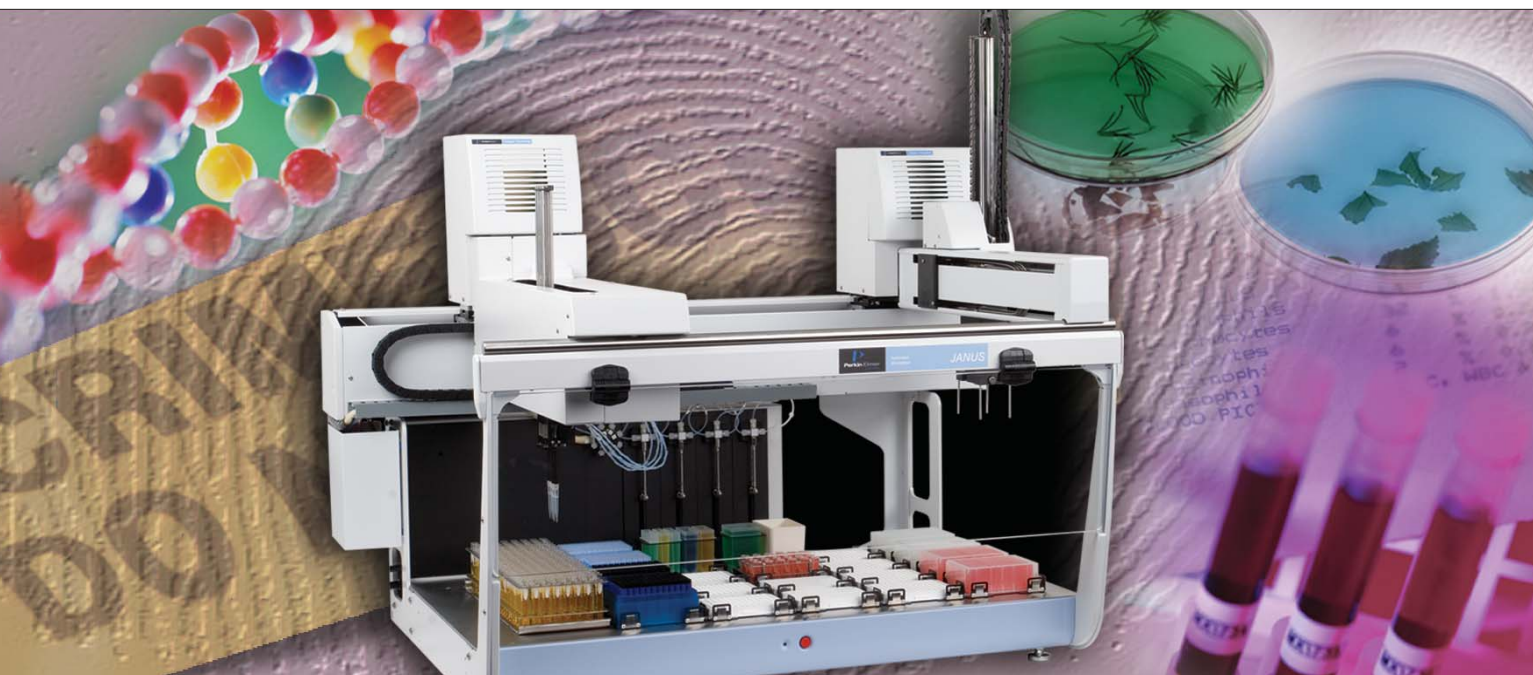


JANUS

Forensic Workstation



for high throughput
DNA sample processing

Now here's an **automation tool** designed for the **unique needs** of the **forensic lab**

The JANUS® Forensic Workstation from PerkinElmer streamlines DNA isolation, quantitation and preparation procedures to increase productivity, eliminate human error, reduce contamination, and increase data integrity and traceability. You'll realize higher throughput sample processing along with reliable and high quality Short Tandem Repeat (STR) typing analyses.

The Forensic Workstation comes with built-in automated DNA protocols that let you:

- Save hours on forensic DNA isolation – extracts DNA from virtually any type of casework or database forensic sample.
- Speed up real-time PCR (QPCR) – quickly determine sample suitability for forensic DNA typing.
- Ensure successful downstream STR typing analysis – automate DNA normalization for widely varying concentrations of DNA (10,000-fold range) and PCR setup.

