



Innovating Epigenetics Solutions



WELCOME TO DIAGENODE

ChIP Workshop

Jessica Apulei & Juri Kazakevych

16-17 February 2021



Jessica Apulei



WHO AM I?

- Inside sales representative and technical support specialist at Diagenode
- PhD at Collège de France, Paris , France 2015-2019
- Postdoc at Harvard University, Cambridge, Boston MA 2019-2020



Juri Kazakevych



WHO AM I?

- Inside sales representative and technical support specialist at Diagenode since 2019
- PhD in Mainz, Germany 2012-2016
- Postdoc in Cambridge, UK 2016-2019

OBJECTIVES

Day 1

- ChIP overview
 - Chip-qPCR vs. ChIP-Seq
- Chromatin preparation
 - Fixation, Cell lysis and Chromatin shearing
- Setting up IP
 - Antibodies, Replicates, inputs, controls

Day 2

- ChIP-qPCR
- ChIP-seq: library prep & sequencing
- ChIP-seq: analysis
- Overview of Alternate methods

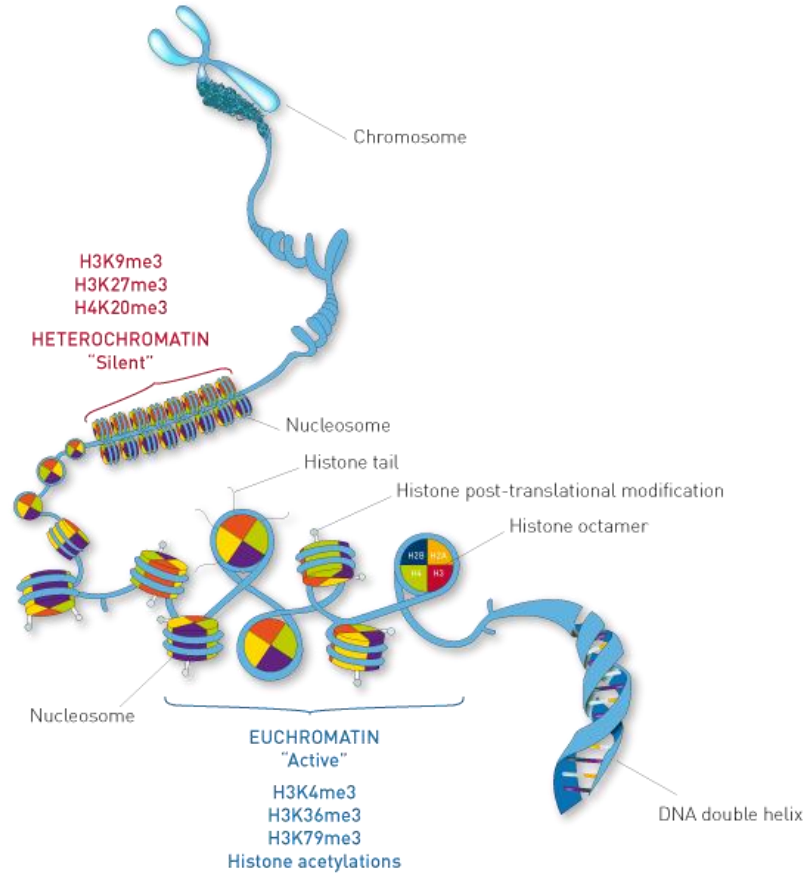




INTRODUCTION

What is Chromatin?

A complex of DNA and proteins found in eukaryotic cells

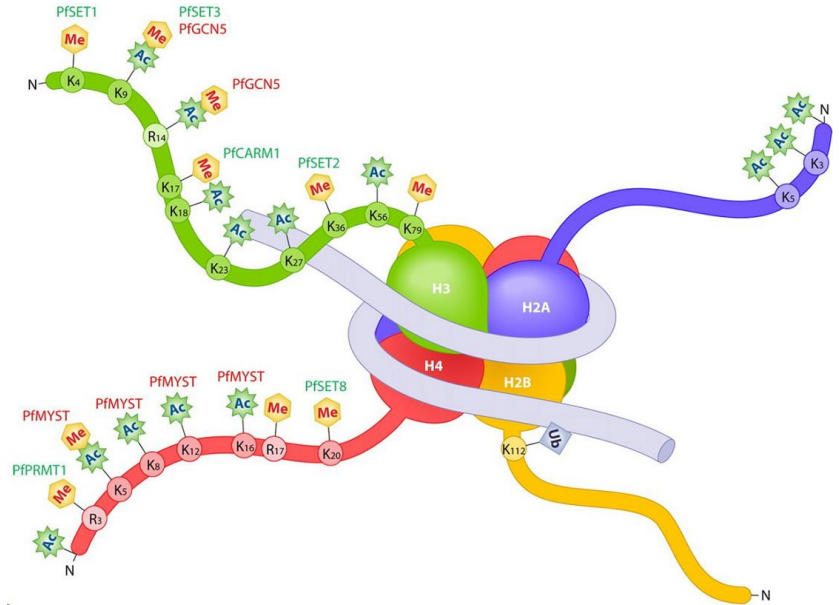




INTRODUCTION

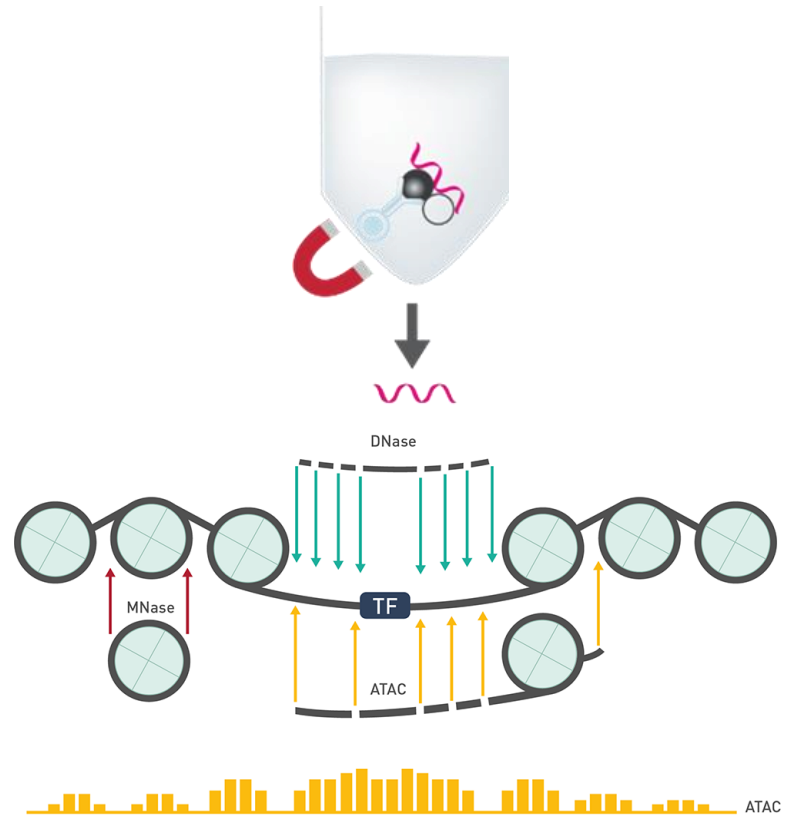
What is Chromatin?

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



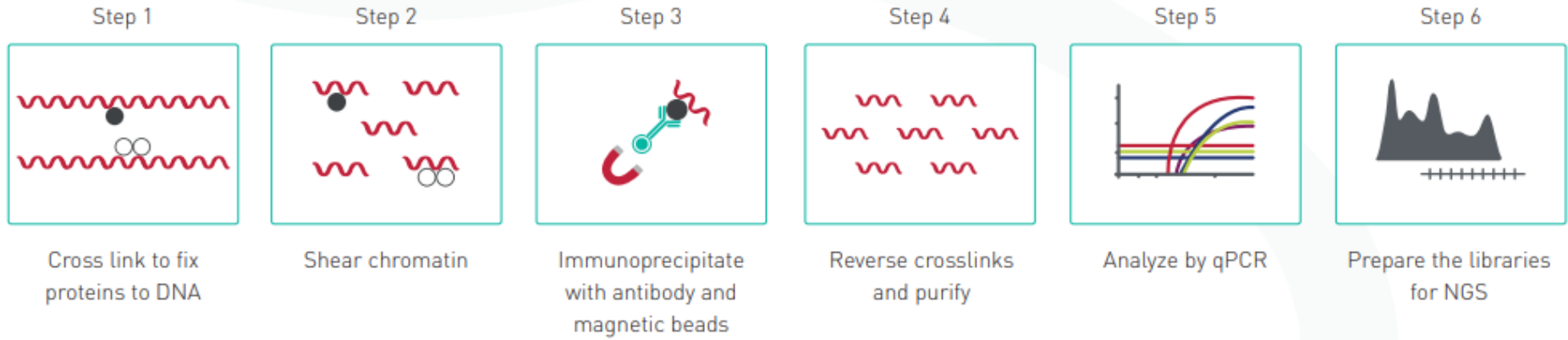
CHROMATIN ANALYSIS

- 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):



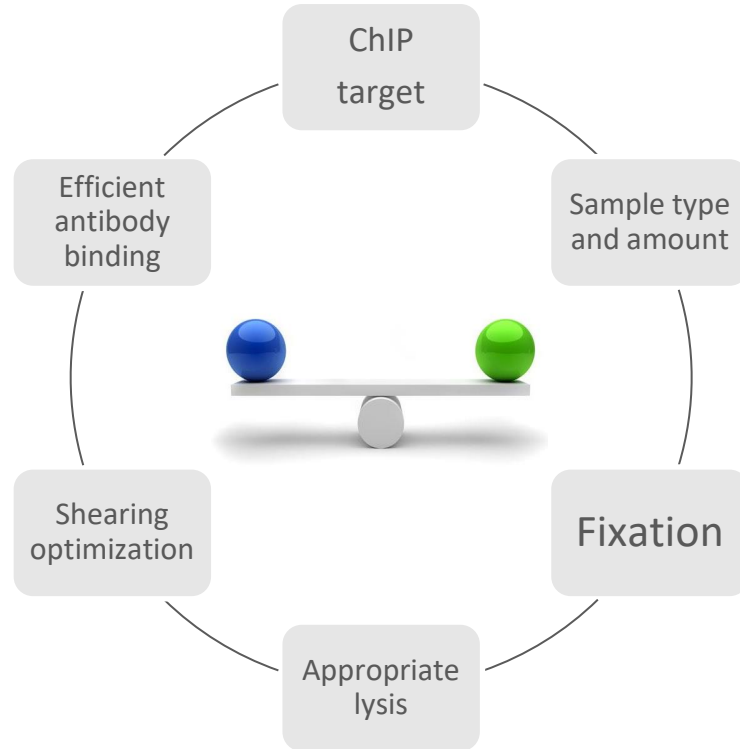


ChIP-qPCR or ChIP-Seq?

