

BOOKING STATION DNA ANALYSIS AND SEARCHING OF CODIS DATABASES

T.F. Callaghan¹, J. Gerst², D. Hares¹

¹FBI Laboratory Division

²FBI Criminal Justice Information Services Division

Commercial Rapid DNA (RDNA) instruments have been available for three years. These platforms were developed for automated typing and STR allele calling of buccal swabs for database enrollment and searching. Results are produced in approximately 90 minutes with minimal operator intervention. The presentation will focus on the FBI's two phase plan for booking suite integration of RDNA instruments into the United States National DNA Database. Electronic linking of arrestee information to the buccal swab for database enrollment will be discussed. Enhancements to CODIS for searching and hit notification to the booking station and investigating agency will also be presented. The concept of prioritizing the searching of "crimes of special concern" for immediate lights out hit notification will be introduced. With CODIS in use by more than 50 countries, rapid enrollment enhancements for CODIS have considered possible use by other countries. The use of an Electronic Biometric Transaction Standard (EBTS) and Arrestee Enrollment Format (AEF) with fingerprint based arrest records may allow for adoption by CODIS counties using electronic fingerprint systems for point of collection DNA registration.