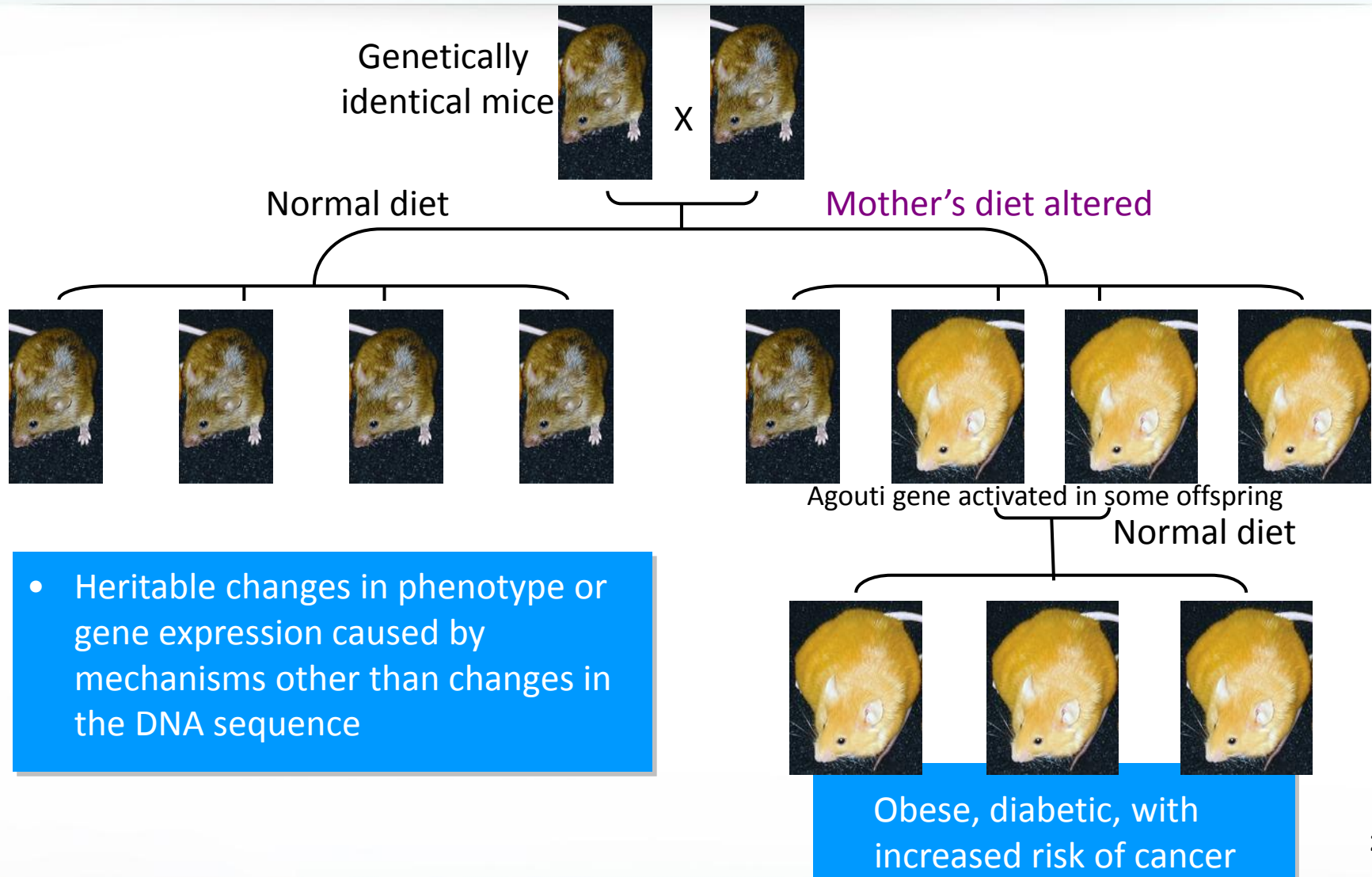


DNA Methylation Mechanisms and Analysis Methods to Study this Key Epigenetic Control

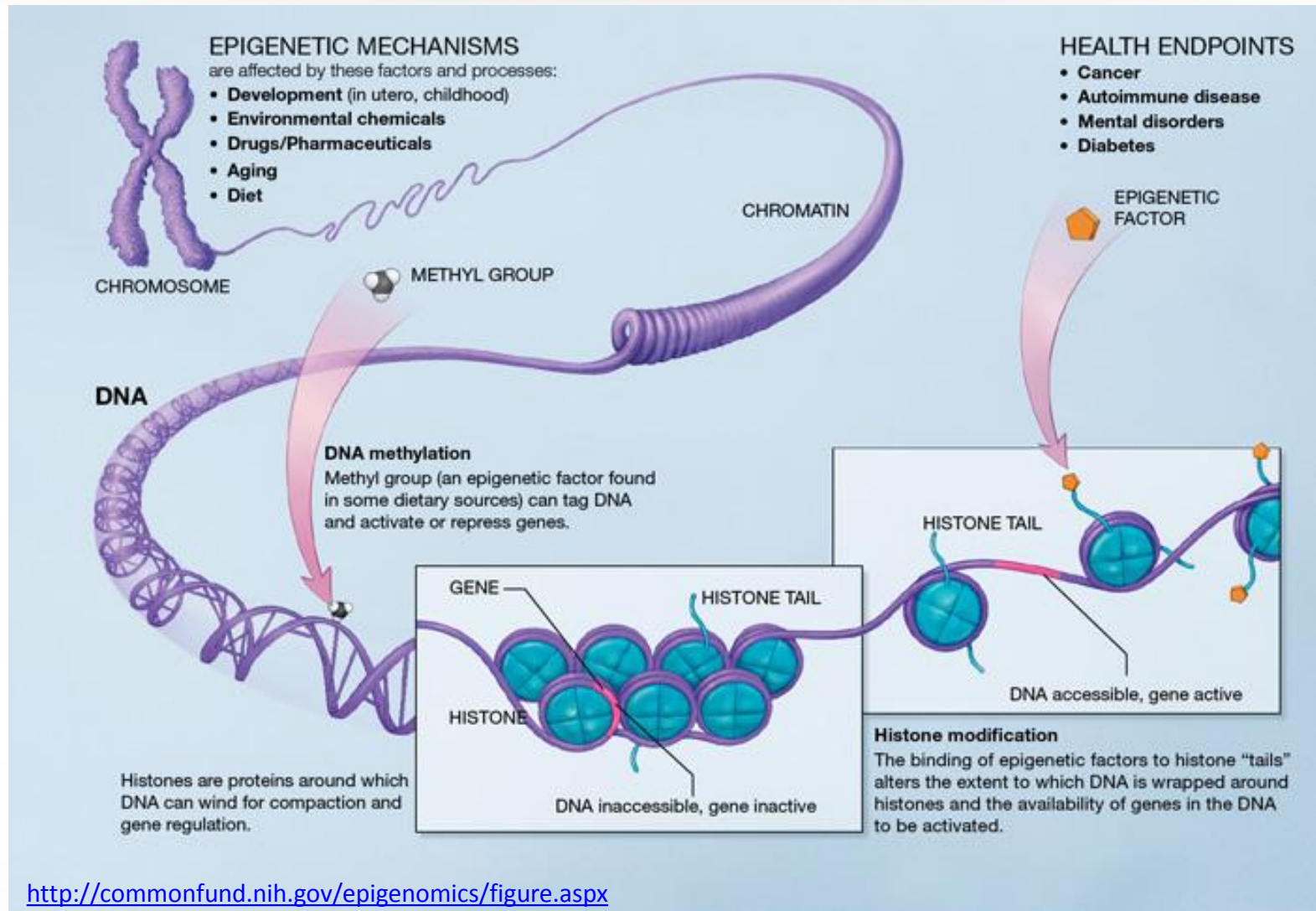
Karen Reece , Ph.D.
September 2012



Epigenetics in Action



Epigenetics Overview



Overview



1. DNA methylation
2. Introduction of bisulfite conversion chemistry
3. Downstream analysis methods
4. Considerations for obtaining quality data

Overview

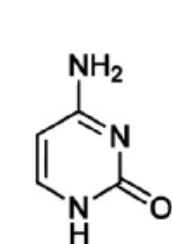


- 1. DNA methylation**
2. Introduction of bisulfite conversion chemistry
3. Downstream analysis methods
4. Considerations for obtaining quality data

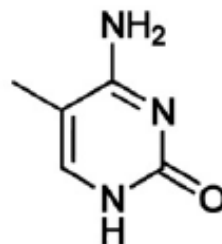
DNA Methylation

Primary Epigenetic Modification of DNA

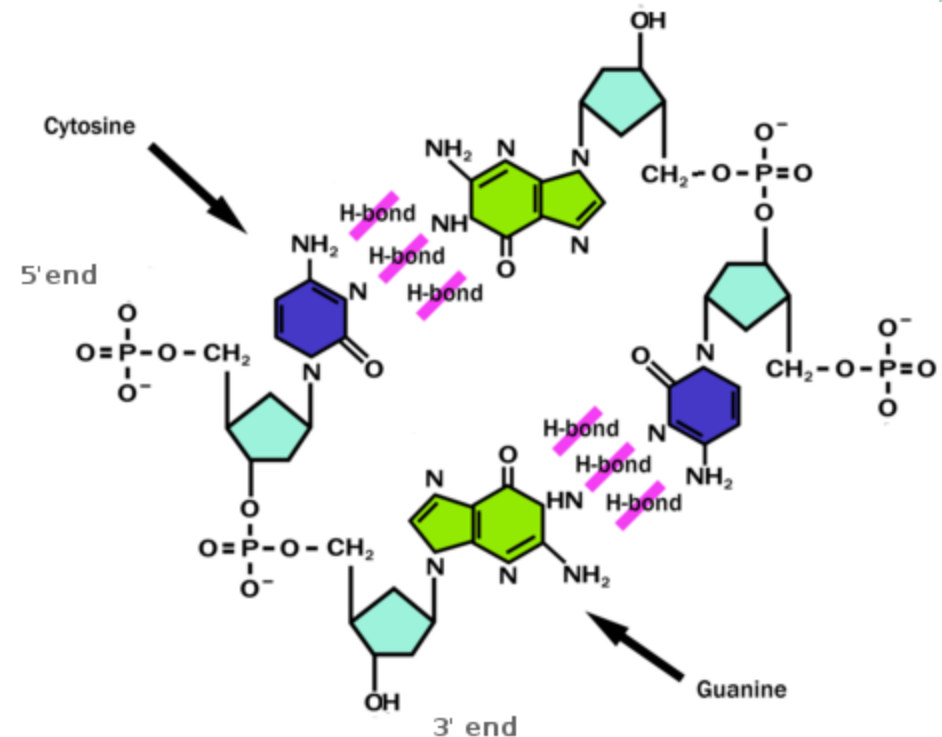
- The predominant epigenetic modification of DNA in mammalian genomes is methylation of cytosine nucleotides (5-MeC).
- The primary target sequence for DNA methylation in mammals is 5'-CpG-3' dinucleotides



