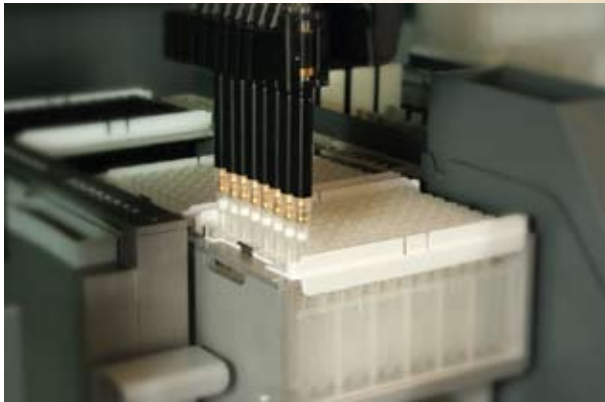


Promega/Tecan Solution



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.

The Tecan Freedom EVO® series offers versatile liquid-handling robotic workstations serving a full range of applications. Tecan presents the Freedom EVO® platform in several different sizes and offers an optional enclosure to minimize contamination.

Low Operating Costs

- Reduce disposable tip usage through reracking.
- Minimize reagent waste by processing anywhere from 1 to 96 samples.

Security

- Verify presence and absence of liquid volumes with liquid-sensing capability.
- Confirm labware positions with optional PosID™ Positive Identification System.

Flexibility

- Process samples from a deep-well plate or individual tubes thanks to adjustable tip spacing.
- Select the lower cost 4-tip instrument or a higher throughput 8-tip instrument.

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 aliquoting

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 1.5 hours.
- 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:

- Heranz, J.A., Poveda, E.G. and Fernandez, F.A. (2008) DNA extraction using the Tecan Freedom EVO® 200 and DNA IQ™ System. *Profiles in DNA*, **11(1)**, 7–8.
- Cowan, C. (2006) The DNA IQ™ System on the Tecan Freedom EVO® 100. *Profiles in DNA*, **9(1)**, 8–10.
- 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.

Freedom EVO and PosID are trademarks of Tecan AG Corporation.

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.