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



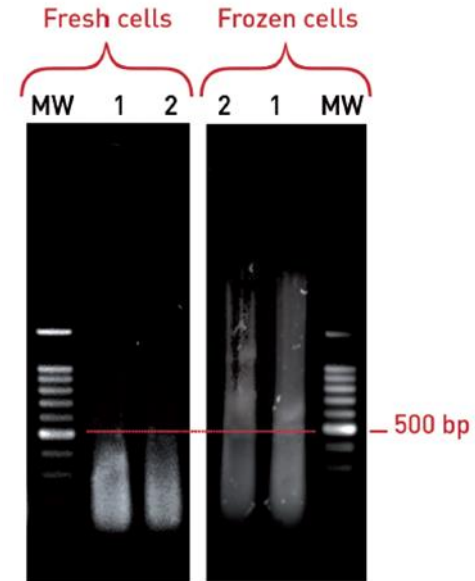
Fixation

- 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

- One step lysis – for low cell numbers
 - Lyse cells directly with an SDS-containing buffer
- Two step lysis – standard protocol, difficult cells
 - Remove soluble cytosolic proteins first
 - Improves sonication efficiency
 - Reduces background
- 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



Optimal: Perform CHIP workflow directly, without freezing

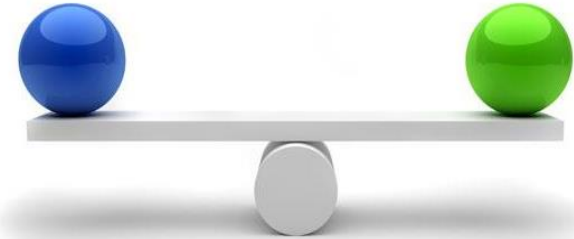


Secrets of ChIP Success

- 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

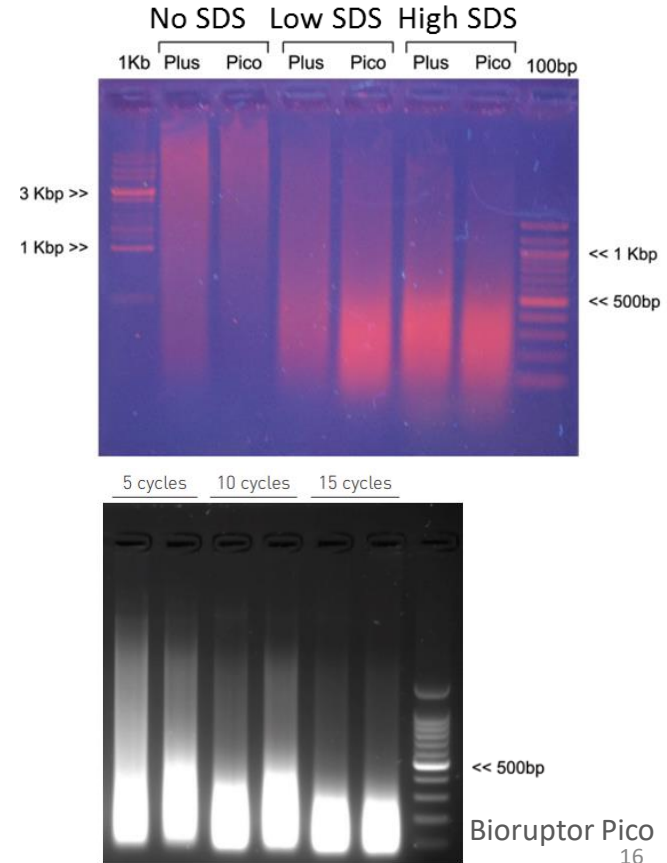
Extracted and
fragmented chromatin

Protein integrity



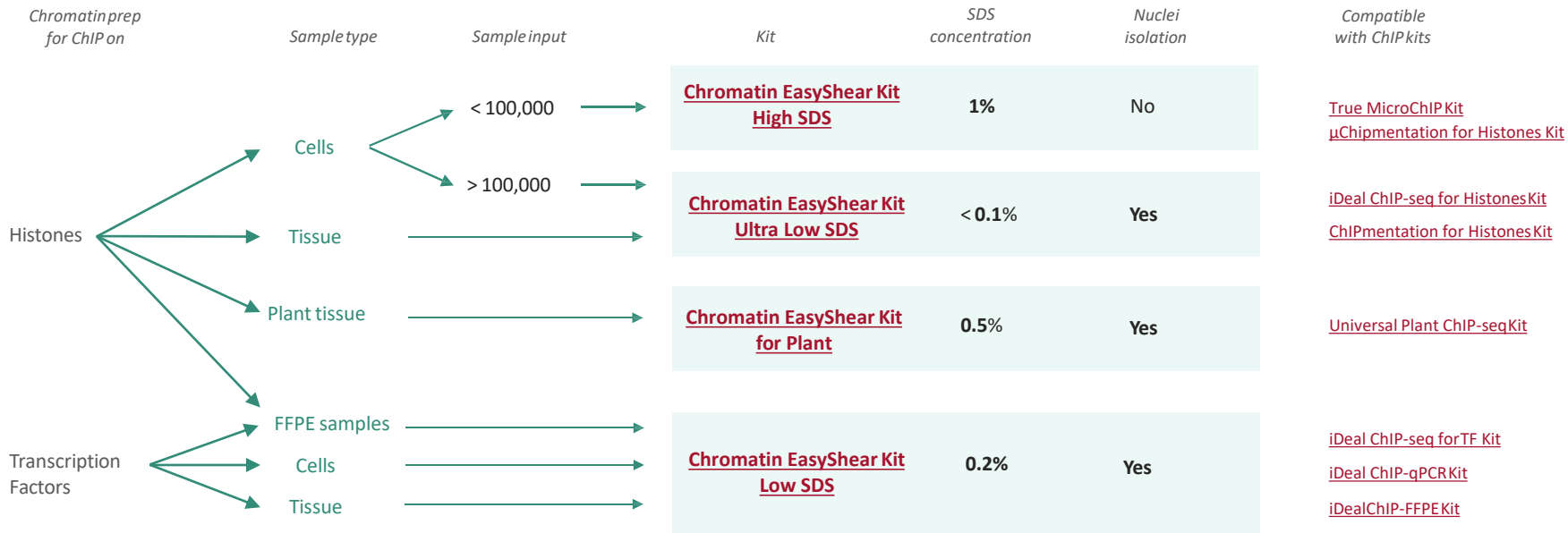
Chromatin Shearing

- 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





Chromatin Shearing



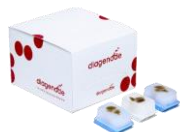


Chromatin Shearing

Chromatin EasyShear Kits

Kit of choice for:

- Chromatin prep with Bioruptor



Chromatin EasyShear Kit

+



Bioruptor

- Chromatin optimization prior to ChIP kits



Chromatin EasyShear Kit

+



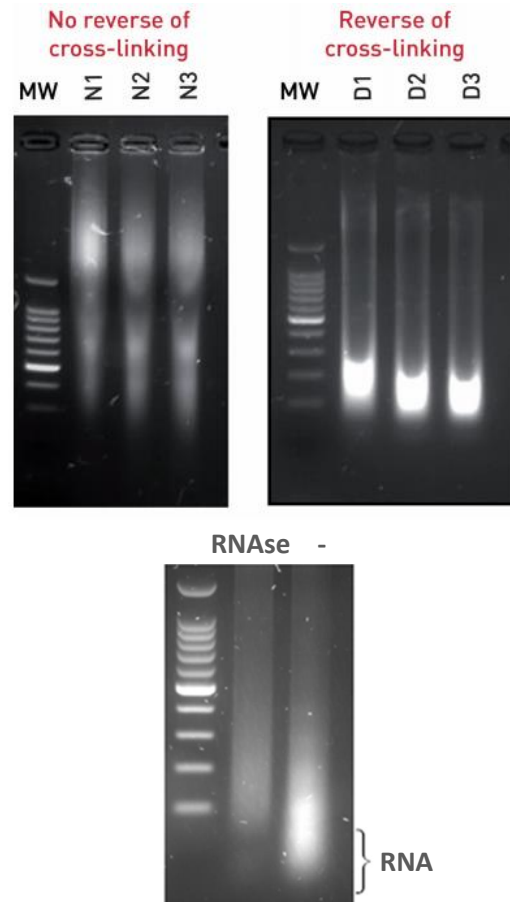
Diagenode's ChIP Kit

Features & Benefits

- Highly optimized for chromatin preparation
- Preserves epitope integrity
- Recommended for the optimization of the chromatin shearing of a new cell line/new sample type prior to ChIP
- Validated: Kit performance has been validated in ChIP-seq

