

## PROTEIN-BASED HUMAN IDENTIFICATION: PROOF OF PRINCIPLE USING THE HAIR SHAFT PROTEOME

Deon Anex<sup>1</sup>, Tami Leppert<sup>2</sup>, Lisa Baird<sup>2</sup>, Nori Matsunami<sup>2</sup>, Mark Leppert<sup>2</sup>, Brad Hart<sup>1</sup>,  
Glendon Parker<sup>1,3</sup>

<sup>1</sup>Forensic Science Center, Lawrence Livermore National Laboratory

<sup>2</sup>Department of Human Genetics, University of Utah

<sup>3</sup>Department of Biology, Utah Valley University

We have developed methodology to extract identifying genetic information from proteomic datasets. DNA-typing has revolutionized forensic practice and jurisprudence, however DNA often is degraded due to biological, chemical or environmental factors. Protein is considerably more stable and more abundant than DNA and persists in the environment for a longer period. Protein also contains genetic information in its primary structure, the result of non-synonymous SNPs (nsSNPs) that manifest as Single Amino-Acid Polymorphisms (SAPs). These SAPs-containing peptides are accessible to shotgun tandem mass spectrometry. We have identified nsSNP-containing peptides from 35 alleles in 26 genes expressed in the forensically informative hair shaft proteome. We obtained complex proteomic datasets from trypsin digests of the hair shafts of 54 validated European American individuals. Peptides corresponding to nsSNPs expressed in this protein population were identified and collated for each individual. The combined probability of each individual nsSNP profile was calculated using genotypic frequencies of each allelic combination in the European population (1000 Genomes Project) and the “product-rule”. The power of genetic discrimination ranged from 1 in 1,002 to 1 in 9,000. The average power of discrimination was 1 in 280. The power of discrimination increased as a function of proteomic dataset quality ( $r^2 = 0.624$ ,  $n = 58$ ,  $p < 0.0001$ ). When the power of discrimination is calculated using genotypic frequencies from the African population increased powers of discrimination are achieved. This is consistent with a decreased likelihood that the samples originate from an African origin. Relative likelihood measurements of European compared to African genetic origin range from 1 to 780 with an average of 50, a median of 18, and a standard deviation of 116. ( $n = 64$ ). Direct validation of the imputed status of each nsSNP allele was achieved with Sanger sequencing. A total of 430 genotype determinations were made from the proteomic data and 426 assignments were confirmed (specificity = 99.1%, FPR = 0.93%). The overall sensitivity was 31%. We have established a framework for the use of proteomic datasets as a source of identifying genetic information, allowing measures of identity and biogeographic background to be made from forensic or anthropological protein sources, including bone, teeth, preserved soft tissue, and trace evidence such as fingerprints.