

Upcoming Meetings

18th International Association of Forensic Sciences Triennial Meeting

New Orleans, Louisiana, USA
July 21–28, 2008
www.iafs2008.com

Canadian Society of Forensic Science Conference

Halifax, Nova Scotia, Canada
September 15–18, 2008
www2.csfs.ca

MAFS 37th Annual Meeting

Des Moines, Iowa, USA
September 28–October 3, 2008
www.mafs.net

19th International Symposium on the Forensic Sciences

Melbourne, Australia
October 6–9, 2008
www.anzfs2008.org.au

19th International Symposium on Human Identification

Hollywood, California, USA
October 13–16, 2008
www.promega.com/hollywood/

Dear Readers,

This issue of *Profiles in DNA* has articles covering an assortment of topics, so there is bound to be something of interest for everyone. In the first article, Chris Maguire and Michael Woodward from The Forensic Science Service present information about DNA-based kinship analysis and include examples of cases where kinship analysis contributed key information to an investigation. They also introduce the FSS-ibd software, which is designed to simplify these types of analyses.

Several articles in this issue deal with DNA extraction from a variety of substrates, including a summary by José Andradás Heranz, Emilio García Poveda and Francisco Alvarez Fernández at the Scientific Police Headquarters in Madrid of their experience isolating DNA from reference and casework samples using the Tecan Freedom EVO® liquid handler. Also, Anthony Tambasco and Rita Simons from the Mansfield Police Department share their protocol for successfully isolating DNA from drug paraphernalia and generating CODIS-eligible DNA profiles in the majority of cases. Finally we announce the availability of the Maxwell® 16 Applications Database, an online tool to obtain and share information about DNA isolation from various sample types using the Maxwell® 16 Instrument.

In this issue's interesting case report, Han Myunsoo from the DNA Identification Center in Seoul, Korea, describes his quest to identify the mother and father of two infants found in a freezer in Seo-rae Village.

Finally we wrap up the issue with a few announcements. Lyn Fatolitis and Jack Ballantyne describe a recent collaborative effort to consolidate several U.S.-based Y-STR databases into a single database that includes Y-STR haplotype data from 13,000 males. We also introduce you to three new Promega Genetic Identity team members.

Didn't see anything of interest to you? Please contact the editor at the address given on page 1 or by e-mail at: profilesindna@promega.com to contribute an article or offer suggestions for content of future issues of *Profiles in DNA*.



Terri Sundquist
Editor,
Profiles in DNA

⁽¹⁾U.S. Pat. Nos. 6,027,945, 6,368,800 and 6,673,631, Australian Pat. No. 732756, European Pat. No. 1 204 741 and Mexican Pat. No. 209436 have been issued to Promega Corporation for methods of isolating biological target materials using silica magnetic particles and simultaneous isolation and quantitation of DNA. Other patents are pending.

⁽²⁾U.S. Pat. Nos. 7,112,552, 7,304,015 and 7,304,016 have been issued to Promega Corporation for compositions for separating molecules. Other patents are pending.

⁽³⁾U.S. Pat. No. 7,319,021 has been issued to Promega Corporation for cell lysis composition, methods of use, apparatus and kit. Other patents are pending.

⁽⁴⁾The purchase of this product does not convey a license to use AmpliTaq Gold® DNA polymerase. You should purchase AmpliTaq Gold® DNA polymerase licensed for the forensic and human identity field directly from your authorized enzyme supplier.

⁽⁵⁾U.S. Pat. Nos. 6,238,863 and 6,767,703 and Korean Pat. No. 691195 have been issued to Promega Corporation for materials and methods for identifying and analyzing intermediate tandem repeat DNA markers. Other patents are pending.

⁽⁶⁾U.S. Pat. Nos. 5,843,660, 6,479,235, 6,221,598 and 7,008,771, Australian Pat. No. 724531, Canadian Pat. No. 2,118,048, Korean Pat. No. 290332, Singapore Pat. No. 57050 and Japanese Pat. No. 3602142 have been issued to Promega Corporation for multiplex amplification of STR loci. Other patents are pending.

⁽⁷⁾STR loci are the subject of U.S. Pat. No. RE 37,984, German Pat. No. DE 38 34 636 C2 and other patents issued to the Max-Planck-Gesellschaft zur Förderung der Wissenschaften, e.V., Germany. The development and use of STR loci are covered by U.S. Pat. No. 5,364,759, Australian Pat. No. 670231 and other pending patents assigned to Baylor College of Medicine, Houston, Texas.

Patents for the foundational PCR process, European Pat. Nos. 201,184 and 200,362, expired on March 28, 2006. In the U.S., the patents covering the foundational PCR process expired on March 29, 2005.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

Maxwell and PowerPlex are registered trademarks of Promega Corporation. Differex, DNA IQ and PowerType are trademarks of Promega Corporation.

ABI PRISM, GeneMapper, GeneScan, Genotyper, Identifier and Quantifier are registered trademarks of Applied Biosystems Corporation. Biomek is a registered trademark of Beckman Coulter, Inc. Freedom EVO is a registered trademark of Tecan AG Corporation. FSS-IP is a trademark of The Secretary of State for the Home Department. GeneAmp is a registered trademark of Roche Molecular Systems, Inc. Hi-Di is a trademark of Applied Biosystems Corporation. Microcon is a registered trademark of Millipore Corporation. UltraPure is a trademark of Invitrogen Corporation.