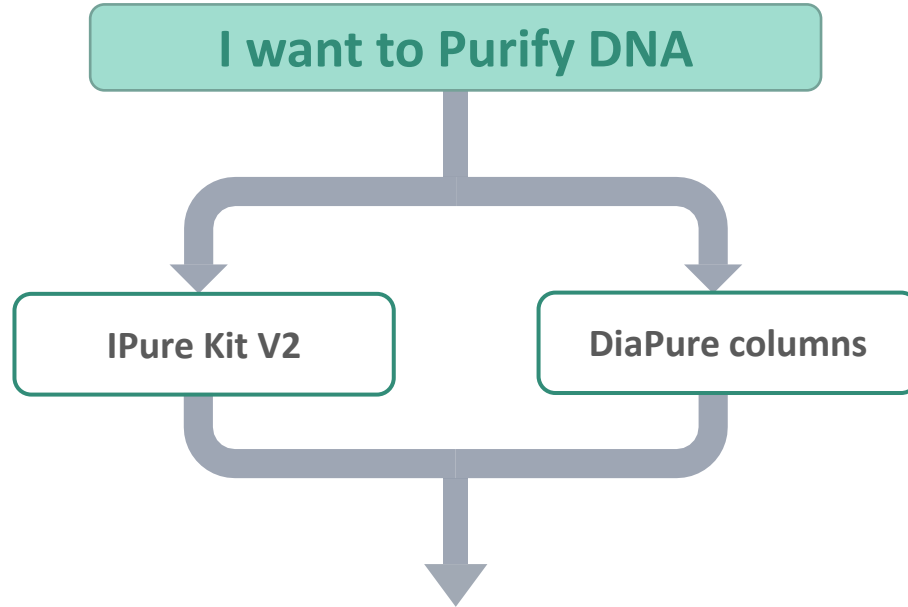
Analyzing Fragment Size

- **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





DNA Purification



Provides pure DNA for any downstream application (e. g. NGS)



DNA Purification

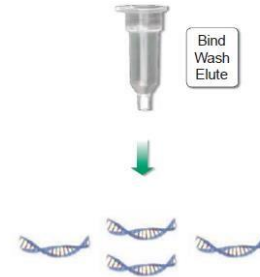
IPure Kit v2

- Best yields
- Recovery of small amounts of DNA
- No toxic reagents (e.g. phenol/chloroform)
- Compatible with automation



MicroChIP DiaPure columns

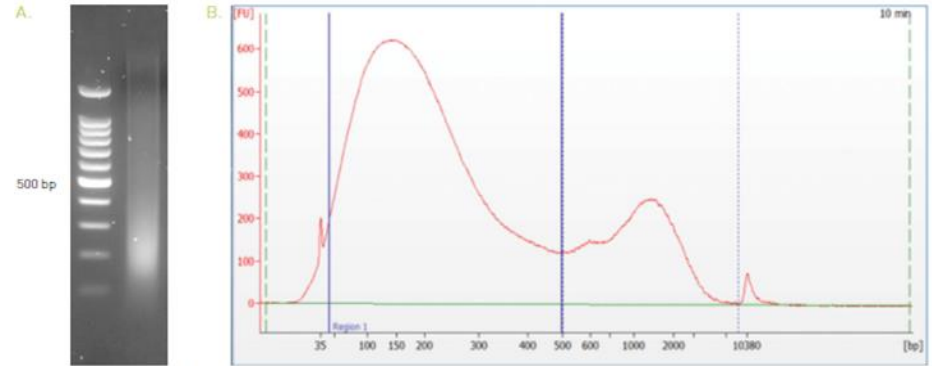
- Perfect for low concentrated samples (elution from 6 μ l)
- DNA recovery 70-90% (50bp - 10kB)
- No toxic reagents (e.g. phenol/chloroform)





Analyzing fragment size

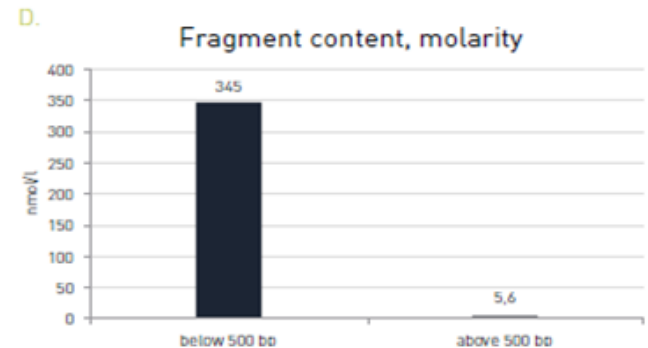
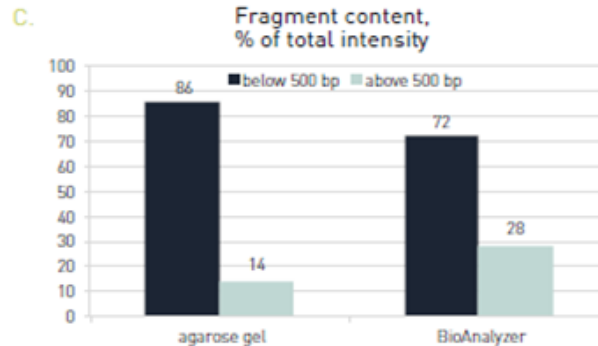
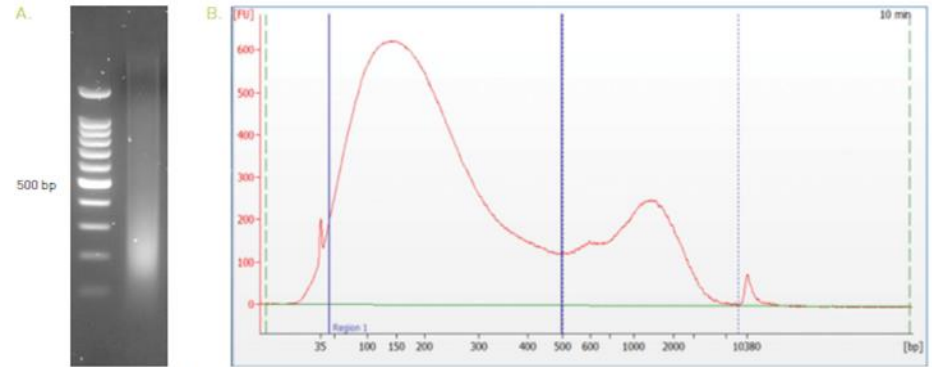
- Use agarose gel or fragment analyzer
- Bioanalyzer or Tapestation:
 - Over-representation of HMW fragments
 - Log-based -> visual misinterpretation of fragment distribution
 - More sensitive to overloading, incomplete reverse crosslinking and residual contaminants





Analyzing fragment size

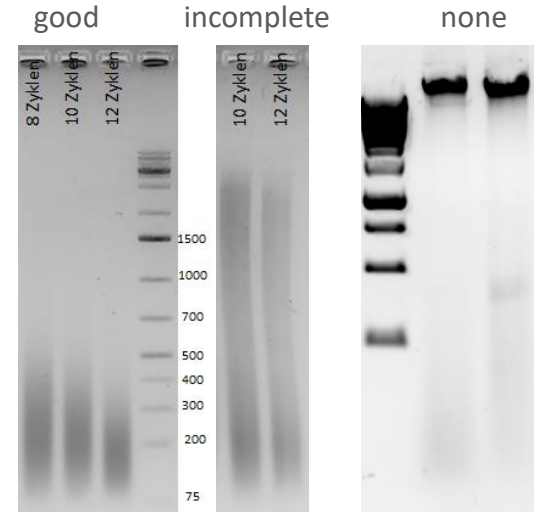
- Use agarose gel or fragment analyzer
- Bioanalyzer or Tapestation:
 - Over-representation of HMW fragments
 - Log-based -> visual misinterpretation of fragment distribution
 - More sensitive to overloading, incomplete reverse crosslinking and residual contaminants





