

STUDY OF DOMINICAN POPULATION SAMPLES CHARACTERIZE WITH 29 AUTOSOMAL STR LOCI AVAILABLE IN COMMERCIAL STR MULTIPLEX KITS

Vilma Diaz, Laboratorio Clinico Patria Rivas

One of most effective tools for individual and population genetic characterization that we will use today are STRs Markers. This work present the results of a population study with 29 autosomal STR loci available in commercial STR multiplex kits that has the aim of creating a local database consider our anterior study with 15 autosomal STR loci included in Powerplex 16 (Promega Corporation). Buccal swabs were obtained from 500 hundred unrelated individuals were born in Dominican Republic and lived there at least two generations.