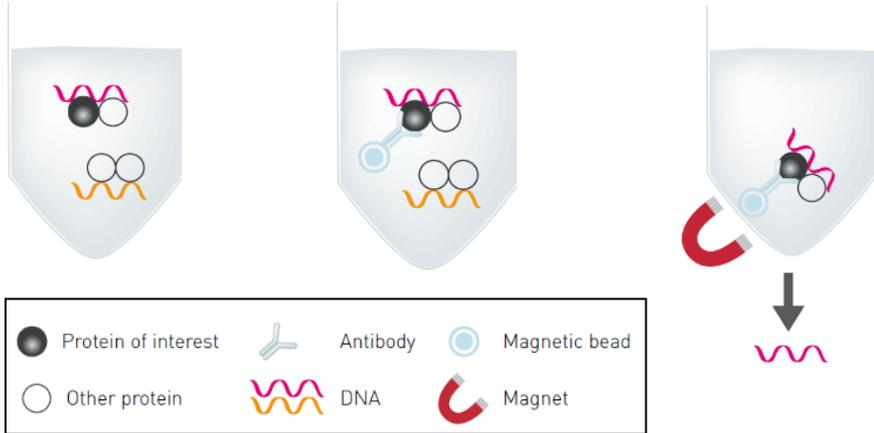
Polyclonal H3K36me3 antibody titration (Diagenode C15410192)

1µg IgG as negative IP control

Chromatin from 100.000 cells



Setting up IP



Components:

- Sheared chromatin
- **ChIP grade antibodies**
-> optimized quantity
- ProteinA/G magnetic beads
- ChIP buffer
- Protease inhibitor cocktail





Antibodies for ChIP – What beads?

Agarose beads

Required: centrifuge

Sensitive to handling

High background

Risk of carry-over

Magnetic beads

Required: magnetic rack

Robust

Low background

Easy separation

Limit antibody amounts to bead capacity!



Protein G or A beads

- Both bind to IgG antibodies and are structurally similar
- Slightly different affinities for IgG subclasses across different species.
- Use appropriate depending on the IgG subtype you are using:

Protein A

Rabbit
Pig
Dog
Cat

Protein G

Mouse
Rat
Human



Setting up IP: Input Sample

- Fraction of sheared chromatin is kept aside as INPUT
 - Processed in parallel with IP-samples from reversed crosslinking
 - Include one input for each chromatin sample
- Key reference for ChIP-qPCR and ChIP-seq analysis
- ChIP-qPCR: used to calculate the recovery (% of input)
- ChIP-seq: mandatory for bio-informatics analysis
 - Normalization for mappability of a region, avoid duplication bias etc.
 - Input pooling can be considered for ChIP-seq on very similar samples

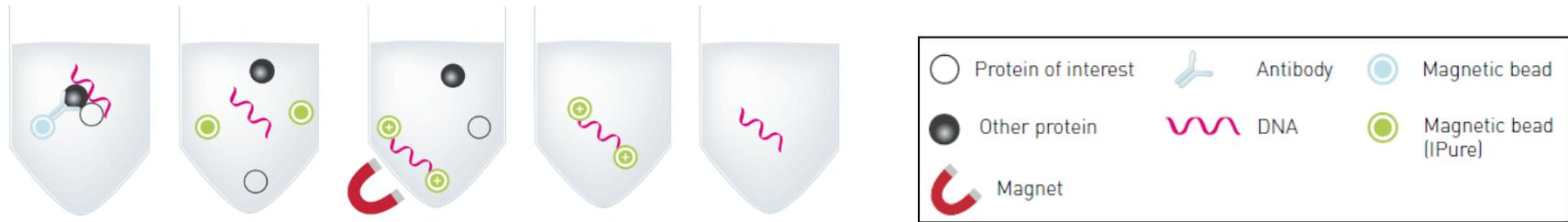


Setting Up IP: Additional Controls

- Positive control (H3K4me3, CTCF):
 - Confirm overall efficiency of ChIP workflow
 - ChIP optimization for new target
- Negative Control (IgG)
 - Measure of non specific IP background
 - Include one negative IgG control in each series of ChIP reactions
- Not necessary to sequence these but good control for qPCR
- Biological Replicates
 - ChIP-qPCR ≥ 3
 - ChIP-seq ≥ 2



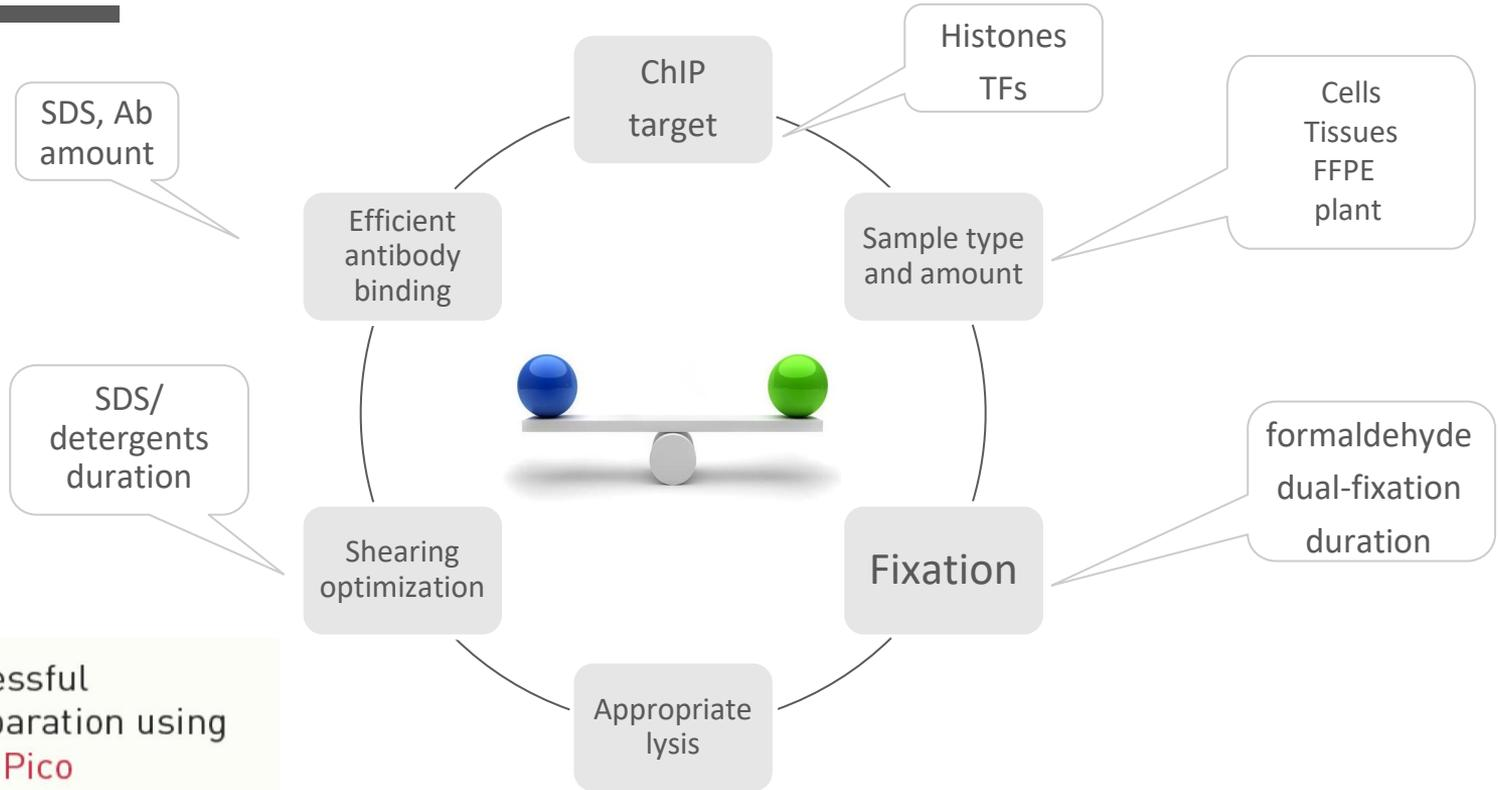
ChIP Protocol – Elution, de-crosslinking and DNA isolation



