

Abstracts

Typing of Samples for DQA1, LDLR, GYPA, HBGG, D7S8, GC, D5S818, D13S317, D7S820, D16S539, vWA, TH01, TPOX, and CSF1PO Loci From a Body Exhumed after 27 Years

Mohammad A. Tahir, Ph.D.¹ and Sudhir Sinha, Ph.D.²

¹ IMC Forensic Services Agency, 40 S. Alabama Street, Indianapolis, IN 46204

² ReliaGene Technologies, Inc., 5525 Mounes Street, New Orleans, LA 70123

A male body was buried in 1970 and it was exhumed in September of 1997. Samples for DNA extraction were taken from various parts of the body including fingernails, teeth and bone. Extracted DNA was quantitated and typed for the above mentioned loci without any problem.

ପୁନଃ