

GENETIC POLYMORPHISM ANALYSIS OF 20 STR LOCI IN THE HAN POPULATION OF GUANGDONG PROVINCE (SOUTHERN CHINA)

Chang-hui Liu, Hong Liu, Xing-yi Yang, Chao Liu, Guangdong Institute of Criminal Science and Technology

To study the genetic polymorphism and evaluate the 20 loci for forensic purpose, allele frequencies for 20 short tandem repeats (STR) loci included in the Powerplex 21 System (Promega, Madison, WI, USA) were obtained from 580 individuals of Guangdong Han. The accumulated powers of discrimination and power of exclusion were $1-3.58 \times 10^{-25}$ and 0.999999996, respectively. It proved the Powerplex 21 System is a powerful tool for individual identification and paternity testing in Guangdong Han population.