- Elution of the chromatin complexes from protein A/G-bound magnetic beads:
elution buffer 30 min at RT
- Reversal of cross-links:
Incubation for at least 4h at 65°C
- Isolation of the ChIP'd DNA:
IPure magnetic beads
Column purification (DiaPure columns for low elution volumes >6μl)



Summary – Tips to Prepare Good Chromatin



Guide for successful chromatin preparation using the **Bioruptor® Pico**

ChIP-qPCR

- Target & Primer selection is key for ChIP-qPCR
- Predict qPCR-targets from ChIP-seq data
- qPCR as QC prior ChIP-seq
- If no ChIP-seq data:
 - estimate binding from similar data, biological function etc.
 - use multiple regions
- PCR program depends on Master Mix, qPCR system and primer pairs

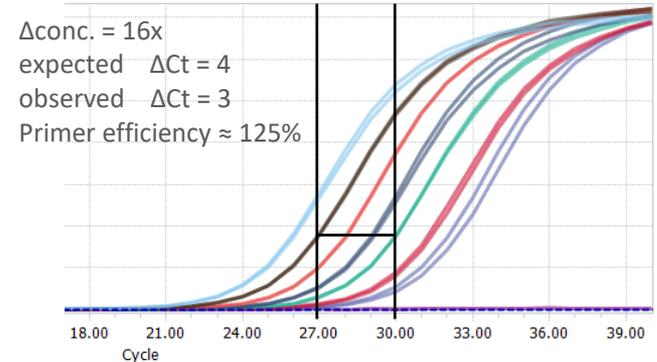
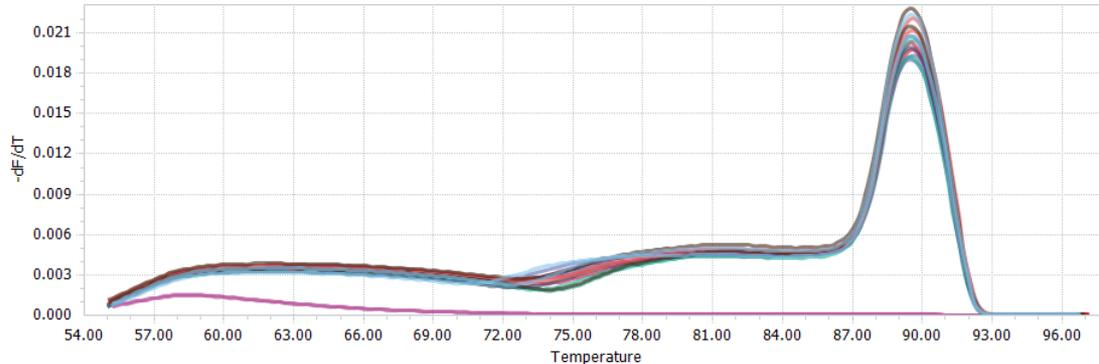




Setting up ChIP-qPCR

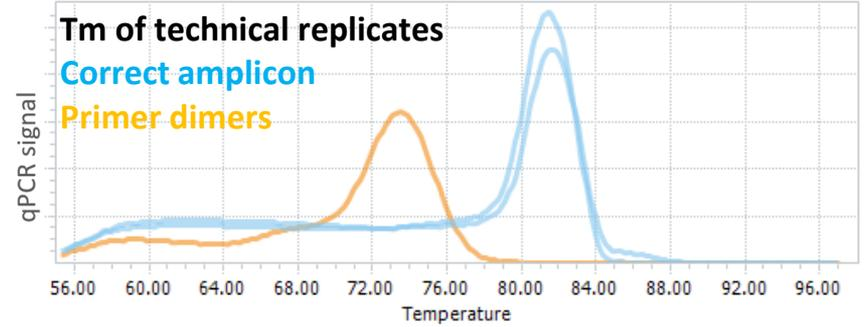
- **Target primer design:**
 - Place primers around binding site
 - 50-150bp amplicons
 - 20-30 bp primers with a similar T_m between 55° and 60°C

- **Primer pair validation:**
 - Check on gDNA/input for T_m profile
 - Check efficiency (95-105% acceptable)



Setting up ChIP-qPCR

- QC
 - T_m : no second peaks, no primer dimers
 - technical replicates within 0.3 Ct
 - Ct values $\gg 30$ are often not reliable