Streamline sample processing with proven JANUS performance

The JANUS system is known for its reliability, flexibility and user-friendly software, and has become a hallmark robotic system for genomic, life science and pharmaceutical laboratories around the world. The powerful Forensic Workstation with the JANUS Automated Workstation is a fully automated sample preparation and liquid handling robotic system that will seamlessly streamline the processing of casework and database samples in forensic laboratories.

- Provides automated protocols and the necessary hardware to conduct the fundamental processes of DNA analysis.
- Reliably performs tedious and repetitive tasks often associated with these applications.
- Uses a single protocol to extract DNA from various sources to deliver reliable, consistent, automated DNA isolation in microplate or single tube format for rapid throughput with decreased human error.

- Isolates DNA from various sample types in the same plate.
- Automates all DNA sample processing including DNA isolation, quantitation, normalization and PCR setup for STR DNA typing analysis.

Forensic applications-ready workstation

The following walk-away, automated protocols are available for all the procedures necessary to prepare samples for forensic DNA typing:

- DNA isolation
- DNA quantification
- DNA normalization and dilution
- PCR setup for STR DNA typing analysis

Save hours on forensic DNA isolation

Automating DNA isolation using the Forensic Workstation and a Promega® DNA IQ™ System protocol increases throughput at least 10-fold compared to manual processing. The system automates the pipet-

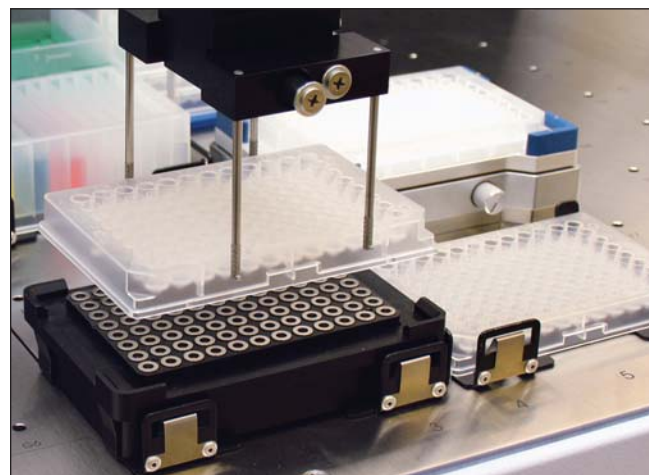


Figure 1. The JANUS gripper provides full automation of applications by providing on and off deck labware movement between a variety of peripheral devices. The toroid magnet, designed specifically with automation in mind, effectively captures paramagnetic particles for efficient washing and elution steps within the DNA IQ application.

ting, magnetic separation, heating, and shaking/mixing steps of the DNA IQ protocol, improving performance and increasing productivity. DNA is extracted from 96 casework samples in less than two hours. DNA contained in blood-stained Whatman® FTA® card punches can also be processed with high sensitivity using the Forensic Workstation.

Speed up QPCR

PCR (RT-QPCR) analysis using the ABI Quantifiler® Kits. Both human DNA and Y-specific DNA sequences can be quantified in the same plate. Additionally, we have automated the Centre for Forensic Sciences (CFS, Toronto, ON) QPCR method using only 5 µl of DNA sample. Eighty samples (plus sixteen standards) can be processed in approximately thirty minutes, providing rapid quantitation of DNA in samples and determination of suitability for forensic DNA typing. Once QPCR DNA concentrations are obtained, JANUS can then provide down-stream processing such as DNA normalization protocols.

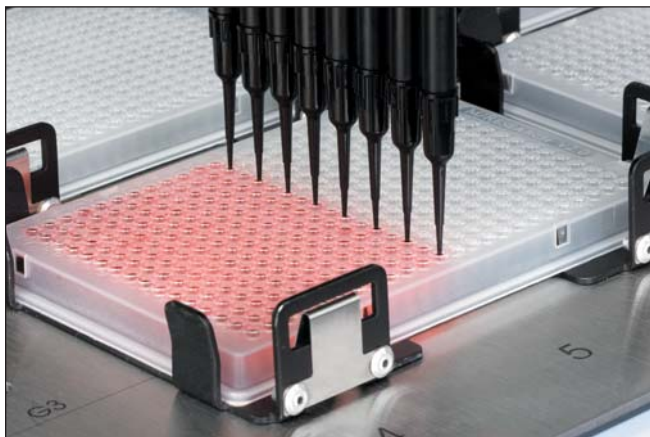


Figure 2. The JANUS Varispan arm (8-tip shown), provides the flexibility and pipetting precision needed for full forensic application processing from DNA purification and normalization to PCR master mix creation and PCR setup.

