

## Abstracts

## Distribution of Polymarker, DQA1 and D5S818, D13S317, D7S820, D16S539, vWA, TH01, TPOX, and CSF1PO loci in Bengali and Punjabi Populations

Sudhir Sinha, Ph.D.<sup>1</sup>, Muhammad Amjad, Ph.D.<sup>2</sup>, Bruce Budowle, Ph.D.<sup>3</sup>, and Mohammad A. Tahir, Ph.D.<sup>1</sup>

<sup>1</sup> ReliaGene Technologies, Inc. 5525 Mounes Street, New Orleans, LA 70123

<sup>2</sup> IMC Forensic Services Agency, 40 S. Alabama Street, Indianapolis, IN 46204

<sup>3</sup> FBI Academy, Quantico, VA



PCR-based typing of polymarker, DQA1 and D5S818, D13S317, D7S820, D16S539, vWA, TH01, TPOX and CSF1PO loci from blood DNA samples collected from Bengali and Punjabi populations of Pakistan and India was used to determine allele frequency and to construct a database. Tests of departure from Hardy-Weinberg equilibrium in these samples showed no significant deviation.

