

Upcoming Meetings

**Forensic Science Society
48th Annual General Meeting**
Wyboston, United Kingdom
November 1, 2008
www.forensic-science-society.org.uk

**Northwest Association of
Forensic Scientists Meeting**
Boise, Idaho, USA
November 3–7, 2008
www.nwafs.org

**American Academy of Forensic
Science 61st Annual Scientific
Meeting**
Denver, Colorado, USA
February 16–21, 2009
www.aafs.org

**29. Spurenworkshop der
Deutschen Gesellschaft fuer
Rechtsmedizin**
Muenster, Germany
February 27–28, 2009
www.r-km.de/spurenworkshop2009/

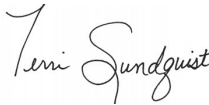
Dear Readers,

Successful DNA typing requires that each step in the process, including DNA isolation, quantitation and amplification, separation of amplified products and data analysis, be performed accurately, consistently and with sufficient throughput to avoid a backlog of samples. The articles in this issue describe processes that can improve workflow and reduce a scientist's hands-on time without sacrificing the quality of the final result.

As one of the first steps toward consistency and high-quality results, DNA-typing laboratories record and validate their processes and often undergo accreditation. One option is ISO accreditation. In the first article of this issue, Patricia Wojtowicz from Forensic Quality Services—International addresses misconceptions about the ISO/IEC 17025 accreditation process.

Three articles in this issue describe tools to reduce a scientist's hands-on time when processing samples, which can result in increased consistency. First Jindřich Novotný and Thorsten Hadrys describe the coupling of laser microdissection and pressure catapulting of single cells from a slide with on-chip PCR, allowing rapid amplification and interpretable DNA profiles from as few as one cell. Next Heike Felske-Zech, Kai Steffen Edler, Igor Martytschan, Christin Gruber and Frank Heidorn relate their experiences with automated DNA isolation using the DNA IQ™ Casework Sample Kit for Maxwell® 16. Finally, Sarah Shultz and Benjamin Krenke present some frequently asked questions about the new automated Plexor® HY System setup methods, which direct robot-mediated quantitative PCR assembly.

In the Tech Tips article, Robert McLaren, Marty Ensenberger, Cindy Sprecher, Dawn Rabbach, Patricia Fulmer, Joe Bessetti and Doug Storts describe a novel solution to eliminate double-stranded DNA artifacts that can appear at the vWA locus. Finally, we present a list of useful citations for laboratories that are validating or considering validating Promega genetic identity products.



Terri Sundquist
Editor,
Profiles in DNA

¹The 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.

²STR 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.

³Allele sequences for one or more of the loci vWA, FGA, D8S1179, D21S11 and D18S51 in allelic ladder mixtures is licensed under U.S. Pat. No. 7,087,380, Australia Pat. No. 2003200444 and corresponding patent claims outside the US.

⁴CAL Fluor® technology is the subject of pending patents and is licensed and sold under agreement with Biosearch Technologies, Inc., for research and development and forensic and paternity testing. These products are sold for use by the end-user only and may not be resold, distributed or repackaged.

⁵This product is sold under licensing arrangements with the USB Corporation for Forensic and Genetic Identity Applications Fields specifically excluding tissue typing related to transplantation or other medical procedures. Further licensing information may be obtained by contacting the USB Corporation, 26111 Miles Road, Cleveland, OH 44128.

⁶This product is sold under licensing arrangements with Stratagene for Forensic and Genetic Identity Applications Fields specifically excluding tissue typing related to transplantation or other medical procedures. Further licensing information may be obtained by contacting the Business Development Department, Stratagene California, 11011 North Torrey Pines Road, La Jolla, CA 92037.

⁷Patents 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.

⁸The purchase of this product conveys to the buyer the limited, nonexclusive, nontransferable right (without the right to resell, repackage, or further sublicense) under U.S. Published Patent Appln. 20020150900 and U.S. Pat. Nos. 5,432,272, 6,617,106 and 6,140,496 to use the product. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of this product does not include or carry any right or license to sell this product. For information on purchasing a license for other uses, please contact Promega Corporation, Business Development, 2800 Woods Hollow Road, Madison, WI 53711, or EraGen Biosciences, Corporate Licensing, 918 Deming Way, Suite 201, Madison, WI 53717. Phone (608) 662-9000; Fax (608) 662-9003.

⁹Use of this product for basic PCR is outside of any valid US or European patents assigned to Hoffman La-Roche or Applera. This product can be used for basic PCR in research, commercial or diagnostic applications without any license or royalty fees.

¹⁰Use of the SE33 locus is licensed under U.S. Pat. Nos. 5,468,610 and 5,721,100.

¹¹PowerPlex® 16 BIO System incorporates dye conjugates made with the Rhodamine Red™-X and Texas Red®-X fluorescent reactive dyes, which are licensed from Molecular Probes, Inc., under U.S. Pat. Nos. 5,798,276 and 5,846,737 for DNA analysis. Rhodamine Red is a trademark of Molecular Probes, Inc., and Texas Red is a registered trademark of Molecular Probes, Inc.

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

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

ABI PRISM, AmpFISTR, Cofiler and SGM Plus are registered trademarks of Applera Corporation. AmpliGrid is a registered trademark of Olympus Corporation. AmpliTaq Gold and GeneAmp are registered trademarks of Roche Molecular Systems, Inc. Biomek is a registered trademark of Beckman Coulter, Inc. CAL Fluor is a registered trademark of Biosearch Technologies, Inc. Chelex is a registered trademark of Bio-Rad Laboratories, Inc. Freedom EVO is a registered trademark of Tecan AG Corporation. FSS-i³ is a trademark of The Secretary of State for the Home Department. HemaColor is a registered trademark of Merck. LMPC is a trademark of P.A.L.M. MicroLaser Technologies. Mastercycler is a registered trademark of Eppendorf AG. PALM is a registered trademark of P.A.L.M. MicroLaser Technologies. POP-4 and Profiler Plus are trademarks of Applera Corporation.