

GENOTYPING OF CHALLENGING SAMPLES USING EIGHT-DYE POWERPLEX® STR AND SPECTRUM CE SYSTEMS

Michael Lauck, Yolanda Stypula, Kristy A. Lenz, Margaret M. Ewing, Cynthia J. Sprecher and Douglas R. Storts, Promega Corporation

Traditional capillary electrophoresis (CE) is still widely used for forensic DNA typing, mainly due to its time- and cost-effectiveness. Currently available CE technology limits multiplex STR systems to 5 or 6 color channels. The new Spectrum CE system offers increased spectral capacity, thus allowing 8 color multiplex STR systems while also supporting existing 4-, 5- and 6-color systems. The Promega PowerPlex® 35GY System simultaneously amplifies 35 different loci, including the 20 CODIS core loci, Amelogenin and DYS391 for gender identification, as well as a Quality Indicator (QI). Ten additional Y-STR loci are included to enable familial searching and to assist with forensic casework on sexual assault evidence. The availability of 8 colors allows the inclusion of smaller, more numerous loci, thus increasing a laboratory's chance of success with challenging samples.

Data will be presented demonstrating the performance of the PowerPlex® 35GY System, focusing in particular on inhibited and degraded samples as well as on mixtures. Our results highlight the advantages of the new 8-color system in comparison to existing 4-, 5- and 6-color systems while also demonstrating the capabilities of the Spectrum CE System.