

## CONVERGE™ SOFTWARE - AN ENTERPRISE SOFTWARE SOLUTION FOR FORENSIC DNA LABS

**J Deng<sup>1</sup>, R Gupta<sup>1</sup>, M Karpagavinayagam<sup>1</sup>, N Rajagopalan<sup>1</sup>, S Vijaychander<sup>1</sup>, S Goli<sup>1</sup>, P Sabadra<sup>1</sup>, J Ge<sup>1</sup>, H Kim<sup>1</sup>, S Purankar<sup>1</sup>**

<sup>1</sup> Thermo Fisher Scientific, 180 Oyster Point Blvd, South San Francisco, CA 94080

Management and analysis of DNA workflow data in forensic labs is a daunting task. The forensic DNA workflow requires numerous intricate steps, including complex analysis, procedures, and decisions. With the development of next generation sequencing (NGS) the need of managing tsunami of NGS data and the complexity of operating traditional DNA workflow with NGS workflow interchangeably makes this even more challenging. In this presentation, we will discuss an integrated software platform that helps solve the above-mentioned challenges, integrates the CE and NGS workflows presenting a unified operational view to provide confidence in the results.

Converge is built on the philosophy to provide best of the breed sample to justice solution focusing on the deep needs of validated DNA labs with functionality that can be configured and customized.

As of now, Converge consists of:

- Platform Module, which addresses the fundamental needs of validated DNA labs, such as security, auditing, e-signature, lab data management, high performance computing (HPC), integration with LIMS and backup/restore.
- Case Management Module, which presents a case centric view in your lab to manage important artifacts like samples, workflows, analysis, DNA profiles (CE and NGS) and reports. It also provides many case related operations such as document management, instrument data management, export/import, and archiving.
- Kinship Module, which can simplify and streamline your laboratory operations and reduce the complexity of the laboratory workflows and statistical analyses. This module integrates all points in the relationship testing workflow, from sample to report, which can significantly increase your lab's efficiency.
- NGS Module, which enables integration of NGS workflows, such as sequencing, data analysis, results review, persist NGS DNA analysis results into local profile database, ancestry inference, phenotypic inference, and mixture analysis. Further the Converge platform enables CE and NGS data interoperability to ease labs transition to NGS.