

HIGH-RESOLUTION MELT ANALYSIS OF DNA METHYLATION PATTERNS CAN DISCRIMINATE BODY FLUID OF ORIGIN IN CRIME SCENE SAMPLES

Joana Antunes^a, Vanessa Aguiar-Pulido^b, Kuppareddi Balamurugan^c, George Duncan^d, Giri Narasimhan^b and Bruce McCord^a

^aDepartment of Chemistry & Biochemistry, Florida International University

^bBioinformatics Research Group (BioRG), School of Computing and Information Sciences, Florida International University

^cSchool of Criminal Justice, University of Southern Mississippi

^dDNA Laboratory, Broward County Sheriff's Office

Body fluid discrimination can aid the forensic investigation by determining the relevance of the DNA samples collected at a crime scene. We have had success in discriminating body fluids by studying epigenetic DNA methylation (1-4), which is characterized by the presence of a methyl group in the 5' carbon of a cytosine followed by a guanine (CpG) (5).

To expand the number of CpGs for forensics, a recently published methylation study involving 450,000 individual CpGs in the human genome was performed (6). The subsequent results were next examined in detail using pyrosequencing (2, 6) or SNaPshot (7); however, not all of those CpGs showed the expected outcome.

As a result, we have applied a novel, *in silico* analysis of array data in order to define useful loci for discrimination of body fluids.

Our *in silico* analysis gives a higher weight to raw array data and uses a combination of Kruskal Wallis and ANOVA followed by Dunn's and Tukey post-hoc tests, respectively. In our initial study, CpGs with the potential to discriminate vaginal epithelia were prioritized and 8 specific CpGs were identified. The next step involved developing primers for these regions using Methprimer (8) or BiSearch (9) according to the guidelines of Wodjacz (10).

Blood, saliva, vaginal epithelia and semen were collected from volunteers according to the approved IRB protocols. DNA was extracted using the BioRobot® EZ1 automated purification workstation (Qiagen, CA) and 100 ng of DNA were bisulfite-modified using the EpiTect® Fast DNA Bisulfite Kit (Qiagen, CA). Amplification and melt-analysis reactions were performed using the Epitect® HRM kit on a Rotor Gene 6000 real time instrument (Qiagen, CA).

The results obtained show that of the initial list of 8 CpGs identified from the array, 4 produced the expected differences in T_m between vaginal epithelia and other body fluids.

Of the remaining 4 CpGs, 3 were not amplified, a result which probably indicated deviations in annealing temperature between forward and reverse primers. The 4th locus showed a similar melt temperature between semen and vaginal epithelia. The array experiment did not include DNA from semen, which may explain why our *in silico* analysis was unable to predict this similarity.

In conclusion, our combination of *in silico* and HRM analysis results in a fast PCR-based method to identify new CpGs for body fluid discrimination. Application of this process should help to accelerate discovery of additional epigenetic loci for forensic analysis.

References

1. Madi T, Balamurugan K, Bombardi R, Duncan G, McCord B. The determination of tissue-specific DNA methylation patterns in forensic biofluids using bisulfite modification and pyrosequencing. *Electrophoresis*. 2012 JUL;33(12):1736-45.
2. Silva DSBS, Antunes J, Balamurugan K, Duncan G, Alho CS, McCord B. Developmental validation studies of epigenetic DNA methylation markers for the detection of blood, semen and saliva samples. *Forensic Science International: Genetics* 2016/04;23:55-63.
3. Antunes,J.,Silva,DSBS.,Kuppareddi,B.,Duncan,G.,Alho,CS.,McCord,B. High Resolution Melt analysis of DNA methylation to discriminate semen in biological stains. *Analytical Biochemistry*. 2016;494:40-45.
4. Antunes J, Silva DSBS, Balamurugan K, Duncan G, Alho CS, McCord B. Forensic discrimination of vaginal epithelia by DNA methylation analysis through pyrosequencing. *Electrophoresis*. 2016:n/a,n/a.
5. Bird A. DNA methylation patterns and epigenetic memory. *Genes Dev*. 2002 JAN 1;16(1):6-21.
6. Park J, Kwon O, Kim JH, Yoo H, Lee H, Woo K, et al. Identification of body fluid-specific DNA methylation markers for use in forensic science. *Forensic Science International-Genetics*. 2014 NOV;13:147-53.
7. Lee HY, An JH, Jung S, Oh YN, Lee EY, Choi A, et al. Genome-wide methylation profiling and a multiplex construction for the identification of body fluids using epigenetic markers. *Forensic Science International: Genetics*. 2015 7;17:17-24.
8. Li LC, Dahiya R. MethPrimer: designing primers for methylation PCRs. *Bioinformatics*. 2002;18.
9. Tusnady GE, Simon I, Varadi A, Aranyi T. BiSearch: primer-design and search tool for PCR on bisulfite-treated genomes. *Nucleic Acids Res*. 2005;33.
10. Wojdacz Tomasz, Hansen Lise, Dobrovic Alexander. A new approach to primer design for the control of PCR bias in methylation studies. *BMC Research Notes*. 2008;1(1):54.