

Newest Members of the Promega Genetic Identity Team

CRISTIAN CAPELLI, PH.D.



We are pleased to announce the addition of a new member to the Genetic Identity team in Europe. Cristian Capelli joins Promega as a Technical Services Scientist supporting the needs of our customers in Europe. Cristian grew up in central Italy and received his degree in Molecular Biology at the University “La Sapienza” in Rome, where he studied the role of small nucleolar RNAs in ribosome biosynthesis. He then earned his Ph.D. in Forensic Genetics at the Institute of Legal Medicine, Catholic University of Rome, where he performed mtDNA sequencing and STR analysis for casework and database setup. As part of his Ph.D. program, he spent one year at the Institute of Zoology of the Ludwig Maximilian University in Munich and one year at the Galton Laboratory of the University College of London, where he focused on ancient DNA and human population genetics. He has extensive experience in STR analysis, with particular focus on Y-chromosome markers. As part of his studies, he developed PCR-RFLP protocols for multiplex genotyping of Y-chromosome SNPs that were used for studies focused on populations in South East Asia, the British Isles and Italy.

In his new role at Promega, Cristian 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.

Cristian enjoys hiking, motorbikes and science fiction. He lives with his wife in Borgo Montello, near Rome, Italy. Please join us in welcoming him to Promega, and feel free to contact him at: cristian.capelli@promega.com.

YASSER DAOUDI, B.S.



We are pleased to introduce Yasser Daoudi, one of the newest members of our field staff. Yasser joins Promega as a Forensic Regional Account Manager. He comes to us from Applied Biosystems, where he was a Forensic Product Specialist. Before joining Applied Biosystems, Yasser worked at the Armed Forces Institute of Pathology (AFIP) performing criminal casework using mtDNA and STR analysis. His casework experience includes being one of two lead analysts on the project to identify the remains interred in the Vietnam Tomb of the Unknown Soldier.

Yasser will be responsible for supporting laboratories in the western United States, including California, Washington, Oregon, Nevada, Idaho, Montana, Wyoming, Utah, Arizona, Colorado, New Mexico, Alaska and Hawaii.

On a personal note, Yasser resides in beautiful San Francisco, where he enjoys playing soccer, golfing and hiking with his golden retriever. He is an avid Cleveland sports fan. Please join us in welcoming him to Promega, and feel free to contact him at: 800-356-9526, extension 1870, or by e-mail at: yasser.daoudi@promega.com to discuss your laboratory's needs.