

PfuUltra II Fusion HS DNA Polymerase

Catalog #600670, 600672, and 600674

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

Technical Services

US and Canada: Call (800) 227-9770 (option 3, 4, 3).

Or send an e-mail to techservices@agilent.com

Agilent's world-wide Sales and Support contact details can be obtained at

www.agilent.com/en/contact-us/page.

MATERIALS PROVIDED

Materials provided	Quantity		
	Catalog #600670	Catalog #600672	Catalog #600674
PfuUltra II Fusion HS DNA Polymerase	40 reactions	200 reactions	400 reactions
10× PfuUltra II Reaction Buffer	1 ml	2 × 1 ml	4 × 1 ml

Storage: Store at –20°C upon receipt.

INTRODUCTION

In Agilent's PfuUltra II fusion HS DNA polymerase, we couple the fusion polymerase technology with our engineered PfuUltra DNA polymerase, hotstart antibody, 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
PfuUltra II fusion HS DNA polymerase	1 µl	1 µl	1 µl
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)
Primers (each)	0.2 µM each primer	0.4 µM each primer	0.2 µM each primer
dNTP concentration	250 µM each dNTP (1 mM total)	500 µM each dNTP (2 mM total)	250 µM each dNTP (1 mM total)
Final reaction buffer conc	1.0×	1.0×	1.0×
Denaturing temperature	95°C	92°C	95°C
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 ≤10 kb. See the Optimization Parameters section for guidelines on amplifying longer targets. The reaction conditions are for one reaction and must be adjusted for multiple samples. The final volume of each sample reaction is 50 µl. Add the components in order into sterile thin-walled PCR tubes while mixing gently.

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

Component	Amount per reaction
Distilled water (dH ₂ O)	40.5 µl
10× PfuUltra II reaction buffer ^a	5.0 µl
dNTP mix (25 mM each dNTP)	0.5 µl
DNA template (100 ng/µl)	1.0 µl ^b
Primer #1 (10 µM)	1.0 µl
Primer #2 (10 µM)	1.0 µl
PfuUltra II fusion HS DNA polymerase	1.0 µl
Total reaction volume	50 µl

^a The 10× buffer provides a final 1× Mg²⁺ concentration of 2 mM.

^b 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 vector (5–30 ng) PCR targets.

Perform PCR using optimized cycling conditions. Suggested cycling parameters are indicated in the table below. The PCR cycling parameters have been tested on the following: the MJ Research DNA Engine PTC-200, the Applied Biosystems GeneAmp PCR system 9700, the Applied Biosystems GeneAmp PCR system 9600, 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.

PCR Cycling Parameters for PfuUltra II fusion HS DNA Polymerase^a

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

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

Targets >10 kb (vector or genomic DNA)

Segment	Number of cycles	Temperature	Duration
1	1	92°C	2 minutes
2	30	92°C	10 seconds
		Primer $T_m - 5^\circ\text{C}$ ^c	20 seconds
		68°C	30 seconds per kb
3	1	68°C	5 minutes

^a Thin-wall PCR tubes are highly recommended. These PCR tubes are optimized to ensure more efficient heat transfer and to maximize thermal-cycling performance.

^b Denaturing temperatures above 95°C are recommended only for GC-rich templates.

^c The annealing temperature may require optimization. Typically annealing temperatures will range between 55° and 72°C.

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.

NOTICES TO PURCHASER

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).

NOTICE TO PURCHASER: LIMITED LICENSE

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 6,258,569, 6,171,785, 6,127,155, 6,030,787, 5,994,056, 5,876,930, 5,804,375, 5,789,224, 5,773,258 (claims 1 and 6 only), 5,723,591, 5,677,152 (claims 1 to 23 only), 5,618,711, 5,538,848, and claims outside the US corresponding to expired US Patent No. 5,079,352. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

LIMITED LABEL LICENSE FOR PFU-CONTAINING DNA POLYMERASE PRODUCTS

Purchase of this product conveys to the purchaser the non-transferable right to use the product for research use only. 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.

MSDS INFORMATION

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