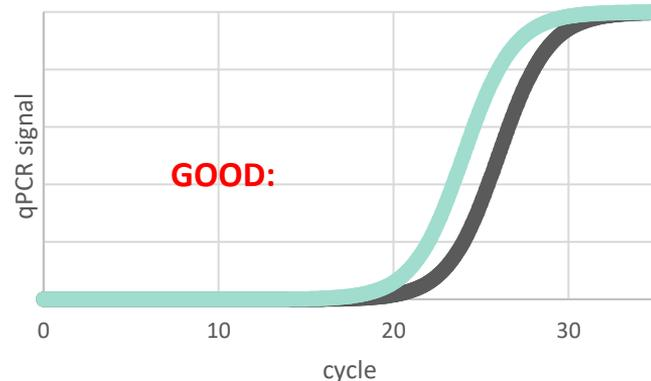
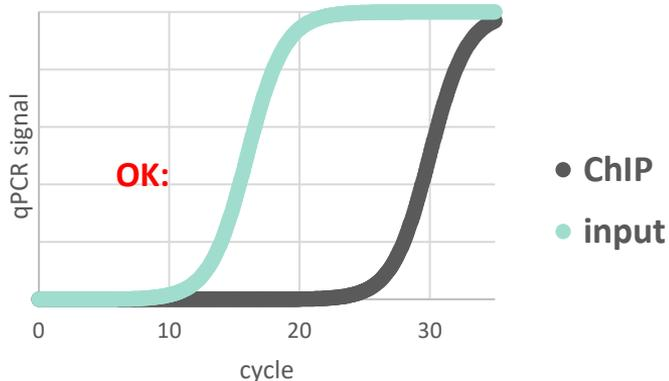
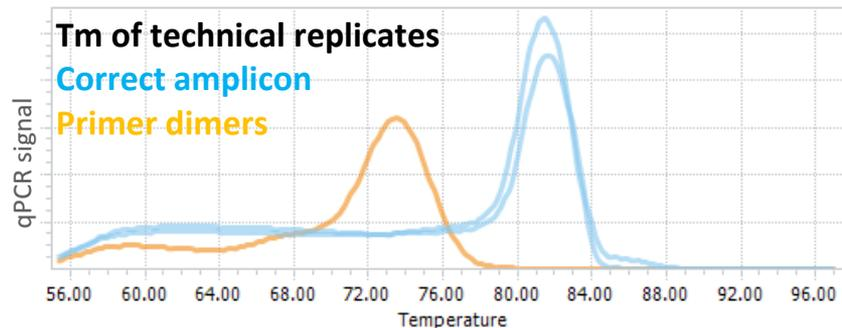
Setting up ChIP-qPCR

QC

- T_m : no second peaks, no primer dimers
- technical replicates within 0.3 Ct
- Ct values $\gg 30$ are often not reliable

ChIP and input samples

- Adjust amount of ChIP-sample/input to obtain comparable Ct values
- Consider primer efficiency for high ΔCt





ChIP-qPCR analysis

ChIP recovery R :

- chromatin recovery as % of Input
- R should be minimal for the IgG control and high for the epitope of interest

$$R = \frac{2^{Ct(input) - Ct(ChIP)}}{100 (input\ fraction)}$$

-> for each target separately

Input fraction is often corrected with a logarithmic compensatory factor, e.g. -6.64 Ct for 1% input

ChIP fold-enrichment F with $\Delta\Delta Ct$ method:

- fold-enrichment of bound vs. epitope-“free” regions
- S varies depending on regions analyzed

$$F = \frac{R(positive\ region)}{R(negative\ region)}$$

-> main success parameter

Successful ChIP?

- If wrong regions targeted – risk of false negative result
- $F > 2$ for ChIP-qPCR analysis
- $F > 4$ for ChIP-seq



ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

Ct values

A

	IgG		H3K4me3		Input (1% of sample)	
P	34.0	36.0	26.0	26.1	27.0	27.1
G1	35.0	35.0	27.0	27.2	26.0	25.7
G2	-	37.0	33.0	34.0	29.0	29.6
N	34.0	35.0	33.0	33.5	28.0	28.1

B

	IgG		H3K4me3		Input (1% of sample)	
	-	36.4	25.4	25.5	26.5	26.6
	35.4	36.4	26.4	26.6	25.4	25.1
	34.4	33.4	28.7	28.5	28.6	28.7
	35.4	-	32.4	33.4	27.4	27.5



ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

Ct values

	A					
	IgG		H3K4me3		Input (1% of sample)	
P	34.0	36.0	26.0	26.1	27.0	27.1
G1	35.0	35.0	27.0	27.2	26.0	25.7
G2	-	37.0	33.0	34.0	29.0	29.6
N	34.0	35.0	33.0	33.5	28.0	28.1

	B					
	IgG		H3K4me3		Input (1% of sample)	
	-	36.4	25.4	25.5	26.5	26.6
	35.4	36.4	26.4	26.6	25.4	25.1
	34.4	33.4	28.7	28.5	28.6	28.7
	35.4	-	32.4	33.4	27.4	27.5

1. Technical sanity check

- values out of range
- high Ct-variation (>0.3) between technical replicates
- A/B inputs shifted
- $Ct(H3K4me3) \approx Ct(input)$



ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

2. Averaging Technical replicates

	A			B		
	IgG	H3K4me3	Input (1% of sample)	IgG	H3K4me3	Input
P	35.0	26.1	27.1	36.4	25.5	26.6
G1	35.0	27.1	25.9	35.9	26.5	25.3
G2	37.0	33.5	29.3	33.9	28.6	28.7
N	34.5	33.3	28.1	35.4	32.9	27.5

3. Biological Sanity Check

- Ct(P) < Ct(N) for H3K4me3
- Ct(H3K4me3) << Ct(IgG)



ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

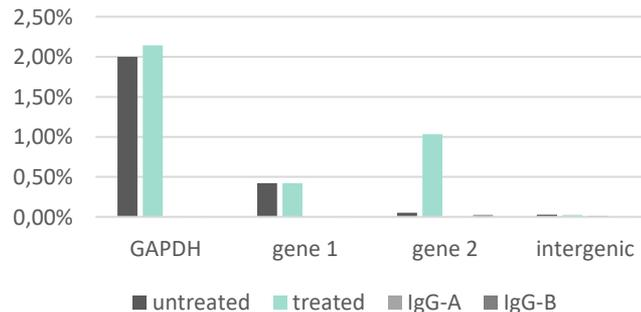
	A			B		
	IgG	H3K4me3	Input (1% of sample)	IgG	H3K4me3	Input
P	35.0	26.1	27.1	36.4	25.5	26.6
G1	35.0	27.1	25.9	35.9	26.5	25.3
G2	37.0	33.5	29.3	33.9	28.6	28.7
N	34.5	33.3	28.1	35.4	32.9	27.5

