

## Newest Genetic Identity Team Members



### DAVID JOHNSON

David joined Promega in January as a Technical Services Scientist—Genetic Identity. His new responsibilities include customer training and providing technical support for Genetic Identity products and applications. He will develop and present educational talks to the DNA-typing community and represent Promega at meetings.

David graduated with a B.S. in Microbiology and a B.A. in Chemistry from Arizona State University. After graduating, he worked as a surgical technician at a local hospital. For the past seven years, he worked for the Phoenix Police Department Crime Lab, where he validated the Biomek® FX method for DNA casework extraction, quantitation and normalization and gained experience in STRs, Y-STRs and mitochondrial DNA testing.

In his free time, he enjoys playing volleyball and kickball, rollerblading and biking.



### CAMI GREEN

Cami joined Promega as a Forensic Regional Account Manager. She supports customers in the Rocky Mountain area of the United States, including Texas, Oklahoma, Kansas, Nebraska, South Dakota, North Dakota, Montana, Wyoming, Colorado and New Mexico. Cami graduated from the University of Utah with a Bachelor of Science in Biology. Her background includes more than seven years of laboratory experience, with five and a half of those as a Forensic DNA Analyst in a forensic laboratory. She has worked at Myriad Genetics, Utah Bureau of Forensic Services and Sorenson Forensics.

Cami's interests outside of work include hiking, mountain biking, skiing and cooking.



### MARTINA VOKURKOVÁ CHCOVÁ

Martina joined the European Genetic Identity team in January. She is based in Prague, Czech Republic, as a Forensic Regional Account Manager, covering Central and Eastern Europe. She graduated with a Masters Degree in Microbiology and Bacterial Genetics in 2003 at Charles University, Prague. While studying for her Masters Degree, she worked at the Academy of Sciences in Prague for three years. After graduation, she worked as a genetic specialist at the Institute of Criminalistics in Prague with the Czech Police for almost five years as a genetics expert. During that time she worked on criminal cases but also on databasing and completed many Police Officer training courses and FBI CODIS training.

In her free time, Martina enjoys almost all sports, including hiking, canoeing, skiing, shooting and squash. She also enjoys reading detective stories and studying English and Russian.