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



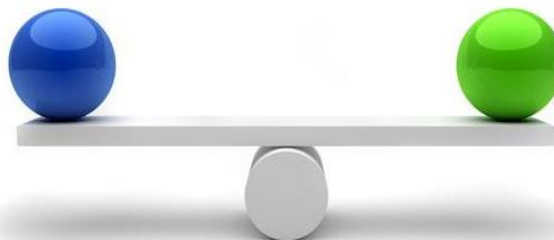


Secrets of ChIP Success

- 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

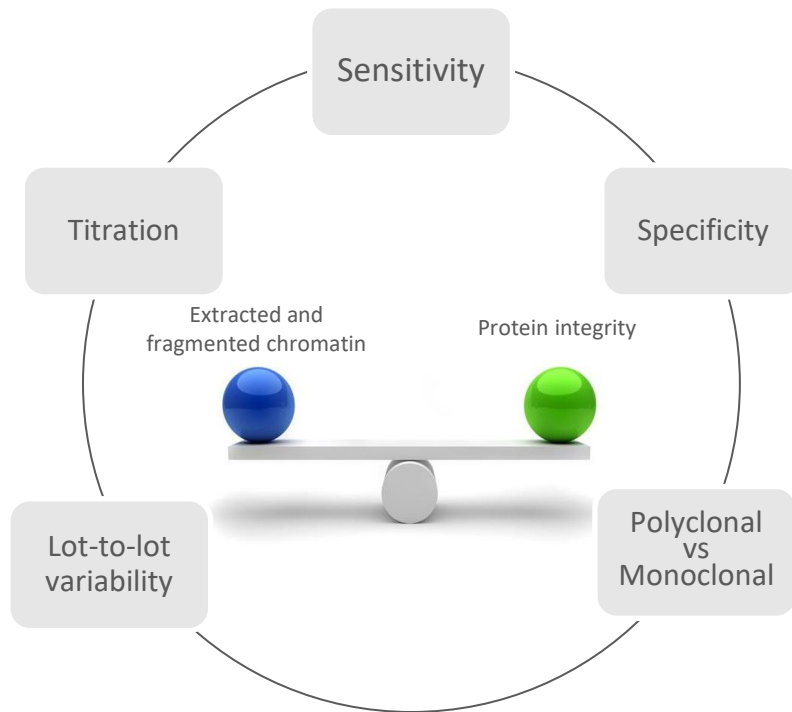
Extracted and
fragmented chromatin

Protein integrity





Antibodies for ChIP



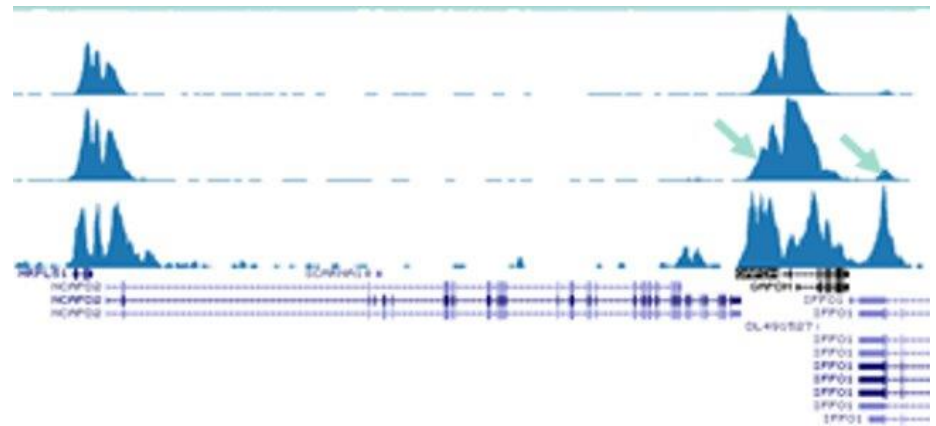
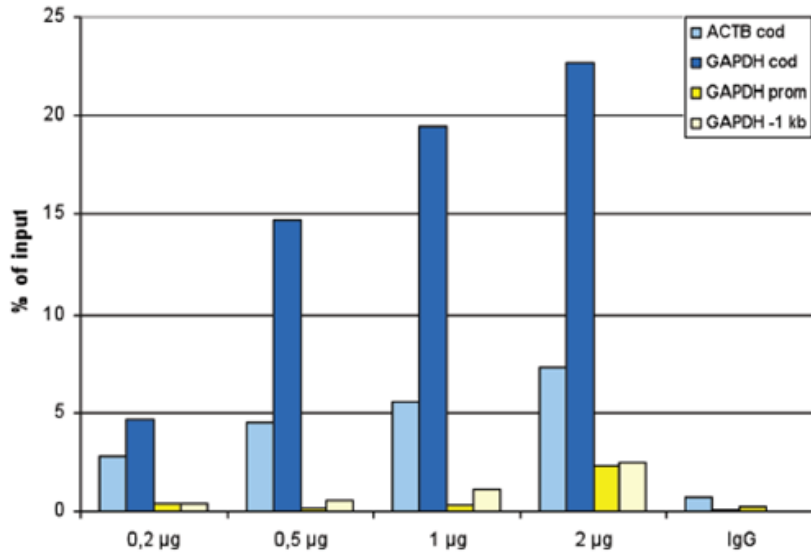


Antibodies for ChIP

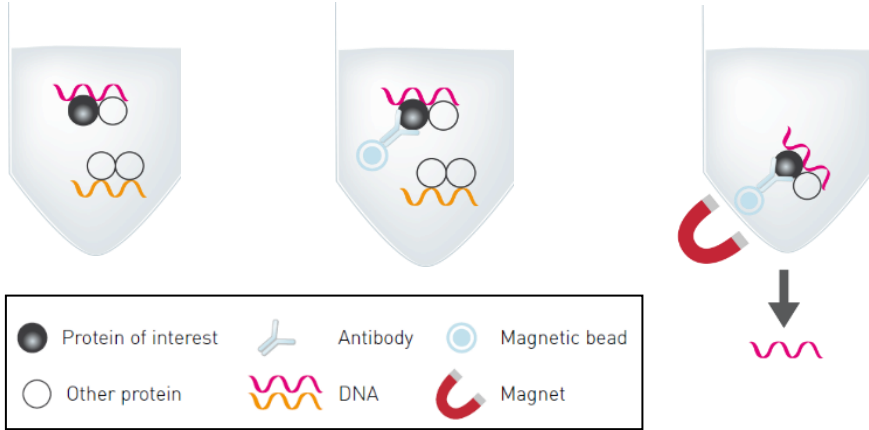
Polyclonal H3K36me3 antibody titration (Diagenode C15410192)

1 μ g 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

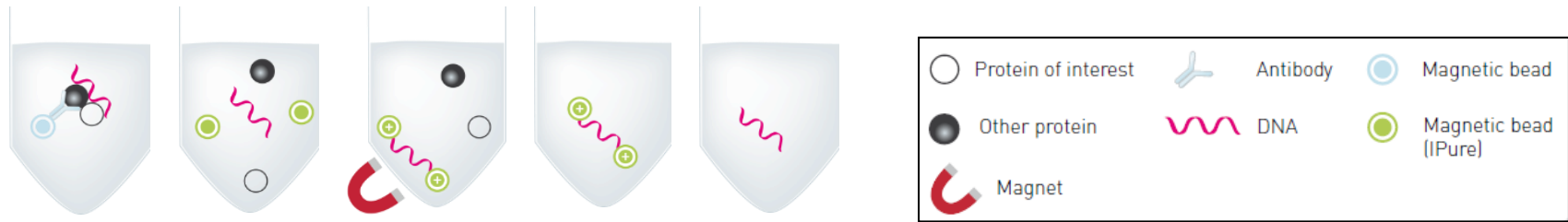


Setting Up IP: Additional Controls

- 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



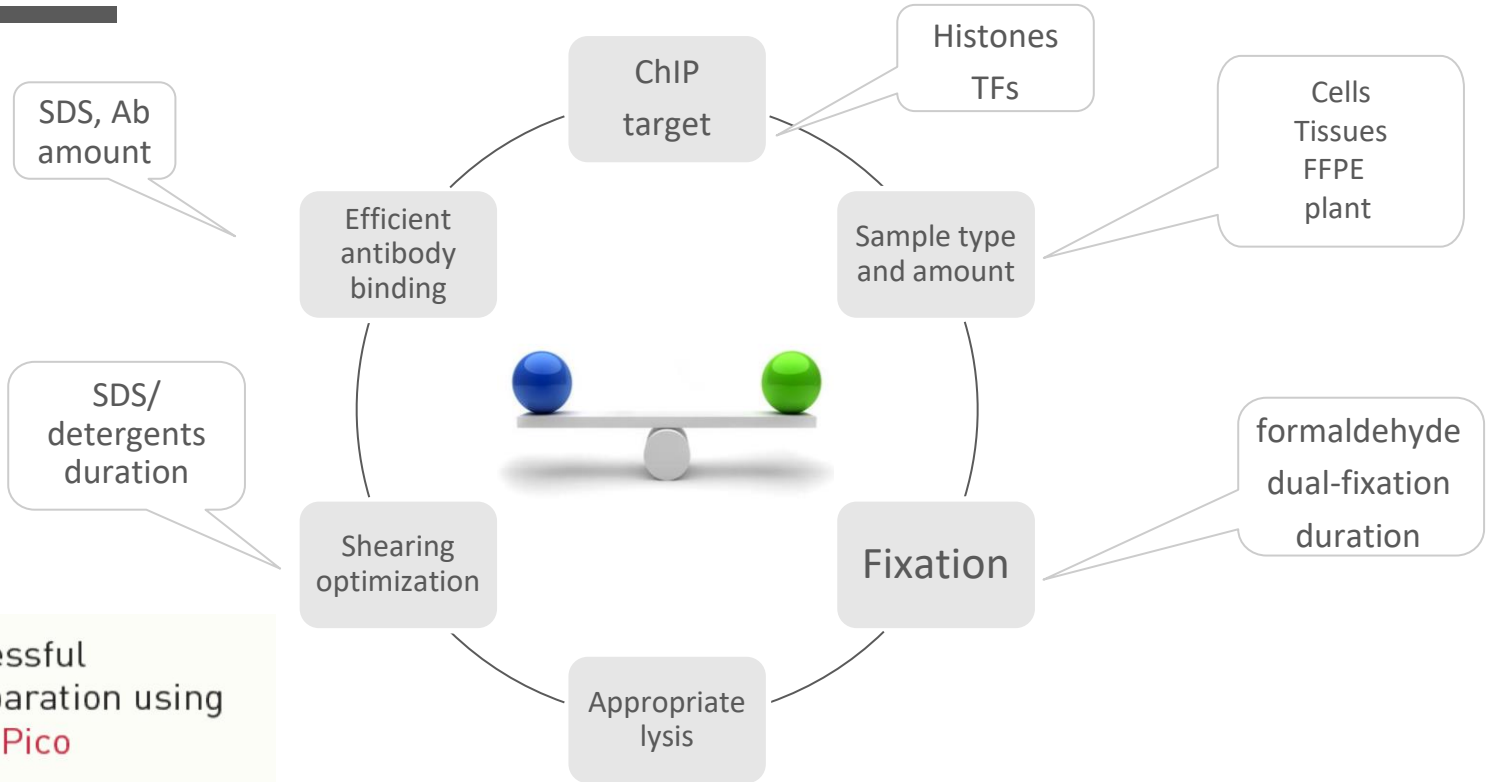
ChIP Protocol – Elution, de-crosslinking and DNA isolation



- 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**



Fastest Fingers Quiz





Fastest Fingers Quiz

Can I use formaldehyde that has been open for 3 months for fixing cells?