Cytidine
C



5-methyl cytidine
5mC



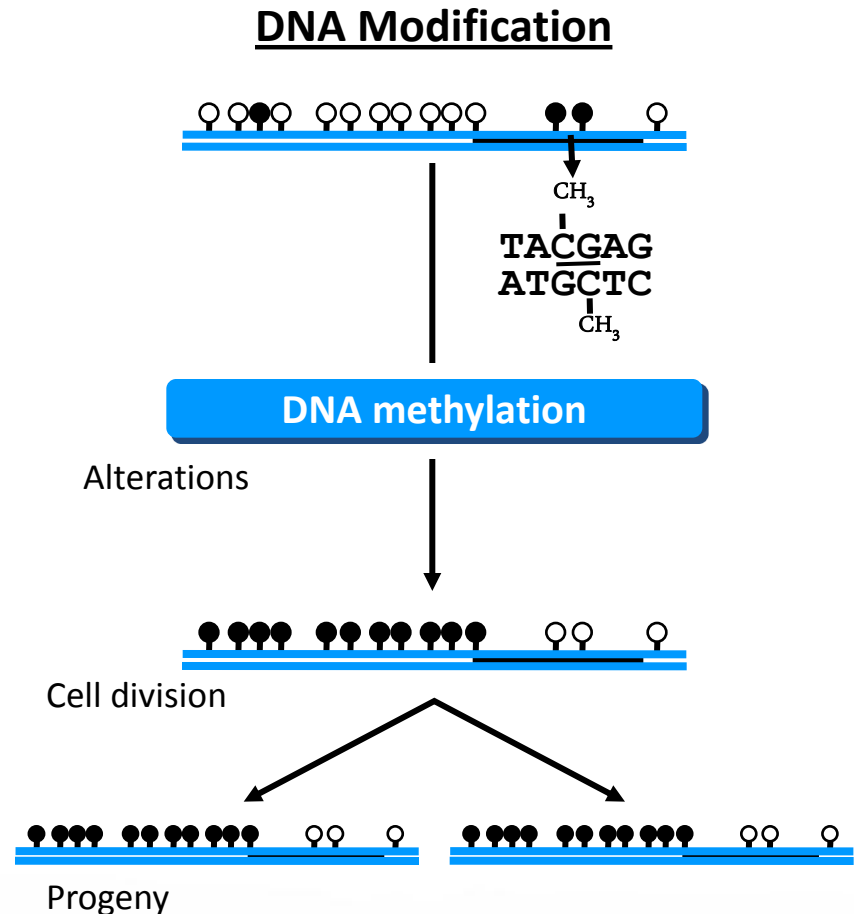
DNA Methylation

Occurs Primarily at CpG Sites in Promoter Regions



CpG islands

- Definition
 - Small stretches of about 300-3000bp
 - >50% GC content, 5' regulatory regions
- Methylation status
 - Generally non-methylated in coding regions allowing gene expression
 - Typically methylated in non-coding regions
- Genomic distribution
 - ~70% of promoter regions contain CpG islands
 - Only 1% of remaining genome contains CpG islands

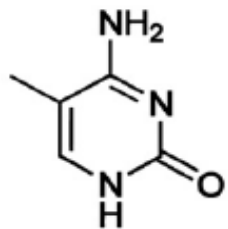


DNA Methylation

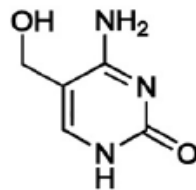
Potential Demethylation Pathway and Intermediates



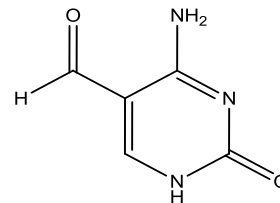
- Covalent addition of $-CH_3$ at the 5' of the cytosine ring by methyltransferases.
- Mechanism for demethylation is unclear, but may involve TET proteins or oxidative pathways leading to intermediates.



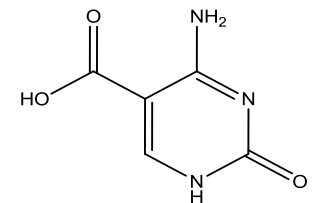
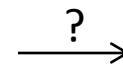
5-methyl cytosine
5mC



5-hydroxymethyl cytosine
5hmC

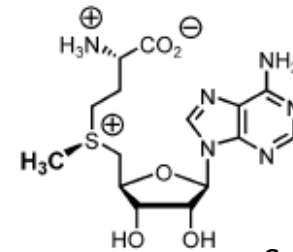
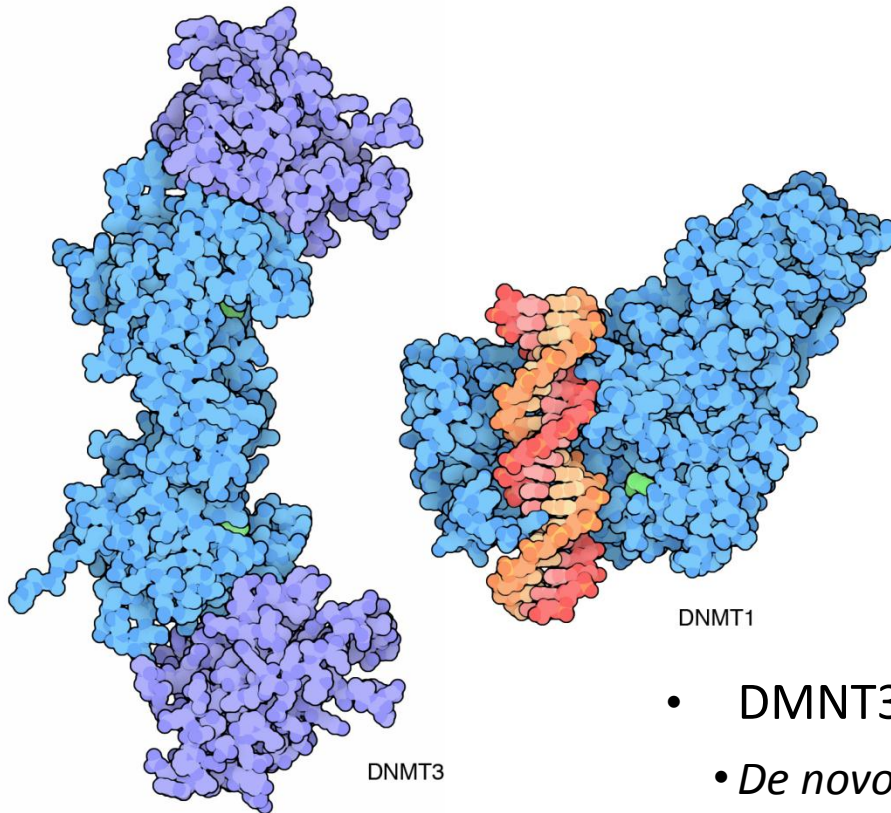


5-formyl cytosine
5fC



5-carboxyl cytosine
5caC

DNA Methylation CpG Methyltransferases



SAM

S-adenosyl-L-
methionine

- DNMT1
 - Maintenance methylation
 - Looks for hemimethylated CpG and maintains methylation pattern following replication
- DNMT3a, DNMT3b
 - *De novo* methylation
 - HCT116 – double knockout cell line
 - DNMT1 and DNMT3b inactive, low level methylation

Overview



1. DNA methylation
- 2. Introduction of bisulfite conversion chemistry**
3. Downstream analysis methods
4. Considerations for obtaining quality data

Studying DNA Methylation

Early Methods Lacked Sensitivity



Methylation-sensitive restriction enzymes combined with Southern blots

- Requires 5 μ g DNA with a low percentage of sites methylated

Restriction digests combined with PCR

- Requires complete digestion of methylated DNA to avoid false positives

Bisulfite modification

- C's are converted to U's
- Methylated C's are NOT converted

Introduction of chemistry:

Shapiro, et al. (1973) JBC vol 248, p4060

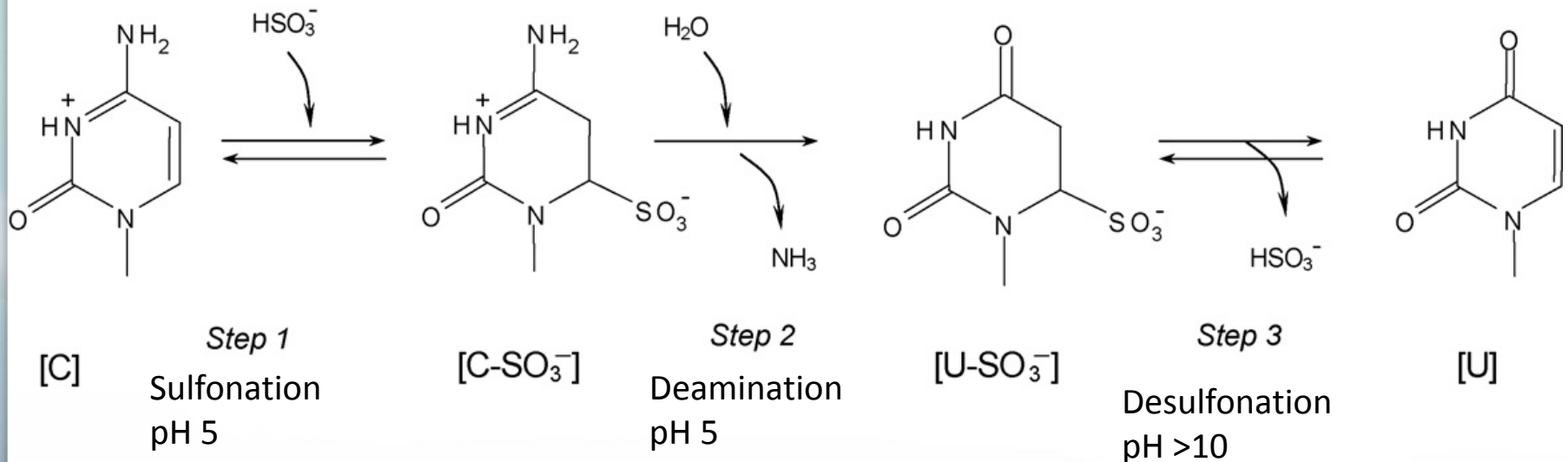
Application to DNA Methylation Detection

