

LATIN-AMERICAN SOCIETY OF FORENSIC GENETICS (SLAGF) RESULTS OF THE INTERLABORATORY QUALITY CONTROL EXERCISE 2010-2011

Cicarelli, R. M. B.¹, De Castro, M.², Penacino, G.³, Vela, M.⁴, Aguirre, V.⁴, Morejon, G.⁴, Burgos, G.⁴, Valenzuela, G.⁴, Builes, J. J.⁵, Gaviria, A.⁴

¹ Laboratório de Investigação de Paternidade, Faculdade de Ciências Farmacêuticas, Universidade Estadual Paulista, Araraquara, SP, Brazil.

² Referencia Laboratorio Clínico, Santo Domingo, República Dominicana.

³ Unidad de Análisis de ADN, Colegio Oficial de Farmacéuticos y Bioquímicos de Capital Federal, Rocamora 4045, Buenos Aires, Argentina.

⁴ Laboratorio de Genética Molecular, Cruz Roja Ecuatoriana, Quito, Ecuador

⁵ Laboratorio Genes Ltda, Medellín, Colombia.

Here are presented the preliminary results of the Interlaboratory Quality Control exercise 2010-2011 of the Latin-American Society of Forensic Genetics (SLAGF). At that time the exercise was composed of a practical part to determine the genetic profiles of 3 blood samples on FTA card, and a theoretical part. The theoretical exercise consisted in two paternity cases for calculations, including an obligatory case with alleged father, mother and child; the other one was a paternity case with a dead alleged father with 3 children and their mother, and an extramarital child with the biological mother. All samples and the CD report formats were sent to 85 participant laboratories on January 14, 2010. The deadline for submission of results was April 15, 2010. To date, 56 reports have been submitted by 13 different countries in Latin America and one from Spain, distributed as follows: 12 from Colombia, 10 from Brazil, 7 from México, 7 from Argentina, 4 from Peru, 3 from Venezuela, Costa Rica and Ecuador, 2 from Bolivia and 1 from Nicaragua, El Salvador, Spain, Uruguay and Dominican Republic. The results will be presented in the poster. With this Interlaboratory Quality Control exercise 2010-2011 of SLAGF, we are going to strengthen the collaboration between all participating countries in Latin America helping to promote scientific exchange, improving the development and knowledge in the Forensic Genetics Science field.