

## Useful Resources for Your Laboratory

Keeping up with the scientific literature can be difficult, considering everything else that requires your time and attention in the laboratory. To make things a little easier, we have compiled a list of useful resources for laboratories that are validating or considering validation of Promega products. This list is not meant to include every useful reference but should provide you with a good starting point in your search for information.

### POWERPLEX® 16 AND 16 BIO<sup>(a-c,k)</sup> SYSTEMS

McLaren, R.S. *et al.* (2008) Post-injection hybridization of complementary DNA strands on capillary electrophoresis platforms: A novel solution for dsDNA artifacts. *Forensic Sci. Int. Genet.* **2**, 257–73.

Krenke, B.E. *et al.* (2002) Validation of a 16-locus fluorescent multiplex system. *J. Forensic Sci.* **47**, 773–85.

Greenspoon, S.A. *et al.* (2004) Validation and implementation of the PowerPlex 16 BIO System STR multiplex for forensic casework. *J. Forensic Sci.* **49**, 71–80.

Levedakou, E. *et al.* (2001) Allele frequencies for fourteen STR loci of the PowerPlex™ 1.1 and 2.1 Multiplex Systems and Penta D locus in Caucasians, African-Americans, Hispanics, and other populations of the United States of America and Brazil. *J. Forensic Sci.* **46**, 736–61.

Budowle, B. *et al.* (2001) STR primer concordance study. *Forensic Sci. Int.* **124**, 47–54.

Budowle, B. and Sprecher, C. (2001) Concordance study on population database samples using the PowerPlex™ 16 kit and AmpF/STR® Profiler Plus™ kit and AmpF/STR® COfiler™ kit. *J. Forensic Sci.* **46**, 637–41.

Masibay, A., Mozer, T.J. and Sprecher, C. (2000) Promega Corporation reveals primer sequences in its testing kits. *J. Forensic Sci.* **45**, 1360–2.

Sprecher, C. *et al.* (2000) The PowerPlex™ 16 System. *Profiles in DNA* **4(1)**, 3–6.

### POWERPLEX® Y SYSTEM<sup>(a,b)</sup>

Krenke, B.E. *et al.* (2005) Validation of a male-specific, 12-locus fluorescent short tandem repeat (STR) multiplex. *Forensic Sci. Int.* **148**, 1–14.

Budowle, B. *et al.* (2005) Twelve short tandem repeat loci Y chromosome haplotypes: Genetic analysis on populations residing in North America. *Forensic Sci. Int.* **150**, 1–15.

Krenke, B. *et al.* (2003) The PowerPlex® Y System. *Profiles in DNA* **6(2)**, 7–10.

Gill, P. *et al.* (2001) DNA Commission of the International Society of Forensic Genetics: Recommendations on forensic analysis using Y-chromosome STRs. *Int. J. Legal Med.* **114**, 305–9.

Roewer, L. *et al.* (2001) Online reference database of European Y-chromosomal short tandem repeat (STR) haplotypes. *Forensic Sci. Int.* **118**, 106–13.

### PLEXOR® HY SYSTEM

Krenke, B.E. *et al.* (2008) Developmental validation of a real-time PCR assay for the simultaneous quantification of total human and male DNA. *Forensic Sci. Int. Genet.*, in press.

Validation Guide for the Plexor® HY System, #GE295, Promega Corporation.

### MAXWELL® 16 SYSTEM

Validation Guide for the DNA IQ™ Reference Sample Kit for Maxwell® 16, #GE181, Promega Corporation.

Krnajski, Z., Geering, S. and Steadman, S. (2007) Performance verification of the Maxwell 16 instrument and DNA IQ reference sample kit for automated DNA extraction of known reference samples. *Forensic Sci. Med. Pathol.* **3**, 264–9.

Promega also offers the Citations Database, a searchable database of peer-reviewed journal articles that describe the use of Promega Genetic Identity products. The Citations Database can be viewed at:

**[www.promega.com/citations/](http://www.promega.com/citations/)**. Simply select “DNA Typing” as the application or enter appropriate keywords to generate a reference list that includes a short description of each article. If you have published an article describing work done with a Promega Genetic Identity product and would like the article included in the database, contact us at: **[profilesindna@promega.com](mailto:profilesindna@promega.com)**