# PfuUltra II Hotstart PCR Master Mix

Catalog #600850 and 600852

## For Research Use Only. Not for use in diagnostic procedures.

## **Technical Services**

#### Telephone:

To find worldwide support telephone numbers visit <u>www.agilent.com/genomics/contactus</u>

Email: techservices@agilent.com

## MATERIALS PROVIDED

	Quantity	
Materials Provided	Catalog #600850	Catalog #600852
PfuUltra II Hotstart 2 $ imes$ Master Mix	100 reactions	400 reactions

**Storage:** Store at  $-20^{\circ}$ C upon receipt. After thawing, store at  $4^{\circ}$ C. Once thawed, full activity is guaranteed for 6 months.

## INTRODUCTION

*PfuUltra* II Hotstart PCR Master Mix\* is a 2× formulation of *PfuUltra* II Fusion HS DNA polymerase,\* an optimized PCR reaction buffer, magnesium, and dNTPs. *PfuUltra* II Fusion HS DNA polymerase couples fusion polymerase technology with our engineered *PfuUltra* DNA polymerase,\* and proprietary ArchaeMaxx PCR enhancing factor to achieve extreme accuracy, high specificity, and long target-length capability while dramatically reducing overall PCR extension times.

## **OPTIMIZATION PARAMETERS (50 µL REACTION VOLUME)**

Parameter	Targets: ≤10 kb (vector or genomic DNA)	Targets: >10 kb (vector or genomic DNA)	cDNA Targets
Extension time	15 seconds for targets ≤1 kb; 15 seconds per kb for targets >1 kb	30 seconds per kb	30 seconds for targets ≤1 kb; 30 seconds per kb for targets >1 kb
Input template	100 ng genomic DNA; 5–30 ng vector DNA	200–250 ng genomic DNA; 5–30 ng vector DNA	1–2 μl cDNA from RT-PCR reaction (50–500 ng starting total RNA template)
Primer concentration	0.2–0.4 µM each primer	0.4 μM each primer	0.2 μM each primer
Denaturing temperature 95°C/20 seconds per cycle and time		92°C/20–40 seconds per cycle	95°C/20 seconds per cycle
Extension temperature	72°C	68°C	72°C

## PCR PROTOCOL

The reaction conditions given here are for amplification of a typical single-copy chromosomal target of  $\leq 10$  kb. See the *Optimization Parameters* section for guidelines on amplifying longer targets, vector DNA targets, or cDNA targets. The reaction conditions are for one reaction and must be adjusted for multiple samples. The final volume of each reaction is 50 µl. Add the components in order into sterile thinwalled PCR tubes while mixing gently.

Important Before dispensing the PfuUltra II Hotstart 2× Master Mix, invert the container several times, then vortex briefly to ensure homogeneity.

Component	Amount per reaction
PfuUltra II Hotstart 2× Master Mix <sup>a</sup>	25.0 μl
Primer #1 (10 μM) <sup>b</sup>	1.0 μl
Primer #2 (10 μM) <sup>b</sup>	1.0 μl
DNA template (100 ng/µl) <sup>c</sup>	1.0 μl
Distilled water (dH <sub>2</sub> O)	22.0 µl
Total reaction volume	50 µl

## Reaction Mixture for a Typical Single-Copy Chromosomal Locus PCR Amplification (≤10 kb)

 $^{\alpha}~$  The final concentration of Mg^{2+} in the 1  $\times$  reaction mixture is 2 mM.

<sup>b</sup> Generally final primer concentrations of 0.2 μM (each primer) are optimal. For longer targets (7–10 kb), increasing the concentration to 0.4 μM (each primer) may substantially increase product yield.

<sup>c</sup> The amount of DNA template required varies depending on the type of DNA being amplified. Generally 100 ng of genomic DNA template is recommended. Less DNA template should be used for amplification of lambda or plasmid (5–30 ng) PCR targets.

Perform PCR using optimized cycling conditions. Suggested cycling parameters for PCR are indicated below. The PCR cycling parameters have been tested on the following: the MJ Research® DNA Engine® PTC-200, the Applied Biosystems® GeneAmp® PCR systems 9600 and 9700, the Bio-Rad iCycler, and the Agilent Mx3000P QPCR system. Optimized cycling parameters are not necessarily transferable between thermal cyclers designed by different manufacturers.

