

INTELLIGENCE LED INVESTIGATIONS: DOES EVERY FORCE ParaDNA?

Tobias Hampshire, LGC

Waiting times for processing of DNA evidence often run into weeks and even months, with delays in the investigative process. The LGC ParaDNA® Intelligence System delivers STR profiles within 75 minutes of sample receipt, with no destruction of sample. Labs can make informed case/resource prioritisation decisions in conjunction with LEAs and district attorneys based on this output; excluding items only containing victim/volunteer DNA, rapidly confirming initial assessments of at-scene blood patterns, and early inclusion of prolific offenders/exclusion of innocent parties, to reference a few examples.

Initial piloting labs were selected based on a plan that utilized experienced DNA scientists to act as gatekeepers to the technology, and as leaders in supporting appropriate development of advisory language to lead investigators. In parallel testing alongside full SGM+ data delivered a positive correlation sufficient for several of these to move to live casework scenarios, with recent developments including the potential for speculative searching against national DNA databases. Early data supports the argument that the system is supporting bringing offenders to justice at the earliest opportunity.

LGC ParaDNA® Instruments provide a simple, fast, robust and reliable way to profile DNA samples in a field-deployable manner, providing rapid intelligence for screening and identification of individuals and biological samples. The ParaDNA® Intelligence System delivers a 5 STR + AMEL profile in 75 minutes.