4. Recovery (% of input)

$$R = \frac{2^{Ct(input)} - Ct(ChIP)}{100(input\ fraction)}$$

	A		B	
	IgG-A	IgG-B	IgG-A	IgG-B
P	2.00%	0.00%	2.14%	0.00%
G1	0.42%	0.00%	0.42%	0.00%
G2	0.05%	0.00%	1.04%	0.03%
N	0.03%	0.01%	0.02%	0.00%

H3K4me3 %input





ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
Antibodies: IgG, H3K4me3

	H3K4me3 % of input	
	A	B
P	2.00%	2.14%
N	0.03%	0.02%

5. ChIP fold-enrichment

$$F_A = \frac{R(P_A)}{R(N_A)} = 74x \text{ enrichment}$$

$$F_B = \frac{R(P_B)}{R(N_B)} = 94x \text{ enrichment}$$



ChIP-qPCR Exercise

Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

	H3K4me3 % of input	
	A	B
P	2.00%	2.14%
N	0.03%	0.02%

5. ChIP fold-enrichment

$$F_A = \frac{R(P_A)}{R(N_A)} = 74x \text{ enrichment}$$

$$F_B = \frac{R(P_B)}{R(N_B)} = 94x \text{ enrichment}$$

7. Assessment

- $F > 2$ ChIP-qPCR qualified
- $F > 4$ ChIP-seq qualified

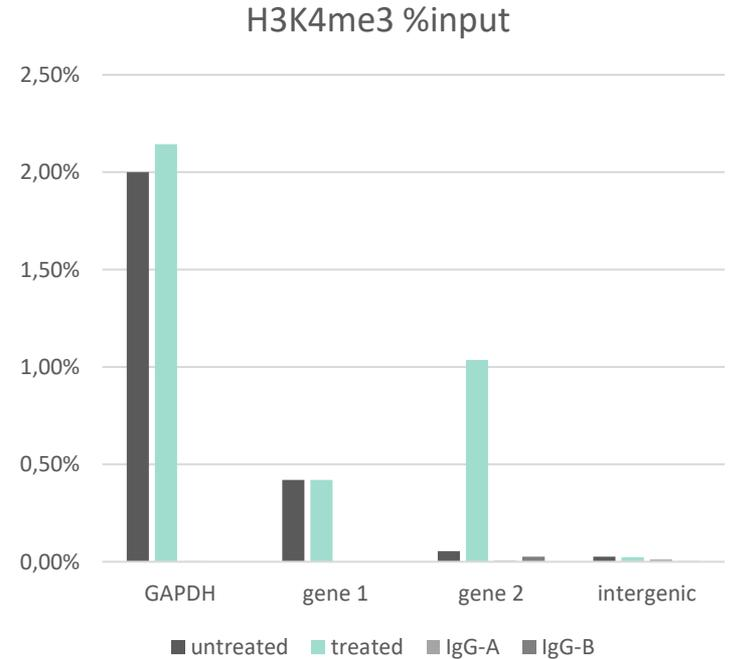
8. Optimization

- if R or F are low, optimize ChIP parameters



ChIP-qPCR Analysis summary

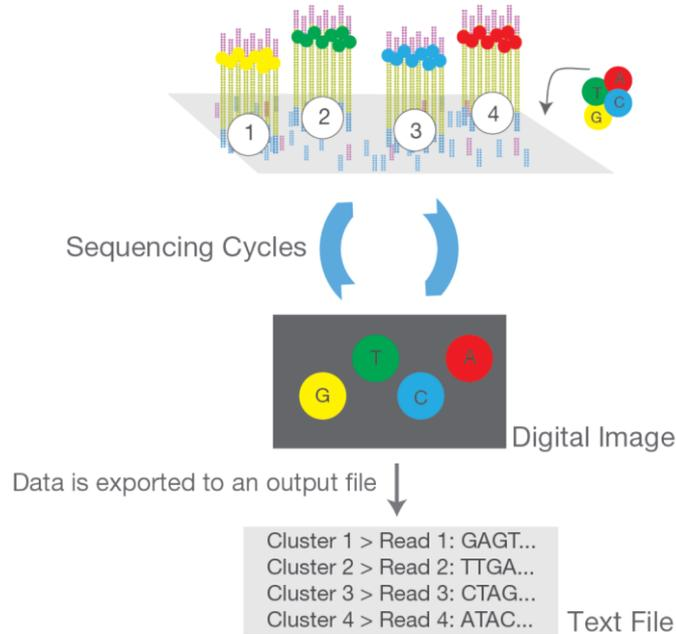
- Determine ChIP-recovery and fold-enrichment
- input used as reference to calculate ChIP-recovery
- Each ChIP-target requires specific control regions
- Suitable control regions can vary among samples





ChIP-seq: library prep

C. Sequencing



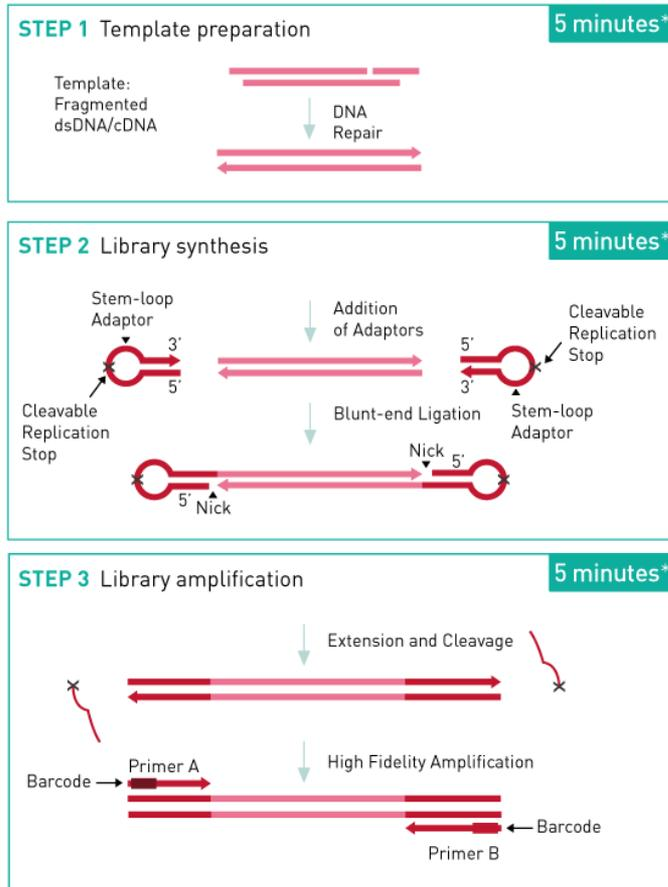
ChIP-seq: Library Prep



- Low input
- Minimal steps
 - To maximize recovery
- Sensitive
 - Minimal PCR amplification
- Suitable for pooling



