

THE VERSAPLEX™ 27PY SYSTEM-A NEW PROMEGA SIX-DYE MULTIPLEX FOR CASEWORK

Dawn Rabbach, Promega Corporation

The VersaPlex™ 27PY System is a 6-color, 27-locus STR system for simultaneously amplifying and genotyping 23 autosomal loci, three Y-STR loci, and Amelogenin. It is intended for forensic casework or kinship STR analysis in regions that prefer the D6S1043 locus. In addition to D6S1043, the autosomal loci are CSF1PO, FGA, TH01, TPOX, vWA, D1S1656, D2S441, D2S1338, D3S1358, D5S818, D7S820, D8S1179, D10S1248, D12S391, D13S317, D16S539, D18S51, D19S433, D21S11, D22S1045, Penta D, and Penta E. This 27-locus STR system will be the only D6S1043-containing multiplex that includes multiple Y-STR loci. These three Y-STR loci allow more confident determination of the number of male contributors in complex mixtures. While designed for an optimal input DNA amount of 1ng, the system is sensitive to low DNA input amounts and is robust against high concentrations of PCR inhibitors. Additionally, nine loci are under 200bp, making it useful for degraded samples. This poster includes preliminary data highlighting the kit's capabilities.