



**Promega**

## Technical Bulletin

---

# PowerPlex® 16 and PowerPlex® ES Monoplex Systems

INSTRUCTIONS FOR USE OF PRODUCTS DC6551, DC6561, DC6571,  
DC6581, DC6591, DC6601, DC6611, DC6621, DC6631, DC6641, DC6651,  
DC6661, DC6671, DC6681, DC6691, AND DC6751.



[www.promega.com](http://www.promega.com)

PRINTED IN USA.  
Revised 1/07



AF9TBD017 0107TBD017

Part# TBD017

# PowerPlex® 16 and PowerPlex® ES Monoplex Systems

All technical literature is available on the Internet at: [www.promega.com/tbs/](http://www.promega.com/tbs/)  
Please visit the web site to verify that you are using the most current version of this  
Technical Bulletin. Please contact Promega Technical Services if you have questions on use  
of this system. E-mail: [techserv@promega.com](mailto:techserv@promega.com)

I. Description.....	1
II. Product Components and Storage Conditions .....	2
III. Amplification Protocols.....	3
IV. Detection Methods .....	4
A. Detection Using the ABI PRISM® 310, 3100 or 3100- <i>Avant</i> Genetic Analyzer and the Applied Biosystems 3130 or 3130 <i>xl</i> Genetic Analyzer ...	4
B. Detection Using the Hitachi FMBIO® II Fluorescence Imaging System. ....	5

## I. Description

The PowerPlex® 16 and PowerPlex® ES Monoplex Systems contain primer pairs that have the same sequence as those used in the PowerPlex® 16 (Cat.# DC6531), PowerPlex® 2.1 (Cat.# DC6471), PowerPlex® 16 BIO (Cat.# DC6541) and PowerPlex® ES Systems (Cat.# DC6731). Refer to the *PowerPlex® 16 System Technical Manual #TMD012*, for locus-specific information, allele size range, allelic ladder components and genotypes of commonly used standard DNA templates. Refer to the *PowerPlex® ES System Technical Manual #TMD017* for SE33 locus-specific information. Technical Manuals are available upon request or at: [www.promega.com/tbs/](http://www.promega.com/tbs/)

## II. Product Components and Storage Conditions

Product	Size	Cat.#
PowerPlex® 16 Monoplex System, Penta E (Fluorescein) <sup>(a,b,c)</sup>	100 reactions	DC6591
PowerPlex® 16 Monoplex System, Penta D (JOE) <sup>(a,b,c)</sup>	100 reactions	DC6651
PowerPlex® ES Monoplex System, SE33 (JOE) <sup>(a,b,d)</sup>	100 reactions	DC6751

Not For Medical Diagnostic Use. Each system includes:

- 150µl Internal Lane Standard
- 300µl Gold ST★R Buffer
- 250µl 10X Primer Pair Mix
- 70µl Allelic Ladder
- 2 Protocols

Product*	Size	Cat.#
PowerPlex® 16 Monoplex System, D3S1358 (Fluorescein) <sup>(a,b)</sup>	100 reactions	DC6551
PowerPlex® 16 Monoplex System, TH01 (Fluorescein) <sup>(a,b)</sup>	100 reactions	DC6561
PowerPlex® 16 Monoplex System, D21S11 (Fluorescein) <sup>(a,b)</sup>	100 reactions	DC6571
PowerPlex® 16 Monoplex System, D18S51 (Fluorescein) <sup>(a,b)</sup>	100 reactions	DC6581
PowerPlex® 16 Monoplex System, D5S818 (JOE) <sup>(a,b)</sup>	100 reactions	DC6601
PowerPlex® 16 Monoplex System, D13S317 (JOE) <sup>(a,b)</sup>	100 reactions	DC6611
PowerPlex® 16 Monoplex System, D7S820 (JOE) <sup>(a,b)</sup>	100 reactions	DC6621
PowerPlex® 16 Monoplex System, D16S539 (JOE) <sup>(a,b)</sup>	100 reactions	DC6631
PowerPlex® 16 Monoplex System, CSF1PO (JOE) <sup>(a,b)</sup>	100 reactions	DC6641
PowerPlex® 16 Monoplex System, vWA (TMR) <sup>(a,b)</sup>	100 reactions	DC6661
PowerPlex® 16 Monoplex System, D8S1179 (TMR) <sup>(a,b)</sup>	100 reactions	DC6671
PowerPlex® 16 Monoplex System, TPOX (TMR) <sup>(a,b)</sup>	100 reactions	DC6681
PowerPlex® 16 Monoplex System, FGA (TMR) <sup>(a,b)</sup>	100 reactions	DC6691

Not For Medical Diagnostic Use. Each system includes:

- 150µl Internal Lane Standard
- 300µl Gold ST★R Buffer
- 250µl 10X Primer Pair Mix
- 2 Protocols

**Storage Conditions:** Store all components at -20°C. The fluorescent 10X Primer Pair Mix is light-sensitive; therefore minimize light exposure and store in the dark.

