

## UPDATED AUTOMATE EXPRESS™ EXTRACTION SYSTEM PROTOCOL CARD WITH VARIABLE ELUTION VOLUMES

Keith Eskinazi, Jason Liu, Harmandeep Gill, Patrick Gray, Jennifer Elliott, Jarrett Roth, Jianye Ge

As a forensic scientist, you face several hurdles in extracting DNA from challenging samples. To help increase your chances of obtaining successful STR profiles, you'll often need to obtain either more or less concentrated DNA. For samples with low amounts of DNA, lower elution volumes provide higher DNA concentrations to help obtain the maximum information from STR analysis. In samples with high amounts of DNA, such as sexual assault case samples, higher elution volumes can automatically dilute samples for downstream analysis.

We're pleased to introduce an automated solution offering a high level of sensitivity and flexibility: the new Applied Biosystems™ AutoMate™ Express™ Forensic DNA Extraction System IC Protocol Card with updated protocol scripts (V1.1), enabling a total of 7 elution volumes on our Applied Biosystems™ AutoMate™ Express™ Forensic DNA Extraction System. Using the new protocol with blood and buccal samples, we've obtained excellent and consistent STR profile quality at different elution volumes.

### Key features

- Maximizes the chance of obtaining an interpretable STR profile from challenging samples
- The greatest range of elution volumes on the market: with 7 elution volumes from 20, 30, 40, 50 100, 200, and 250  $\mu$ L
- Multiple elution volume options allow modulation of final DNA concentration
- Less time-consuming and more consistent than manual extraction
- Works for both Applied Biosystems™ PrepFiler™ Express standard and BTA protocols
- Improved STR peak heights with lower elution volumes and higher concentration

For Research, Forensics and Paternity only. Not for use in diagnostic procedures.