

Newest Members of the Genetic Identity Team



LEN GOREN

Len joined Promega in August as Director of Genetic Identity. His role includes global responsibility for marketing, sales and technical support. Len has an extensive background in marketing, business development, sales, engineering and operations management. He spent the last 16 years with Sartorius, most recently as a Senior Vice President. At Sartorius, Len's focus was the design, manufacturing, marketing and support of instruments and equipment.

Len recently completed his MBA at IMD (University of Lausanne) in Switzerland. His graduate studies include Masters work in molecular biology at Wayne State University School of Medicine and a B.S. in Microbiology from Michigan State University. He is a member of the Biopharm International editorial board, served on the board of the State of Pennsylvania Bioprocess Resource Center, and was an instructor at Pennsylvania State University where he led biotech workshops.

Len enjoys spending time with his wife and children, home-remodeling projects, skiing, biking, racquetball, theatre and travel.



MATTHIAS LINDNER, DR. RER. NAT.

Matthias joined Promega Germany in 2003 as a Technical Services Scientist. Over the years, he assumed more responsibilities for supporting the DNA-typing community by providing PowerPlex® Systems support. In March, Matthias officially joins the European Genetic Identity team as a Technical Services Scientist.

Matthias received his Diploma in Biology from the University of Heidelberg, Germany, where he also earned his Ph.D. at the German Cancer Research Center. In his new role, Matthias will be involved in customer training and technical support for Genetic Identity products and applications. He will also develop educational talks and represent Promega at meetings.

Matthias enjoys parks and gardens, hiking and graphical art. He will be based in Paris, France. Please join us in welcoming him to the GI team, and feel free to contact him at: matthias.lindner@promega.com



MELISSA SCHWANDT

Melissa joined Promega in February as a Technical Services Scientist—Genetic Identity. Melissa received her B.S. in Biology from the University of Wisconsin—Stevens Point and her Ph.D. in Experimental Pathology from the University of Utah. Before joining Promega, she worked at the Wisconsin State Crime Laboratory for 5.5 years and the New York City Office of Chief Medical Examiner for 3 years. Most recently she was involved in validating the Tecan Freedom EVO® 100 for DNA casework at the Wisconsin State Crime Laboratory. She has experience extracting, quantifying and amplifying forensic DNA samples, interpreting results and presenting those results to the legal community.

At Promega, Melissa will be responsible for customer training and support for Genetic Identity products and applications. She will represent Promega at meetings and develop and present educational talks to the DNA-typing community.

Melissa and her husband live in Roxbury, Wisconsin, where she enjoys reading the classics, cooking and spending time with her nieces and nephews.