- A. Yes, it is not an issue as long as you make a fresh dilution
- B. No, open fresh bottle





Fastest Fingers Quiz

I am working with tissues, but I do not have time to shear chromatin today. What should I do?

- A. Fix tissues, lyse and freeze chromatin - shear later
- B. Fix tissues and freeze - lyse and shear chromatin later
- C. Freeze fresh tissue - fix, lyse and shear chromatin later

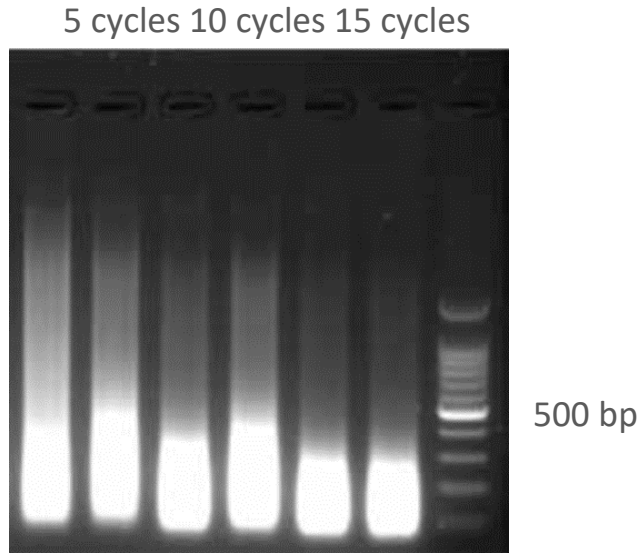




Fastest Fingers Quiz

Which cycle in the image would you use for chromatin shearing?

- A. 5 cycles
- B. 10 cycles
- C. 15 cycles





Innovating Epigenetics Solutions



WELCOME TO DIAGENODE

ChIP Workshop

Jessica Apulei & Juri Kazakevych

16-17 February 2021

OBJECTIVES

Day 1

- CHIP overview
 - Chip-qPCR vs. ChIP-Seq
 - Chromatin preparation
 - Fixation, Cell lysis and Chromatin shearing
- Setting up IP
 - Antibodies, Replicates, inputs, controls

Day 2

- ChIP-qPCR: set-up
- ChIP-qPCR: analysis
- ChIP-seq: library prep & sequencing
- ChIP-seq: analysis
- Overview of Alternate methods



ChIP-qPCR

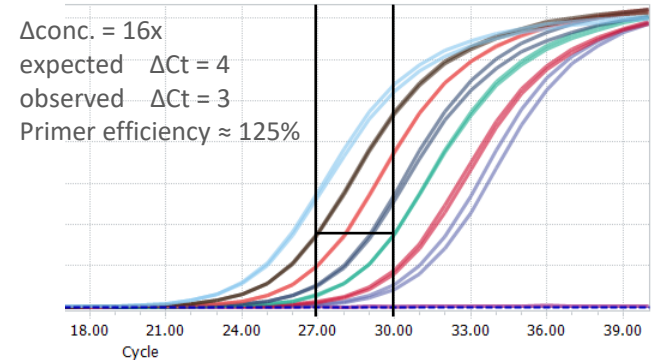
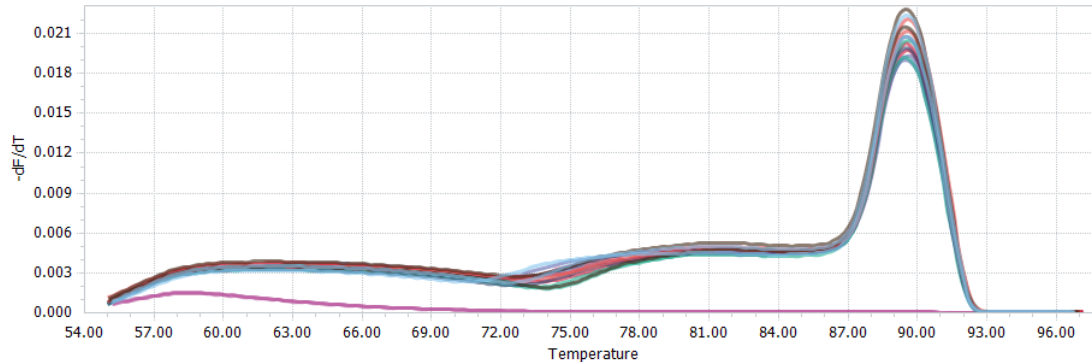
- Target & Primer selection is key for ChIP-qPCR
- Predict qPCR-targets from ChIP-seq data
- qPCR as QC prior ChIP-seq
- If no ChIP-seq data:
 - estimate binding from similar data, biological function etc.
 - use multiple regions
- PCR program depends on Master Mix, qPCR system and primer pairs





Setting up ChIP-qPCR

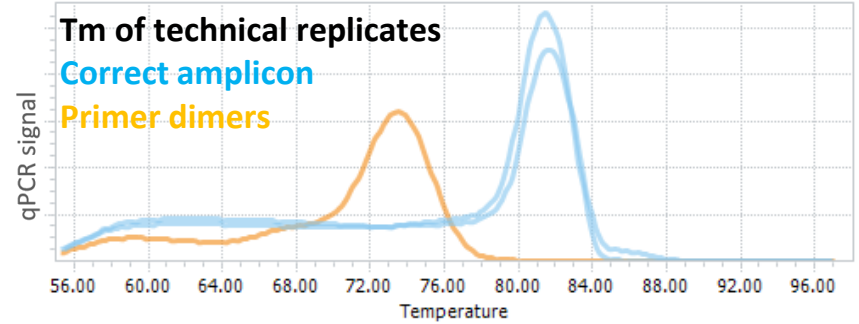
- **Target primer design:**
 - Place primers around binding site
 - 50-150bp amplicons
 - 20-30 bp primers with a similar T_m between 55° and 60°C
- **Primer pair validation:**
 - Check on gDNA/input for T_m profile
 - Check efficiency (95-105% acceptable)



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

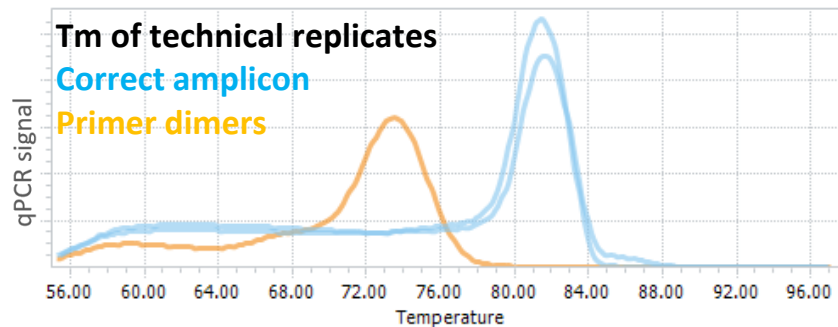




Setting up ChIP-qPCR

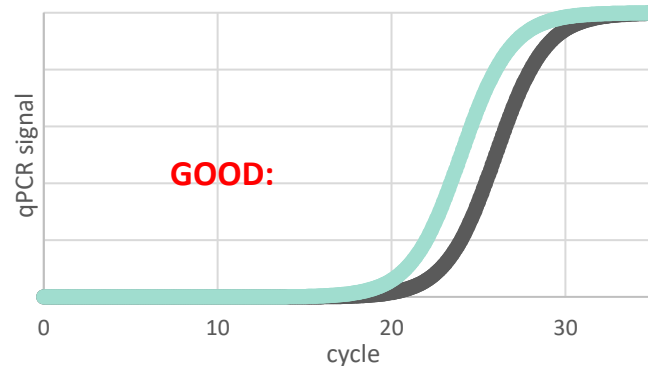
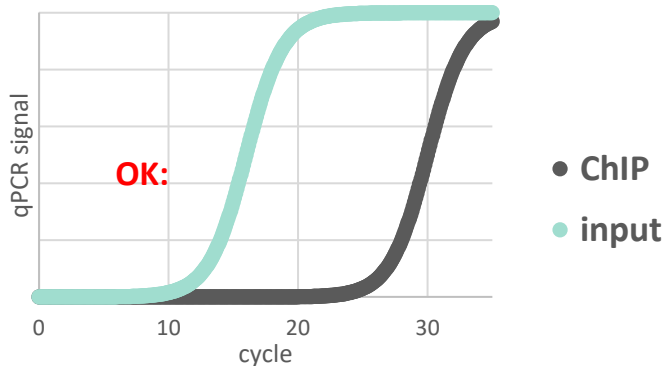
QC

- 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
- $Ct(H3K4me3) \ll Ct(IgG)$



ChIP-qPCR Exercise

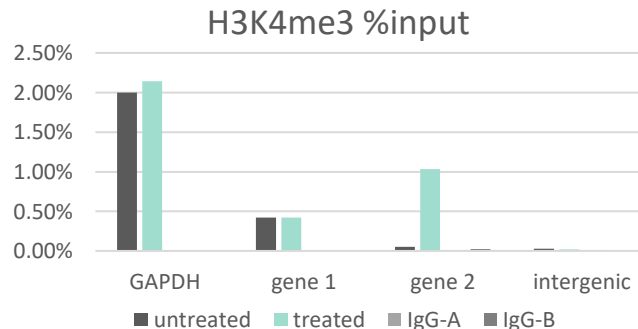
Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

