

Promega/PerkinElmer Solution



Promega products are compatible with the most popular instrument platforms and come complete with optimized methods and the best technical support in the industry. Our products allow you to increase throughput and efficiency in ways that fit your laboratory's needs, whether you want to automate individual processes or your full workflow.

PerkinElmer® has developed a suite of methods for DNA extraction, qPCR setup, normalization and STR setup using Promega chemistries. The JANUS™ Forensic Workstation increases lab efficiency and comes close to complete workflow automation.

Speed

- Increase throughput tenfold compared to manual DNA isolation.
- Set up 80 samples and two DNA standards in less than 40 minutes.

Confidence

- Ensure security by 21 CFR Part 11 compatibility.
- Minimize malfunction with liquid-level sensing.

Flexibility

- Choose between disposable or fixed washable tips.
- Select the throughput you need by choosing either the 4-tip or 8-tip configuration.

Automation will allow you to

- Increase consistency and reduce human errors
- Focus on data analysis instead of sample preparation
- Improve overall laboratory efficiency by automating additional processes such as dispensing

Automated applications

- DNA extraction using the trusted DNA IQ™ System
- Quantitative PCR setup with the Plexor® HY System for human and male DNA quantification
- Normalization and PCR setup for PowerPlex® Systems

To find out more, please call us at: 1-800-356-9526 (U.S.) or at: 608-274-4330 (International) or send us an e-mail at: genetic@promega.com

References:

Pizzamiglio M. *et al.* (2006) Robotic DNA extraction system as a new way to process sweat traces rapidly and efficiently. *International Congress Series*. **1288**, 598–600

Biondi M. *et al.* (2003) Forensic sample processing using a robotic workstation: Automated paper-based spotting of whole blood convicted offender samples and high throughput DNA isolation for STR analysis. *Proceedings from 14th International Symposium on Human Identification*.

Tack L. and Grunst T. (2002) Automated forensic DNA isolation from blood using Promega's DNA IQ System and the MultiPROBE II Forensics DNA Workstation. Poster: SAFS Fall Annual Meeting

Tack L. (2003) PerkinElmer's MultiPROBE II Forensic Workstation. Application Note: AAFS 55th Annual Meeting

DNA IQ is a trademark, and Plexor and PowerPlex are registered trademarks of Promega Corporation.

JANUS is a trademark of Perkin Elmer, Inc.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

