

## **FBI Vision for Booking Station Rapid DNA Analysis and Searching of CODIS**

Thomas Callaghan<sup>1</sup>, James Gerst<sup>2</sup>, Douglas Hares<sup>1</sup>

<sup>1</sup>FBI Laboratory Division, Quantico, Virginia

<sup>2</sup>FBI Criminal Justice Information Services Division, Clarksburg, West Virginia

Commercial Rapid DNA (RDNA) instruments have been available for five years and have been used to submit reference profiles to CODIS from accredited laboratories since 2015. 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 planned integration of Rapid DNA following a change in Federal Law allowing submission of booking station generated arrestee CODIS profiles to 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. Requirements and milestones for arrestee Rapid DNA implementation and pilot studies will be introduced. The concept of prioritizing the searching of "crimes of special concern" for immediate lights out hit notification will be explained.