	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%





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

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

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

6. Assessment

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

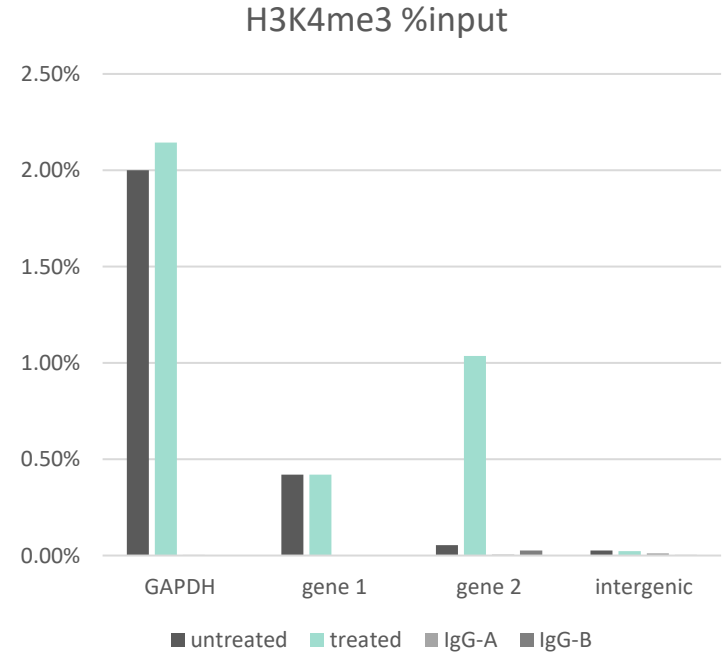
7. Optimization

- 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





Fastest Fingers Quiz





Fastest Fingers Quiz

You did a histone ChIP-qPCR and got 1.5-fold enrichment. Should you go forward with seq?

A. Yes

B. No





Fastest Fingers Quiz

You performed a CHIP on a novel Transcription factor. You do not see an enrichment of positive control regions over negative (intergenic) ones. Which could be the reasons? (select all applicable)

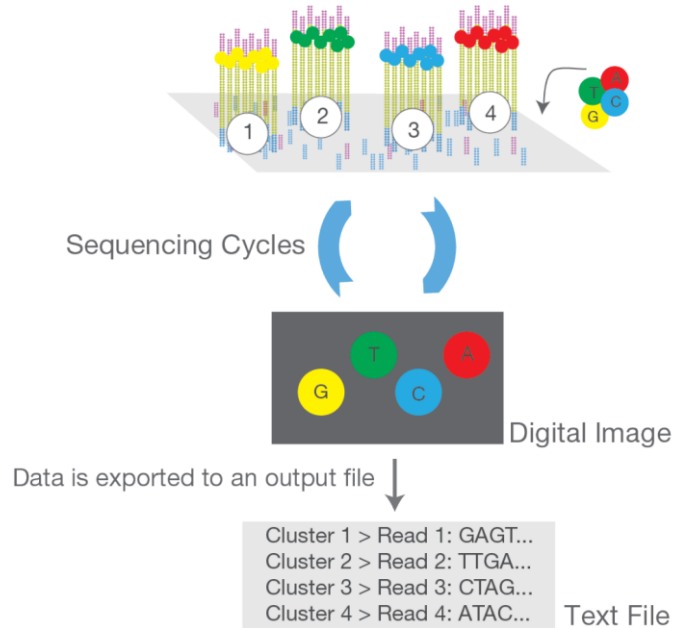
- A. antibody used is not suitable for CHIP
- B. CHIP ok but positive control regions are not bound by the TF
- C. fixation protocol not adequate for targeted TF
- D. Covid19





