



Innovating Epigenetics Solutions



WELCOME TO DIAGENODE

ChIP Workshop

Jessica Apulei, Ekaterina Gracheva & Juri Kazakevych

June 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

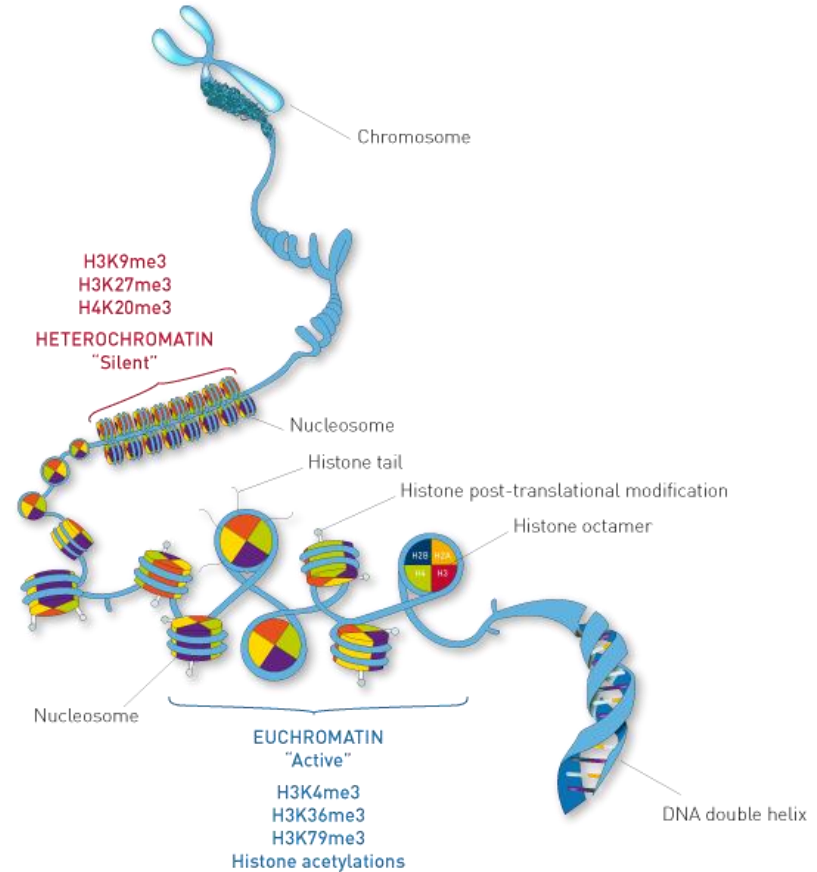
Day 2

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



What is Chromatin?

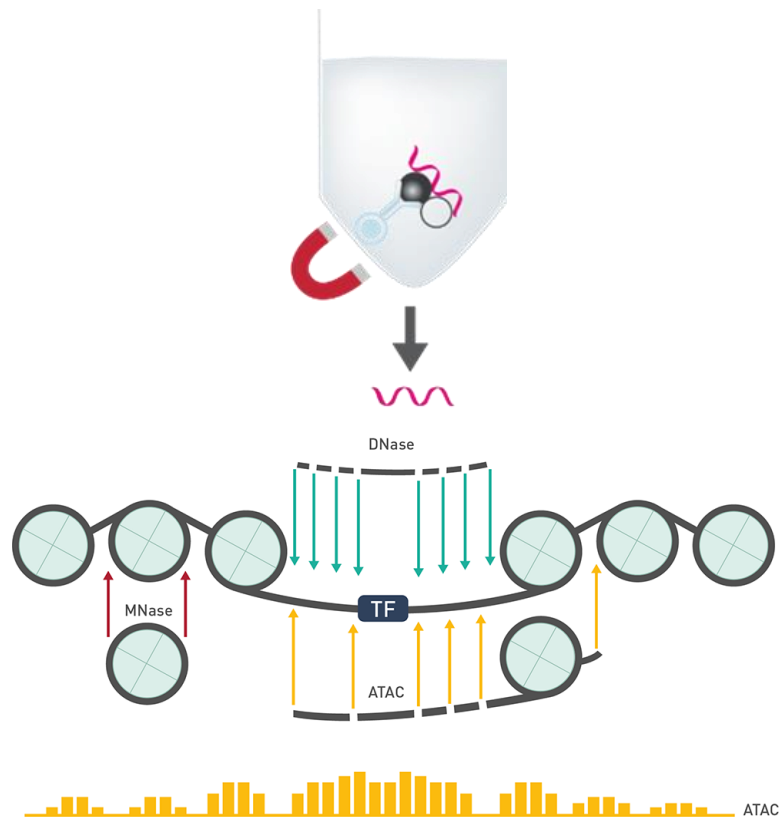
A complex of DNA and proteins found in eukaryotic cells



-

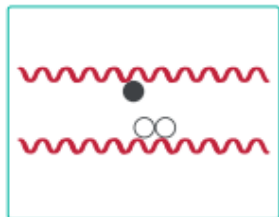
CHROMATIN ANALYSIS

- 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



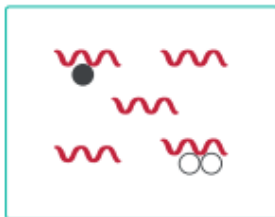
Workflow: Chromatin ImmunoPrecipitation (ChIP):

Step 1



Cross link to fix
proteins to DNA

Step 2



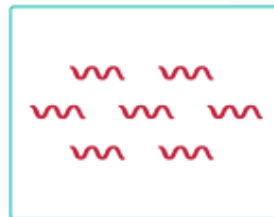
Shear chromatin

Step 3



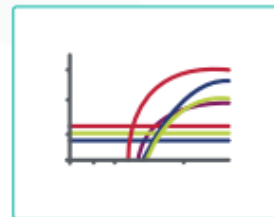
Immunoprecipitate
with antibody and
magnetic beads

Step 4



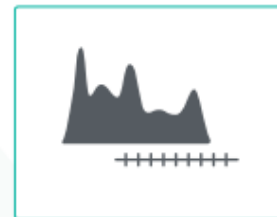
Reverse crosslinks
and purify

Step 5



Analyze by qPCR

Step 6



Prepare the libraries
for NGS

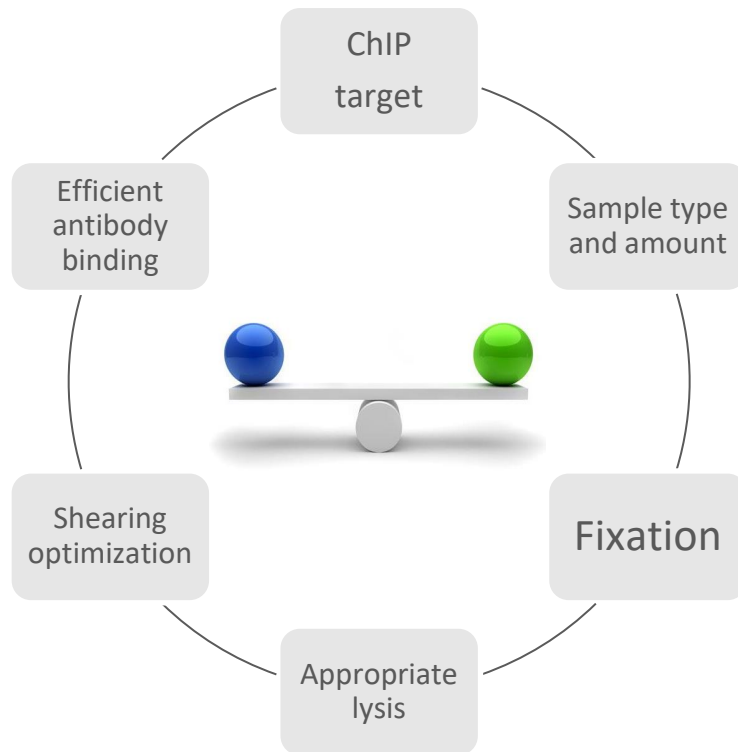


ChIP-qPCR or ChIP-Seq?

