

## New Members of the Genetic Identity Team



### LAURA BIMBASHI

Laura Bimbashi joined Promega in November as a Forensic Regional Account Manager for Europe. Laura is based in the United Kingdom and will be covering the United Kingdom, Scandinavia and other parts of Europe. She will provide technical and sales support for Genetic Identity customers.

Laura graduated with BSc (Hons) in Pharmacology from University College London and completed her postgraduate studies in Forensic Medicine at the University of Glasgow. Her background includes eight years of laboratory experience as a Forensic DNA Analyst for LGC Teddington, Sacramento County Forensic Lab and Kosovo Police Services Forensic Lab. For the past two years, she also worked as an account manager for Starlab UK.

In her free time, Laura enjoys skiing, swimming, playing basketball and spending time with her family.



### BENJAMIN KRENKE

Many of you know Ben from his years in the Promega Research and Development group. Ben joined Promega in 1999 and has been instrumental in developing multiple PowerPlex® and Plexor® products.

In his new role as Product Associate, Ben will be managing products from a different perspective as he works to identify new products and applications that will benefit the forensic scientist. Ben graduated from the University of Wisconsin with a Bachelor of Science in Biology and a Masters of Science in Molecular Biology.

While not at work, Ben enjoys biking, racing, camping and spending time with his wife and daughter.



### ERIN McCOMBS

Erin joined Promega as a Product Manager for Genetic Identity in October. In her new role, Erin will be responsible for managing overall product life cycle, which includes collecting and communicating requirements for new product development, as well as developing marketing and commercialization plans for various Genetic Identity products.

Erin graduated with a Bachelor of Science degree in Biology from Marquette University. She worked for five years as a Transplant Services Technologist at the Blood Center of Wisconsin. She has worked most recently at Invitrogen, with two years as a Technical Services Supervisor and seven years as a Product Manager for various transplant-related technologies.

In her spare time, Erin enjoys spending time with her family, gardening and hiking.