Frommer, et al. (1992) PNAS vol 89, p 1827

Studying DNA Methylation

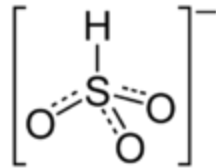
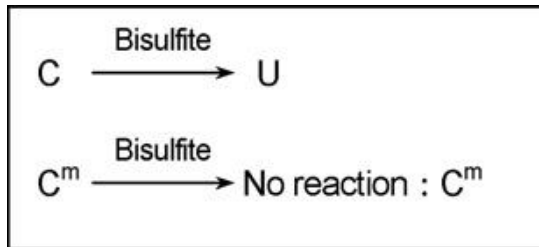
Bisulfite Conversion Converts Cytosines to Uracils

- Bisulfite Conversion is the most widely used technique for studying DNA methylation
 - Converts non-methylated cytosines to uracil
 - No distinction between 5-methyl cytosine (5mC) and 5-hydroxymethylcytosine (5hmC)

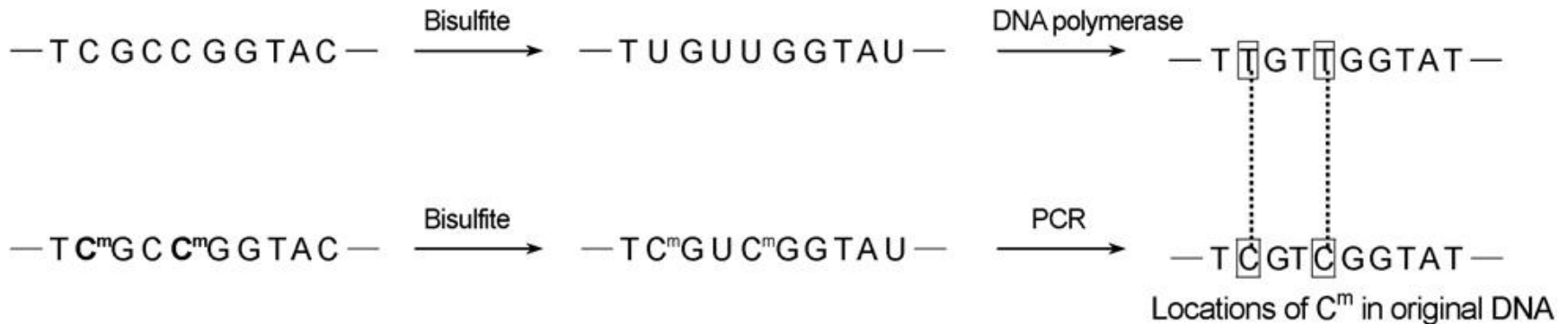


Studying DNA Methylation

Bisulfite Conversion Converts Cytosines to Uracils



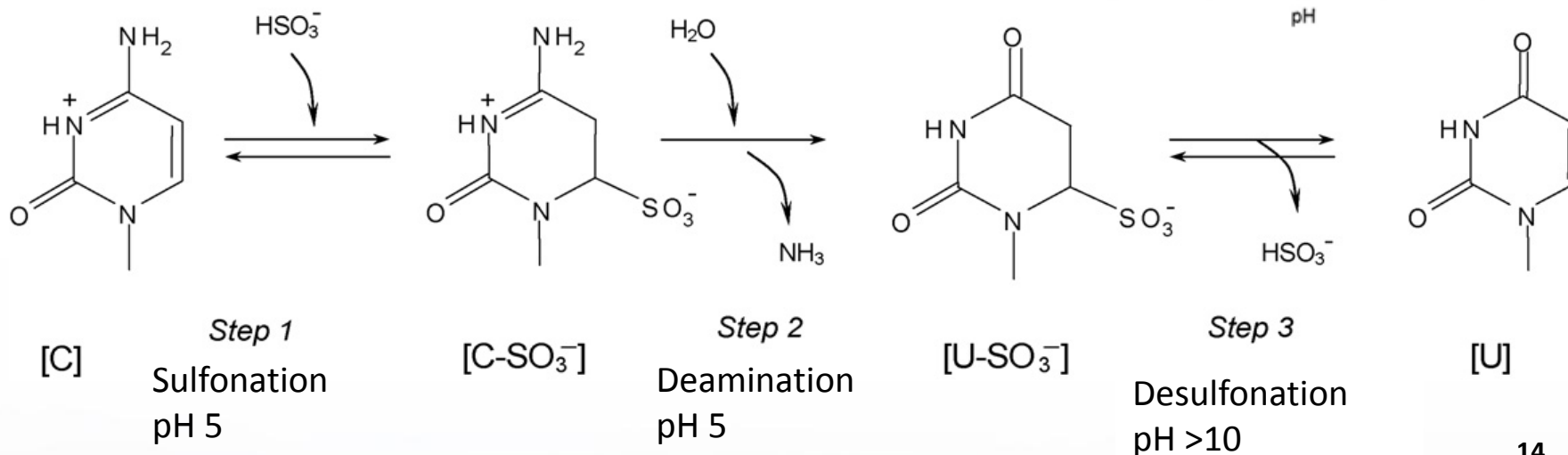
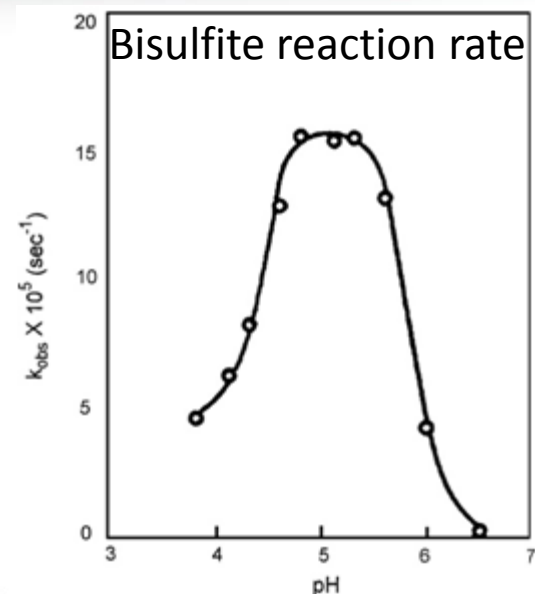
DNA :



Bisulfite Conversion Chemistry

Harsh Reaction Conditions Degrade DNA

- Original conditions:
 - Denature with NaOH or high heat
 - Incubate at pH 5 for 16-20 hours at $\sim 55^\circ\text{C}$
 - Causes depurination, fragmentation
- Early protocols led to loss of >90 % starting material

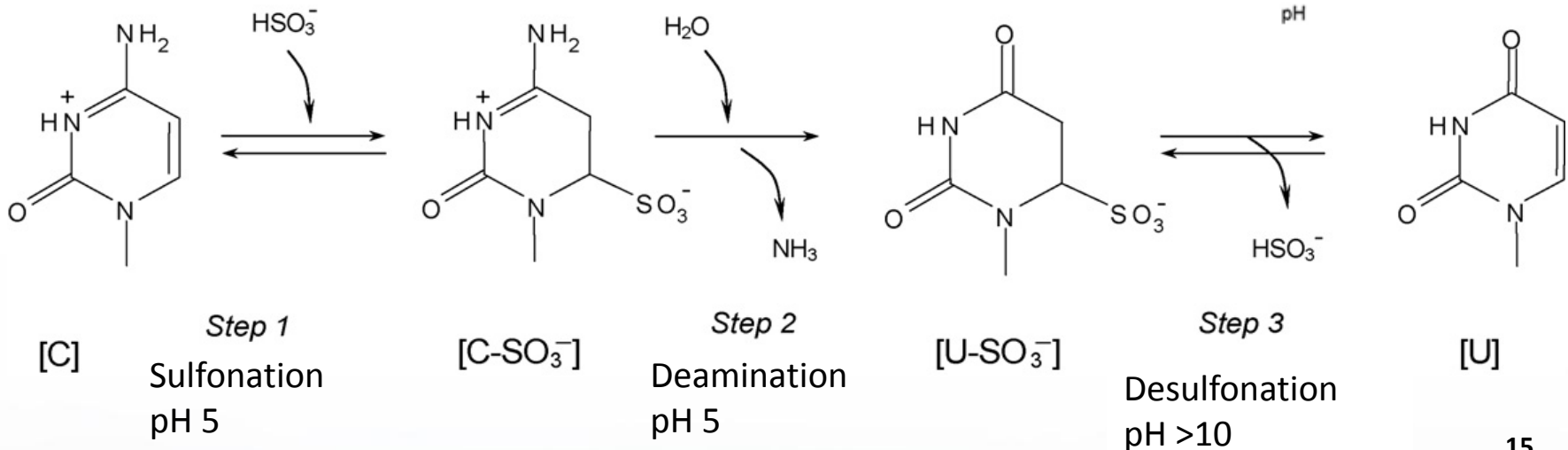
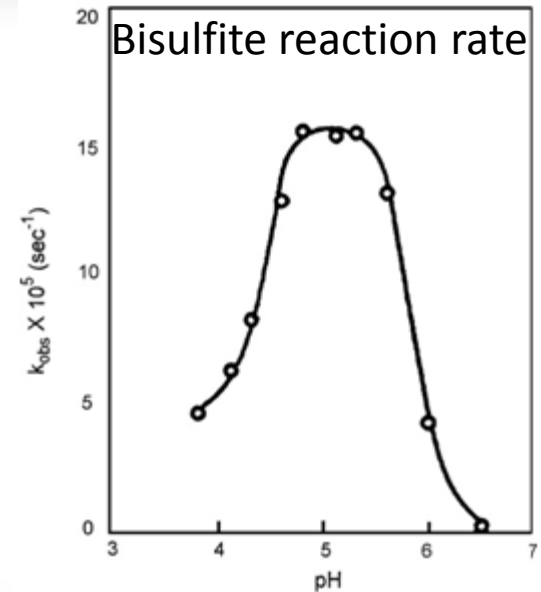