ChIP-seq: library prep

C. Sequencing



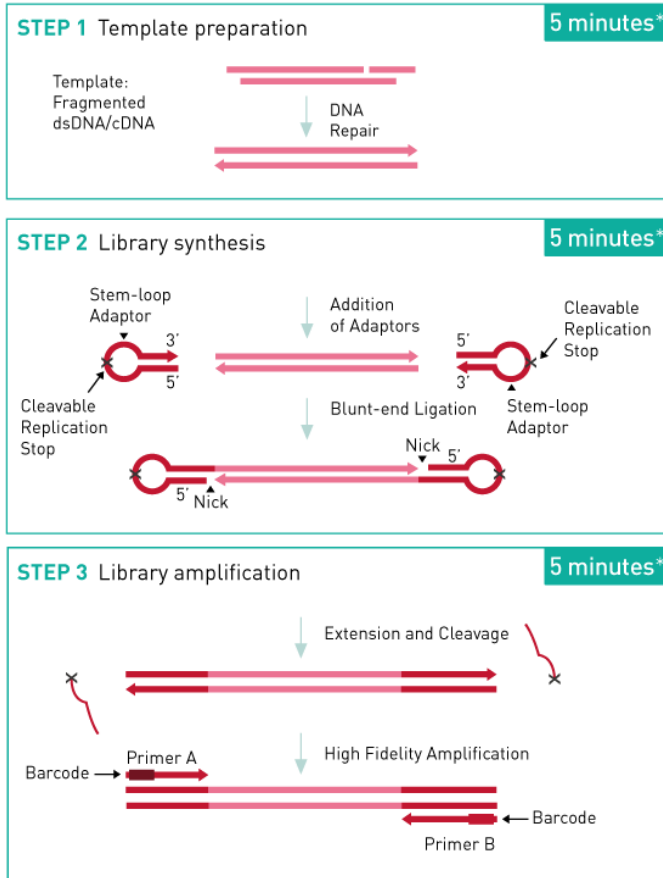
ChIP-seq: Library Prep



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



MicroPlex kit workflow



* hands-on-time



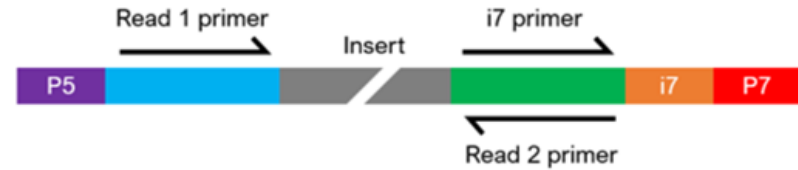
Benefits of Multiplexing

- **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

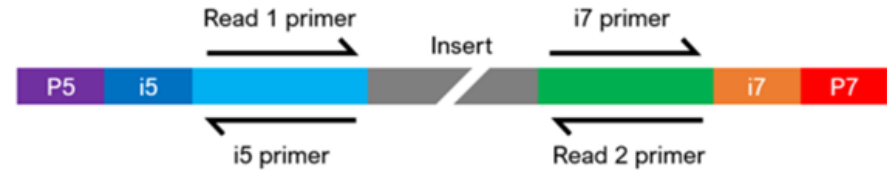


Single and Dual-indexed Libraries

- **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



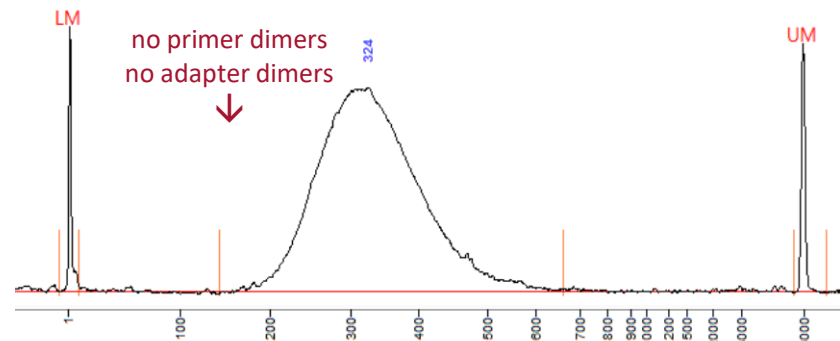
Dual-Indexed Sequencing





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





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

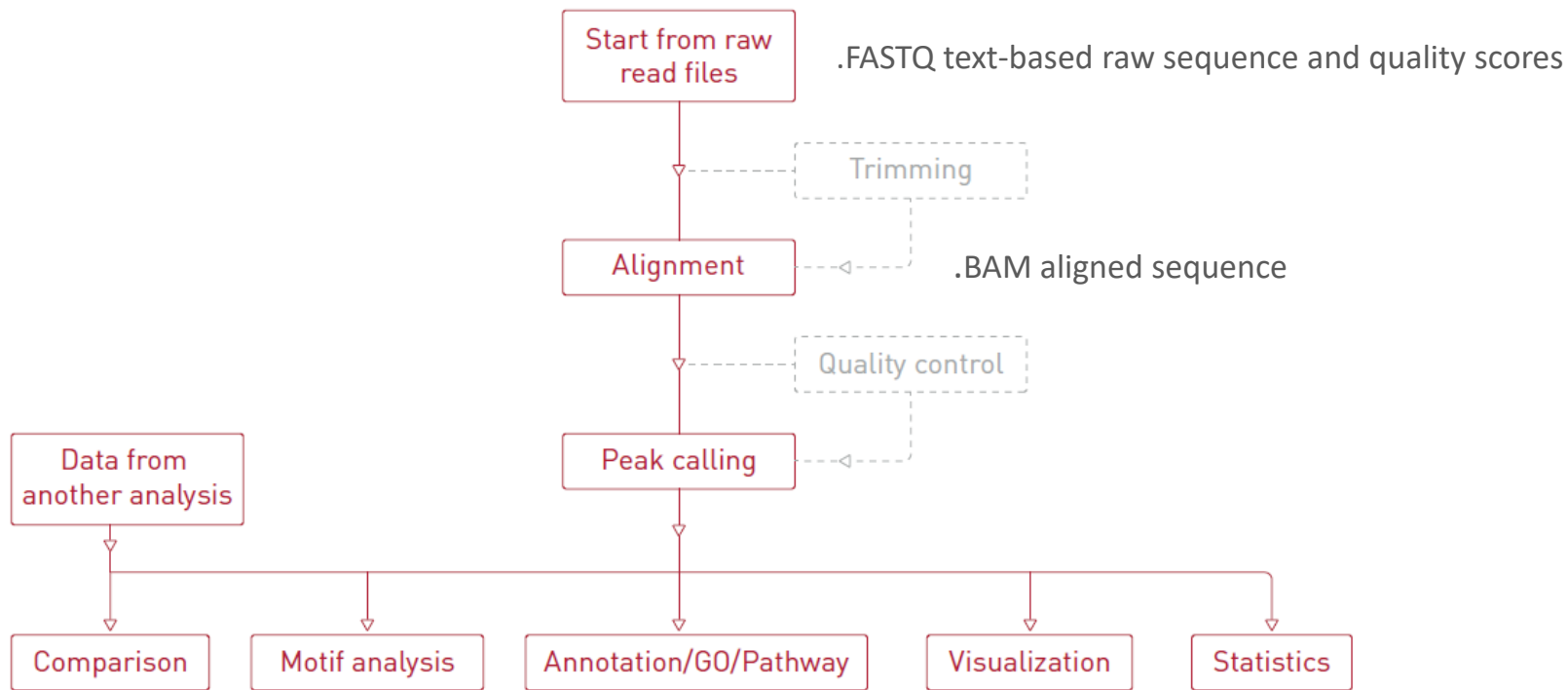


ChIP-seq: Analysis

Bioinformatician	R	https://www.r-project.org/
	free-ware and online tool kits	www.bioinformatics.babraham.ac.uk/projects/seqmonk/ biit.cs.ut.ee/gprofiler/gost
Wet-lab expert with free time	standard bio-informatic services	https://www.diagenode.com/en/categories/Services Comprehensive Multi-Omic and bio-info services
	free-ware and online tool kits	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



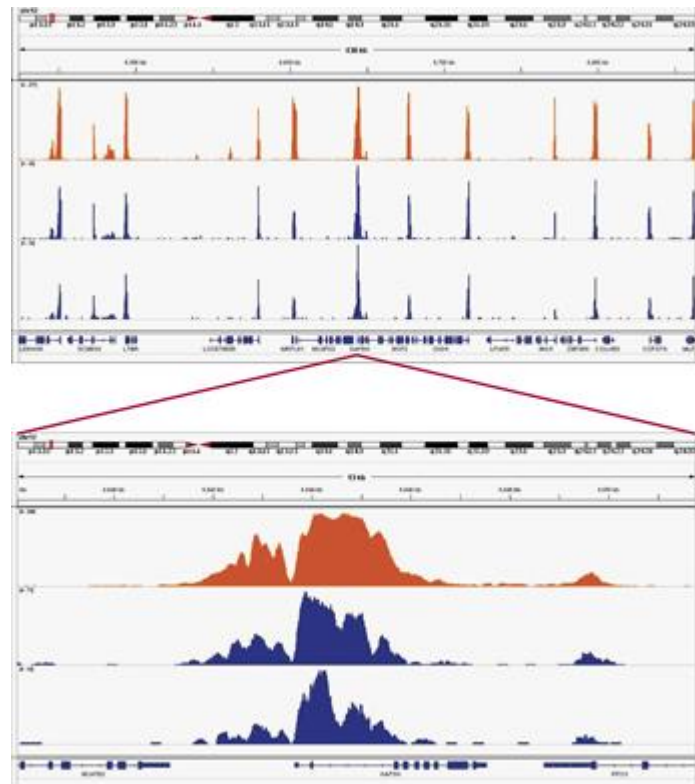
ChIP-seq: Analysis

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





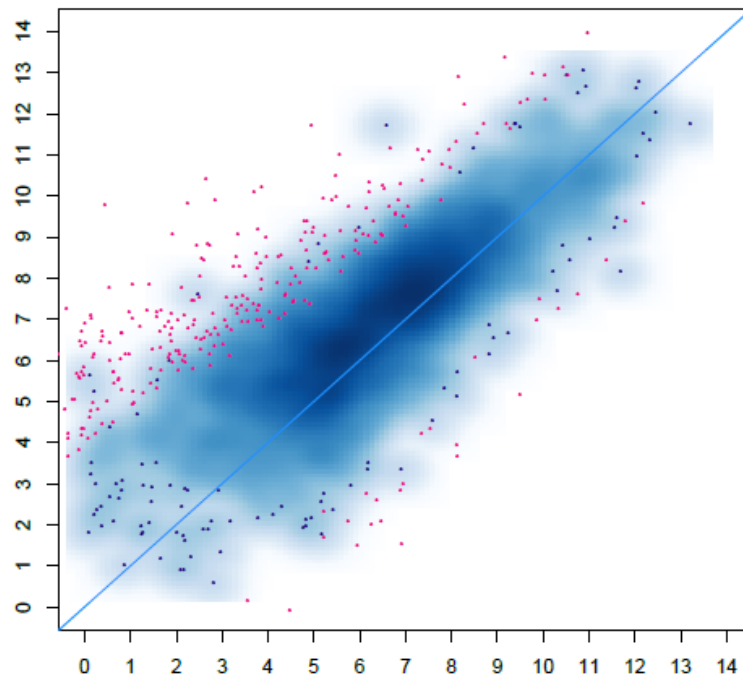
ChIP-seq: Analysis

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





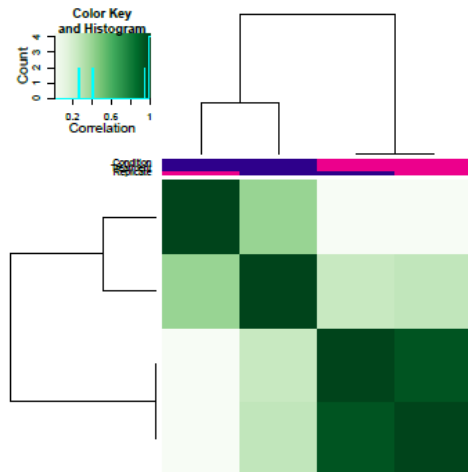
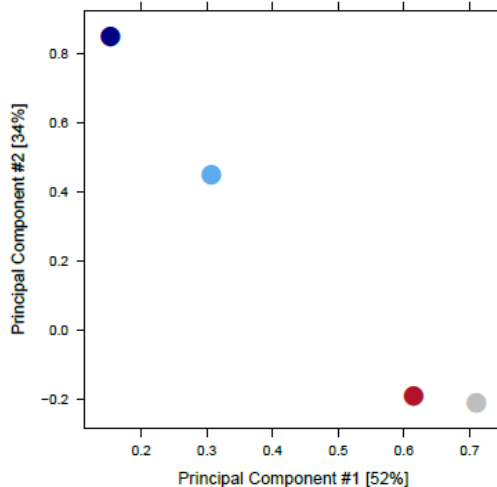
ChIP-seq: Analysis

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



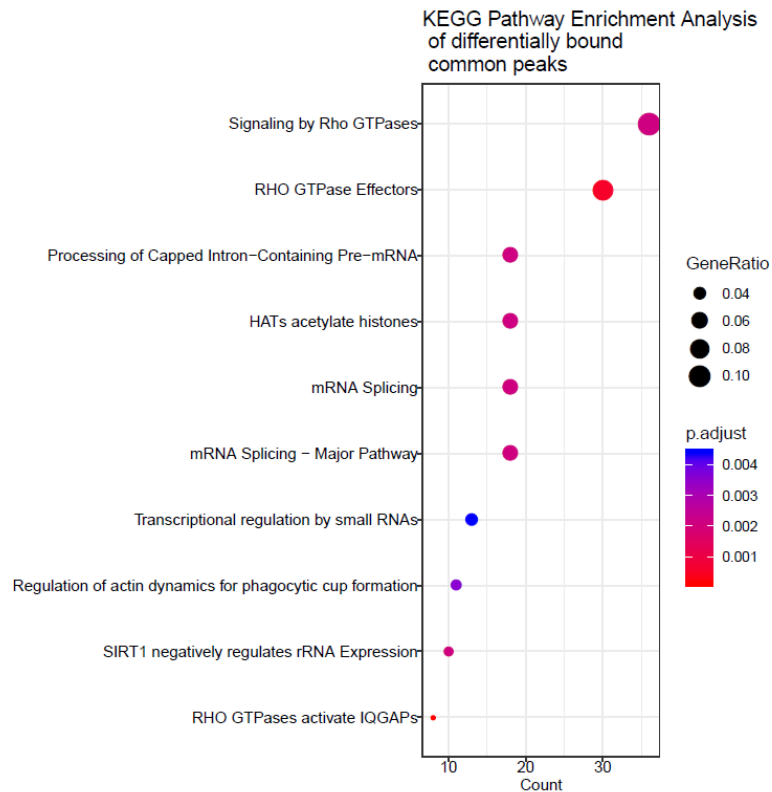
ChIP-seq: Analysis

Standard bioinformatic analysis:

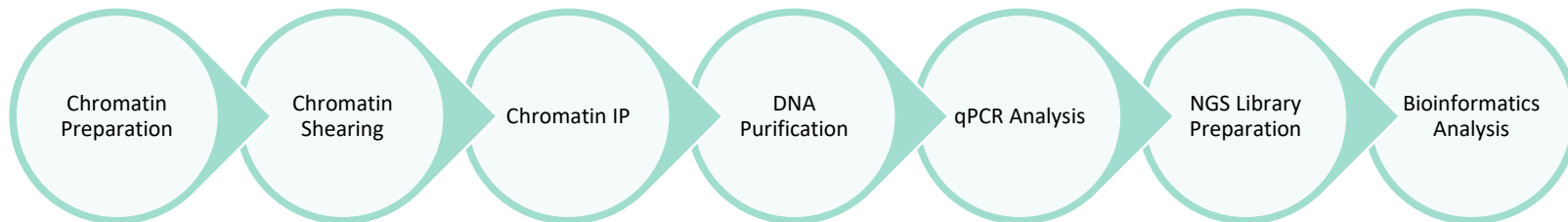
- alignment to reference genome
- peak calling

