

CASE STUDY



"With Promega's automated stocking program, we can focus on the most important task—making the correct diagnosis. We don't have to worry about running out of reagent or tracking down lot numbers. With Promega, it's all hands-free, automatic and documented."

*- John Fulmer, Operations Director,
Avero Diagnostics*

The Company -- Private Reference Diagnostics and Research Laboratory

Avero Diagnostics is a private reference diagnostics and research laboratory. The company, founded in 2004, conducts diagnostic testing for clients across the United States. Avero's research and development arm is committed to improving patient diagnosis and outcome for a variety of cancers by advancing diagnostics.

The Need -- Streamline Workflow and Increase Efficiency, Reproducibility

Avero's diagnostics laboratory conducts approximately 15,000 HPV tests annually. The growing laboratory's 31 technicians manage a huge clinical workload. In addition to performing diagnostic tests, they are responsible for tracking lot-specific information for all of the reagents used during sample processing and testing.

The laboratory director wanted to automate as much of the testing process as possible to streamline workflow and increase efficiency and reproducibility. He visited the Promega web site and discovered a potential solution--the Maxwell® 16 Instrument and the Maxwell Total Viral Nucleic Acid Purification System. The lab became a beta test site for the bench top automation system.

The Solution -- Automated Stocking System for Ordering, Tracking

Based on workload, the lab calculated they would use 16 Maxwell reagent systems per month. By using Promega Helix™ intelligent on-site stocking, Avero technicians were free from tracking lot-specific reagent information on the nucleic acid purification systems. Helix automatically managed ordering, inventory and tracking which saved countless hours of administrative time.

The Helix system offers additional benefits:

- Wireless units can be moved to where they are most needed, providing flexibility as the laboratory grows and changes.
- The stocking system can expand as the laboratory process new sample types.

Avero Diagnostics is considering automating processing of additional sample types, including tissue and blood. The company may add a new Maxwell Instrument and third Helix unit to manage the reagents for these new sample types.