Bisulfite Conversion Chemistry

Protocol Modifications Preserve DNA Quality



- Sodium bisulfite is most common, but formulations include K and NH_3 bisulfite
- Heat and incubation time are important considerations for preserving DNA quality which impact downstream analysis methods
- New commercial kits incubate for 30-60min

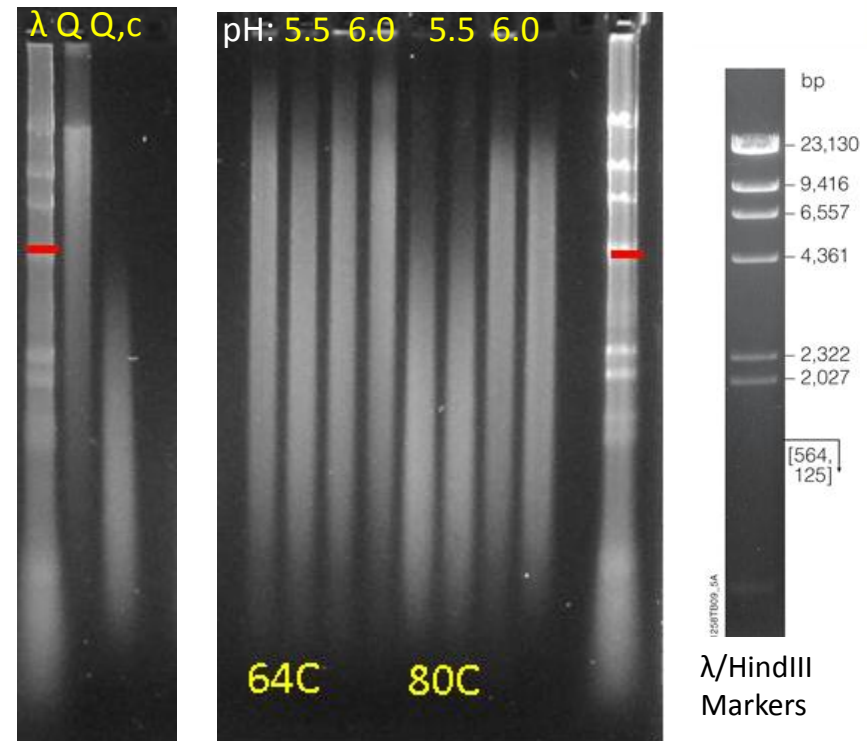


Bisulfite Conversion Chemistry

Protocol modifications Preserve DNA Quality



- λ /HindIII markers with commercially available DNA from Qiagen (Q = unconverted, Q,c = converted)
- Samples converted at either 64°C or 80°C at pH 5.5 or pH 6.0
 - Higher temperature improves denaturation, but increases fragmentation
 - Higher pH preserves DNA, but lowers conversion efficiency.



Bisulfite Conversion Chemistry

Summary



- Bisulfite Conversion is the most widely used technique for studying DNA methylation
- Harsh conditions (low pH/high temperature)
 - Variation in these parameters affect reaction rate and extent of DNA fragmentation
- Commercial kits and new protocols take only a few hours to complete and often yield less fragmented DNA compared to earlier methods

Overview



1. DNA methylation
2. Introduction of bisulfite conversion chemistry
- 3. Downstream analysis methods**
4. Considerations for obtaining quality data

Studying DNA Methylation

Downstream Analysis Methods

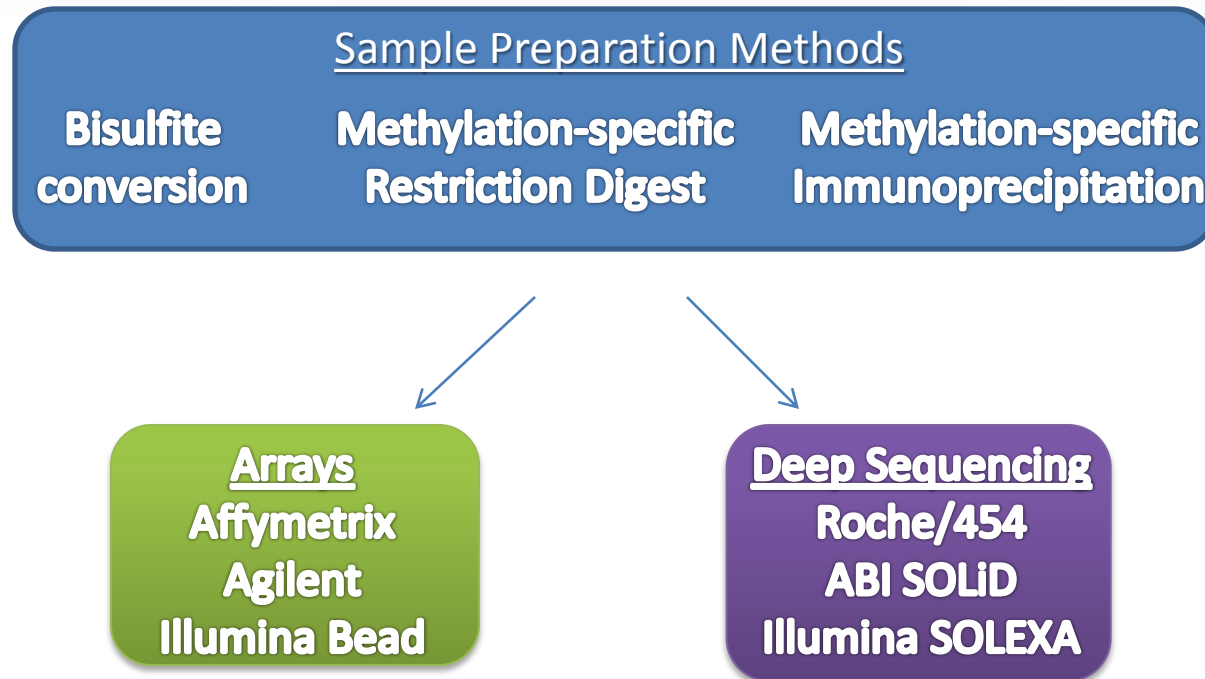


Wide range of techniques used to study DNA post-bisulfite conversion

- Methylation Specific Restriction Enzymes
- Sanger Sequencing or Pyrosequencing
- Microarrays
- PCR Techniques
 - Bisulfite Specific PCR (BSP)
 - COBRA—determination of methylation at specific RE sites within PCR amplicon

Studying DNA Methylation

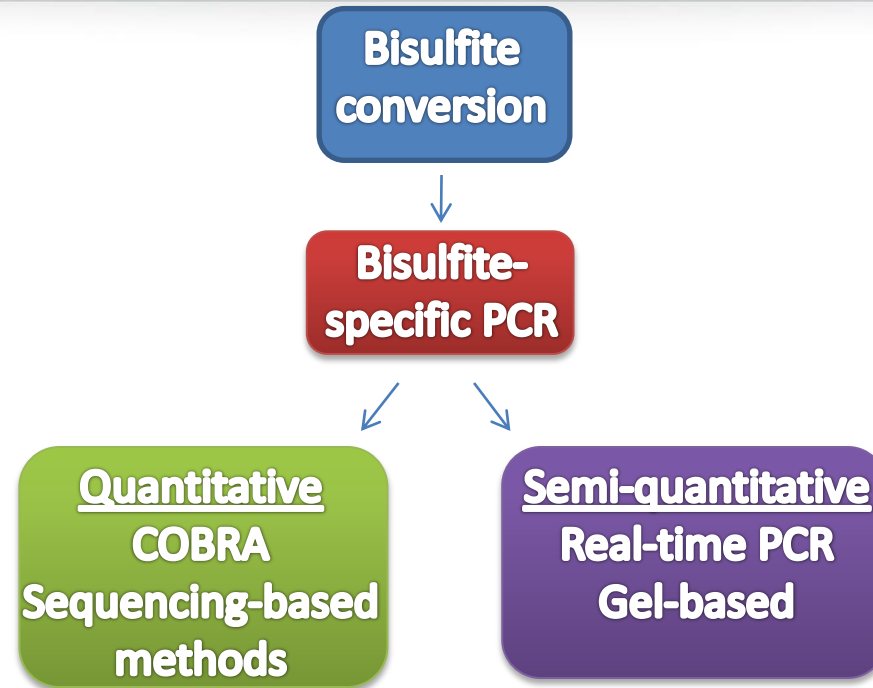
Genome-wide Approaches



- Arrays: High-throughput, Hybridization/probe-based, Lower specificity
- Sequencing: High-throughput, polymerase or ligase-based, very complicated data analysis

Studying DNA Methylation

Gene-specific Approaches



- COBRA – Combined Bisulfite Restriction Analysis
- Sequencing-based methods: pyrosequencing, CE-based sequencing, cloning and Sanger sequencing
- Semi-quantitative: Uses real-time or gel-based visualization to estimate percentages and general location of methylation

Bisulfite-specific PCR

Primers Must Be Specific to Converted Sequence



- Bisulfite-specific primer pair (blue): amplifies both bisulfite-converted methylated and bisulfite converted unmethylated DNA
- Wildtype-specific primer pair (orange): amplifies unconverted DNA, both methylated or unmethylated sequences
- Primers do not contain any CpG sites, but do contain non-CpG cytosines

NON-CONVERTED FORWARD SEQUENCE (77 Cs, T_m = 98C)

```
AAAATGGGCTAGACACAAAGGACTCGGTGCTTGTCCCAGCCAGGCGCCCTCGGCGACCGGGCAGCTGGG
AGGGGAATGGGCGCCCGGACCCAGCTGGGAACCCCGGGTGCAGACTCCACCTACCTAGTCCGGCGCCAGGC
CGGGTCCGACAGCTCCGGCAGCCAGCCCGCCCGCCGTGTCCAGATGTCGCGTCAGAGGCGTGACCGGTT
TAGTTTAATTTCCGCTTGTTTTTCCAAATCTAGAAGAGG
```

BISULFITE CONVERTED FORWARD STRAND (UNMETHYLATED) (0 Cs, T_m = 84C)

```
AAAATGGGTTAGATATAAAGGATTGGTGTGTTTGTGTTTAGTTAGGTGTTTTTGGTGATGTGGGTAGTTGGG
AGGGGAATGGGTGTTTGGATTTAGTTGGGATTTTTGGGTGTGATTTTATTTATTTAGTTTGGTGTAGGT
TGGGTTGATAGTTTTGGTAGTGTTAGTGTTGTGTTGTGTTTAGATGTTGTGTTAGAGGTGTGTAGTGGTT
TAGTTTAATTTGGTTTGTGTTTTTAAATTTAGAAGAGG
```

