

## **RAPID DNA WORKS IN THE BOOKING ENVIRONMENT**

Gray Amick, Ph.D.<sup>1</sup>, Capt. Washava Moyer<sup>2</sup>, Sgt. Tormesha Haynes<sup>2</sup>, Sgt. Samuel Jackson<sup>2</sup>,  
Roy Swiger, Ph. D.<sup>3</sup>, Shanin Lodhi, M.S.<sup>3</sup>

<sup>1</sup> Richland County Sheriff's Department

<sup>2</sup> Alvin S. Glenn Detention Center

<sup>3</sup> IntegenX, Inc.

In collaboration with the Alvin S. Glenn Detention Center, the Richland County Crime Lab (RCSD) investigated the use of the RapidHIT ID system in the booking environment. The RAPID Trial (Richland Assesses Potential of near-Instant DNA) is a proof-of-concept study that compares time to results of arrestee DNA testing of the in-jail system relative to the current program, demonstrating the ability to generate high quality DNA profiles that meet QAS requirements for CODIS upload by non-scientist operators. In this talk, principal investigator Gray Amick discusses the study design, operator training, sample inclusion criteria, and other gritty details behind the first Rapid DNA jail pilot study in the U.S. By the end of this presentation, forensic scientists considering implementation of a Rapid DNA program will have a framework for working with their local agencies to set up their own jail pilot.