

Abstracts

A Genetic Population Study of Chromosome Polymorphisms Amplified by Means of PCR in South-West Spain

J.J. Gamero², J.L. Romero², L. Souto¹, M.I. Arufe², D.N. Viera¹, M.C. Vide¹

¹Institute of Legal Medicine of Coimbra, Coimbra 3000, Portugal

²Department of Legal Medicine, Faculty of Medicine, 11003 Cadiz, Spain



The recent discovery of a series of polymorphisms in the Y chromosome that are only inherited paternally is of great interest within the field of forensic genetics. Such polymorphisms are now considered to be an interesting tool in individual diagnosis by a large number of authors in, for example, those cases derived from forensic practice which involve crimes of a sexual nature and a fraction of male DNA belonging to the assailant is mixed with the victim's DNA in the analysed stain.

This study presents the results obtained from the genetic population study for two polymorphisms recently identified in the Y chromosome, specifically DYS385 and DYS390.