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



Summary – Tips to Prepare Good Chromatin



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



Starting material: Cells and Tissues

	Standard	Low-input
Cells	Histones: 1 million/IP TF: 4 million/IP for TF (from 100.000 cells depending on target) iDeal ChIP kits	Histones only 10-100k/IP True MicroChIP-seq Kit
Tissue	soft tissues: Dounce homogenization (liver, brain) hard/frozen/fibrous: bead beater like TissueLyser Histones: 5-7 mg/IP TF: 30mg/IP (from 1.5mg/IP depending on target) iDeal ChIP kits	Histones only amounts depend on tissue type
FFPE-tissue	Challenging due to extensive crosslinking de-paraffinization with Heptane instead of xylene -> easier, non-toxic workflow from 300ng DNA per IP iDeal ChIP-FFPE kit	
Plant tissue	65mg - 2g / chromatin preparation , depending on sample Universal Plant ChIP-seq kit	

Low-input ChIP

True MicroChIP-seq kit

10k-100k cells/histone-IP

10k-700k cells /chromatin prep

Suitable for FACS-sorted samples

Single step lysis & minimal handling

LEGEND



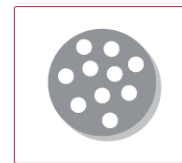
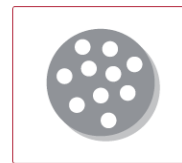
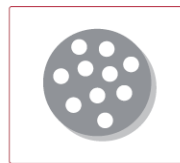
Protocol for:

Batch

Individual
samples

FACS-sorted
cells

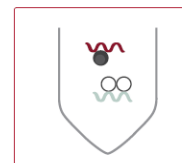
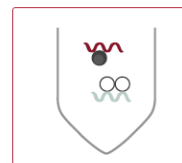
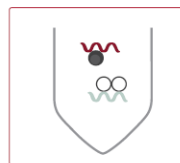
STEP
1



Cell or tissue collection
and DNA-protein
cross-linking

30 minutes
to 1 hour

STEP
2



Cell lysis and
chromatin shearing

1 to 2
hours

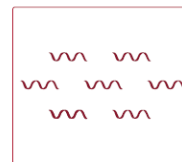
STEP
3



Magnetic
immunoprecipitation

Overnight

STEP
4



Elution, decross-linking
and DNA purification

5 hours



Fixation

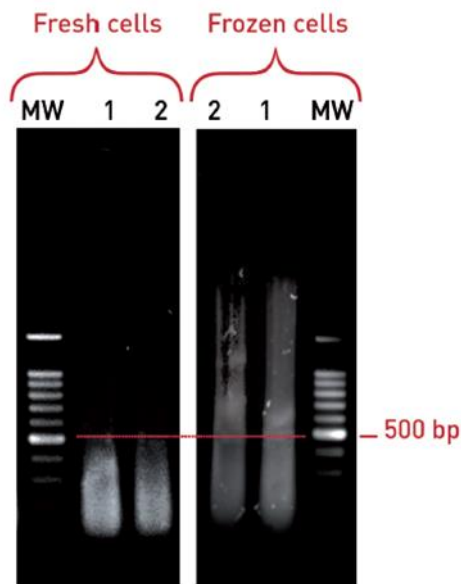
- Covalent stabilization of protein-DNA interactions; Reversible
Directly in medium for weak or rare protein-DNA interaction
For histone marks, cells can be resuspended by trypsinization before fixation
- Common fixative: Formaldehyde
Fresh
Methanol-free not mandatory

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



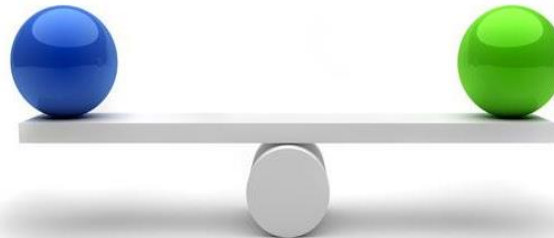
Optimal: Perform ChIP workflow directly, without freezing

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

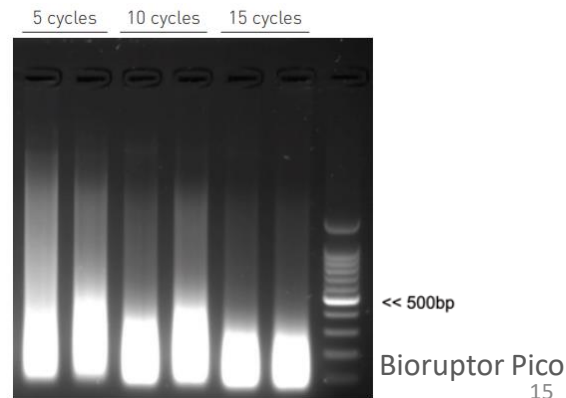
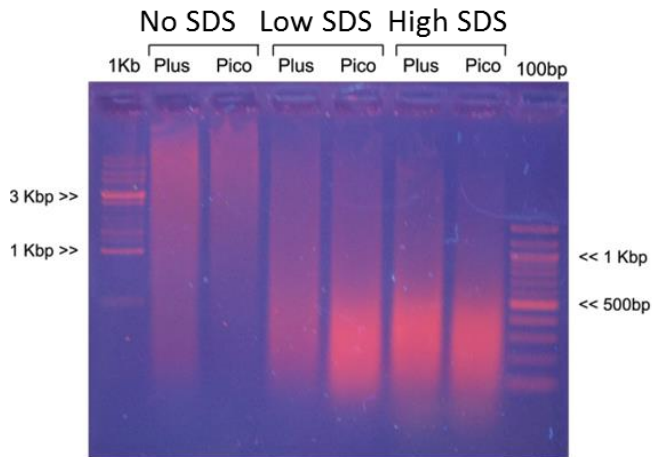
Extracted and
fragmented chromatin

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 prep for ChIP on	Sample type	Sample input per IP	Kit	SDS concentration	Nuclei isolation	Compatible with ChIP kits
Histones	Cells	10-100k	Chromatin EasyShear Kit High SDS	1%	No	True MicroChIP-seq Kit μChipmentation for Histones Kit
		> 100k				
	Tissue		Chromatin EasyShear Kit Ultra Low SDS	< 0.1%	Yes	iDeal ChIP-seq for Histones Kit ChIPmentation for Histones Kit
	Plant tissue		Chromatin EasyShear Kit for Plant	0.5%	Yes	Universal Plant ChIP-seq Kit
Transcription Factors	FFPE					
	Cells		Chromatin EasyShear Kit Low SDS	0.2%	Yes	iDeal ChIP-seq for TF Kit iDeal ChIP-qPCR Kit
	Tissue					iDealChIP-FFPE Kit

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

+



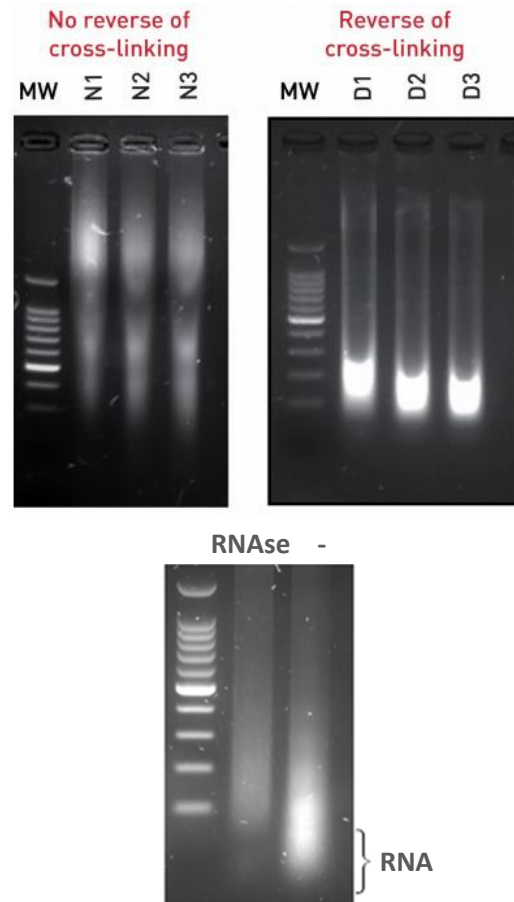
Diagenode's ChIP Kit

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

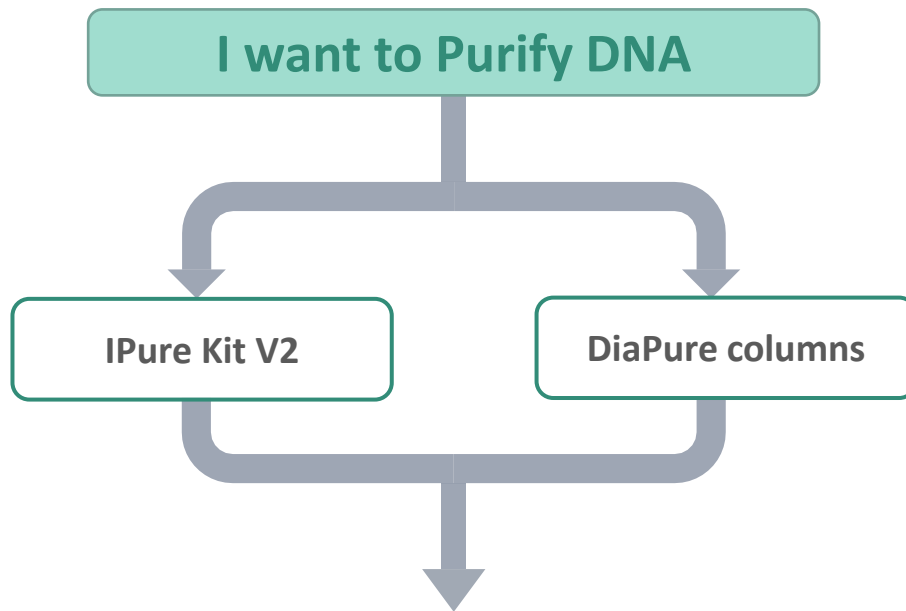
Analyzing Fragment Size

- **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

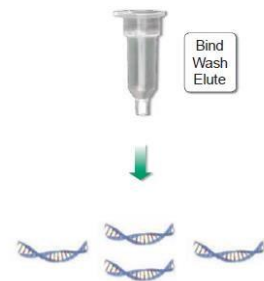
IPure Kit v2

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



MicroChIP DiaPure columns

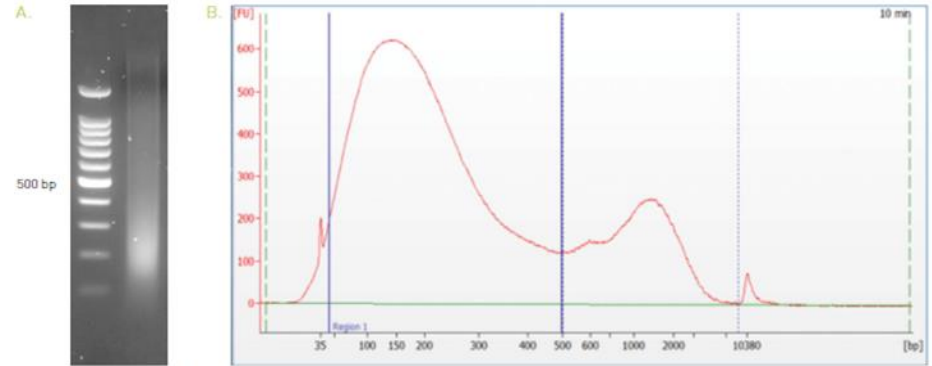
- Perfect for low concentrated samples (elution from 6 μ l)
- DNA recovery 70-90% (50bp - 10kB)
- No toxic reagents (e.g. phenol/chloroform)





