

Newest Members of the Genetic Identity Team



DORA GRANSDEN

Dora joined the European Genetic Identity team at Promega in January as a European Account Manager based in the UK. Her role involves providing both sales and technical support for Genetic Identity customers.

Dora has over 14 years of experience in laboratory work with 10 years at the Institute for Animal Health, where she worked on a project mapping developmental and drug resistance genes in *Eimeria tenella*. Before joining Promega, she spent more than four years in a private forensic DNA laboratory in the UK, extracting, profiling and analyzing forensic casework samples.

Dora enjoys socializing, shopping, reading and drinking the occasional glass of red wine.



LOTTE DOWNEY

Lotte joined Promega in June as the Genetic Identity Marketing Manager to develop and implement worldwide marketing strategies for Genetic Identity products and technologies.

She comes to us from Applied Imaging where she was the Marketing Manager for Cytogenetics. There she led efforts to develop automated systems for cytogenetic laboratories. Lotte has also worked at MiraiBio, Inc., and Genomic Solutions. Originally from the Netherlands, she has a Masters Degree in biology and business administration.

When she is not working, Lotte loves spending time with her family and scrapbooking. She recently relocated from California to the Midwest and is looking forward to a "White Christmas".



MATT SELIGA

We are pleased to introduce Matt as the new Field Regional Account Manager for the western United States. Matt's responsibilities include providing product information and customer training, managing sales and representing Promega at various human identity meetings throughout the year.

Before joining Promega, Matt worked at Applied Biosystems, where he was a Forensic Technical Application Specialist. His experience includes supporting gel-based and capillary-based instruments and associated software programs, including GeneScan®, Genotyper® and GeneMapper®.



BRET LIGHT

Promega welcomes Bret as the new Manager, Global Sales and Tech Support.

Bret has a bachelor's degree in Zoology and master's degrees in Molecular Pathology and Business Administration. Bret has spent the last three years as the Eastern Regional Manager in the Promega Life Science area. Before that, he worked at Hitachi Genetic Systems/MiraiBio as a sales manager and at Oncor in Research and Development and Technical Support for molecular diagnostic products.

He enjoys the outdoors, hiking, biking, kayaking, running, flying and walking his dog in the woods. He loves to spend time with his wife and two children, playing games or watching movies. He also enjoys making sawdust in his workshop on a Saturday afternoon.