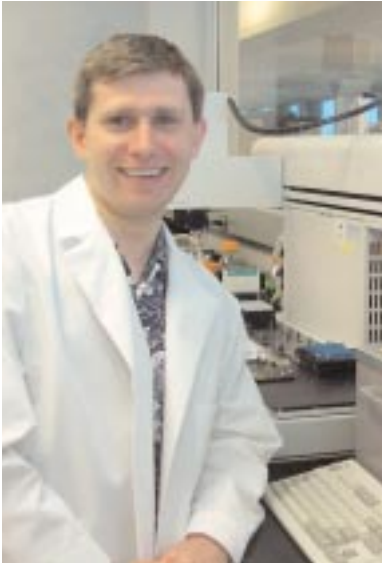


Newest Member of the Promega Genetic Identity Team



BOB MCLAREN, PH.D.

Please welcome the newest member of the Genetic Identity team: Bob McLaren.

Bob joined Promega in 1998 as a Technical Services Scientist. Over the years, he took on more responsibilities supporting the DNA-typing community, including the automation of the DNA IQ™ System on many instrument platforms and PowerPlex® System support. In September of 2005, Bob officially joined the Genetic Identity team as a Technical Services Scientist, Genetic Identity. Many of you have already met Bob during the setup of the DNA IQ™ System on a Biomek® 2000 or Biomek® 3000 workstation in your laboratory or during FSS-i3™ Expert Systems Software training.

In his new role, Bob will be responsible for training customers and providing support for Genetic Identity products and applications. He will also develop educational talks and represent Promega at meetings.

Bob received his B.S. degree in Microbiology from University College London, England, and his Ph.D. in Biochemistry from the University of Dundee in Scotland. He and his wife live in Oregon, Wisconsin, where he enjoys sailing, golfing and shoveling snow.

PROFILES IN DNA READERS SURVEY

We would like to thank all of you who took the time to complete the *Profiles in DNA* Readers Survey. We received a lot of valuable feedback about past issues and many ideas for new articles in future issues. This feedback will help us to ensure that *Profiles in DNA* articles continue to be of interest to our readers. We would like to congratulate the recipients of the *Profiles in DNA* Readers Survey prizes:

Grand Prize: Nancy Hill Jones, Alabama Department of Forensic Science

Second Prize: Trevor Seifert, University of Rochester, Center for Oral Biology

Third Prize: David Fisher, Office of Chief Medical Examiner, New York