

## **Nine STR Loci Co-Amplified: Portuguese Population Study**

Souto, L.; Carvalho, M.; Anjos, M.J., Vieira, D.N.; Vide, M.C.  
Instituto de Medicina Legal, Largo da Sé Nova 3000 Coimbra Portugal



The recent development of new STR multiplex kits with co-amplification of increasing number of STRs, provides both higher discrimination information and productivity for forensic labs dealing with large number of samples.

In this communication a first study for a Portuguese population (Central Area) is presented for the multiplex STRs: D3S1358; vWA; FGA; D8S1179; D21S11; D18S51; D5S818; D13S317; D7S820.

A total of 100 unrelated samples were chelex extracted and amplified with the PE Applied Biosystems AmpFISTR Profiler Plus kit and detected on an ABI Automated Sequencer 377.