

**A NINE STR LOCUS MATCH BETWEEN TWO APPARENTLY UNRELATED
INDIVIDUALS USING AmpF/STR® PROFILER PLUS™ AND COFILER™**

Kathryn Troyer, Theresa Gilboy, and Brian Koeneman
Arizona Department of Public Safety, Phoenix, AZ



The Arizona Department of Public Safety Crime Laboratory is reporting a nine STR locus match between two apparently unrelated individuals at vWA, D21S11, D5S818, D13S317, D7S820, D16S539, TH01, TPOX, and CSF1PO. The samples were analysed by Myriad Genetic Laboratories, using Applied Biosystem's AmpF/STR® Profiler Plus™ and COfiler™ kits. The nine STR locus match was discovered using a duplicate search by Myriad Genetic Laboratories. Further investigation by the Arizona Department of Public Safety revealed that the two offenders varied by age and race. In addition, it was noted that the two offenders shared one allele at three of the remaining four loci, D3S1358, D8S1179, and D18S51. The match is in the process of being confirmed with Applied Biosystem's AmpF/STR® Identifier™ kit.