

## Promega/Beckman Coulter 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.

Beckman Coulter offers a range of automation platforms. The Biomek<sup>®</sup> 3000, Biomek<sup>®</sup> NX<sup>P</sup> and Biomek<sup>®</sup> FX<sup>P</sup> workstations allow labs to choose the automation capabilities and throughput level appropriate to their needs.

### Speed

- Process up to 40 differential extractions from sample to eluted DNA in under 5 hours with less than 1 hour of hands-on time.
- Isolate DNA from 96 samples in 2 hours.

### Efficiency

- Eliminate the need for pipette-based mixing with the integrated shaker.
- Count on methods developed and supported by chemistry automation experts.

### Flexibility

- Process a variety of casework and database sample types.
- Choose the platform that best fits the needs of your laboratory.

## Beckman Coulter Systems

### Biomek® 3000

- Interchangeable tools system with gripper, single channel and 8-channel pipetting capability
- Tube-to-plate pipetting capability
- Liquid-level sensing with single channel tools
- Cost effective entry-level instrument

### Biomek® NX<sup>P</sup>

- Single head with Span-8 independent pipetting and gripper
- Tube-to-plate pipetting capability
- Liquid-level sensing
- Independent well access
- Optional integrated barcode reader

### Biomek® FX<sup>P</sup>

- Dual-head system incorporating 96-well pipetting and gripper on one head with Span-8 independent pipetting on second head
- Tube-to-plate pipetting capability
- Liquid-level sensing with Span-8 tips
- Independent well access
- Optional integrated barcode reader
- Increased sample throughput

### 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

- Automated differential extractions featuring the Differex™ System.  
*Process up to 40 differential extractions in under 5 hours with less than 1 hour of hands-on time.*
- DNA extraction using the trusted DNA IQ™ System  
*Isolate DNA from 96 samples in 2 hours or less depending on the platform used.*
- Quantitative PCR setup with the Plexor® HY System for human and male DNA quantification
- Normalization and PCR setup for PowerPlex® Systems for the NX<sup>P</sup> and FX<sup>P</sup> robotic platforms
- Full workflow automation with complete sample tracking

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](mailto:genetic@promega.com)

### References:

- McLaren, B., Bjerke, M. and Tereba, A. (2006) Automating the DNA IQ™ System on the Biomek® 3000 laboratory automation workstation. *Profiles in DNA*, **9(1)** 11–13.
- Olson, R. and Cowan, C. (2007) Automating the Differex™ System. *Profiles in DNA*, **10(2)** 19–20.
- McLaren, B. (2006) Automation in a forensic laboratory: An update. *Profiles in DNA*, **9(2)** 19–20.

Biomek is a registered trademark of Beckman Coulter, Inc.  
Differex and DNA IQ are trademarks, and Plexor and PowerPlex are registered trademarks 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.