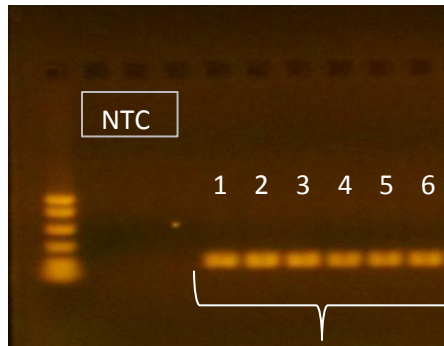
BISULFITE CONVERTED FORWARD STRAND (METHYLATED) (25 Cs, T_m = 88C)

```
AAAATGGGTTAGATATAAAGGATTGGTGTGTTTGTGTTTAGTTAGGCGTTTTCGGCGACCGGGTAGTTGGG
AGGGGAATGGGCGTTCGGATTTAGTTGGGATTTTCGGGTGCGATTTTATTTATTTAGTTCCGGCGTTAGGT
CGGGTCCGATAGTTTCGGTAGCGTTAGCGTCCGCGTCCGTGTTTAGATGTCGCGTTAGAGGCGTGTAGCGGTT
TAGTTTAATTTGGTTTGTGTTTTTAAATTTAGAAGAGG
```

Bisulfite-specific PCR

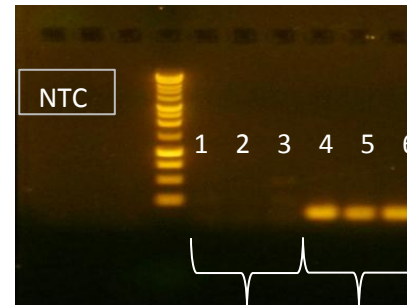
Detecting Underconverted DNA

Bisulfite-specific primers: a band should be visible if DNA is converted



Band indicate bisulfite-converted DNA

Wildtype-specific primers : No bands should be visible if conversion is complete



No bands suggest complete conversion

Bands here shows unconverted DNA is present

- Primers do not contain any CpG sites, but do contain non-CpG cytosines
- **Bisulfite-specific primer pair:** amplifies both bisulfite-converted methylated and bisulfite converted unmethylated DNA
- **Wildtype-specific primer pair:** amplifies unconverted DNA, both methylated or unmethylated sequences

Bisulfite-specific PCR

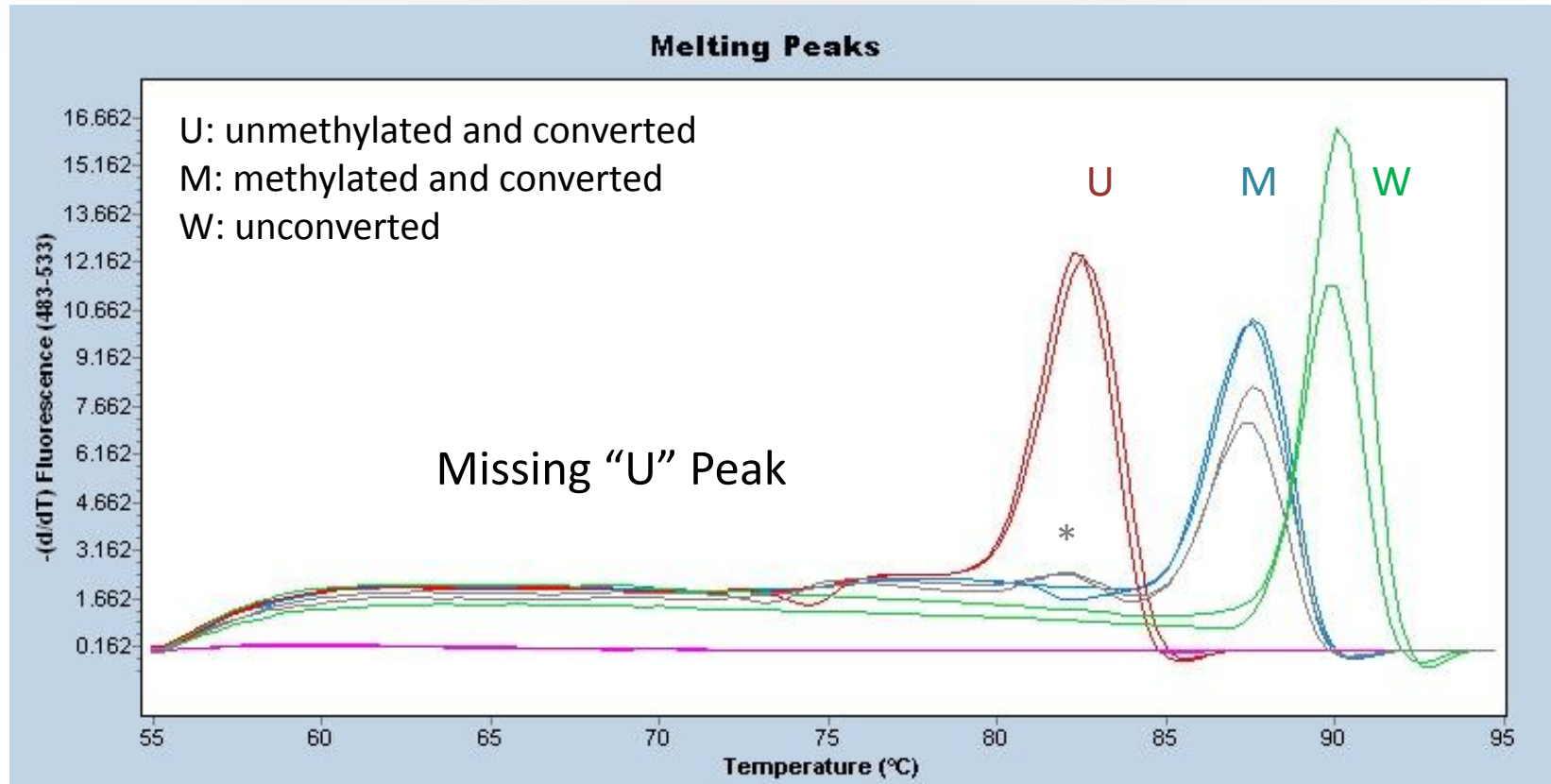
Evidence of PCR Bias



- Most literature examples show preferential amplification of bisulfite-converted unmethylated DNA over bisulfite-converted methylated DNA
- Common techniques to overcome bias are based on destabilizing GC rich regions and secondary structure of methylated DNA
- No universal approach has been reported to overcome problem
- Correction of PCR bias by means of cubic polynomial regression¹
 - Involves running control samples varying in % methylation and calculating a regression curve
 - Equation of the best-fitting curve is then used for correction of the data obtained from the samples of interest

Bisulfite-specific PCR

Preferential Amplification of Methylated DNA



- Template DNA contains a mixture of fully methylated and converted, fully unmethylated and converted DNA.
- In this assay, methylated DNA is preferentially amplified over unmethylated DNA

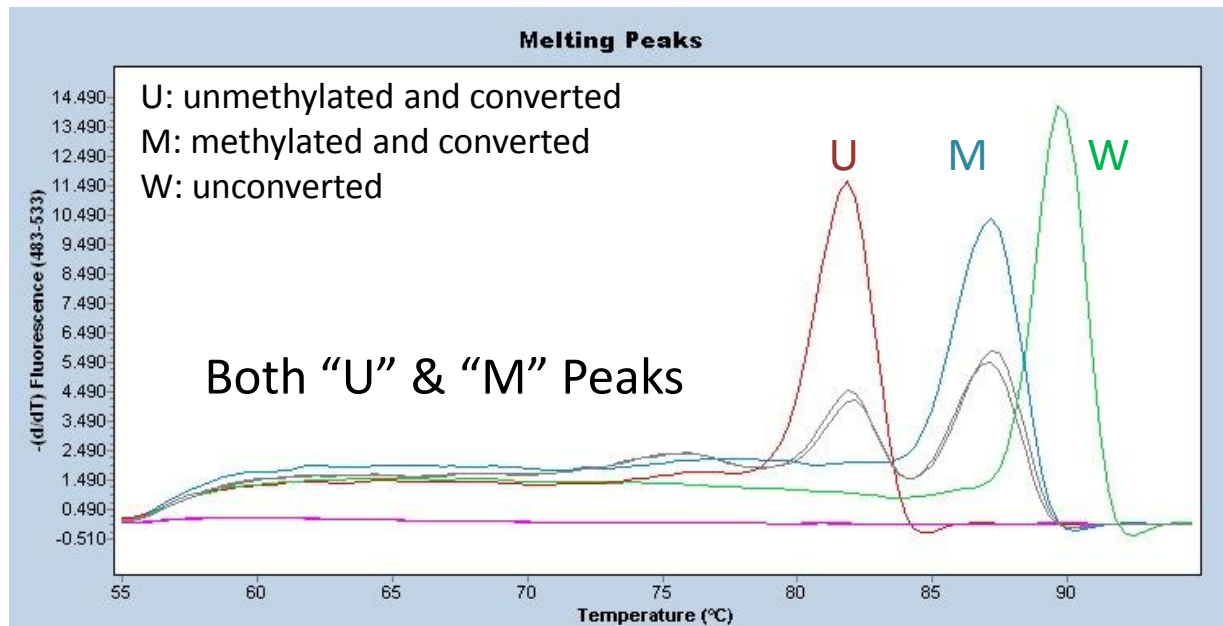
Bisulfite-specific PCR

Correction of PCR Bias by Modifying Cycling Protocol



Rand, et al Epigenetics. 2006 Apr-Jun;1(2):94-100 used a differential denaturation in real time PCR to amplify unmethylated DNA in a methylated DNA background, after treatment with bisulfite.

