## EARLY CUSTOMER DATA USING PROMEGA CHEMISTRIES ON THE RAPIDHIT™ 200

Stevan Jovanovich¹, Rolfe Anderson¹, Rich Belzinski¹, Keith Elliott¹, Helen Franklin¹, Stephanie Gagnano¹, Dennis Harris¹, Yuan Li¹, Roger McIntosh¹, Bill Nielsen¹, Charles Park¹, Alex Kindwall¹, Charles Troup¹, Ezra Van Gelder¹, Mattias Vangbo¹, Jason Werking¹, Steve Williams¹, David Wyrick¹, Douglas Storts², Robert McLaren², Jonelle Thompson², Brandy Cook³, Lesley Ives⁴

<sup>1</sup>IntegenX Inc, Pleasanton, CA, USA

<sup>2</sup>Promega Corporation, WI, USA <sup>3</sup>Palm Beach County Sheriff's Office, FL, USA

<sup>4</sup>Kev Forensic Services, UK

IntegenX has developed an automated DNA analyzer, capable of producing database quality STR profiles, sample in, answer out, in less than 90 minutes.

In partnership with Promega, IntegenX has designed and manufactured cartridges and assays using Promega's DNA IQ™ System for DNA purification and PowerPlex® 16HS System multiplex PCR chemistries.

In July 2012 early customers from forensic laboratories were invited for the first time to the IntegenX laboratories in Pleasanton, CA for training, followed by a technical evaluation of the RapidHIT in their own hands. The objective was to enable customers already skilled in analyzing DNA results to experience the operation of the RapidHIT first-hand, and to obtain their early feedback on the quality of results generated using the Promega DNA IQ™ System and PowerPlex® 16HS chemistries on the RapidHIT 200. Data generated by customers on several parameters such as success rates, ease of use and sensitivity were generated, back-to-back against standard manual processing of the same samples using the same Promega reagents, and are presented here.

These results demonstrate the ability of the RapidHIT to work with existing commercially available and accepted chemistries, in addition to demonstrating the flexibility of Promega's DNA IQ™ System and PowerPlex® 16HS to be applied to emerging rapid DNA platforms.