USE OF THE PROMEGA POWERPLEX® 18D SYSTEM AND SWAB-2 SOLUTION FOR HIGH-THROUGHPUT DNA DATABASING OF OMNI SWABS

John Teske Texas Department of Public Safety, Austin, TX

As the benefits of DNA for solving crimes continues to grow so does the need to increase the speed at which DNA profiles are generated. Changes in technology over the years have allowed DNA databasing labs to not only decrease the time necessary to generate a DNA profile, but also to decrease the number of steps involved. The PowerPlex 18D System and Swab-2 solution is one of those recent changes in technology that allows analysts to go from swab to profile in as little as one day with only five steps. This study by the Texas Department of Public Safety CODIS Lab, presented here, shows the decrease time and steps required for omni swab processing.