



# ***PfuTurbo* Hotstart DNA Polymerase**

## **Instruction Manual**

**Catalog #600320, #600322, and #600324**

Revision G0

**Laboratory Reagent.**

600320-12



## 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.

## ORDERING INFORMATION

Please visit [www.agilent.com](http://www.agilent.com).

## TECHNICAL SERVICES

### United States and Canada

Email: [techservices@agilent.com](mailto:techservices@agilent.com)

Telephone: 800 227 9770 (option 3,4,3)

### All Other Locations

Please visit [www.agilent.com/en/contact-us/page](http://www.agilent.com/en/contact-us/page).

# **PfuTurbo Hotstart DNA Polymerase**

## **CONTENTS**

<b>Materials Provided</b> .....	<b>1</b>
<b>Storage Conditions</b> .....	<b>1</b>
<b>Notices to Purchaser</b> .....	<b>1</b>
<b>Additional Materials Required</b> .....	<b>1</b>
<b>Introduction</b> .....	<b>2</b>
<b>Critical Optimization Parameters for <i>PfuTurbo</i> Hotstart DNA Polymerase-Based PCR</b> .....	<b>4</b>
Extension Time.....	4
Enzyme Concentration .....	4
Primer-Template Purity and Concentration.....	5
Primer Design.....	5
Deoxynucleoside Triphosphates.....	6
Reaction Buffer .....	6
Magnesium Ion Concentration .....	6
PCR Cycling Parameters .....	6
Amplification of Genomic Targets >6 kb .....	7
<b>Application Notes</b> .....	<b>7</b>
Thermostability .....	7
Terminal Transferase Activity.....	7
Reverse Transcriptase Activity.....	7
<b>PCR Protocol Using <i>PfuTurbo</i> Hotstart DNA Polymerase</b> .....	<b>8</b>
<b>Troubleshooting</b> .....	<b>10</b>
<b>Preparation of Media and Reagents</b> .....	<b>10</b>
<b>References</b> .....	<b>11</b>
<b>MSDS Information</b> .....	<b>11</b>

# PfuTurbo Hotstart DNA Polymerase

## MATERIALS PROVIDED

Materials provided	Quantity		
	Catalog #600320	Catalog #600322	Catalog #600324
PfuTurbo hotstart DNA polymerase (2.5 U/μl)	100 U	500 U	1000 U
10× Cloned Pfu DNA polymerase reaction buffer <sup>a</sup>	1 ml	2 × 1 ml	4 × 1 ml

<sup>a</sup> See Preparation of Media and Reagents.

## STORAGE CONDITIONS

All components: -20°C

## NOTICES TO PURCHASER

### 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.

## ADDITIONAL MATERIALS REQUIRED

Temperature cycler  
PCR tubes  
PCR primers  
Deoxynucleoside triphosphates (dNTP's)

## INTRODUCTION

---

*PfuTurbo* hotstart DNA polymerase is a special formulation of Agilent's high fidelity *PfuTurbo* DNA polymerase with hot start activity. (See Table I for a comparison of Agilent hot start PCR enzymes.) *PfuTurbo* hotstart DNA polymerase is formulated with heat labile monoclonal antibodies that, at room temperature, effectively neutralize DNA polymerase and 3'-5' exonuclease (proofreading) activities. Full enzyme activity is regained upon denaturation of the antibody during the initial denaturation step. *PfuTurbo* hotstart DNA polymerase retains the high fidelity, sensitivity and yield of *PfuTurbo* DNA polymerase, while facilitating room temperature PCR assembly. *PfuTurbo* hotstart DNA polymerase provides reduced background and improved detection sensitivity by preventing priming until stringent primer annealing temperatures are reached. Amplification systems that are most likely to benefit from the hot start capability of *PfuTurbo* hotstart DNA polymerase are those (1) designed to detect low-copy-number targets in complex DNA backgrounds, (2) prone to primer-dimer formation, and (3) assembled robotically during high throughput PCR procedures.

