

Introducing the Newest Member of the Promega Genetic Identity Team

At Promega, we know that providing excellent technical support enables customer success. We are pleased to announce the newest member of our technical support team.



CURTIS KNOX, B.S.

Curtis Knox joined the Genetic Identity team in January of 2004 as a Technical Services Scientist. Curtis received his B.S. degree in Genetics from Iowa State University. Before joining Promega, he worked at the Wisconsin State Crime Laboratory in Madison, Wisconsin, for 7 years and the North Louisiana Crime Laboratory in Shreveport, Louisiana, for 3 years.

In his new role at Promega, Curtis will be responsible for training customers, providing support for Genetic Identity products and applications. He will also develop educational talks and represent Promega at meetings. His extensive experience includes extracting and amplifying DNA samples, interpreting and reviewing DNA evidence, and presenting this data to law enforcement officials and attorneys. His 10 years of experience as a forensic scientist makes him ideally suited to be a technical resource for our customers.

Curtis and his wife live in Madison, where he enjoys fly fishing and playing soccer.



Career Opportunities

TECHNICAL SERVICES—GENETIC IDENTITY, EUROPE

We have an opening in Europe for an experienced scientist to provide technical support for Promega Genetic Identity products and related applications, coordinate and participate in projects, and assume a leadership role in supporting our customers. Successful candidates will have a B.S., M.S. or Ph.D. degree, a minimum of 3 years of lab experience in molecular biology, cell biology, forensic biology or a related area, and experience with the ABI PRISM® 310 or 3100 capillary electrophoresis platform with the ability to troubleshoot its use. Candidates should be able to travel approximately 40% of the time.

FORENSIC REGIONAL ACCOUNT MANAGER

We are looking for an experienced account manager to plan, execute and monitor sales, marketing and technical activities for Promega Genetic Identity products in the United States. Job responsibilities include direct sales efforts, attending key meetings, and organizing and conducting workshops, focus groups, and product alpha test or validation activities. Successful candidates will have a B.S., M.S. or Ph.D. degree in molecular biology or forensic biology, 2–3 years of lab experience with DNA-typing protocols and, preferably, at least 3 years of sales and marketing experience. Candidates should be able to travel domestically and internationally approximately 50% of the time. For more information, see: www.promega.com/hr/