- T_m differences of 2.3 - 5°C allowed selective amplification of unmethylated amplicons over corresponding methylated amplicons



Bisulfite-specific PCR

Designing Primers to Avoid PCR Bias



Suggestions for bisulfite-specific primer design:

1. The fewer CpG sites, the better
2. When CpG sites are present, keep them away from the 3' end of the primer
3. Salt-adjusted T_m of primer should be around 65°C to run PCR annealing step at 60°C
4. Include one or more T's near 3' end of primer (not originating from CpG site)
5. Evaluate primers for secondary structure or primer dimer complications

Wojdacz, et al., 2008, A new approach to primer design for the control of PCR bias in methylation studies, BMC Research Notes, 1:54

Bisulfite-specific PCR

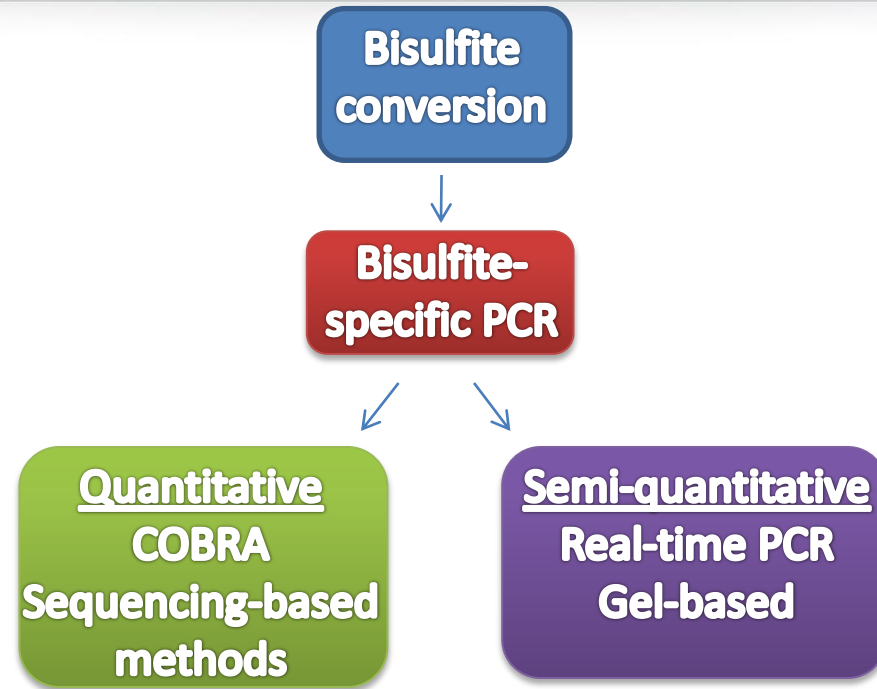
Summary



- Important step for any gene-specific DNA methylation study
- Uses two primer pairs
 - To detect DNA (methylated or unmethylated) that has been bisulfite converted
 - To detect DNA that has not be converted
- PCR bias is a common problem
 - Follow published primer design recommendations
 - Qualify primers up front using validated control DNA sources
 - Can test bias by amplifying mixtures of fully methylated and unmethylated DNA in various ratios
 - If bias is still present, consider additives or experimental redesign

Studying DNA Methylation

Gene-specific Approaches



- COBRA – Combined Bisulfite Restriction Analysis
- Sequencing-based methods: pyrosequencing, CE-based sequencing, cloning and Sanger sequencing
- Semi-quantitative: Uses real-time or gel-based visualization to estimate percentages and general location of methylation

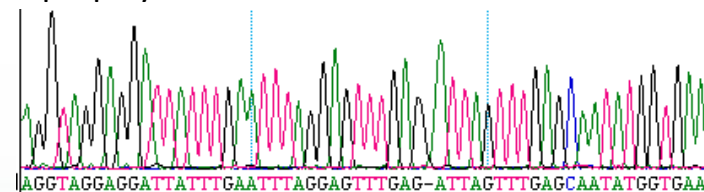
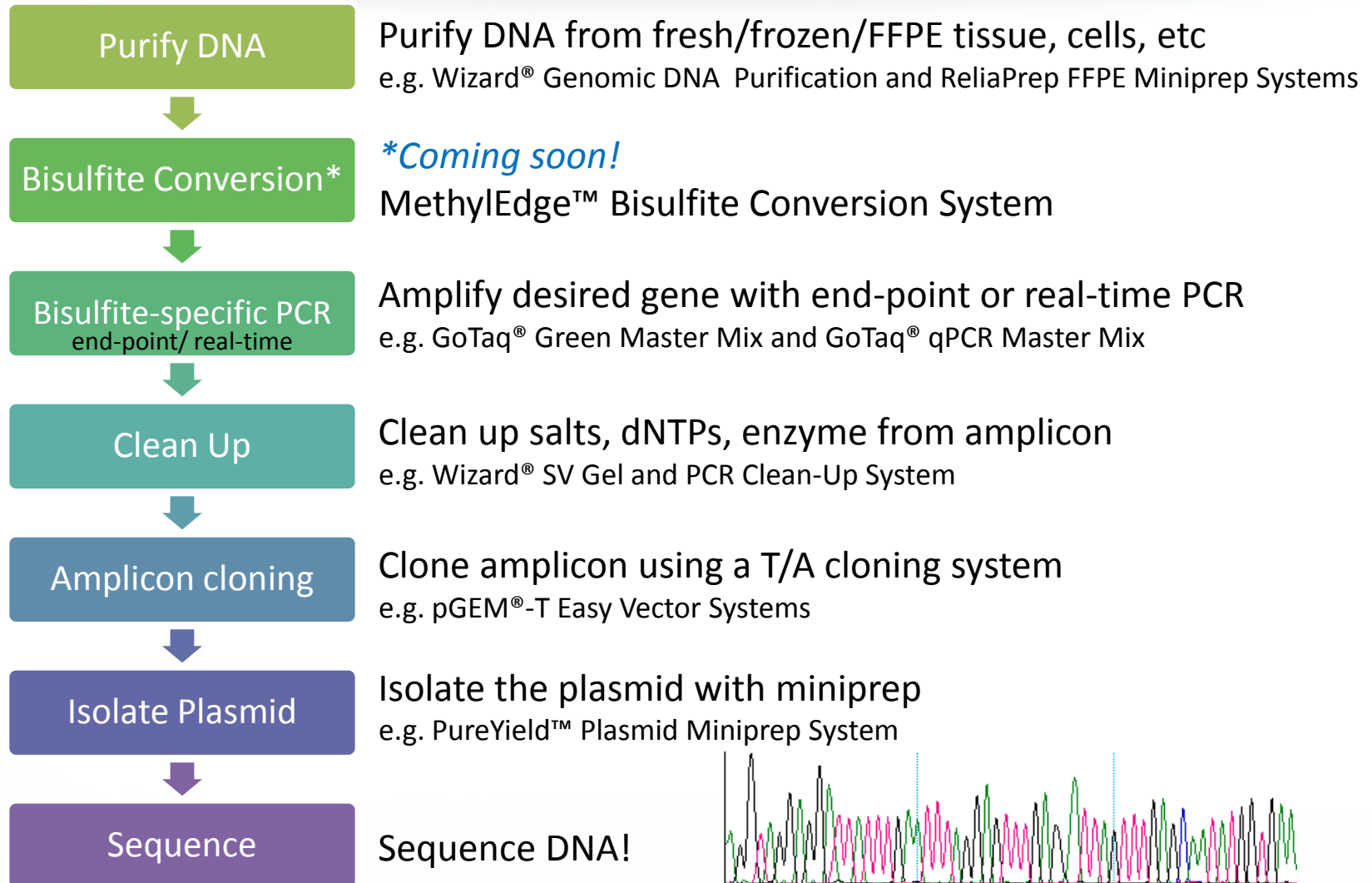
Studying DNA Methylation

Sequencing Methods for Bisulfite-Converted DNA



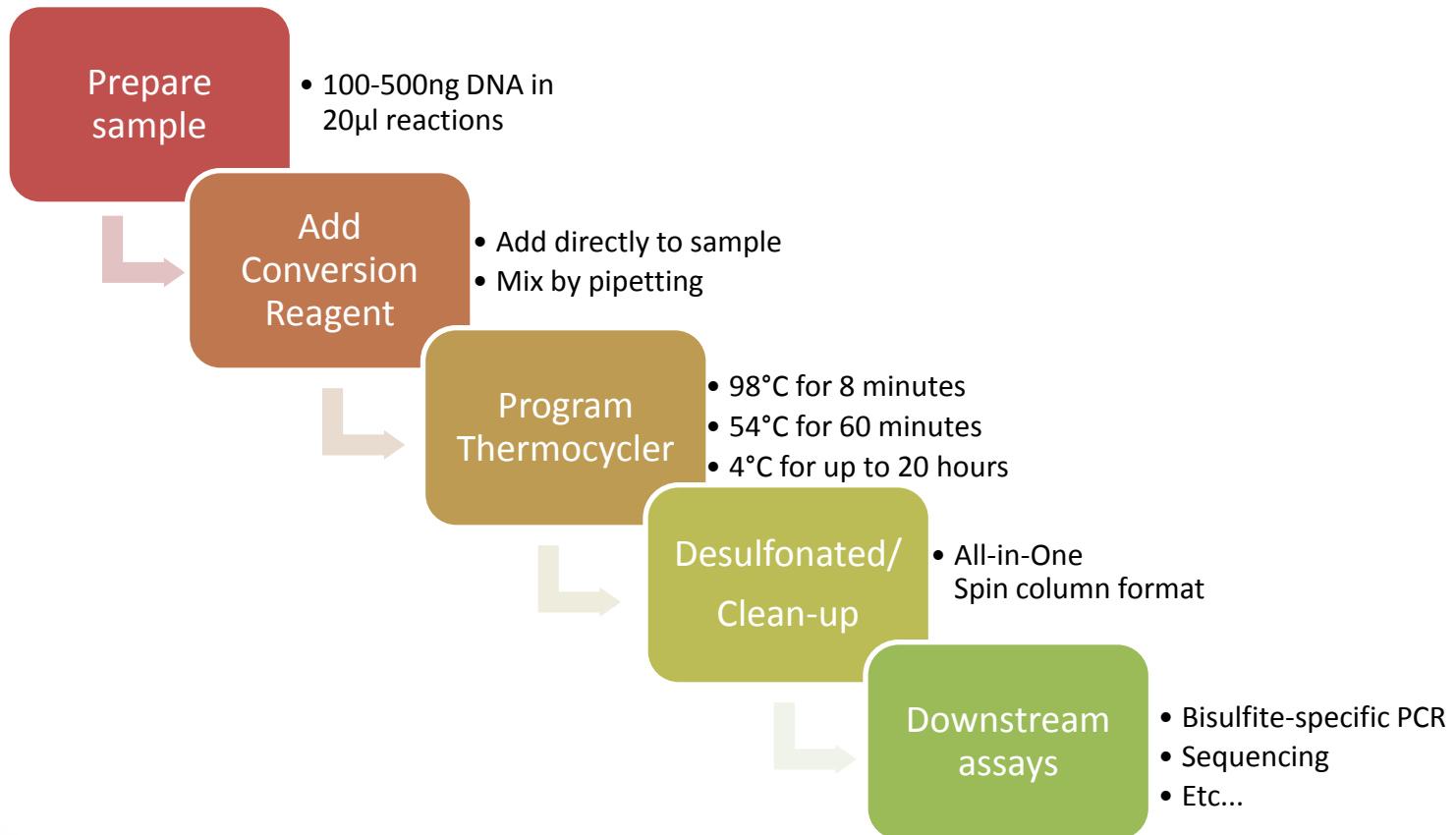
- **Pyrosequencing**
 - Detection of pyrophosphate upon nucleotide incorporation using ATP-coupled luciferase reaction
 - Light produced and measured only when complementary base is added to strand
 - No PCR reaction necessary
 - High-throughput, very expensive
- **Dye-terminator/Sanger Sequencing**
 - Each nucleotide is labeled with a different fluorophore
 - Sequence is read by chromatogram following capillary electrophoresis
 - Cloning and sequencing necessary
 - Low-throughput, more affordable