MicroPlex kit workflow

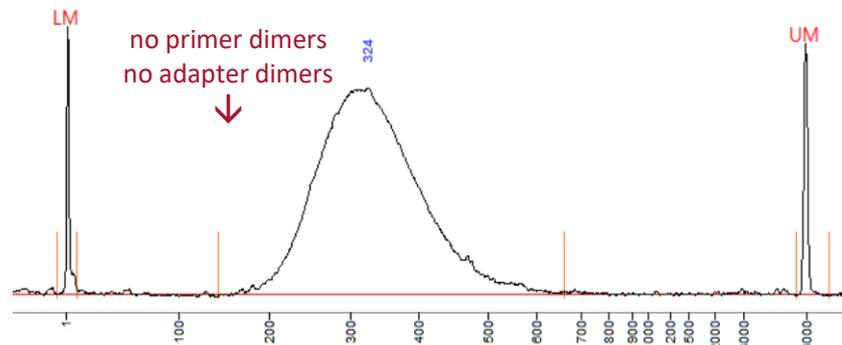


* hands-on-time



Library Pooling

- **Determine library size**
 - Bioanalyzer or Fragment Analyzer
 - Identify adapter dimers or unexpected library sizes
- **Quantify**
 - Qubit
 - qPCR – quantify sequencable library
 - Convert from ng/μl to nM using average library size
- **Dilute and Pool normalized libraries**
 - Same size for best clustering





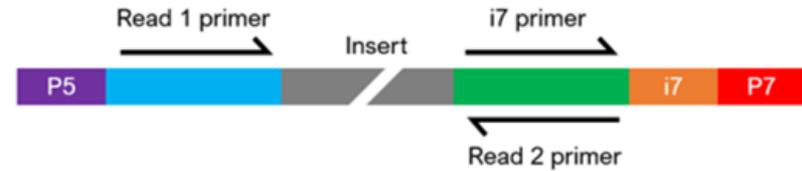
Benefits of Multiplexing

- **Fast High-Throughput Strategy:**
 - Large sample numbers can be simultaneously sequenced
- **Cost-Effective Method:**
 - Reduces time and reagent use
 - Cluster detection more efficient with different bases in beginning of read
- **Simplified Analysis:**
 - Automatic sample identification with "**barcodes**" using Illumina software

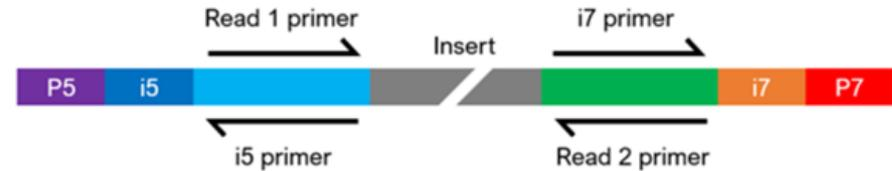


Single and Dual-indexed Libraries

- **Single-index sequencing**
 - Low level of multiplexing
- **Dual indexing**
 - Higher multiplexing - more samples per lane possible
 - Higher accuracy of sample identification
- **Unique dual indexing (UDI)**
 - Allows filtering of index-hopping events



Dual-Indexed Sequencing





ChIP-seq: Sequencing Settings

Read length

- 50 bp sufficient for most ChIPs
- adjust fragment-size to read length

Sequencing depth

- mainly set by samples/flow cell and flow cell type
- 30 M reads for sharp peaking targets e.g. H3K4me3, H3K27ac
- 50 M for broadly distributed and abundant targets e.g. H3K27me3
- use same depth for input

Replicates

- \geq duplicates
- increased replicate number will improve sensitivity of the downstream analysis

Input sequencing

- one input per sample is gold standard
- pooling inputs from replicates can often be considered