\*Items listed are available from Promega as “special order” items. A minimum order is not required, but items may not be immediately available for shipment. Contact Promega Customer Service for more information.

### III. Amplification Protocols

Follow the protocols in the *PowerPlex® 16 System Technical Manual #TMD012* for the PowerPlex® 16 Monoplex Systems and in the *PowerPlex® ES System Technical Manual #TMD017* for the PowerPlex® ES Monoplex System. However, follow the directions below for the amount of AmpliTaq Gold® DNA polymerase to use per 25µl reaction.

1. Prepare the amplification mix as directed in Tables 1 and 2.
2. Add template DNA as directed in Tables 1 and 2. For optimal amplification results, we recommend 0.5–2ng of DNA per reaction.

**Table 1. Amplification Mix for a Single Sample.**

PCR Component	Volume Per Sample
Nuclease-Free Water	17.4µl
Gold ST★R 10X Buffer	2.5µl
10X Primer Pair	2.5µl
AmpliTaq Gold® DNA polymerase <sup>1</sup>	0.1µl (0.5 units)
<b>Total amplification mix volume</b>	<b>22.5µl</b>
Template DNA <sup>2</sup> (to be added)	2.5µl
<b>Total Reaction Volume</b>	<b>25.0µl</b>

<sup>1</sup>Assumes the AmpliTaq Gold® DNA polymerase is 5u/µl. If the enzyme concentration is different, adjust the volume of enzyme used accordingly.

<sup>2</sup>Assumes the template DNA volume is 2.5µl.

**Table 2. Preparation of Amplification Mix for Multiple Samples.**

PCR Component	Volume Per Sample	× Number of Reactions	= Final Volume (µl)
Nuclease-Free Water	µl		
Gold ST★R 10X Buffer	2.5µl		
10X Primer Pair	2.5µl		
AmpliTaq Gold® DNA polymerase <sup>1</sup>	0.1µl (0.5 units)		
<b>Total amplification mix volume<sup>2</sup></b>	<b>µl</b>		
Template DNA (to be added)	µl		
<b>Total Reaction Volume</b>	<b>25.0µl</b>		

<sup>1</sup>Assumes the AmpliTaq Gold® DNA polymerase is 5u/µl. If the enzyme concentration is different, adjust the volume of enzyme used accordingly.

<sup>2</sup>The volume of template DNA plus the volume of amplification mix should equal 25.0µl.

#### IV. Detection Methods

**Note:** Allelic ladders are only included with the PowerPlex® 16 Monoplex System, Penta E (Fluorescein), PowerPlex® 16 Monoplex System, Penta D (JOE) and PowerPlex® ES Monoplex System, SE33 (JOE). Allelic ladders for the other systems are available by custom order. Allelic ladder options are included in Tables 3 and 4.

##### IV.A. Detection Using the ABI PRISM® 310, 3100 or 3100-Avant Genetic Analyzer and the Applied Biosystems 3130 or 3130xl Genetic Analyzer

Specific directions for detection using the ABI PRISM® 310, 3100 or 3100-Avant Genetic Analyzers and the Applied Biosystems 3130 or 3130xl are given in the *PowerPlex® 16 System Technical Manual #TMD012* and the *PowerPlex® ES System Technical Manual #TMD017*. Please follow these protocols when using the PowerPlex® Monoplex Systems. For additional questions, contact Promega Technical Services.

**Table 3. Allelic Ladder Options for the ABI PRISM® 310, 3100 and 3100-Avant Genetic Analyzers and the Applied Biosystems 3130 or 3130xl Genetic Analyzer.**

<b>PowerPlex® 16/ES Monoplex Loci</b>	<b>Allelic Ladder Options<sup>1,2</sup></b>
D18S51, D21S11, TH01, D3S1358,	
FGA, D8S1179, vWA	PowerPlex® 16 or PowerPlex® ES
Penta E	Penta E or PowerPlex® 16
Penta D	Penta D or PowerPlex® 16
CSF1PO, D16S539, D7S820,	
D13S317, D5S818, TPOX	PowerPlex® 16
SE33	SE33 or PowerPlex® ES

<sup>1</sup>The PowerTyper™ Macros (Cat.# DG3470) can be used for data analysis.

