

## **STACS DNA DELIVERS THE DNA LAB ESSENTIALS**

Malena B. Jimenez, B.S., Hays Young, Ph.D., STACS DNA

Let's cut straight to the point. You have more work than available resources. You have rush cases. You have trainees. Someone used an expired reagent and no one ordered additional consumables to replace the expiring ones. You forgot your thermal cycler's annual certification was due. Management wants your stats for last month and the NIJ wants your grant reports. New legislation is being proposed for sexual assault kit turnaround within 30 days. You have an audit next week. And, you just spilled coffee down your shirt.

DNA life is hard. Let us make it easy. STACS DNA delivers the DNA lab essentials: Lab Management, Sample Processing, Compliance, and Integration. Current customers report a 100% time savings with reagent and instrument tracking, an 80% time savings with managing samples and instrumentation and resolving troubleshooting issues, and a 40-100% time savings with documenting cases, performing peer review, preparing documents for legal proceedings and preparing for and conducting audits. This allows users to process 43% more samples and complete 30% more cases. STACS DNA improves documentation quality, detail, and standardization. About the only thing we can't do for you is remove that coffee stain.