

Upcoming Meetings

2009 Mid-Atlantic Association of Forensic Scientists Annual Meeting

Hunt Valley, Maryland, USA
May 4–8, 2009
www.maafs.org

5th European Academy of Forensic Science Conference

Glasgow, Scotland
September 8–11, 2009
www.eafs2009.com

23rd World Congress International Society for Forensic Genetics

Buenos Aires, Argentina
September 23–26, 2009
www.isfg2009.org

20th International Symposium on Human Identification

Summerlin, Nevada, USA
October 12–15, 2009
www.promega.com/geneticsymp20/

IVth Mediterranean Academy of Forensic Sciences Meeting

Antalya, Turkey
October 14–18, 2009
www.mafs2009.org

Dear Readers,

It has been said that we are moving from the industrial age to the information age. To keep pace with our workload and life in general, it is increasingly important to find specific information quickly and easily. Computers and other digital devices are ideally suited to this purpose. Rather than sorting through stacks of paper and visually scanning page after page to find specific information, you can perform a simple search and get the desired result much faster. In addition, unlike a static printed page, electronic media allow us the opportunity to customize the interface and interact with scientists around the globe. In this way, electronic media have changed the way that we gather information forever.

In light of this, *Profiles in DNA* is evolving: shedding its paper constraints and going completely digital. The web-based version of *Profiles in DNA* will include all of your favorite articles, as well as informative new articles, useful resources, opportunities for interaction and even some fun activities. Don't miss your opportunity to stay connected to this valuable resource. Visit the Promega Web site at: www.promega.com/stayconnected/ and update your profile.

Keeping with the theme, we are ushering in a new era of information-gathering methods by presenting you with information about a new era of Promega products, including the new PowerPlex® European Systems, which amplify 15 ENFSI-recommended STR loci plus Amelogenin with or without SE33, and the new PowerPlex® 16 HS System, which amplifies the same 15 STR loci found in the PowerPlex® 16 System and includes hot-start *Taq* DNA polymerase in a convenient master mix. First, however, we start this issue with an article about the discovery of a 100- to 150-year-old mummy in Chile and how a multidisciplinary group of scientists performed anthropological and genetic analyses to learn more about the origin of this individual and his modern descendants.



Terri Sundquist
Editor,
Profiles in DNA

^(a)U.S. Pat. Nos. 6,027,945, 6,368,800 and 6,673,631, Australian Pat. No. 732756 and other patents and patents pending.

^(b)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.

^(c)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.

^(d)U.S. Pat. Nos. 6,238,863 and 6,767,703 and other patents and patents pending.

^(e)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 and other patents and patents pending.

^(f)Allele 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.

^(g)Licensed under U.S. Pat. Nos. 5,338,671 and 5,587,287 and corresponding patents in other countries.

^(h)CAL 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.

⁽ⁱ⁾This 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.

^(j)This 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.

^(k)U.S. Pat. No. 6,242,235, Australian Pat. No. 761757, Canadian Pat. No. 2,335,153 and other patents and patents pending.

^(l)Use 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.

^(m)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.

⁽ⁿ⁾The 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.

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

Maxwell, Plexor, PowerPlex and Wizard are registered trademarks of Promega Corporation. DNA IQ is a trademark of Promega Corporation.

ABI PRISM, AmpFISTR, Identifier and Quantifiler are registered trademarks of Applied Biosystems Corporation. AmpliTaq Gold is a registered trademark of Roche Molecular Systems, Inc. CAL Fluor is a registered trademark of Biosearch Technologies, Inc. MegaBACE is a registered trademark of GE Healthcare Bio-Sciences. MiniFiler is a trademark of Applied Biosystems Corporation.