GENETIC DATA OF 10 X-CHROMOSOMAL LOCI IN VITÓRIA CITY POPULATION (ESPÍRITO SANTO STATE, BRAZIL)

Martins JA^{1*}, Costa JC¹, Paneto GG¹, Gusmão L², Sánchez-Diz P³, Carracedo A³, Cicarelli RMB¹

X-chromosome STR analysis has become of increasing interest in forensic practice, especially in complex cases of kinship. Brazilian genetic data on autosomal and Y-chromosome STRs are already present in the literature, but regarding X-STRs, few studies have been published in Brazil to date. The present work investigated 10 X-STRs (DXS8378, DXS9898, DXS7133, GATA31E08, GATA172D05, DXS7423, DXS6809, DXS7132, DXS9902 and DXS6789) in a sample of 224 unrelated individuals belonging to Vitória (Espírito Santo State, Brazil). Peripheral blood was collected in FTA cards and DNA was extracted using FTA protocols. Amplification was performed in a single PCR multiplex reaction as proposed by Gusmão et al. in the GEP-ISFG collaborative study. The Hardy-Weinberg equilibrium (HWE) was examined using the exact test with Arlequin software; the mean exclusion chance of trios involving daughters (MEC_T) as well as in father/daughter duos (MEC_D) and power of discrimination in females (PD_F) and in males (PD_M) were computed as proposed by Desmarais et al. Allele frequencies between men and women samples were not significantly different. For a significance level of 0.005 (after Bonferroni's correction) no deviations from the Hardy-Weinberg equilibrium were observed. DXS7423 was the least polymorphic marker whereas the most diverse was DXS6809. Power of discrimination in females ranged from 0.952 (DXS6809) to 0.837 (DXS8378/DXS7423) and in males ranged from 0.831 (DXS6809) to 0.676 (DXS7423). The combined MEC_D, MEC_T, PD_F and PD_M were 0.99988233, 0.99999778, 0.9999999996 and 0.9999995, respectively. The 10 X-STRs analyzed constitute a powerful tool for human identification purposes in Vitória city population, Brazil.

*JAM supported by CAPES.

Financial support: PADC/FCF-UNESP and FAPESP.

¹UNESP - São Paulo State University, School of Pharmaceutical Sciences, Laboratório de Investigação de Paternidade, Araraquara, Brazil

²IPATIMUP - Institute of Pathology and Molecular Immunology of University of Porto, Portugal ³Institute of Legal Medicine. Genomics Medicine Group. University of Santiago de Compostela. A Coruña. Spain