Ensure successful downstream STR typing analysis

To ensure successful downstream STR typing analysis, the Forensic Workstation automates DNA normalization and PCR setup. The protocol carries out all steps for forensic DNA normalization in the 40-367,000 pg/µL range, and PCR setup using the ABI AmpFLSTR® Profiler Plus® Amplification Kit.

Using the results from the QPCR analysis, the workstation dilutes and normalizes all DNA samples to a fixed concentration (1 ng/5 µL). It creates the PCR master mix and assembles 15 µL - 25 µL reactions in a 96-well PCR plate by using automated protocols. Processing time is less than forty minutes for eighty samples and two DNA control standards. The process is easily amenable for use with any commercial forensic STR, Y-STR and mtDNA analysis kit such as Promega® PowerPlex® and ABI COfiler®.

Move your forensic lab forward with JANUS automation

The Forensic Workstation is the answer to your DNA sample processing challenges. You'll eliminate backlogs and improve quality and consistency with the proven user-friendly automation this instrument provides. It's everything you need when you cannot do without:

Enhanced reliability for error-free processing and superior data quality

- Delivers consistent, reliable and proven protocols from DNA isolation to PCR set-up.
- Reduces human handling to help eliminate contamination.
- Ensures sample integrity with no cross-contamination or carry-over from adjacent wells.
- Handles automated spotting of barcode-labeled blood samples onto FTA® cards for database archival of convicted offender whole blood samples.

More flexibility to handle several protocols and higher throughput reducing processing times and backlog

- Analyzes 96 forensic DNA samples in less than two hours – a 10-fold increase in throughput when compared to manual processing.
- Processes numerous sample types, including swabs, fabric, whole blood, blood stains, and gum.
- Varispan™ feature enables reformatting of samples from tubes to microplates.
- Handles magnetic particle or filtration-based protocols for either large database or small case-work samples.
- VersaTip® PLUS technology enables use of disposable or fixed washable tips and changes “on the fly” between or within an assay; several volume options. Disposable filter tips reduce sample cross-contamination.
- Delivers high throughput processing with either 4-tip or 8-tip JANUS configurations.
- Independently programs heating of up to 4 microplates from ambient to 105 °C with the heater option.

Reliable security and safety for preserving the chain of evidence throughout the process

- 21 CFR Part 11 compatibility ensures security of the system throughout a process by providing electronic signature, instrument access security, data security and audit logs.

- Includes Microsoft Access® database software for real-time, user-friendly cataloging capability to store run log files related to the protocol for auditing, post-run analysis, transfer of data to Excel® for manipulation, and to provide input data for other instruments.
- Real-time error feedback with Liquid Level Sensing (LLS) detects and pauses the protocol if a malfunction is detected, saving time and precious reagents.
- Reagent batch ID entry allows for assigning reagent lot numbers to reagent troughs for tracking and report generation. Assign and record batch ID's for all the reagents used in a protocol for easy reproducibility at a later date.

Automation designed to improve forensic laboratory productivity

With ever increasing demand on analysis of DNA evidence for database and casework samples, automation has become a necessity. The Forensic Workstation from PerkinElmer provides a complete automation solution for meeting today's demand for forensic DNA analysis. Not only is the workstation reliable and flexible, it increases throughput, minimizes human error and provides superior data quality to manual processing. It comes with standardized protocols for a variety of procedures that will simplify sample handling, DNA extraction, quantification and PCR set-up in a timely fashion. The Forensic Workstation is the answer to reducing your backlog without compromising quality.

For more information, visit www.perkinelmer.com/forensicworkstation or call us at 1-800-762-4000.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/asooffices

©2007 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. Varispan is a trademark and JANUS, PerkinElmer, PlateStak and VersaTip are registered trademarks of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. DNA IQ is a trademark and MagnaBot, PowerPlex and Promega are registered trademarks of Promega Corporation. FTA and Whatman are registered trademarks of Whatman International Ltd. AmpFLSTR, COfiler, Profiler Plus and Quantifiler are registered trademarks of Applied Biosystems Corporation. Microsoft, Access and Excel are registered trademarks of Microsoft Corporation. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.