

## **STR DATA FOR 15 LOCI FOR THE CHINESE POPULATION IN HONG KONG**

Christina Li, Lai-fan Fung, Kwong-yuk To

Forensic Science Division, Government Laboratory, The HKSAR Government, China

Samples were collected from 309 unrelated Chinese residing in Hong Kong for analysis of allele frequencies at STR loci D3S1358, vWA, D16S539, D2S1338, D8S1179, SE33, D19S433, TH01, FGA, D21S11, D18S51, LPL, F13B, FESFPS and F13A01. No significant deviation from Hardy-Weinberg equilibrium was found in all loci. The observed heterozygosity (OH) ranged from 0.398 (F13B) to 0.927 (SE33) and the power of discrimination (PD) ranged from 0.593 (F13B) to 0.990 (SE33). The power of exclusion (PE) ranged from 0.194 (F13B) to 0.883 (SE33) and the combined power of exclusion was 0.9999973.