THE DISTRIBUTION OF ALLELE FREQUENCIES OF 15 STRS IN DOMINICAN POPULATION

Vilma Díaz¹, Patria Rivas¹, Carmen Olivero¹, Angel Carracedo²
Laboratorio Clinico Lic. Patria Rivas. Santo Domingo, Dominican Republic ²Instituto de Medicina Legal, Facultad de Medicina, Universidad de Santiago de Compostela, Galicia, Spain

The most effective tools for individual and population genetic characterization that we'll use today are STRs. This work presents the results of a population study of 15 autosomal STRs loci included in PowerPlex 16 (Promega Corporation) that has the aim of creating a local database. Buccal swabs and blood samples were obtained from more than two thousands unrelated individuals were born in Dominican Republic and lived there for at least two generations.