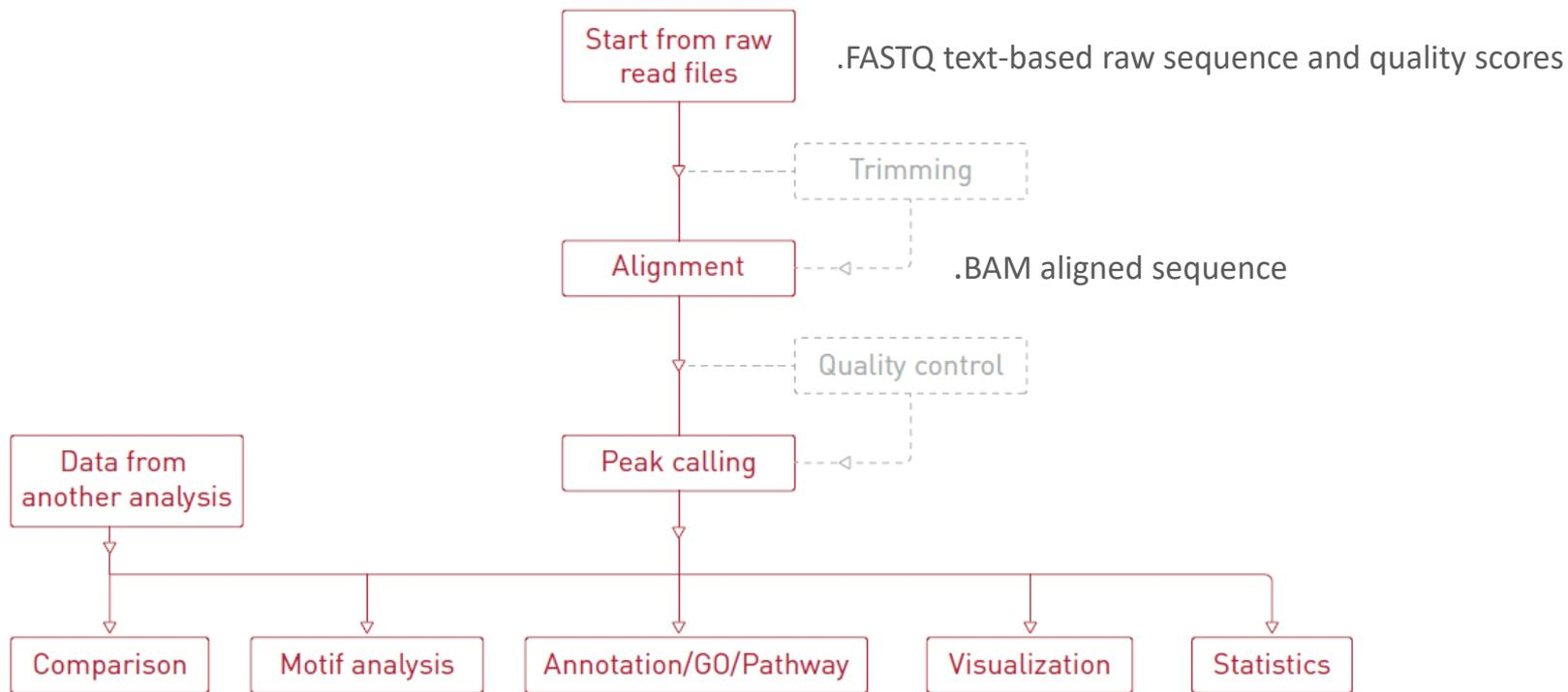


ChIP-seq: Analysis

Bioinformatician	R free-ware and online tool kits	https://www.r-project.org/ www.bioinformatics.babraham.ac.uk/projects/seqmonk/ biit.cs.ut.ee/gprofiler/gost
Wet-lab expert with free time	standard bio-informatic services free-ware and online tool kits	https://www.diagenode.com/en/categories/Services Comprehensive Multi-Omic and bio-info services www.bioinformatics.babraham.ac.uk/projects/seqmonk/ initial & advanced data analysis, genome browser, graphical presentation of data https://biit.cs.ut.ee/gprofiler/gost Functional profiling tool
no expertise or no free time	advanced bio-informatic services	https://www.diagenode.com/en/categories/Services Comprehensive Multi-Omic and bio-info services



ChIP-seq: Analysis



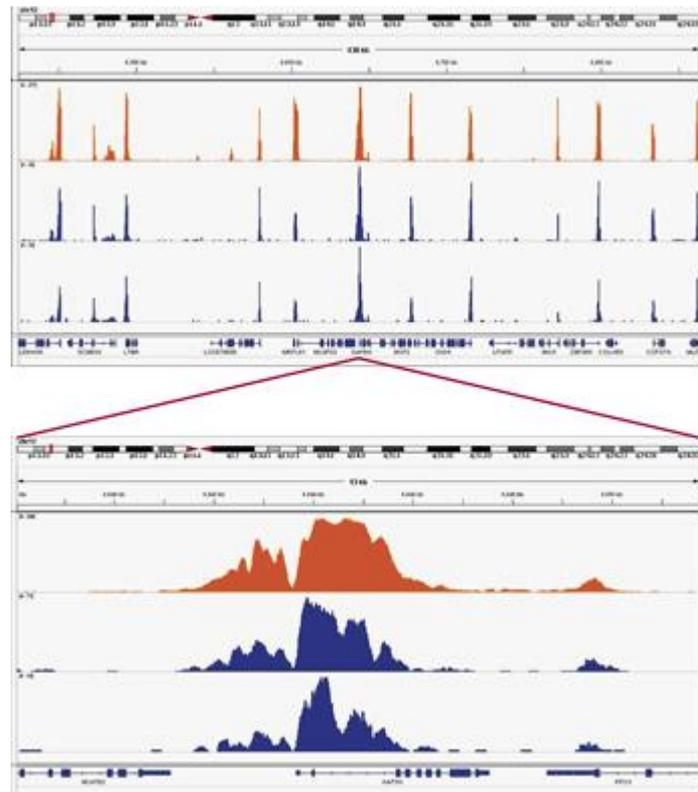
ChIP-seq: Analysis

Standard bioinformatic analysis:

- alignment to reference genome
- peak calling

Advanced bioinformatic analysis:

- annotation of peaks and genes
- differential analysis of peak/gene lists
- unsupervised analysis (PCA, clustering)
- functional enrichment analysis
(e.g. Pathway analysis, Gene ontology)
- Machine learning
- integrative analysis
(RNA-seq, ATAC-seq, more ChIP-seq targets)
- publication-ready Visualization of genomic regions





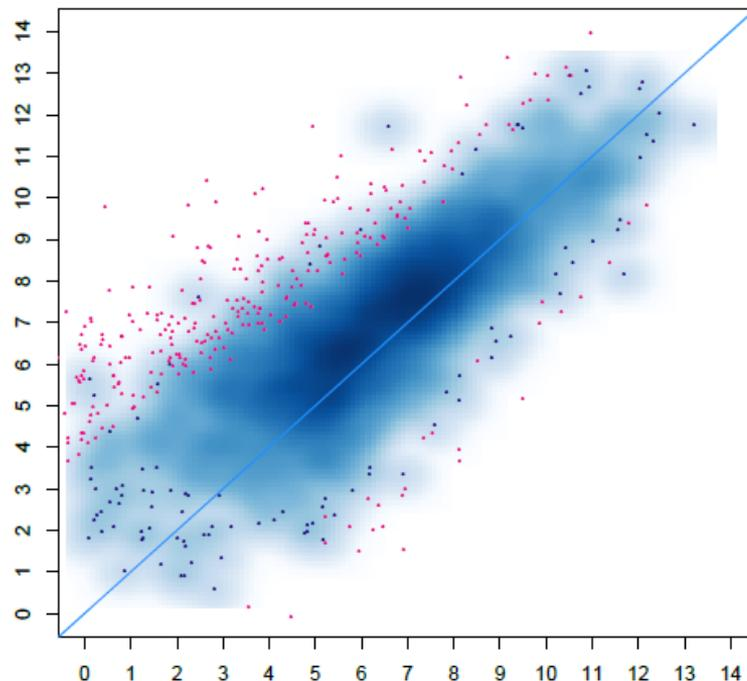
ChIP-seq: Analysis

Standard bioinformatic analysis:

- alignment to reference genome
- peak calling

Advanced bioinformatic analysis:

- annotation of peaks and genes
- **differential analysis of peak/gene lists**
- unsupervised analysis (PCA, clustering)
- functional enrichment analysis
(e.g. Pathway analysis, Gene ontology)
- Machine learning
- integrative analysis
(RNA-seq, ATAC-seq, more ChIP-seq targets)
- publication-ready Visualization of genomic regions





