CHOOSING AND IMPLEMENTING NEW TEST SYSTEMS

Robert C. Giles, Ph.D. GeneScreen, Inc.



An overview of possible DNA testing platforms will be presented with specific emphasis placed on the use of PowerPlex[®] 16 run on the SpectruMedix 9610 capillary electrophoresis instrument. The effect of front-end automation and the 9610 on laboratory efficiency will be discussed.

In addition to the information regarding the use of a 96 capillary electrophoresis instrument, the implementation of single nucleotide polymorphisms (SNPs) in parentage testing will also be discussed. GeneScreen's parent company, Orchid Biosciences, has developed a high-throughput platform for testing SNPs known as the UHT (ultra high throughput). The methodology, known as SNP-IT, involves single base, primer extension of nucleic acid polymorphisms followed by either indirect or direct detection of fluorescently labeled nucleic acid chain terminators. Based on several characteristics, a large panel of selected SNPs has been constructed to examine its usefulness as a potential paternity testing panel. This platform will allow for rapid, high volume testing that, when fully implemented, should provide a completely automated and cost efficient parentage test.