VALIDATION OF CASELYNX A NOVEL LABORATORY INFORMATION MANAGEMENT SYSTEM INCORPORATING RFID TAGS

<u>Shelley Johnson</u>, William Miller, Thomas R. Reynolds Fairfax Identity Laboratories a Division of AlBiotech, 601 Biotech Drive Richmond, VA 23235

This presentation will detail the forensic sample tracking through a new laboratory information management program called CaseLynx. The program tracks evidence from receipt to return and has incorporated Radio Frequency Identification Tags for evidence and reference sample tracking. This system was designed to include all critical aspects of laboratory processes. It provides web based integration, and contains modules for case processing, quality management features to assist with quality assurance, document control, auditing training, validations, equipment maintenance, testimony monitoring, proficiency testing and quality control. This presentation will detail the proof of concept and validation studies.