## UTILITY OF STR GENOTYPING SOFTWARE IN FORENSIC CASEWORK AND DATABANKS

Sean Walsh, Rhonda Roby, Mike Troutman, Candia Brown, Cydne Holt

PE Biosystems, Foster City, CA



DNA Profiling via STR analysis can be accomplished by using an integrated system that includes optimized AmpF/STR® kit reagents, multicolor fluorescent detection instruments, and specialized software for sizing and genotyping. Custom templates are available for Genotyper® software, which designate sample alleles by comparison to alleles in the corresponding allelic ladder. The role of the genotyping software as a tool in casework and convicted offender databanks will be discussed. Updated template features will be explained, including stutter filtering, off-ladder allele designations for CODIS, and automatic comparison of overlap loci when using AmpF/STR® Profiler Plus™and AmpF/STR® Cofiler™ kits. Limitations of genotyping software and the role of the analyst in the process will also be discussed.