Analyzing fragment size

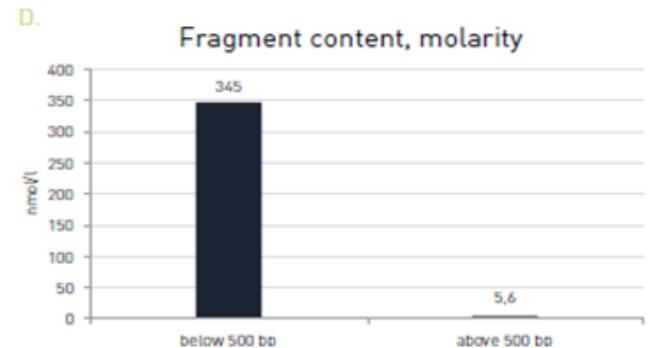
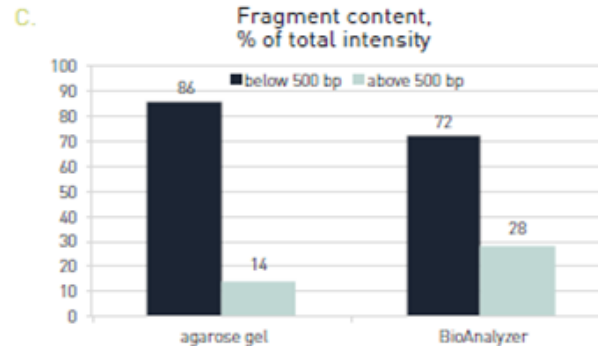
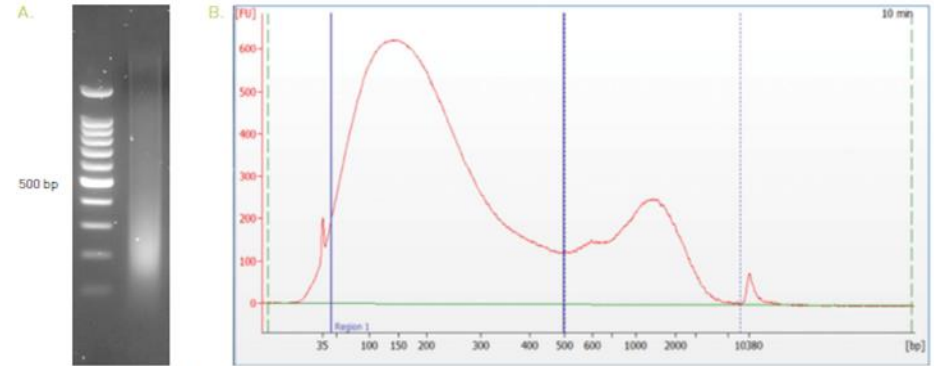
- Use agarose gel or fragment analyzer
- Bioanalyzer or Tapestation:
 - Over-representation of HMW fragments
 - Log-based -> visual misinterpretation of fragment distribution
 - More sensitive to overloading, incomplete reverse crosslinking and residual contaminants





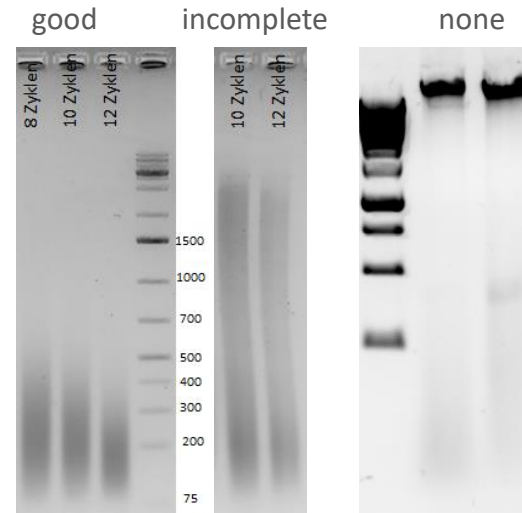
Analyzing fragment size

- 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

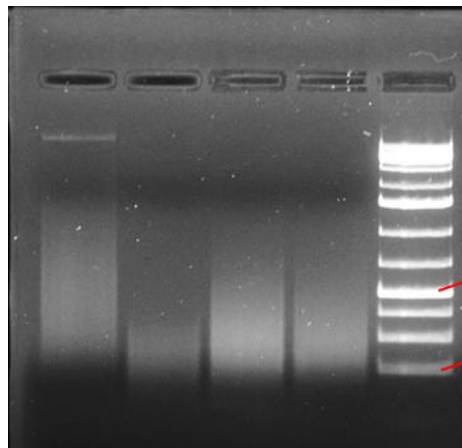


Concentration matters



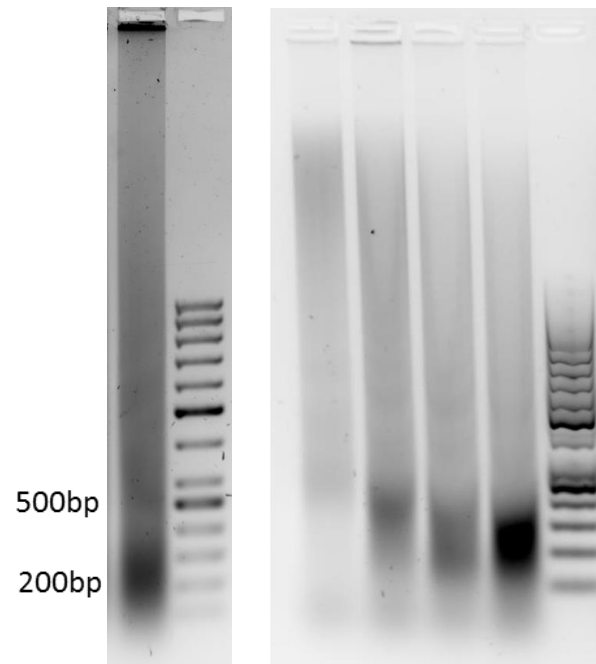
A. Thaliana root

3x-diluted		undiluted (0.5g/600μl)	
cycles: 8	12	12	15



2 Mio Monocytes

volume:	100μl	200μl			
cycles:	12	4	6	8	10

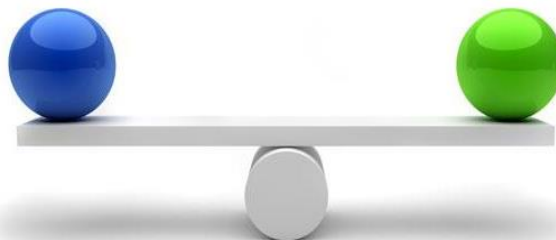


Secrets of ChIP Success

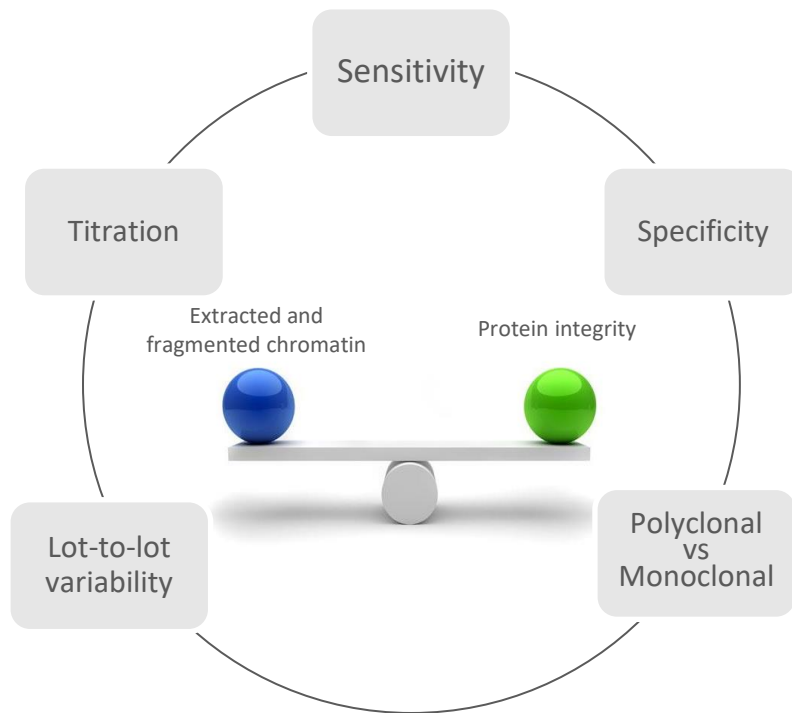
- Prepare “good” chromatin
 - Suitable fragment size and available epitopes
- Use a good antibody at the right concentration
- Optimize for highest specific signal and the lowest background