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



Hardware:	Bioruptor	IP-Star		IP-Star
Reagents:	Cross-link Gold	ChIP-seq grade antibodies	IPure kit MicroChIP DiaPure columns	Primer Pairs
Kits:	EasyShear kits			
	iDeal ChIP kits			MicroPlex kit or TAG kits
	True MicroChIP kit			
	ChIPmentation kits			
Services:	Epigenomic Profiling Services			
				Data Analysis Service



Other Methods to Study Chromatin

- Interaction between proteins and DNA (immuno- assays)
 - ChIPmentation
 - Cut&Run and Cut&TAG

- Methods to study chromatin accessibility (non-immuno assays)
 - ATAC-seq



ChIPmentation™ & μChIPmentation™

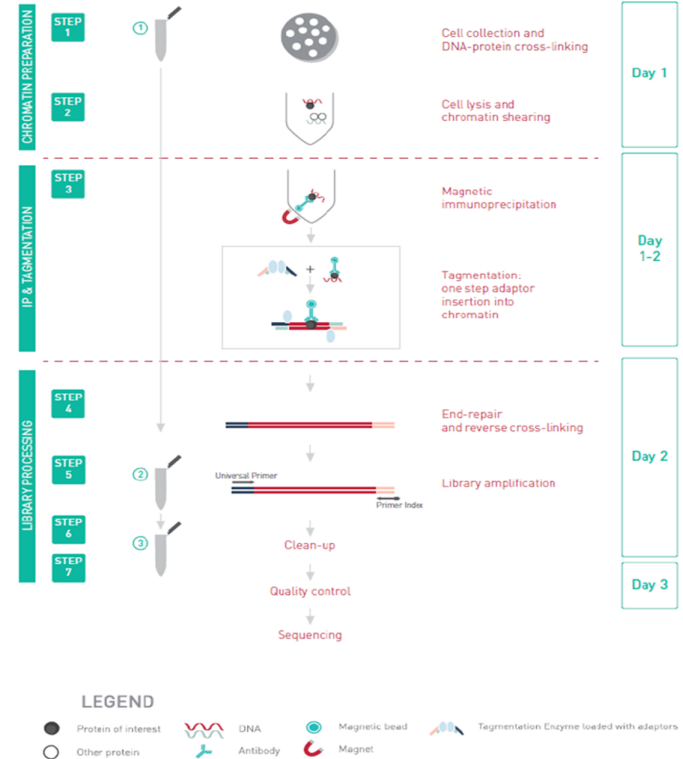
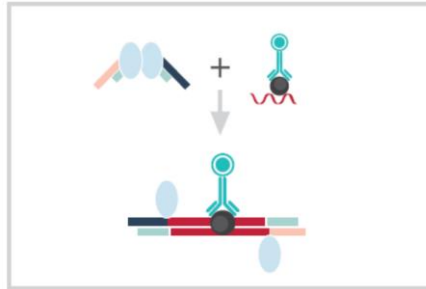
Easier and faster than classical ChIP-seq

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

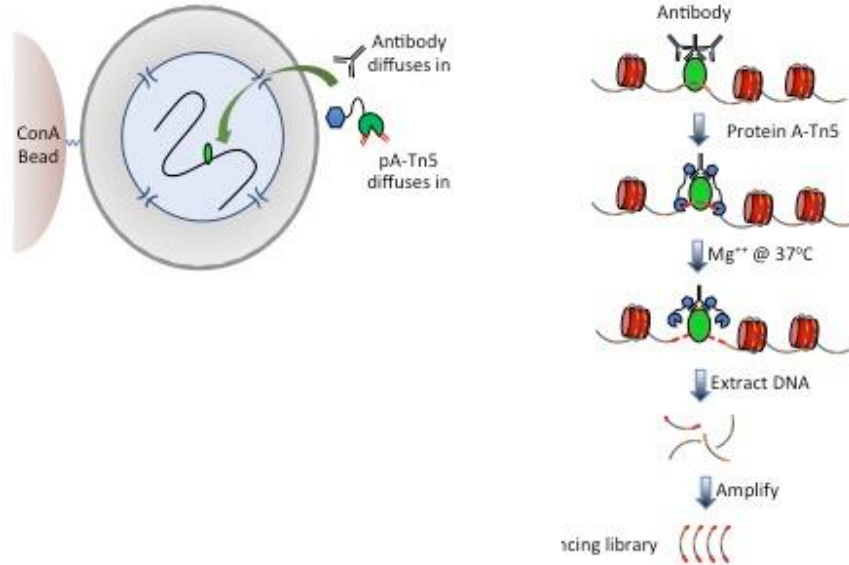
μChIPmentation for 10,000 cells





CUT&Tag: Cleavage Under Targets and Tagmentation

CUT&Tag (Cleavage Under Targets & Tagmentation)



Key features:

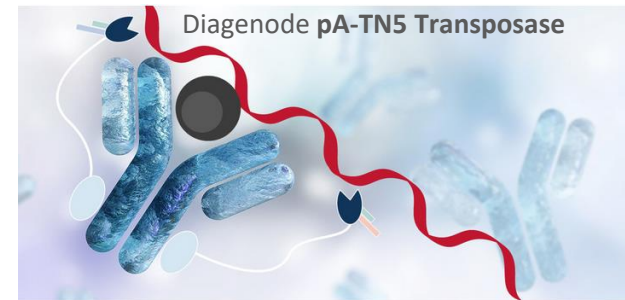
Crucial reagent:

- proteinA-Tn5

Fast and easy protocol:

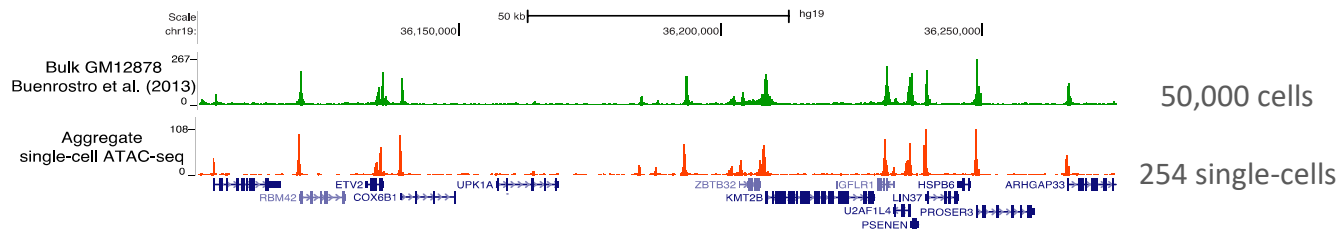
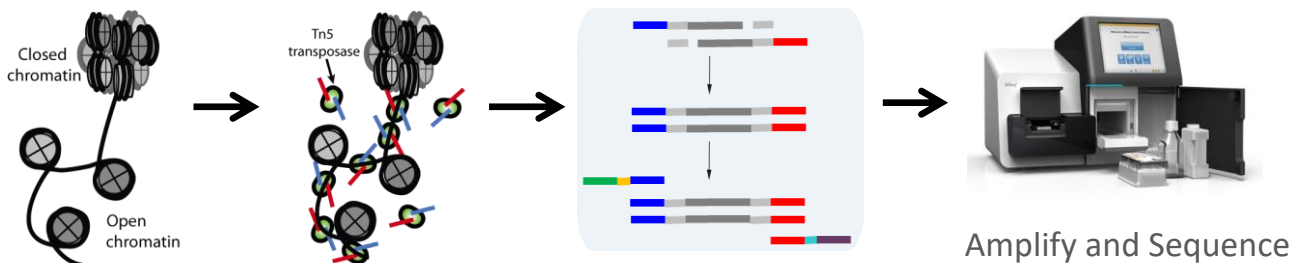
- fast tagmentation-based library prep
- No chromatin prep

Suitable for low cell numbers





Assay for Transposase-Accessible Chromatin



Buenrostro et al., Nature, 2013

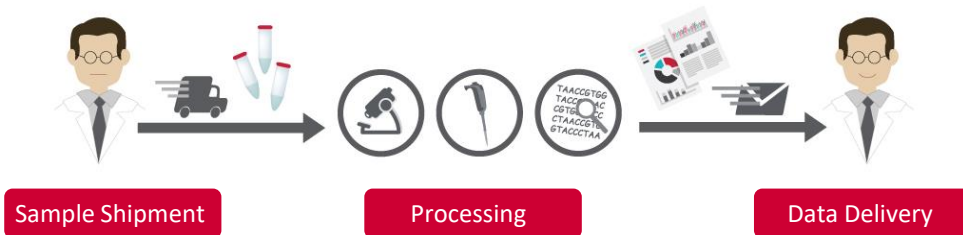
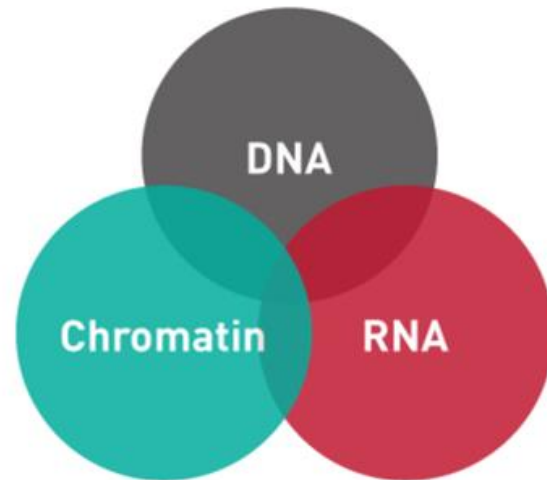




EPIGENOMICS PROFILING SERVICES

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 :-))

www.diagenode.com

