

INNOVATIONS THAT RESHAPE **SAMPLE PREP** FOR BIOANALYSIS



SAME PROCESS, SAME RESULTS, LESS HANDS-ON TIME

Proteomics research presents issues of limited and variable sample material and degradation, vast dynamic range and a plethora of post-translational modifications. Whether analyzing a single protein or an entire proteome, automation can help achieve enhanced reproducibility while also delivering cost- and time-saving benefits to proteomics researchers. Automation solutions are central to these improvements, and also to linking proteomics-based information with related areas of research.

The Biomek automated workstations offered by Beckman Coulter offer a simple, rapid, accurate and reproducible solution for the quantitative analyses that are routinely processed manually in analytical laboratories.



Biomek 4000



Biomek NX[®]

The Biomek liquid handling systems present a dynamic range of pipetting capabilities that can be configured for semi- or complete automation of protein assays and meticulous management of precious protein samples.

By automating the sample preparation prior to analysis with the Biomek automated workstation, you can achieve:

- Higher throughput
- Reduced active bench time, freeing up resources
- Minimized opportunities for human error and user-to-user variability
- Repeatability; obtain crucial samples and answers, detect smaller difference in samples, and reduce re-tests, leading to faster insights

With innovations in automation for sample preparation, Beckman Coulter products are at the forefront of research in biopharma research, clinical research, biomarker discovery and biologics analysis.

For more information, visit beckman.com.



Biomek FX[®]



Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at beckman.com
© 2016 Beckman Coulter, Inc.

AAG-1467FLY02.16-A