Extracted and
fragmented chromatin

Protein integrity



Antibodies for ChIP



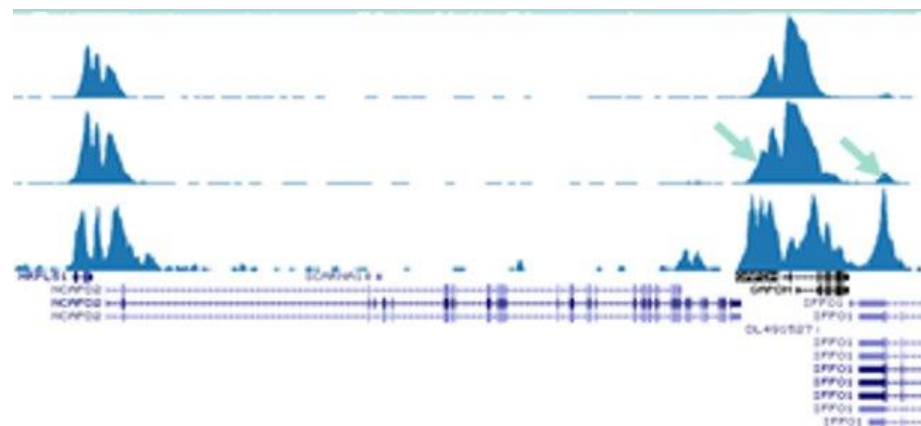
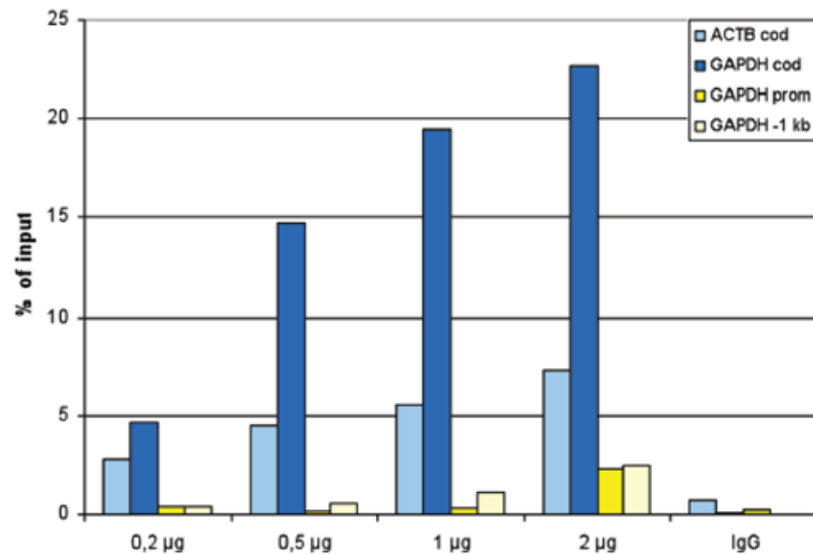


Antibodies for ChIP

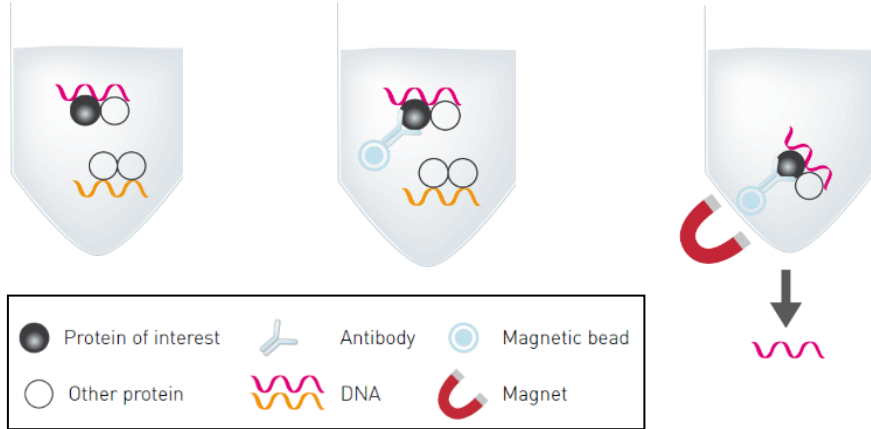
Polyclonal H3K36me3 antibody titration (Diagenode C15410192)

1µg IgG as negative IP control

Chromatin from 100.000 cells



Setting up IP



Components:

- Sheared chromatin
- **ChIP/ChIP-seq 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

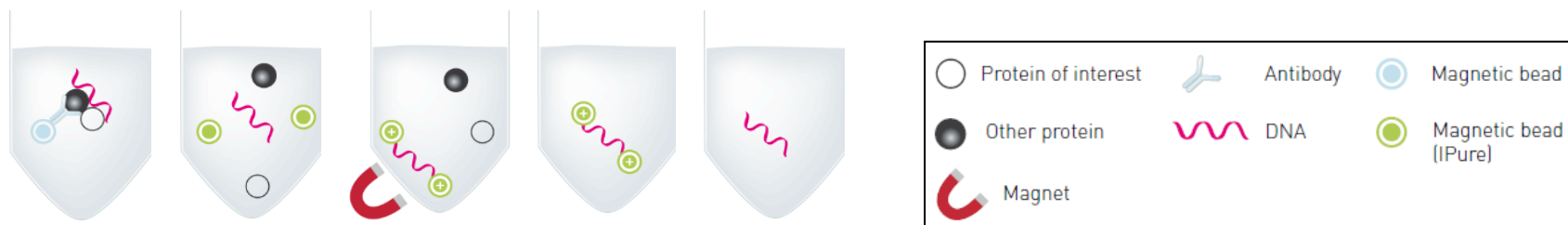
- Fraction of sheared chromatin is kept aside as INPUT
 - Processed in parallel with IP-samples from reversed crosslinking
 - Include one input for each chromatin sample
- Key reference for ChIP-qPCR and ChIP-seq analysis
- ChIP-qPCR: used to calculate the recovery (% of input)
- ChIP-seq: mandatory 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

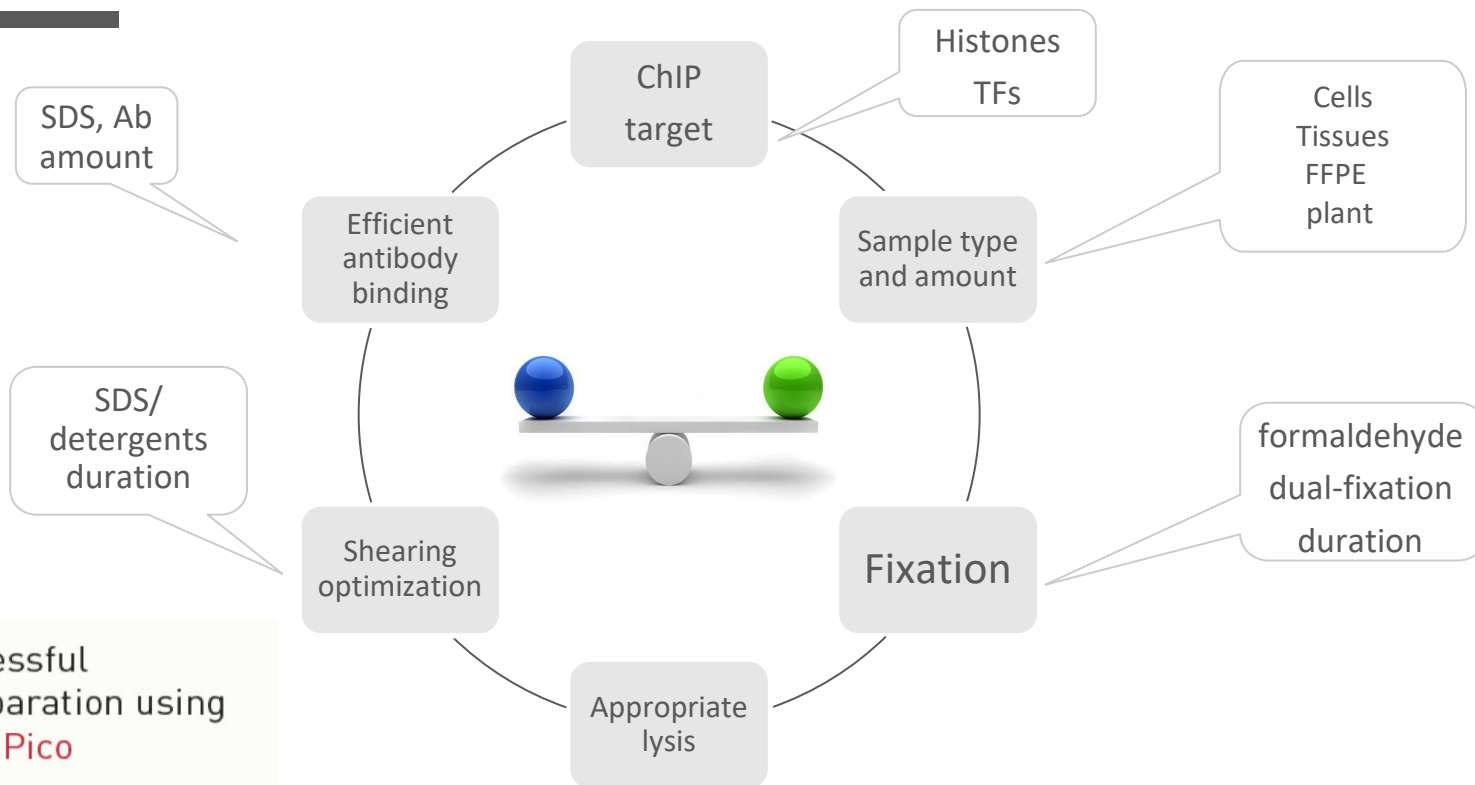
ChIP Protocol – Elution, de-crosslinking and DNA isolation