Bisulfite Sequencing Direct Sequencing Workflow



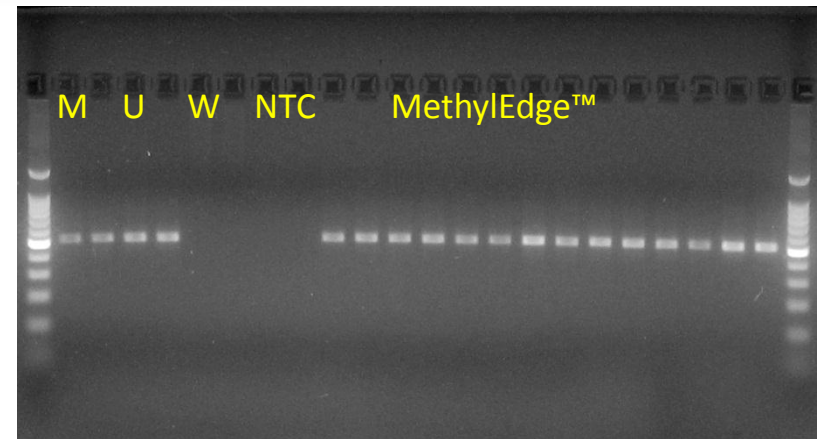
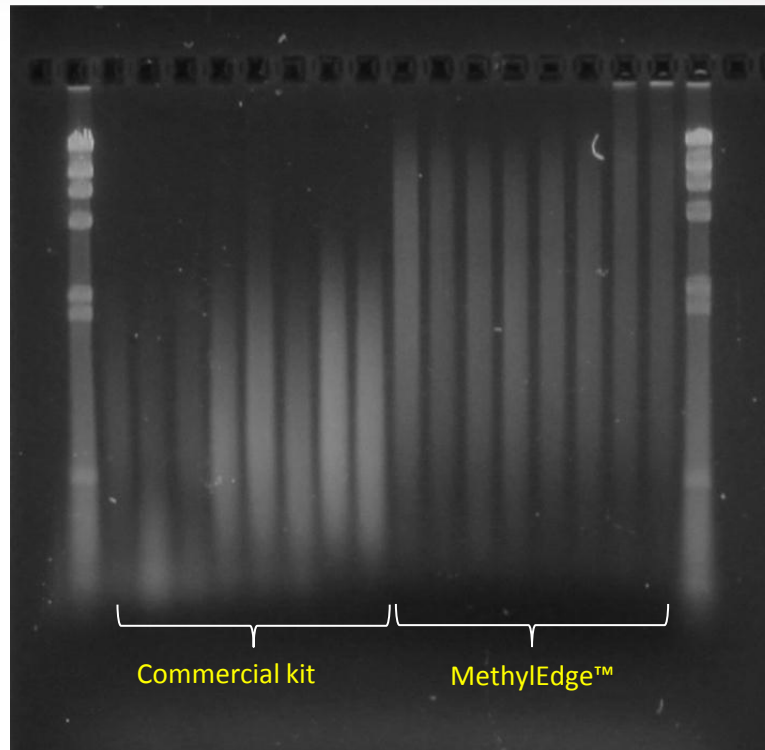
MethylEdge™ Bisulfite Conversion System

Streamlined Protocol Complete in Under Two Hours



MethylEdge™ Bisulfite Conversion System

Maintaining DNA Integrity



U: unmethylated and converted
M: methylated and converted
W: unconverted

- MethylEdge™ Bisulfite Conversion System leaves DNA less fragmented than the leading commercial kit
- Bisulfite-specific PCR shows amplification of a 524bp fragment
- Direct sequencing indicates >99% conversion

Downstream Analysis Methods Summary



Genome-wide approaches

- Arrays and deep sequencing
- Facilitate epigenomic mapping

Gene-specific approaches

- Many rely on bisulfite-specific PCR prior to analysis
- COBRA, Bisulfite Sequencing, real-time and end-point PCR
- Elucidate epigenetic changes in genes important for specific processes and conditions

Overview



1. DNA methylation
2. Introduction of bisulfite conversion chemistry
3. Downstream analysis methods
4. **Considerations for obtaining quality data**

Improving Bisulfite-based Analysis and Data Quality

DNA Concentration Can Affect Conversion Efficiency



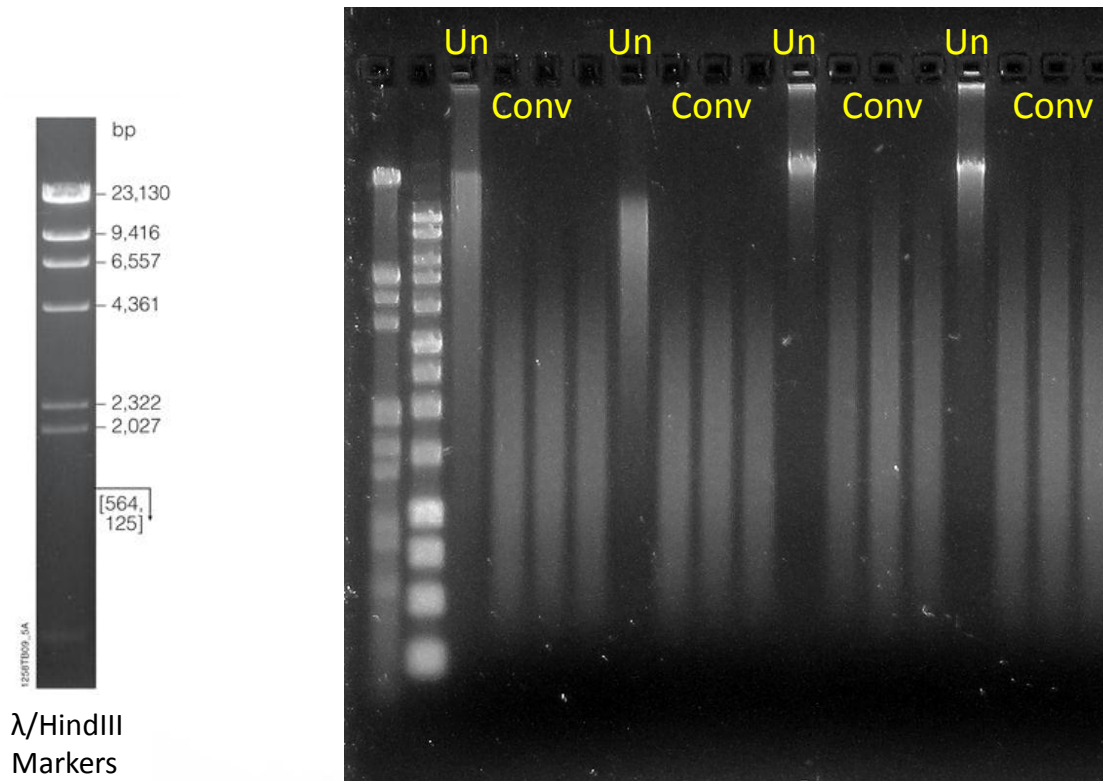
- Most commercial kits recommend converting 100-500ng per 20 μ l reaction for optimal results
 - Low end = 50pg
 - Issues: loss of sample, incomplete conversion due to bisulfite:DNA ratio
 - High end = 2 μ g
 - Issues: Incomplete conversion due to bisulfite:DNA ratio and possibly incomplete denaturation.
- Sheared vs unsheared DNA
 - Early protocols recommended extreme shearing
 - Advantage: better denaturation, avoid incomplete conversion
 - Disadvantage: starting with smaller fragments leads to even smaller fragments following conversion
 - Recommendation: light shearing with 28G needle

Improving Bisulfite-based Analysis and Data Quality

Sheared DNA Results in Smaller Fragment Size



Starting with DNA that is sheared will result in smaller fragments following bisulfite conversion.



