

## GENETIC VARIATION FOR 15 AUTOSOMAL STR LOCI (IDENTIFILER) IN A PUERTO RICO DATABASE POPULATION

**Fernando Mercedes Fernández<sup>1</sup> M.S., Carmen A. Tirado Neris<sup>1</sup> M.S., Jodi Irwin<sup>4</sup>,  
Joseph Bloom<sup>1,2</sup> Ph.D., José Rodríguez Orengo<sup>1,3</sup> Ph.D.**

<sup>1</sup>*Institute of Forensic Science of Puerto Rico, San Juan, PR*

<sup>2</sup>*Department of Pharmaceutical Sciences, School of Pharmacy, San Juan, PR*

<sup>3</sup>*Department of Biochemistry, School of Medicine, University of Puerto Rico, San Juan, PR*

<sup>4</sup>*Armed Forces DNA Identification Laboratory, Armed Forces Institute of Pathology, Rockville, MD*

Amplification of short tandem repeat (STR) loci has become a useful tool for human identification and forensic case analysis because of the large number of STR loci in the human genome and their highly polymorphic nature.

The allele frequencies of 15 autosomal short tandem repeat loci included in the AmpFISTR Identifiler kit (Applied Biosystems) were obtained from a sample set of unrelated individuals living in Puerto Rico (n=211) from different geographic areas of the island taken at random, to determine the allele frequency and genotype of fifteen loci plus Amelogenin contained in the AmpFISTR Identifiler kit. For all loci "D8S1179, D21S11, D7S820, CSF1PO, D3S1358, TH01, D13S317, D16S539, D2S1338, D19S433, vWA, TPOX, D18S51, D5S818, Amelogenin and FGA.", no deviation from Hardy-Weinberg equilibrium was found. The values of heterozygosity, polymorphic information content (PIC), power of discrimination (PD), power of exclusion (PE), paternity index (PI) and matching probability (PM) were calculated.

Using statistical tool, samples were analyzed and interpreted, in order to determine the significance of allele frequency for given locus following chi-square test. In conclusion, in this study was demonstrated that Puerto Rican Data follows Hardy Weinberg equilibrium under assumptions of random mating.