

A NEW VISION OF THE DISTRIBUTUMTION OF Y-CHROMOSOME STRs IN DOMINICAN REPUBLIC USING POWERPLEX® Y23 SYSTEM

Vilma Diaz, Laboratorio Clinico Patria Rivas

We have studied the distribution of Y-chromosome STRs and established a database within population samples from Dominican Republic by analyzing male-specific markers using Powerplex® Y23 System. Laboratorio Clinico Patria Rivas collected in the population samples and this population is truly representative of the entire country. The distribution of Y-chromosome is a powerful tool to study human evolution pathways and to infer about major and local male migration movement or patters. The aim of the present study is to present a new and complete vision of Dominican samples who is reflect the largely heterogeneous population and our work was analyze more samples plus revision of our anterior publication of Dominican Male individuals to evaluate the allelic frequency, thus, the possibility to implement a genetic database.