Un: unconverted DNA

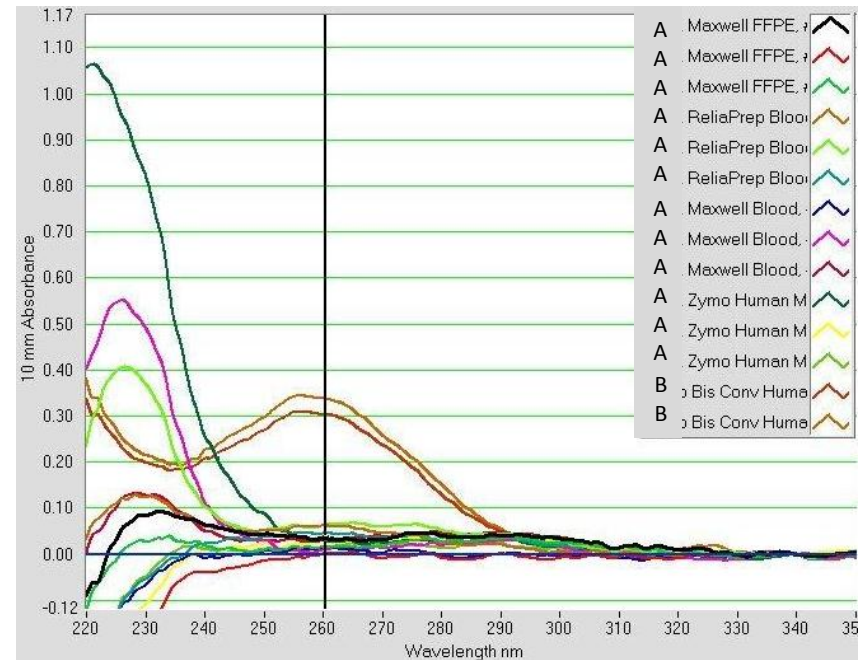
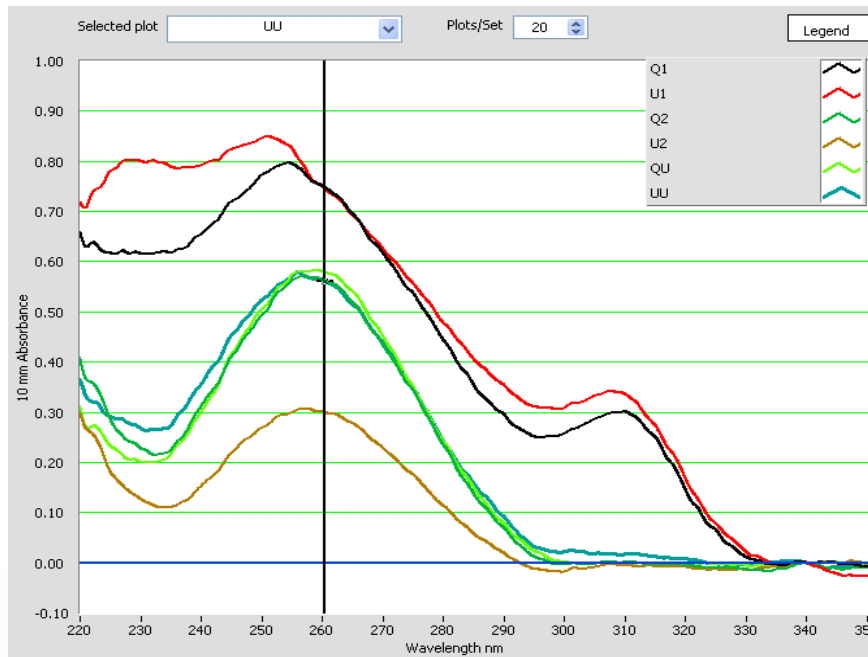
Conv: bisulfite-converted DNA

Improving Bisulfite-based Analysis and Data Quality

Absorbance Scans Can Detect Impurities



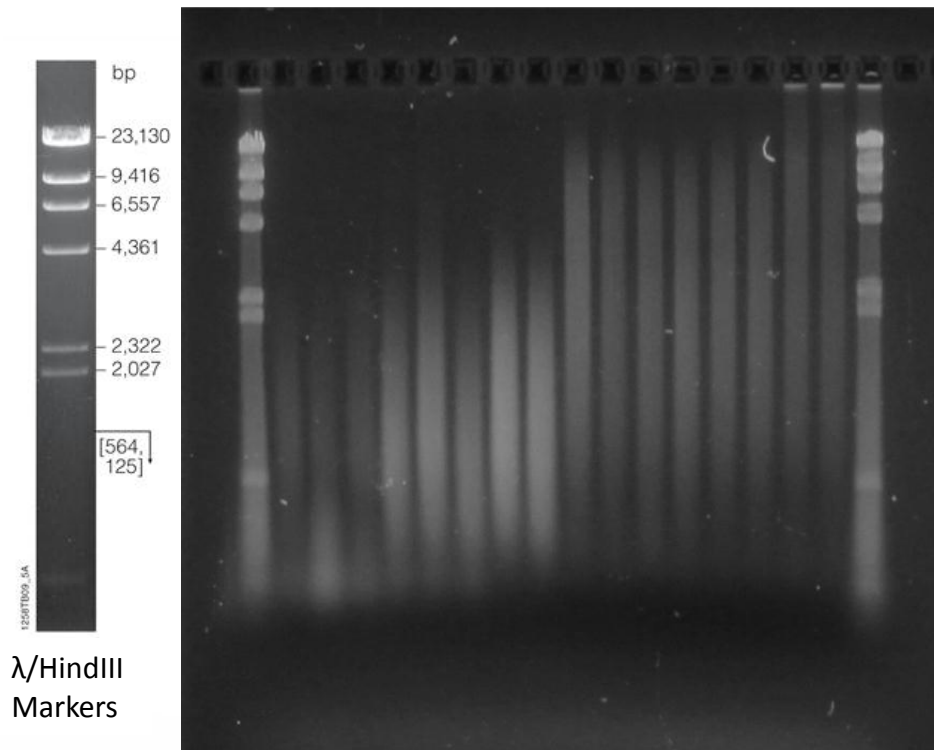
- Assess DNA quality following bisulfite conversion
 - Run absorbance scan to estimate concentration and look for impurities



Improving Bisulfite-based Analysis and Data Quality Visualizing DNA Provides Insight for Future Analysis



- Assess DNA quality following bisulfite conversion
 - Run samples on an agarose gel stained with a fluorescent total nucleic acid stain to visualize fragmentation



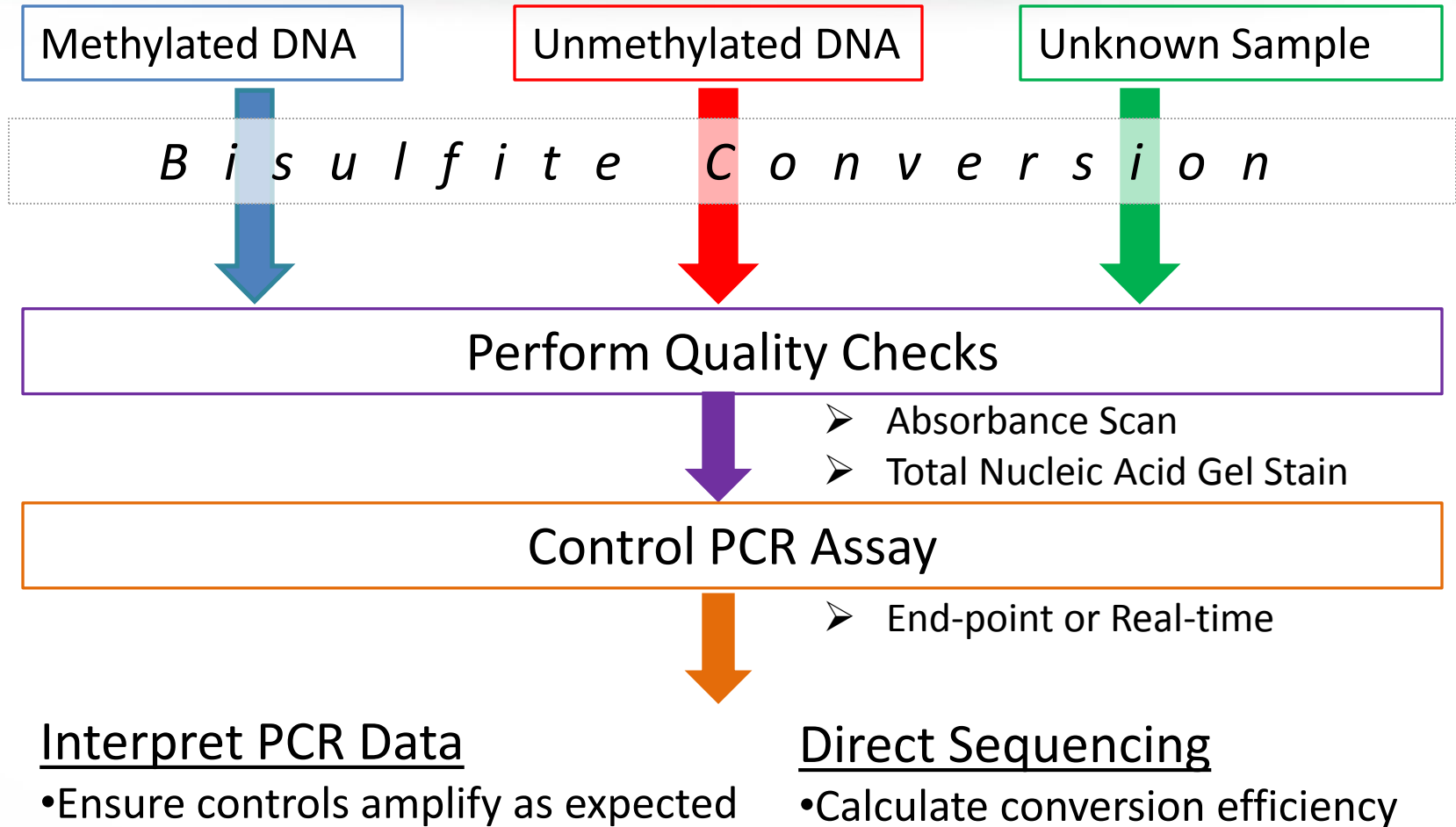
Improving Bisulfite-based Analysis and Data Quality ***Control Assays and DNA Sources Are Critical***



- Control Assays
 - Because the presence of a C following conversion is interpreted as a methylated site, conversion efficiency is critical!
 - Few published studies reference including control samples in bisulfite conversion, but studies looking at error rates indicate there are a number of factors that can influence conversion efficiency
- Always include a qualified DNA source in experiments
 - Qualified fully methylated or unmethylated DNA to run alongside unknowns during conversion
 - Qualified *bisulfite-converted* fully methylated or unmethylated DNA to run alongside *bisulfite-converted* unknowns in downstream assays.

Improving Bisulfite-based Analysis and Data Quality

Bisulfite Control Assay Workflow



Improving Bisulfite-based Analysis and Data Quality

Summary



- Bisulfite:DNA ratio is important for conversion efficiency
- Most protocols are optimal at 100-500ng DNA/20uL reaction
- Denaturation is critical for full conversion
- Light shearing of source DNA is recommended
- Control assays are critical to ensure proper interpretation of data
- If sample is available, always assess purity and quality of bisulfite-converted DNA with an absorbance scan and agarose gel

DNA Methylation Mechanisms and Analysis Methods to Study this Key Epigenetic Control



1. DNA methylation
2. Introduction of bisulfite conversion chemistry
3. Downstream analysis methods
4. Considerations for obtaining quality data

Questions?