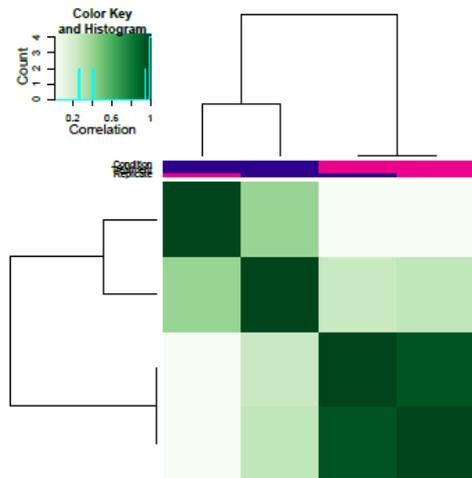
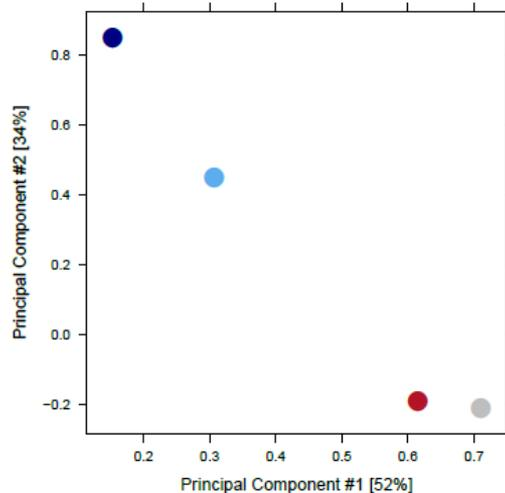
ChIP-seq: Analysis

Standard bioinformatic analysis:

- alignment to reference genome
- peak calling

Advanced bioinformatic analysis:

- annotation of peaks and genes
- differential analysis of peak/gene lists
- **unsupervised analysis (PCA, clustering)**
- functional enrichment analysis
(e.g. Pathway analysis, Gene ontology)
- Machine learning
- integrative analysis
(RNA-seq, ATAC-seq, more ChIP-seq targets)
- publication-ready Visualization of genomic regions



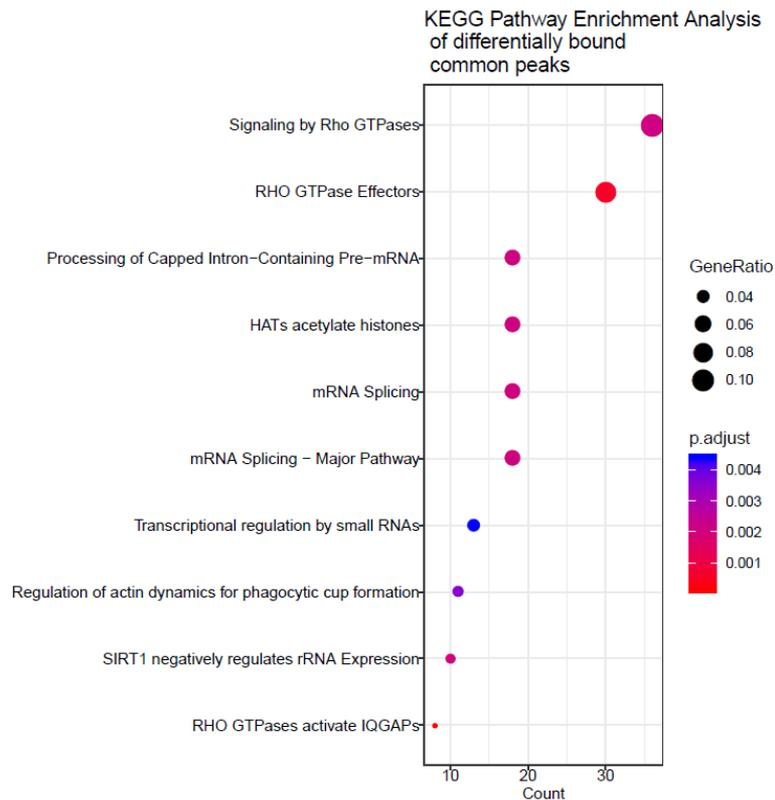
ChIP-seq: Analysis

Standard bioinformatic analysis:

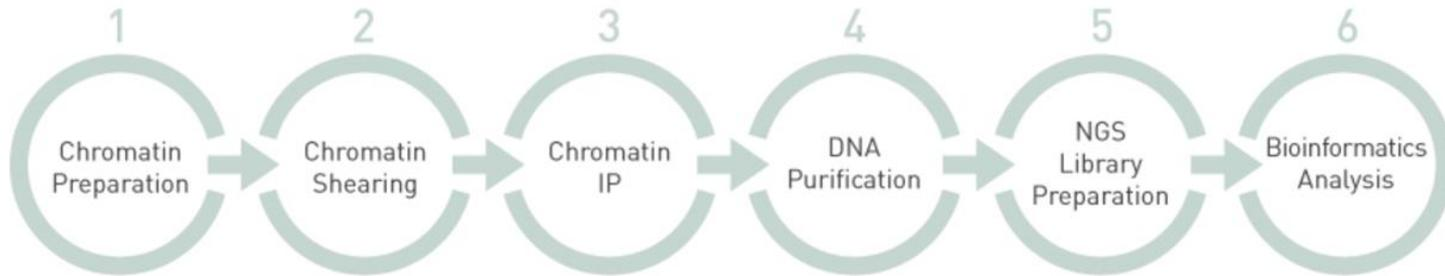
- alignment to reference genome
- peak calling

Advanced bioinformatic analysis:

- annotation of peaks and genes
- differential analysis of peak/gene lists
- unsupervised analysis (PCA, clustering)
- **functional enrichment analysis**
(e.g. Pathway analysis, Gene ontology)
- Machine learning
- integrative analysis
(RNA-seq, ATAC-seq, more ChIP-seq targets)
- publication-ready Visualization of genomic regions



Summary



Solutions for each step of the way : Bioruptor, Kits, Antibodies, NGS library prep and Services

www.diagenode.com



Other Methods to Study Chromatin

- Methods to study the **interaction** between proteins and DNA (immuno- assays):
 - ChIPmentation
 - Cut&TAG

