

Promega Maxwell® 16 System



The Maxwell® 16 instrument from Promega offers forensic laboratories fast, reliable sample processing and provides high-quality automated DNA and RNA purification at a fraction of the cost of high-throughput instruments.

Comprised of a pre-programmed instrument and pre-filled reagent cartridges, the system delivers consistent purification that saves time and labor. At just the touch of a button, you achieve a greater level of confidence in your results. Fast purification of up to 16 samples, we call it Personal Automation™.

Better Productivity

- Achieve consistent results time after time
- Truly “plug-and-play” technology

Efficiency

- Add samples or lysate to pre-filled reagent cartridges
- Obtain purified DNA in 30-45 minutes

Flexibility

- Process 1 to 16 samples at a time
- Easily switch between casework and reference sample setup



Maxwell® 16 System

- Compact, bench-top instrument (12.8 × 17.3 × 12.9 inches)
- Pre-filled reagent cartridges available for both DNA and RNA purification
- Proven DNA IQ™ chemistry
- Integrated system from one supplier
- Affordable automation ideal for use as a stand-alone instrument or to complement a larger robotic platform

Automation will allow you to

- Increase consistency and reduce human errors
- Focus on data analysis instead of sample preparation

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:

- Felske-Zech H., et al. (2008) The Maxwell® 16 Low Elution Volume System for Forensic Casework: Implementation and Routine Use in a Forensic Laboratory. Profiles in DNA, 11(2), 9-11.
- Krnajski Z., Geering S., and Steadman S., (2007) Performance verification of the Maxwell 16 instrument and DNA IQ Reference Sample Kit for automated DNA extraction of known reference samples. Forensic Science, Medicine, and Pathology. Vol. 3. 264-269.
- Bjerke, M. et al. (2006) Forensic applications of the Maxwell® 16 instrument. Profiles in DNA, 9(1), 3-5.

Maxwell is a registered trademark of Promega Corporation.

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

