

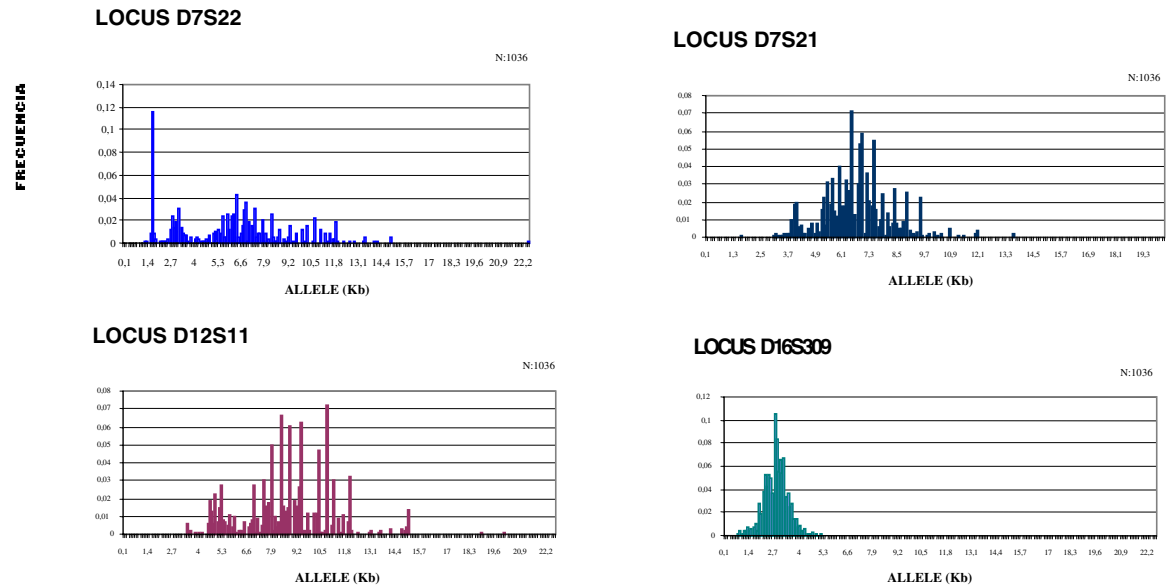
ALLELE FREQUENCIES ANALYSIS OF FOUR VNTRs LOCI IN ARGENTINE POPULATION

Ana María Di Lonardo, Sergio Fernando Valente, Andrea Hilda Szöcs, Carlos Gregorio Echenique, Marcelo Florentino, and Oscar Alberto Santapá

Banco Nacional de Datos Genéticos, Hospital Carlos G. Durand, Buenos Aires, Argentina



We analyzed allele frequencies of four VNTRs loci= D7S21 locus (MS31 probe), D12S11 locus (MS43a probe), D7S22 locus (G3 probe) and D16S309 locus (MS205 probe) by NICE--no isotopic chemiluminiscent (Lifecodes Corporation). We used HINF I restriction enzyme for the digestion of DNA. We studied 1,036 biological unrelated individuals taken from 527 caseworks of paternity testing.



The statistical parameters confirm the usefulness of these markers to exclude or assign biological relationship.