

## **SAMPLE TO ANALYSIS AUTOMATION WITH PROMEGA CHEMISTRIES**

Aaron Bonk, Ph.D., Joseph Bessetti, M.S., Cristopher Cowan, Ph.D.  
Promega Corporation, 2800 Woods Hollow Rd, Madison WI 57311

Processing forensic and reference samples typically involves extracting DNA , quantitating the extracted DNA and preparing an STR amplification reaction for each sample. Manual sample processing is slow, labor-intensive and subject to human error. Automated liquid handlers can process samples quickly with minimal possibility for errors or cross-contamination. Promega offers automated methods for the Biomek 3000 for extracting DNA with the DNA IQ™ System, for quantitating samples with Plexor® Systems, and for STR analysis with PowerPlex® 16 Systems. In summary, Promega offers a complete package of chemistry and methods for automated processing from extraction to analysis.