

Implementation of Next Generation DNA Expert Systems

Martin Bill
Forensic Science Service, Ltd.

The UK Forensic Science Service® is a long-standing leader in the use of expert systems for the assessment of DNA evidence. For more than a decade the FSS has used expert systems to streamline or remove human review stages in the DNA process and the level of automation and efficiency has increased year on year as software and scientific solutions have improved.

In June 2010 the FSS began the implementation of FSS DNA INSIGHT™ a revolutionary new approach to the interpretation and management of DNA data. The project has re-engineered the role of the DNA experts, the forensic process, training, workflow and culture through the introduction of novel software and theory. This presentation discusses the advances made and challenges faced by the project team during the implementation of the technical and process improvements, cultural transformation and large scale changes to 'ways of working'.

We describe the role of FSS DNA INSIGHT™ in delivering improved science and technology to approximately three hundred DNA experts and improved casework services to the criminal justice system.