- 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





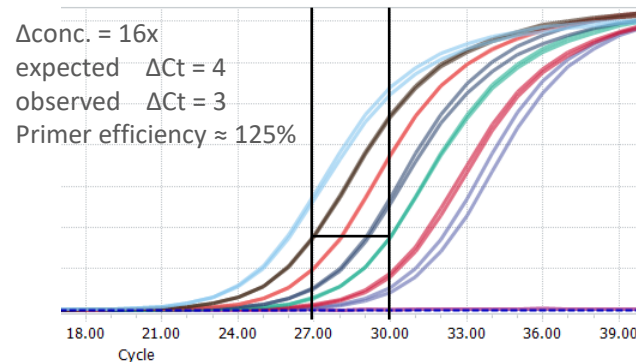
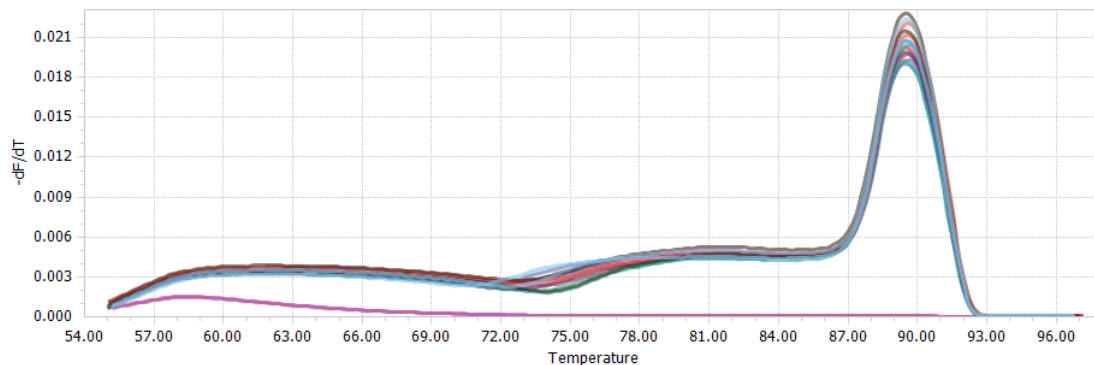
Setting up ChIP-qPCR

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:

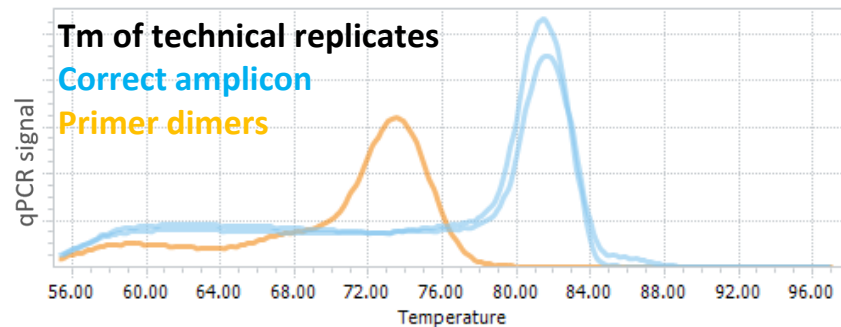
- Check on gDNA/input for T_m profile
- Check efficiency (95-105% acceptable)



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

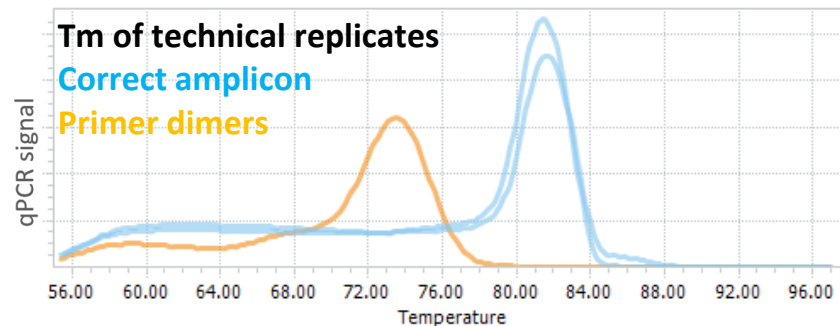




Setting up ChIP-qPCR

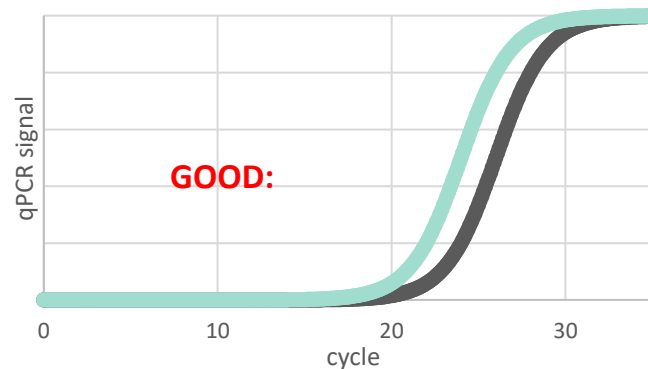
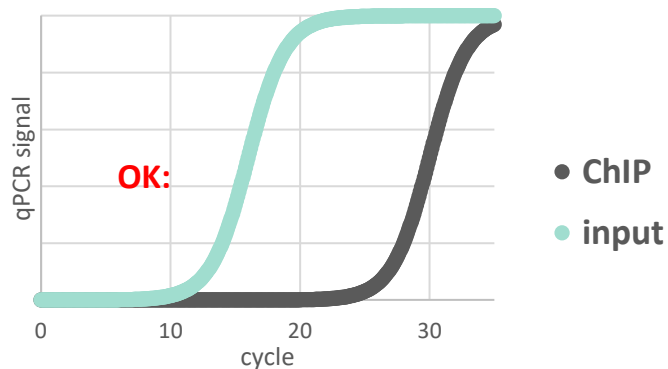
■ QC

- 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
- F 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

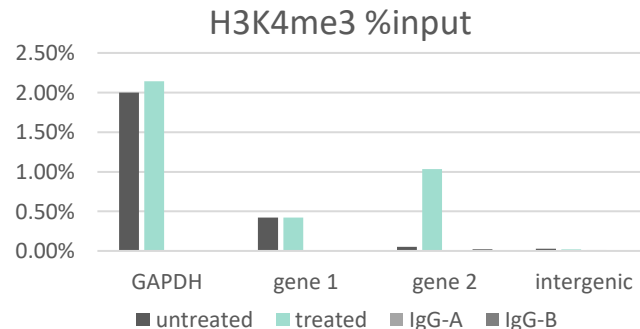
Conditions: untreated (A), treated (B)
 Targets: positive control (e.g. GAPDH) (P), genes of interest (G1, G2) intergenic region (N)
 Antibodies: IgG, H3K4me3

