

# Newest Members of the Promega Genetic Identity Team



## LISA LANE, M.S.

Promega is pleased to introduce Lisa Lane, our newest Regional Technical Specialist. Lisa joined Promega in August of 2000. She supports customers in the Western United States in a variety of sales and technical programs. Most recently she has been working with customers to organize regional Educational Forums, which are programs that discuss issues involved in implementing new technologies into customer laboratories.

Lisa received a Bachelor of Science with an emphasis in microbiology and a minor in chemistry from the University of Oklahoma. She also has a Masters degree in human relations from the same institution.

Prior to joining Promega, Lisa directed operations for the Advanced Center for Genome Technology in Norman, Oklahoma. In this position she established quality control methodologies, scientific protocols and operational procedures. Additionally, she has written and established both scientific and managerial programs for state and private facilities.

Contact information: Tel: 405-364-6214. E-mail: [llane@promega.com](mailto:llane@promega.com).



## PAUL NEWMAN, PH.D.

Paul Newman joined Promega's team in September of 2000 as Genetic Identity's Customer Support and Training Manager. In his position, Paul will develop and manage Promega's Working Groups and Educational Forums. His experience with a variety of instrument platforms, new technologies, and reagent systems makes him ideally suited to help customers and field personnel with a wide variety of technical issues.

Originally from New York, Paul received his B.S. in biology at the State University of New York, Stony Brook. He worked as a research assistant at Cold Spring Harbor in New York and then completed his Ph.D. in molecular genetics at the University of Cincinnati College of Medicine. Afterwards, he was a Postdoctoral Fellow at Tufts University School of Medicine in Boston. More recently, Paul was the director of training for Visible Genetics in Atlanta. In this position he was responsible for setting up their training facility and training curriculum for HIV genotyping.

Contact information: Tel: 608-277-2646. E-mail: [pnewman@promega.com](mailto:pnewman@promega.com).



## STEPHEN LUPTON, B.S.

Stephen Lupton is the new Genetic Identity Key Account Manager in Europe. Stephen manages Genetic Identity product sales and technical support activities from his office in Lyon, France. His responsibilities include providing product information, customer training, sales management and participation as Promega's representative in the various European human identity symposia throughout the year.

Stephen received a fifth year degree in biology from the University of Lille. He has extensive experience in PCR, sequencing, DNA analysis and DNA purification. Stephen previously represented Promega as a Field Applications Specialist serving France's Life Science Research market. Stephen is fluent in English, French and German.

Contact information: Tel: +33-0-437225078. E-mail: [slupton@fr.promega.com](mailto:slupton@fr.promega.com).



## NICOLE ZIEMER, M.S.

Nicole Ziemer joined our team in September of 2000 as the Marketing Assistant and Customer Service Representative for Genetic Identity in Europe. Nicole studied Business Management at the University of Mannheim and European Business Management at the University of Applied Science in Worms. She received her Masters in International Business Management at the University of Lyon.

Nicole resides in Mannheim, Germany and is fluent in German, Spanish, French and English.

Contact information: Tel: 49-621-8501. E-mail: [nziemer@de.promega.com](mailto:nziemer@de.promega.com).