<sup>2</sup>All allelic ladders are available by custom order.

#### IV.B. Detection Using the Hitachi FMBIO® II Fluorescence Imaging System

For detection using the Hitachi FMBIO® II Fluorescence Imaging System, see the *PowerPlex® 2.1 System Technical Manual #TMD011* or *PowerPlex® 16 BIO System Technical Manual #TMD016*. The filter set used for detection of the PowerPlex® 16 and PowerPlex® ES Monoplex loci is shown in Table 4.

**Table 4. Allelic Ladder Options for the Hitachi FMBIO® II Fluorescence Imaging System.** The filter set used for detection of each PowerPlex® 16 or PowerPlex® ES Monoplex locus is shown in bold in the “Filters Needed for Detection” Column.

<b>PowerPlex® 16/ES Monoplex Locus</b>	<b>Allelic Ladder Options<sup>1</sup></b>	<b>Filters Needed for Detection (nm)</b>	<b>PowerPlex® Filter Set</b>
D18S51, D21S11, TH01, D3S1358	PowerPlex® 2.1	<b>505, 585, 650</b>	1.1/2.1
	PowerPlex® 16 BIO <sup>2</sup>	<b>505, 577, 598, 665</b>	16 BIO
CSF1PO, D16S539, D7S820, D13S317, D5S818	PowerPlex® 16 BIO <sup>2</sup>	<b>505, 577, 598, 665</b>	16 BIO
	JOE Ladder Mix	<b>585, 650</b> <b>577, 665</b>	1.1/2.1 16 BIO
FGA, TPOX, D8S1179, vWA	PowerPlex® 2.1	<b>505, 585, 650</b>	1.1/2.1
		<b>505, 577, 665</b>	16 BIO <sup>3</sup>
SE33	SE33 Allelic Ladder	<b>585, 650</b> <b>577, 665</b>	1.1/2.1 16 BIO
Penta E	PowerPlex® 2.1	<b>505, 585, 650</b>	1.1/2.1
	PowerPlex® BIO	<b>505, 577, 598, 665</b>	16 BIO
	Penta E Allelic Ladder	<b>505, 650</b> <b>505, 665</b>	1.1/2.1 16 BIO
Penta D	PowerPlex® BIO	<b>505, 577, 598, 665</b>	16 BIO
	JOE Ladder Mix	<b>585, 650</b>	1.1/2.1
		<b>577, 665</b>	16 BIO
	Penta D Allelic Ladder	<b>585, 650</b> <b>577, 665</b>	1.1/2.1 16 BIO

<sup>1</sup>All allelic ladders are available by custom order.

<sup>2</sup>The Matrix 16 BIO is required for use with the PowerPlex® 16 BIO Allelic Ladder Mix. The Internal Lane Standard 600, labeled with CXR, can be detected with the 665nm filter.

<sup>3</sup>When using the PowerPlex® 16 BIO filter set for TMR-labeled loci, do not include the data from the 598nm filter in the FMBIO® Analysis Project or do not scan using the 598nm filter. Both TMR and CXR will be detected with the 598nm filter making color separation impossible. The TMR-labeled loci will be detected in the 577nm scan (JOE channel).



**Promega**

---

**Notes**

**Notes**

<sup>(a)</sup>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.

<sup>(b)</sup>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.

<sup>(c)</sup>U.S. Pat. Nos. 6,238,863 and 6,767,703 have been issued to Promega Corporation for materials and methods for identifying and analyzing intermediate tandem repeat DNA markers. Other patents are pending.

<sup>(d)</sup>Use of the SE33 locus is licensed under U.S. Pat. Nos. 5,468,610 and 5,721,100.

© 2002-2007 Promega Corporation. All Rights Reserved.

PowerPlex is a registered trademark of Promega Corporation. PowerTyper is a trademark of Promega Corporation.

ABI PRISM is a registered trademark of Applied Biosystems Corporation. AmpliTaq Gold is a registered trademark of Roche Molecular Systems, Inc. FMBIO is a registered trademark of Hitachi Software Engineering Company, Ltd.

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

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.