ChIPmentation™ & μChIPmentation™

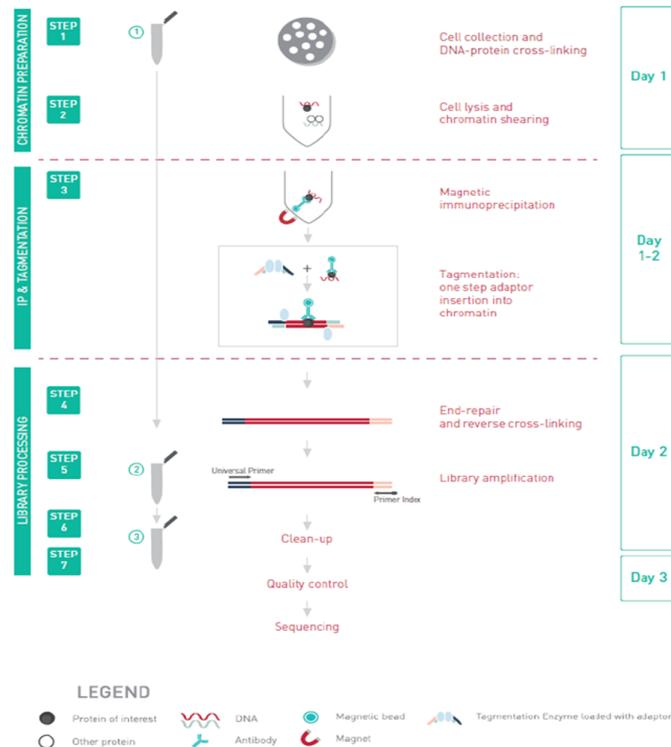
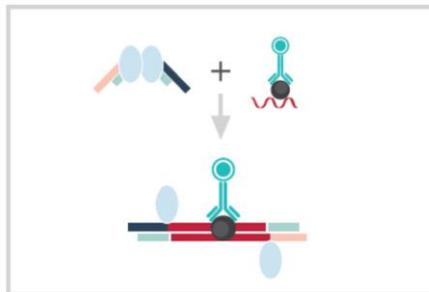
Easier and faster than classical ChIP-seq

Validated for various histone marks

Ideal for analysis of large cohorts of samples (easy and fast)

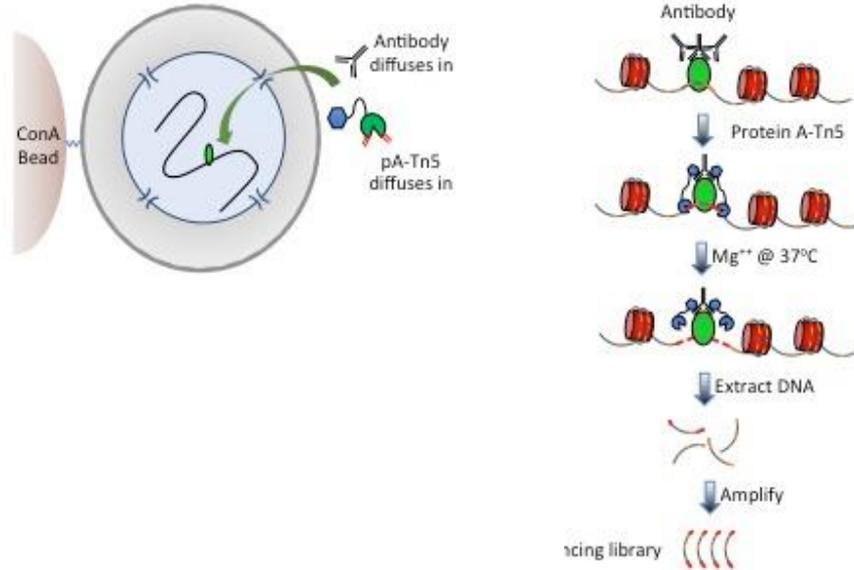
Ideal for analysis of large number of marks on a unique sample

μChIPmentation for 10,000 cells



CUT&Tag: Cleavage Under Targets and Tagmentation

CUT&Tag (Cleavage Under Targets & Tagmentation)



Key features:

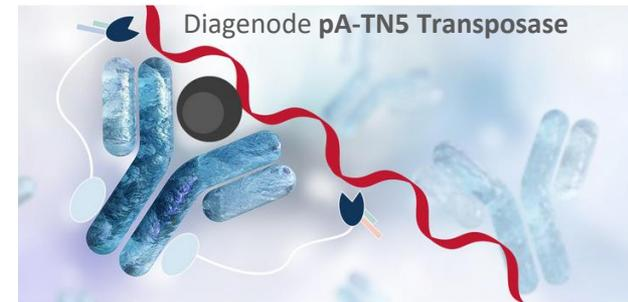
Crucial reagent:

- proteinA-Tn5

Fast and easy protocol:

- fast tagmentation-based library prep
- No chromatin prep

Suitable for low cell numbers

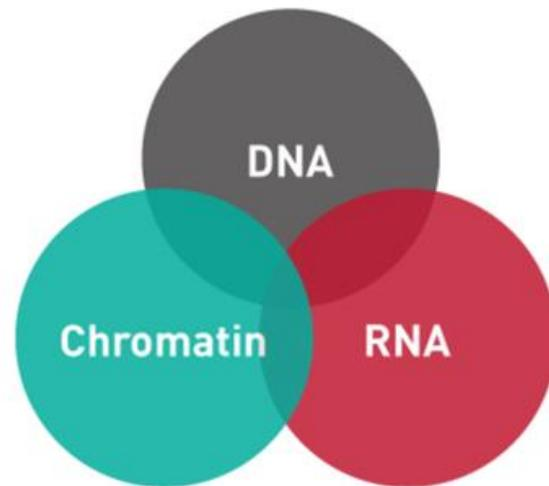




EPIGENOMICS PROFILING SERVICES

15 YEARS OF EXPERTISE IN EPIGENETICS

- End-to-end epigenetic service and analysis
- Collaborative and customized project design
- Dedicated in-house expert for your project
- Presentation-quality data and graphs



CHIP WORKSHOP

THANK YOU!

Thank you for taking part in our ChIP workshops! (more coming)

Presentation will be sent to each participant

Watch for a little survey in your inbox – your feedback is invaluable

Stay in the know

Join our newsletter



@Diagenode

Spread the word
#EpiWorkshopsWithDiagenode
(more gifts for you :-))

customer.support@diagenode.com

www.diagenode.com

