

RAPID DNA PROFILING AT THE SCENE OF THE CRIME: DEVELOPMENT AND VALIDATION OF THE ParaDNA® INTELLIGENCE SYSTEM

Simon Wells, LGC

Implementation of a forensics method always requires validation, but if you're developing a new approach, proving the suitability and performance of the new measurement is a multi-step process. Here we report on the developmental validation of the ParaDNA intelligence test – defining the performance of the analysis; we progress to working collaboratively with external laboratories to ensure transfer of the assay and finally progress to verification in police forces and validation of the particular approach they adopt.

Underpinning this work we use theoretical modelling and experimental measurement to ensure that the systems are fit for purpose – enabling law enforcers to quickly find the most useful evidence, include and exclude individuals and ultimately to compare profiles against a local DNA database.