



PowerPlex[®] 18D System

More Profile. Less Process.

The Power to Solve...*from Sample to Analysis*

Dramatically improve your laboratory efficiency. PowerPlex® 18D streamlines database and paternity testing by reducing manipulations and total processing time.

- **Eliminate Extraction**—Simplify and shorten sample processing with direct amplification from FTA® cards. Simplified protocols for other sample types are available.*
- **Reduce PCR Time**—Amplify in less than 1.5 hours using rapid cycling technology.
- **Upload More Markers**—Amplify D2S1388, D19S433, Penta D, Penta E, Amelogenin and the 13 CODIS loci for more discrimination and more loci in common with neighboring databases.

More Profile.

The PowerPlex® 18D System amplifies 18 of the most common human identification markers (Figure 1). With D2S1338, D19S433, Penta D, Penta E plus the 13 CODIS STR loci and Amelogenin, more discrimination and enhanced database compatibility is possible. PowerPlex® 18D uploaded profiles include all of the loci found in the commonly used PowerPlex® 16 HS System and Identifiler® Kit. Additionally, D2S1338 and D19S433 increase discrimination over the PowerPlex® 16 HS System for difficult paternity testing.

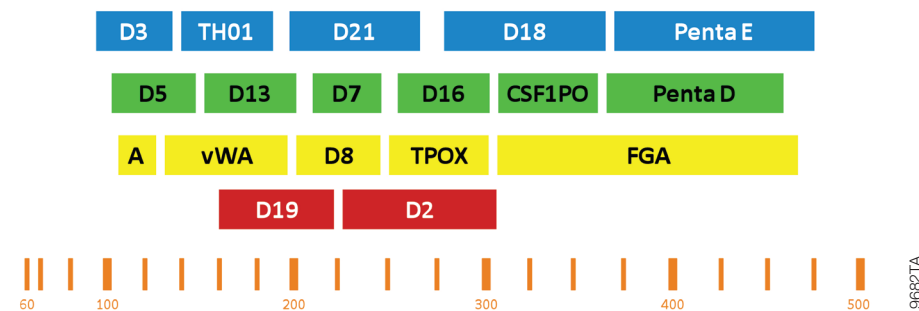


Figure 1. Configuration of the PowerPlex® 18D System. The PowerPlex® 18D System contains all of the 13 CODIS loci: D3S1358, TH01, D21S11, D18S51, D5S818, D13S317, D7S820, D16S539, CSF1PO, vWA, D8S1179, TPOX, and FGA, plus Amelogenin, Penta E, Penta D, D19S433 and D2S1338.

Less Process.

The PowerPlex® 18D System saves labor, time and cost with direct amplification and rapid PCR technology by removing extraction or washing for FTA® cards (or simplified procedures for swabs and other storage cards*) and shortening PCR to <1.5 hours. Less processing means fewer manipulations and savings on purification or wash reagents. Rapid PCR technology significantly shortens a laboratory’s daily workflow and allows a whole plate of samples to move from “punch” to capillary electrophoresis (CE) analysis *in the same day*.

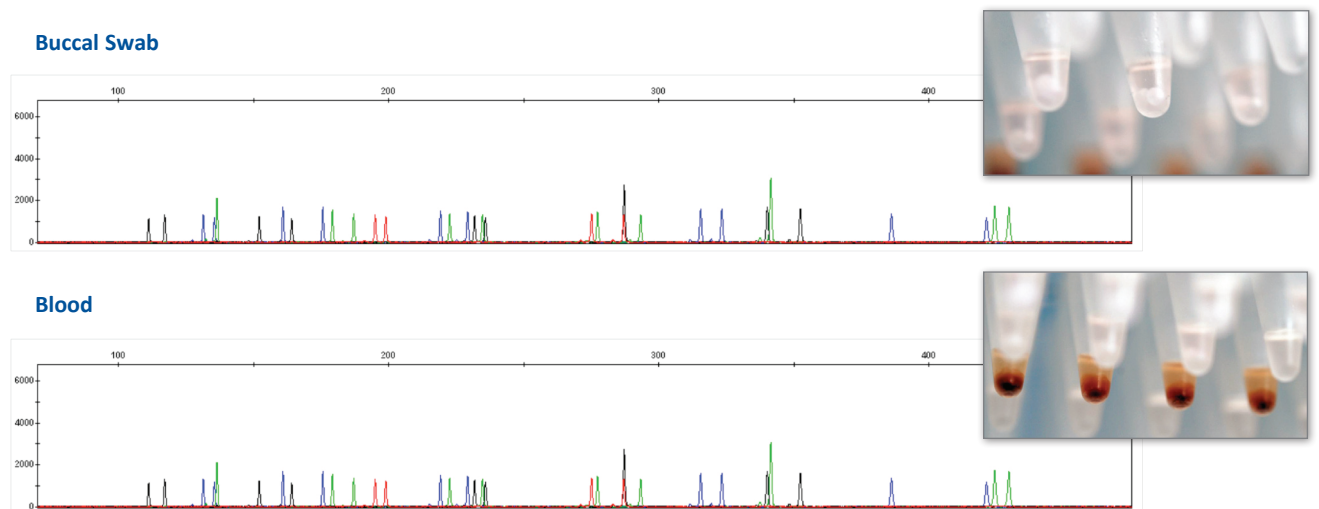


Figure 2. Amplification of samples on FTA® punches. Amplification of two 1.2mm FTA® punches with a buccal sample (top panel) and one 1.2mm FTA® punch containing blood (bottom panel) was performed using 27 cycles on a GeneAmp® PCR System 9700 and analyzed with the Applied Biosystems 3130xl Genetic Analyzer.

For more information visit: www.promega.com/powerplex18D/

“...move from ‘punch’ to CE analysis in the same day”


* Samples collected using *Omni Swabs* and *Bode Buccal DNA Collector™* devices can be quickly processed with simple pretreatment procedures. Contact Promega Technical Services at: genetic@promega.com for details. Analysis of extracted samples possible.

Ordering Information

PRODUCT	SIZE	CAT.#
PowerPlex® 18D System	200 reactions	DC1802
	800 reactions	DC1808
PowerPlex® 5-Dye Matrix Standards, 3100/3130		DG4700
CC5 Internal Lane Standard 500		DG1521

The Power to Solve



 **Identity Automation™:** Fully developed and tested high-throughput solutions with Promega installation, training and support

 Automation  Chemistry  Software

*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, DNA IQ, Identity Automation, Slicprep and STR Normalization Manager are trademarks of Promega Corporation. Bode Buccal DNA Collector is a trademark of the Bode Technology Group, Inc. FSS-i³ is a trademark of The Secretary of State for the Home Department and is only available in select countries. FTA is a registered trademark of Flinders Technologies, Pty, Ltd., and is licensed to Whatman.