	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

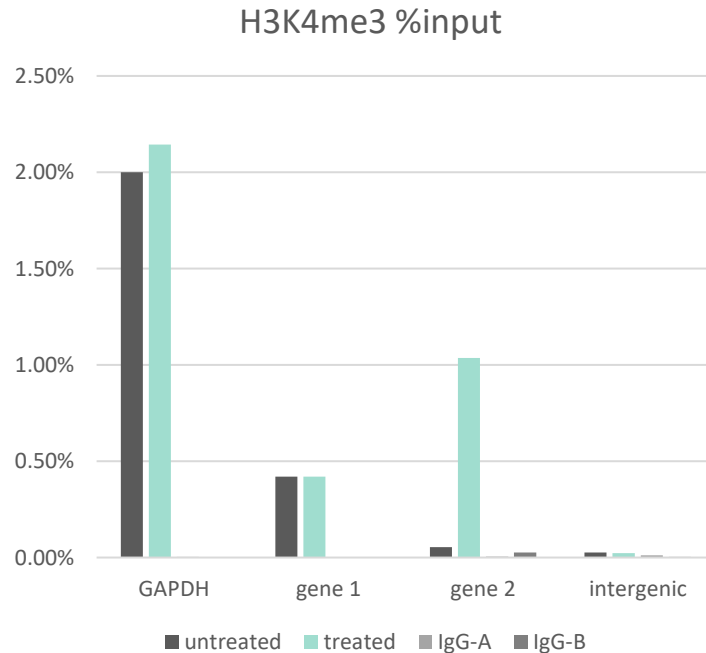
7. Optimization

- if R or F are low, optimize ChIP parameters



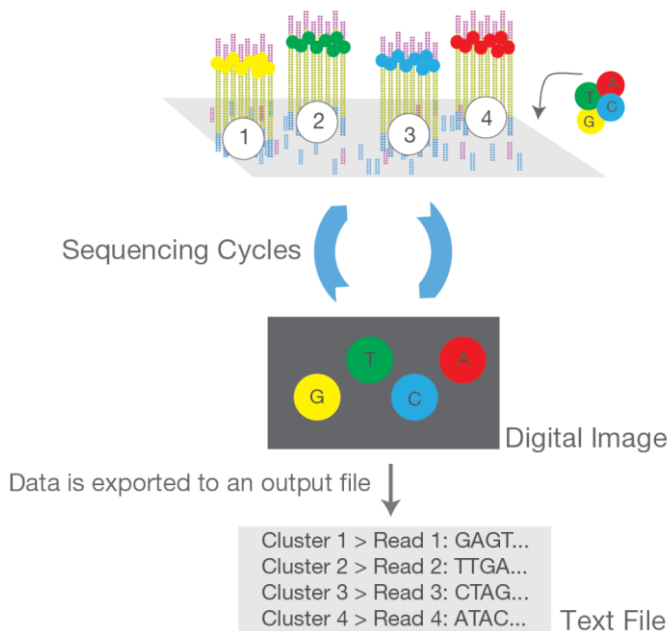
ChIP-qPCR Analysis summary

- 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

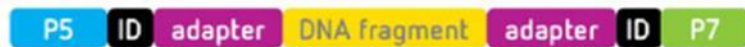


ChIP-seq: library prep

C. Sequencing



ChIP-seq: Library Prep



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

MicroPlex kit workflow

STEP 1 Template preparation

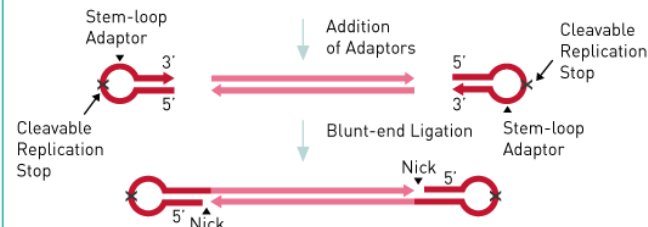
5 minutes*

Template:
Fragmented
dsDNA/cDNA



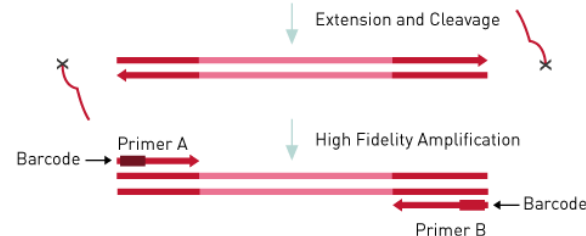
STEP 2 Library synthesis

5 minutes*



STEP 3 Library amplification

5 minutes*



* hands-on-time



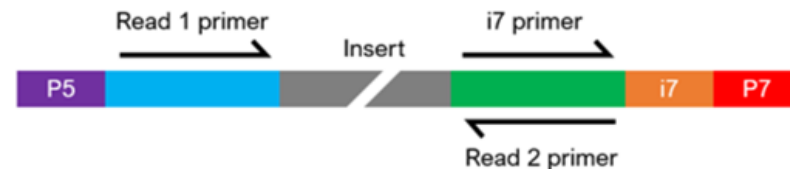
Benefits of Multiplexing

- **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

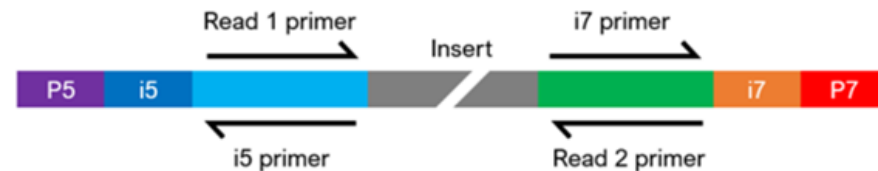


Single and Dual-indexed Libraries

- **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



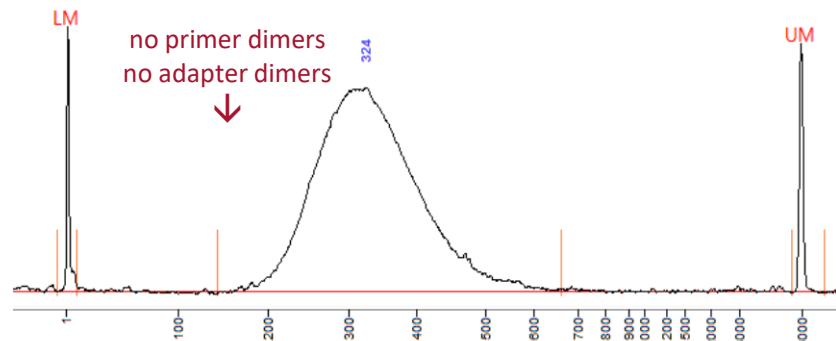
