

A U.S. GOVERNMENT FIELD TRIAL USING THE RAPIDHIT™ 200 HUMAN IDENTIFICATION SYSTEM

Stefanie Gangano, Jason Werking, Roy Swiger, and Keith Elliott
IntegenX Inc., Pleasanton, CA, USA

In spring 2012, IntegenX undertook, for the second year running, a field trial of the RapidHIT™ 200 with interested parties/agencies within the U.S. government. The trial consisted of users evaluating the RapidHIT 200 capability in an operational context.

The objectives of the trial were to demonstrate ease of training and ease of use; demonstrate the ability to use the system to analyze samples and compare these with a database within 90 minutes; test robustness of both the instrument and the consumables during long haul shipment; and to demonstrate the ability of the RapidHIT to be operated in a mobile environment.

Over 70 samples were added to an existing IntegenX database (126 participants were originally added in last year's field trial), generating a database of nearly 200 subjects. Mock "crime" samples were then analyzed on the RapidHIT by the participating field operatives for comparison against this database. These samples consisted of reference cheek (buccal) swabs, other known sample types (such as ear, nose, and blood samples) and unknown high-content (such as blood, hats, and hair follicle) and low-content samples (such as touch, copper wire, and cell phones).

Results from this trial are presented and findings with respect to the trial objectives are discussed. ☘