

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

References

Introduction of chemistry:

- Shapiro, R., Braverman, B., Louis, J. B., & Servis, R. E. (1973). Nucleic acid reactivity and conformation. II. Reaction of cytosine and uracil with sodium bisulfite. *J Biol Chem*, *248*(11), 4060-4064.
- Shapiro, R., & Weisgras, J. M. (1970). Bisulfite-catalyzed transamination of cytosine and cytidine. *Biochem Biophys Res Commun*, *40*(4), 839-843.

Application to DNA Methylation Detection

- Clark, S. J., Statham, A., Stirzaker, C., Molloy, P. L., & Frommer, M. (2006). DNA methylation: bisulphite modification and analysis. *Nat Protoc*, *1*(5), 2353-2364. doi: 10.1038/nprot.2006.324
- Frommer, M., McDonald, L. E., Millar, D. S., Collis, C. M., Watt, F., Grigg, G. W., . . . Paul, C. L. (1992). A genomic sequencing protocol that yields a positive display of 5-methylcytosine residues in individual DNA strands. *Proc Natl Acad Sci U S A*, *89*(5), 1827-1831.

PCR bias:

- Moskalev, E. A., Zavgorodnij, M. G., Majorova, S. P., Vorobjev, I. A., Jandaghi, P., Bure, I. V., & Hoheisel, J. D. (2011). Correction of PCR-bias in quantitative DNA methylation studies by means of cubic polynomial regression. *Nucleic Acids Res*, *39*(11), e77. doi: 10.1093/nar/gkr213
- Rand, K. N., Mitchell, S. M., Clark, S. J., & Molloy, P. L. (2006). Bisulphite differential denaturation PCR for analysis of DNA methylation. *Epigenetics*, *1*(2), 94-100.
- Warnecke, P. M., Stirzaker, C., Melki, J. R., Millar, D. S., Paul, C. L., & Clark, S. J. (1997). Detection and measurement of PCR bias in quantitative methylation analysis of bisulphite-treated DNA. *Nucleic Acids Res*, *25*(21), 4422-4426.
- Wojdacz, T. K., & Hansen, L. L. (2006). Reversal of PCR bias for improved sensitivity of the DNA methylation melting curve assay. *Biotechniques*, *41*(3), 274, 276, 278.
- Wojdacz, T. K., Hansen, L. L., & Dobrovic, A. (2008). A new approach to primer design for the control of PCR bias in methylation studies. *BMC Res Notes*, *1*, 54. doi: 10.1186/1756-0500-1-54

Control Assays and Error

- Genereux, D. P., Johnson, W. C., Burden, A. F., Stoger, R., & Laird, C. D. (2008). Errors in the bisulfite conversion of DNA: modulating inappropriate- and failed-conversion frequencies. *Nucleic Acids Res*, *36*(22), e150. doi: 10.1093/nar/gkn691
- Grunau, C., Clark, S. J., & Rosenthal, A. (2001). Bisulfite genomic sequencing: systematic investigation of critical experimental parameters. *Nucleic Acids Res*, *29*(13), E65-65.
- Reece, K., Mook, J., & Storts, D. (2012). A Real-Time Dye-Based PCR Assay for Assessing Bisulfite Conversion Efficiency. *In Preparation*.
- Warnecke, P. M., Stirzaker, C., Song, J., Grunau, C., Melki, J. R., & Clark, S. J. (2002). Identification and resolution of artifacts in bisulfite sequencing. *Methods*, *27*(2), 101-107.

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

Comprehensive epigenetics resource: <http://epigenie.com/>

Epigenetics overview: <http://www.pbs.org/wgbh/nova/body/epigenetics.html>

NIH epigenomics website: <http://sis.nlm.nih.gov/enviro/epigenomics.html>

NIH Common Fund website: <https://commonfund.nih.gov/epigenomics/>

Agouti gene in mice: <http://www.pbs.org/wgbh/nova/body/epigenetic-mice.html>

DNA Methylation Database: <http://www.methdb.net/>