Dual-Indexed Sequencing





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

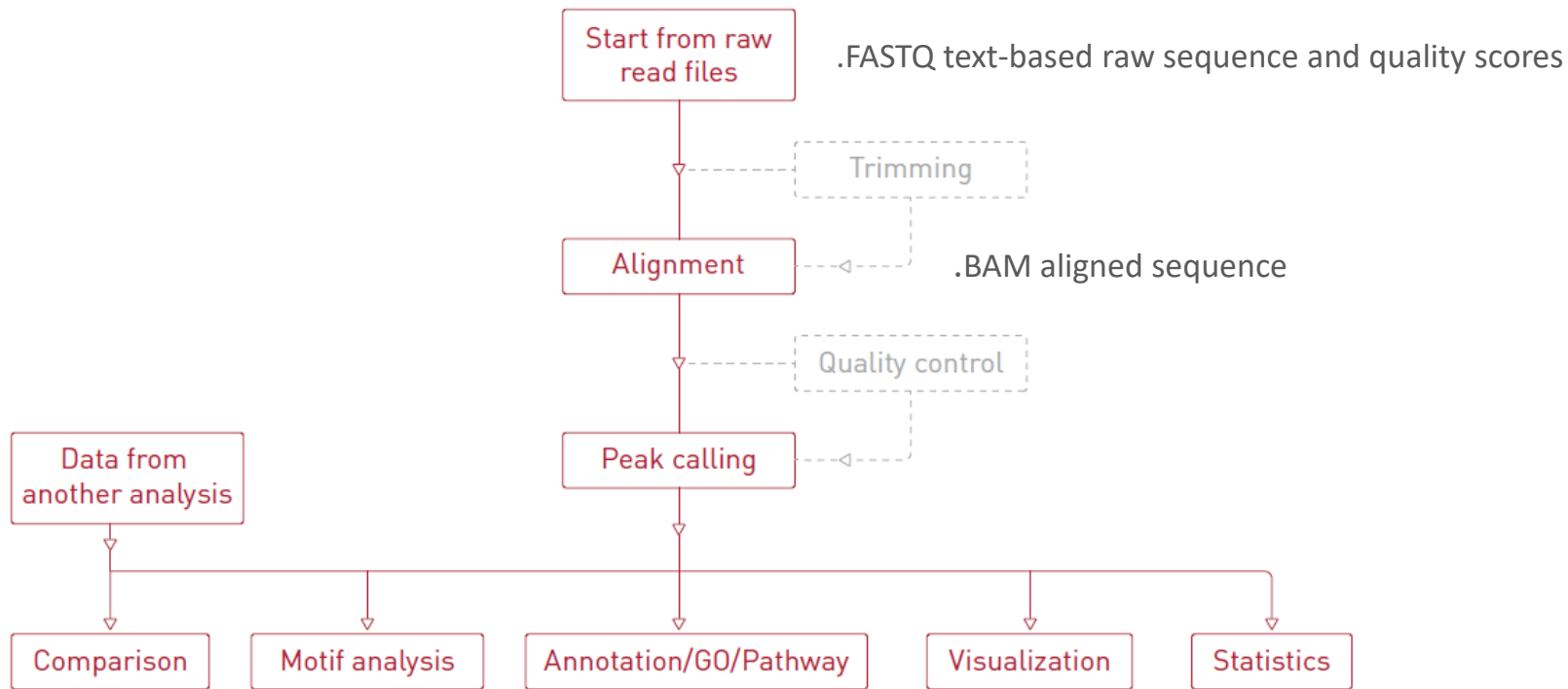


ChIP-seq: Analysis

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



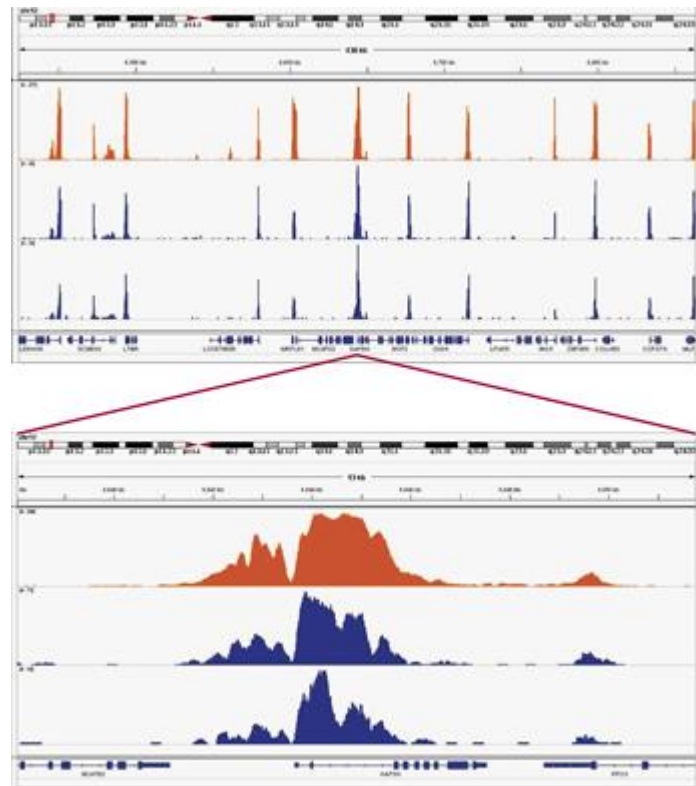
ChIP-seq: Analysis

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





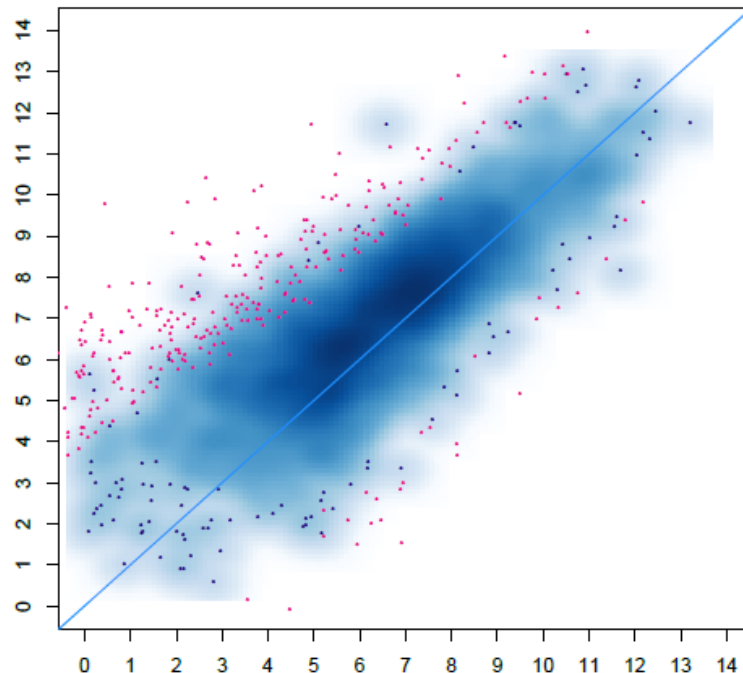
ChIP-seq: Analysis

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



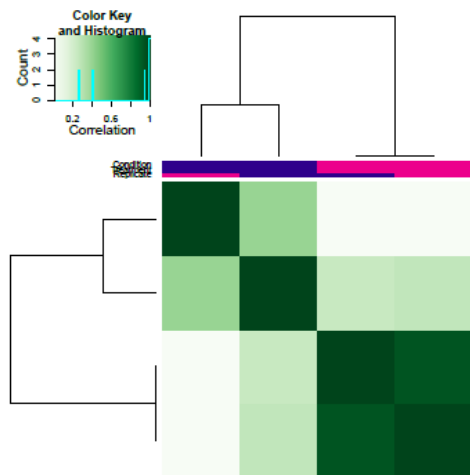
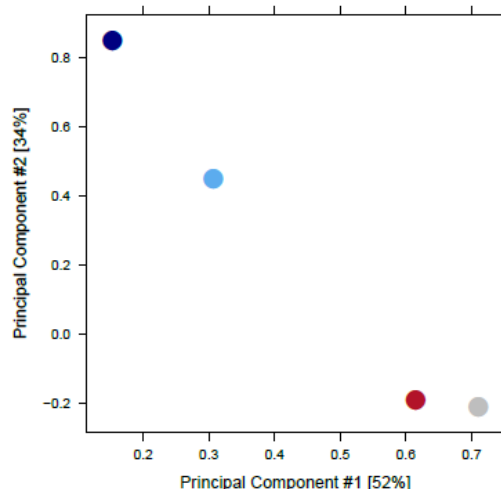
ChIP-seq: Analysis

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



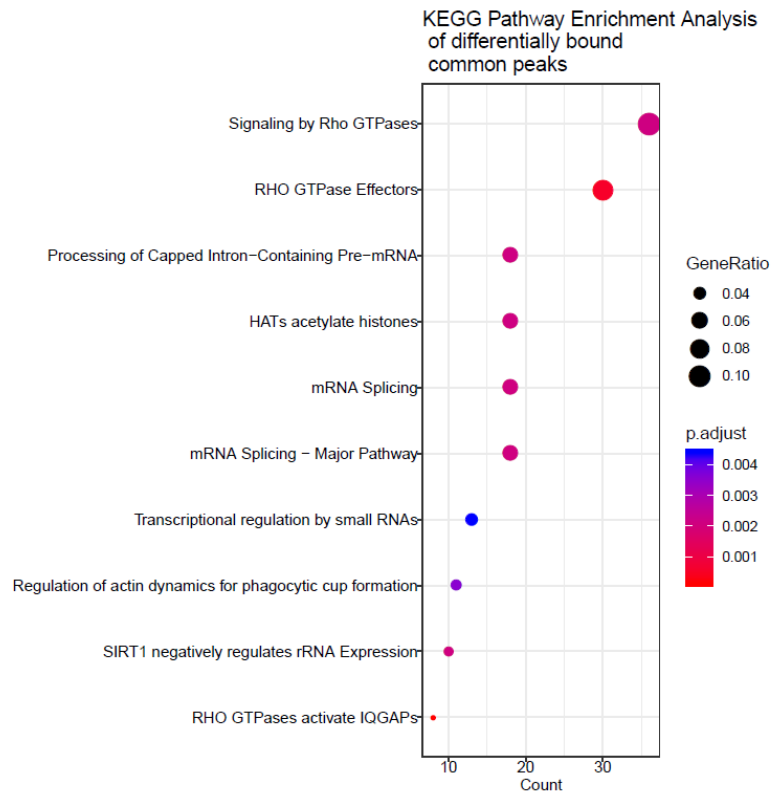
ChIP-seq: Analysis

Standard bioinformatic analysis:

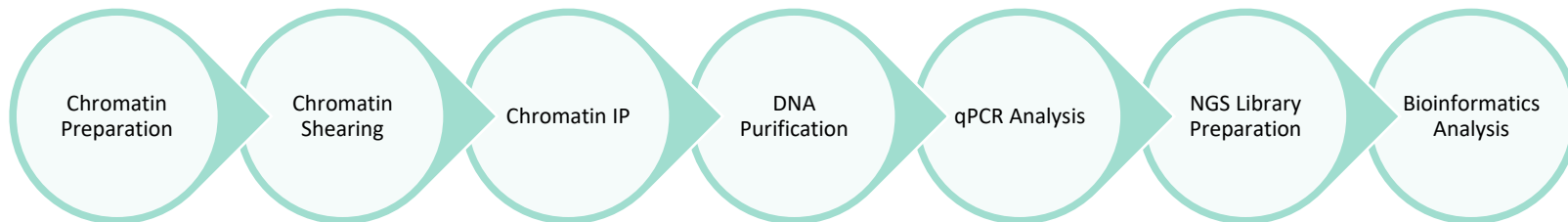
- alignment to reference genome
- peak calling

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-seq 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™

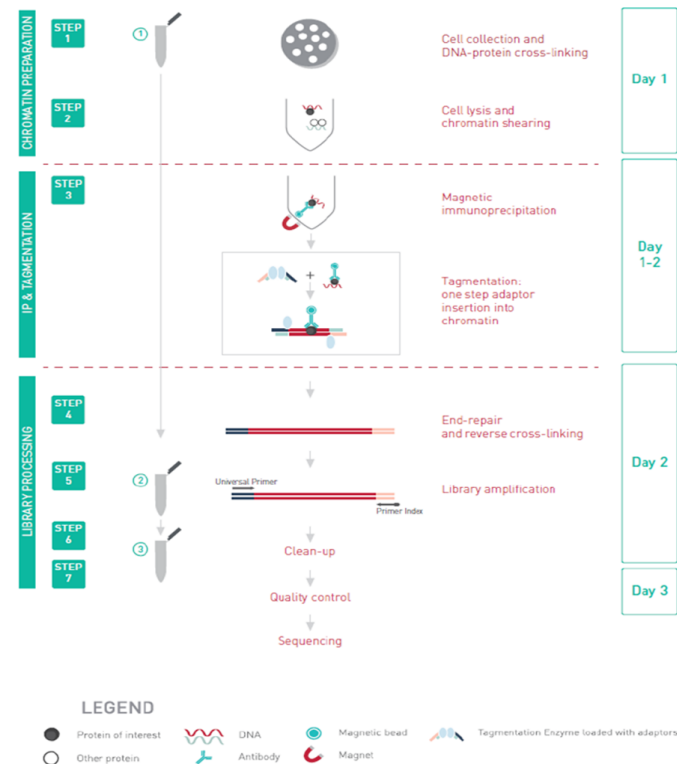
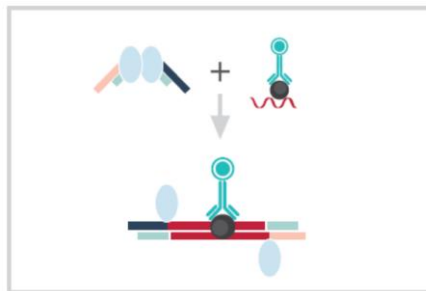
Easier and faster than classical ChIP-seq