Analyze the PCR amplification products on a 0.7–1.0% (w/v) agarose gel.

\* U.S. Patent Nos. 6,734,293, 6,444,428, 6,183,997, and 5,948,663.

## PCR Cycling Parameters for PfuUltra II Hotstart PCR Master Mix<sup>a</sup>

	Number of		Duration	
Segment	cycles	Temperature	(vector or genomic DNA)	Duration (cDNA)
1	1	95°C <sup>♭</sup>	2 minutes	1 minute
2	30 cycles for	95°C	20 seconds	20 seconds
	vector or genomic	Primer $T_m - 5^{\circ}C^{\circ}$	20 seconds	20 seconds
	DNA; 40 cycles for cDNA	72°C	15 seconds for targets ≤1 kb	30 seconds for targets ≤1 kb
			15 seconds per kb for targets >1 kb	30 seconds per kb for targets >1 kb
3	1	72°C	3 minutes	3 minutes

#### Targets ≤10 kb (vector DNA, genomic DNA, and cDNA)

## Targets >10 kb (vector or genomic DNA)

Segment	Number of cycles	Temperature	Duration
1	1	92°C	2 minutes
2	30	92°C	20 seconds <sup>d</sup>
		Primer $T_{\rm m} - 5^{\circ}{\rm C}^{\circ}$	20 seconds
		68°C	30 seconds per kb
3	1	68°C	5 minutes

<sup>a</sup> Thin-wall PCR tubes are highly recommended. These PCR tubes are optimized to ensure more efficient heat transfer and to maximize thermalcycling performance.

<sup>b</sup> Denaturing temperatures above 95°C are recommended only for GC-rich templates.

<sup>c</sup> The annealing temperature may require optimization. Typically annealing temperatures will range between 55° and 72°C.

<sup>d</sup> For long targets (10–19 kb), increasing the duration of the denaturing step to 40 seconds per cycle may increase PCR product yield.

## LIMITED PRODUCT WARRANTY

This warranty limits our liability to replacement of this product. No other warranties of any kind, express or implied, including without limitation, implied warranties of merchantability or fitness for a particular purpose, are provided by Agilent. Agilent shall have no liability for any direct, indirect, consequential, or incidental damages arising out of the use, the results of use, or the inability to use this product.

#### LIMITED LABEL LICENSE FOR PFU-CONTAINING DNA POLYMERASE PRODUCTS

This product is covered by the claims of one or more of the following U.S. Patents: 5,948,663; 6,183,997; 6,444,428. Purchase of this product conveys to the purchaser only the non-transferable right under these patents to use the product for research use only by the purchaser. No rights are granted to the purchaser hereunder to sell, modify for resale or otherwise transfer this product, either alone or as a component of another product, to any third party. Agilent reserves all other rights, and this product may not be used in any manner other than as provided herein. For information on obtaining a license to use this product for purposes other than research, please contact Agilent Technologies, Business Development, 11011 North Torrey Pines Road, La Jolla, California 92037. Phone (858) 535-5400.

Use of this product is licensed under U.S. Patent Nos. 5,587,287 and corresponding foreign patents Europe, Hong Kong, and Japan

This product is provided under an agreement between Bio-Rad Laboratories and Agilent Technologies, Inc., and the manufacture, use, sale or import of this product is subject to US. Pat. No. 6,627,424 and EP Pat. No. 1 283 875 B1, owned by Bio-Rad Laboratories, Inc. Purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in PCR (but not real-time PCR) in the Research Field including all Applied Research Fields (including but not limited to forensics, animal testing, and food testing).

## **ENDNOTES**

Applied Biosystems<sup>®</sup> is a registered trademark of Applera Corporation or its subsidiaries in the US and certain other countries. DNA Engine<sup>®</sup> and MJ Research<sup>®</sup> are registered trademarks of Bio-Rad Laboratories, Inc. GeneAmp<sup>®</sup> is a registered trademark of Roche Molecular Systems, Inc.

## **MSDS INFORMATION**

Material Safety Data Sheets (MSDSs) are provided online at *http://www.genomics.agilent.com*. MSDS documents are not included with product shipments.