

Regional Technical Specialist Joins Promega Genetic Identity Team

Promega's Regional Technical Specialists (RTS) serve as your principal source of information for *GenePrint*® products. Promega's RTS are available to:

- Answer technical questions about *GenePrint*® products.
- Assist in selection of Promega products.
- Provide up-to-date information on new products from Promega.
- Coordinate and present excellent seminars, workshops and training programs.
- Represent Promega at regional, national and international meetings.

ARNI MASIBAY, PH.D.

We are pleased to announce that Arni Masibay has joined the Genetic Identity team as a Regional Technical Specialist (RTS) covering the United States. Arni is responsible for all product sales to *GenePrint*® customers in the United States.

Arni grew up in Chicago, Illinois, and received his bachelor's degree in biology from Illinois Wesleyan University. He then moved to the Washington, D.C., area where he received his master's degree in Forensic Science from George Washington University and his doctorate in Microbiology from Georgetown University. He has since been a staff fellow at the National Institutes of Health and an adjunct professor at Broward Community College, Florida International University and Florida Atlantic University. Arni was a Forensic Scientist with the Broward Sheriff's Office and a member of the Forensic DNA Team in Broward County. He also served as a guest reviewer for the *Journal of Forensic Science* and a CODIS administrator. Arni joins the Promega *GenePrint*® team well qualified to meet the needs of our customers.

Arni lives with his wife, Becky, in Fort Lauderdale, Florida. In his free time, Arni is an avid tennis player and fan. He participates in many outdoor activities such as camping, hiking, diving and other water sports. Arni enjoys traveling, meeting new people and exchanging ideas.

TO CONTACT ARNI

Please contact him by Tel: 800-356-9526, extension 1707, or E-mail: amasibay@promega.com.



Arni Masibay, Ph.D.