

**CODIS Booking Station Rapid DNA Analysis and non-CODIS Crime Scene
Rapid DNA Applications**

Thomas Callaghan

FBI Laboratory Division, Quantico, Virginia

tfcallaghan2@fbi.gov

Rapid DNA instruments were developed for the “fully automated - hands free” typing of reference samples in under two hours for DNA profile submission to the Combined DNA Index System (CODIS). Status of the FBI’s Booking Station implantation of the Rapid DNA Act of 2017 with the Booking Station Rapid DNA Pilot will be reviewed. Rapid DNA has also been used by some agencies in non-CODIS applications to analyze crime scenes in an effort to develop investigative leads and to include or exclude a suspect in near real-time. An overview of non-CODIS crime scene Rapid DNA applications will be presented, along with an update on the on the FBI’s Crime Scene Rapid DNA Task Force that was established in 2018.