

### **Analysis of Y Chromosomal Locus DYS390 in a Chinese Population**

Xiufen Zheng, Weiwei Wu, Zhiping Xie, Jianpo Chen, Jian Ye, Anqian Ji  
Institute of Forensic Science ,Beijing,100038,P.R.China



Y-chromosomal locus DYS390 was studied in 97 unrelated Chinese from the northwest area in China. By using PCR, vertical denatural PAGE and silver stain, six alleles were identified and the frequencies ranged from 0.0103 to 0.3916. The alleles' frequency distribution in this population showed significant difference from that of Holland, not from Germany. The gene diversity (GD) is 0.6875. The analysis of the real father-son pairs showed that there were no mutations occurring. This locus has high specificity to Y chromosomal DNA, even when there exists a high ratio of female's DNA. DYS390 will be widely used in forensic individual identification.