
Abstracts

Y Chromosome Polymorphisms in the Population of Galicia (NW Spain)

C Pestoni; MJ Brión; ML Cal; A González; MV Lareu; A Carracedo.
Institute of Legal Medicine. University of Santiago de Compostela.
San Francisco s/n. 15705 - Santiago de Compostela, Galicia, Spain



Different Y chromosome polymorphisms were investigated in the population of Galicia (NW Spain). The Galician population is a relatively isolated population located at the most westerly European continental edge and is therefore attractive for population studies. The lack of recombination and the mode of inheritance of the Y chromosome non-pseudoautosomal part make the Y chromosome particularly interesting for these studies. Different types of Y chromosome polymorphisms, such as STRs (DYS19, DYS385, DYS389-I, DYS389-II, DYS390, DYS391, DYS392 and DYS393), the haploid minisatellite MSY1, and one biallelic polymorphism (YAP-DYS287) were explored in order to construct highly discriminative Y chromosome haplotypes. These studies contribute to the exploration of the potential of the Y chromosome in both the anthropological and forensic fields.