Validated for various histone marks

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

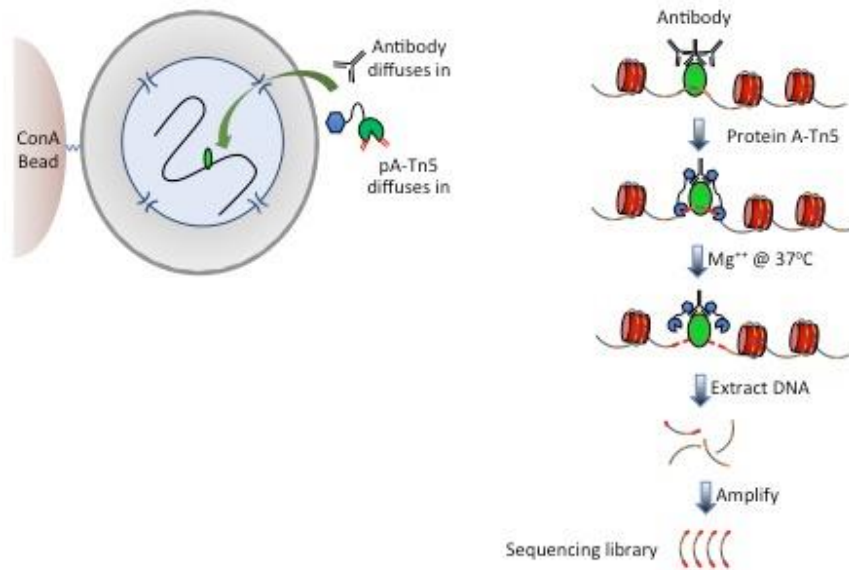
Ideal for analysis of large number of marks on a unique sample

μChIPmentation for 10,000 cells



CUT&Tag: Cleavage Under Targets and Tagmentation

CUT&Tag (Cleavage Under Targets & Tagmentation)



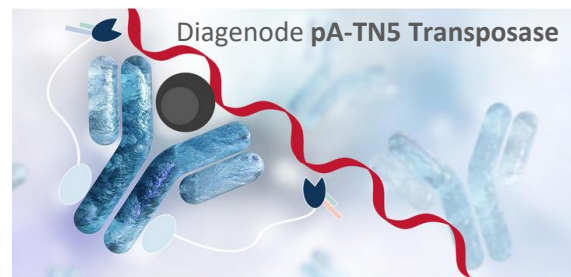
Key features:

Crucial reagent:

- proteinA-Tn5

Fast and easy protocol:

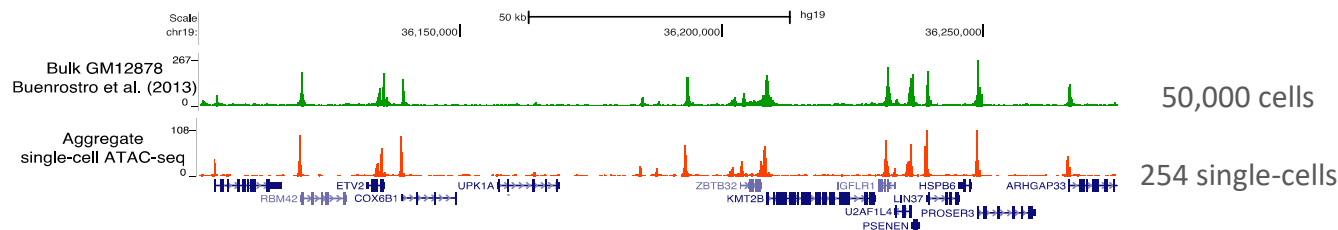
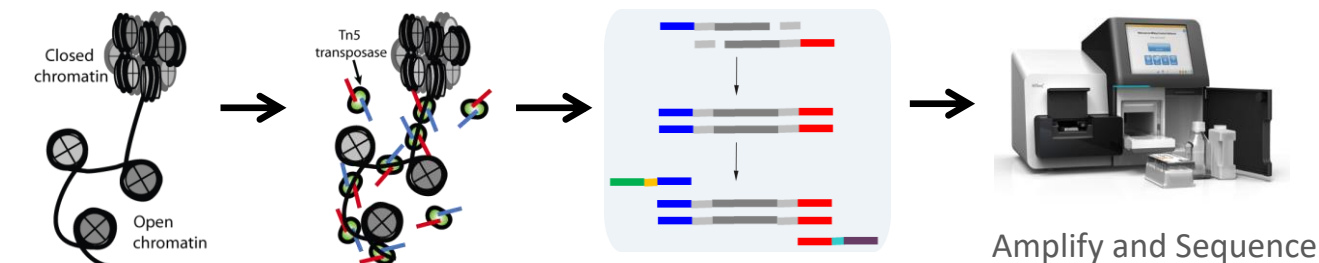
- fast tagmentation-based library prep
- No chromatin prep
- Suitable for low cell numbers



New application note available [here](#)



Assay for Transposase-Accessible Chromatin



Buenrostro et al., Nature, 2013



New! ATAC-seq Kit




Starting material: 200k-500k/sample

Minimal: 20k cells/reaction

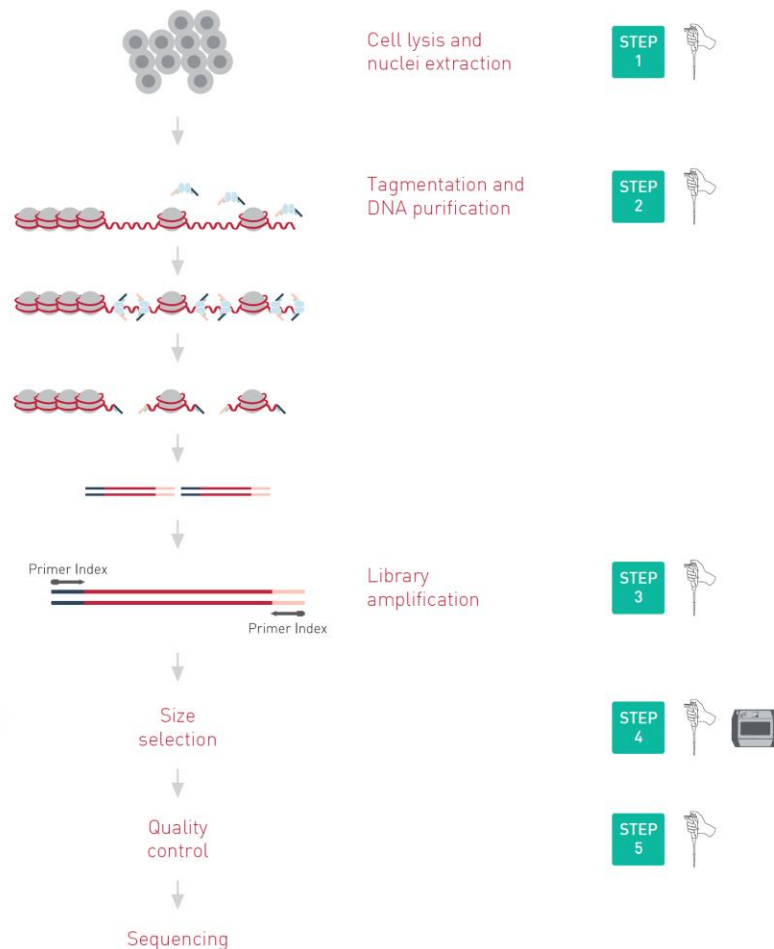
Optimal: 50k cells/reaction

Validated on mammalian cells

LEGEND

 Chromatin  Tagmentase  Fragmented DNA with adaptors

 Manual processing  Auto processing

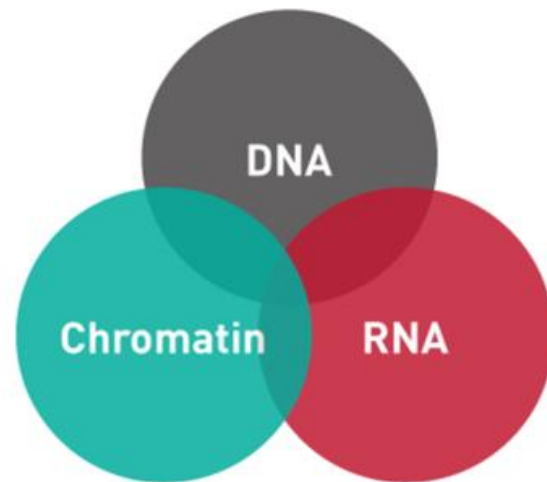




EPIGENOMICS PROFILING SERVICES

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



Sample Shipment

Processing

Data Delivery

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