*PfuTurbo* DNA polymerase has an error rate six-fold lower than *Taq* DNA polymerase, and significantly lower than the error rates of most other proofreading enzymes or DNA polymerase mixtures.<sup>1-3</sup> (See Table II for a comparison of the fidelity characteristics of many of the commercially available DNA polymerases.) Finally, the enhanced performance of *PfuTurbo* DNA polymerase allows the use of shorter extension times, fewer PCR cycles, and lower DNA template concentrations than are required for *Pfu* DNA polymerase, making it ideally suited for high-performance PCR applications.

*PfuTurbo* hotstart DNA polymerase provides robust amplification of long, complex genomic targets. A key component of *PfuTurbo* hotstart DNA polymerase is the exclusive thermostable ArchaeMaxx polymerase-enhancing factor. The ArchaeMaxx factor eliminates a PCR inhibitor and promotes shorter extension times, higher yield, and greater target length capabilities. The ArchaeMaxx factor improves the yield of products by overcoming dUTP poisoning, which is caused by dUTP accumulation during PCR through dCTP deamination.<sup>4</sup> Once incorporated, dU-containing DNA inhibits *Pfu* and most archaeal proofreading DNA polymerases, such as Vent and Deep Vent DNA polymerases, limiting their efficiency.<sup>4</sup> The ArchaeMaxx factor functions as a dUTPase, converting poisonous dUTP to harmless dUMP and inorganic pyrophosphate, resulting in improved overall PCR performance.

**TABLE I**

**Properties of Agilent Hot Start PCR Enzymes**

Hot Start PCR enzyme	Hot Start Method	Activities Neutralized	Activation Procedure <sup>a</sup>	Enzyme Applications
<i>PfuTurbo</i> hotstart DNA polymerase	Antibody	DNA polymerase, 3'-5' exonuclease	PCR Activation 30 cycles	<ul style="list-style-type: none"> <li>highest fidelity</li> <li>genomic DNA templates up to 19 kb</li> </ul>
Herculase hotstart DNA polymerase	Antibody	DNA polymerase, 3'-5' exonuclease	PCR Activation 30 cycles	<ul style="list-style-type: none"> <li>challenging cloning targets</li> <li>long and/or GC-rich targets</li> <li>higher fidelity than <i>Taq</i> DNA polymerase</li> </ul>
SureStart <i>Taq</i> DNA polymerase	Chemical	DNA polymerase, 5'-3' exonuclease	Pre-PCR Activation (9–12 minutes @ 95°C) 30 cycles or PCR Activation 40 cycles	<ul style="list-style-type: none"> <li>routine PCR up to 3 kb</li> </ul>

<sup>a</sup> PCR activation means that full enzyme activity is recovered during temperature cycling, either during the initial denaturation step (antibody-based formulations) or within the first 5–15 cycles (chemical hot start). For SureStart *Taq*, slow enzyme activation during temperature cycling typically necessitates the use of additional PCR cycles to achieve desired product yield (35–45 cycles). In the Pre-PCR activation method, the enzyme is activated prior to temperature cycling, and no additional cycles are necessary.

**TABLE II**

**Comparison of Thermostable DNA Polymerases Using a *lacIOZ $\alpha$* -Based Fidelity Assay<sup>a</sup>**

Thermostable DNA polymerase	Error rate <sup>b</sup>	Percentage (%) of mutated 1-kb PCR products <sup>c</sup>
<i>PfuUltra</i> high-fidelity DNA polymerase	$4.3 \times 10^{-7}$	0.9
<i>PfuTurbo</i> DNA polymerase	$1.3 \times 10^{-6}$	2.6
<i>Pfu</i> DNA polymerase	$1.3 \times 10^{-6}$	2.6
Deep Vent <sub>R</sub> DNA polymerase	$2.7 \times 10^{-6}$	5.4
Vent <sub>R</sub> DNA polymerase	$2.8 \times 10^{-6}$	5.6
PLATINUM <i>Pfx</i>	$3.5 \times 10^{-6}$	5.6
KOD DNA polymerase	$3.5 \times 10^{-6}$	5.6
<i>Taq</i> DNA polymerase	$8.0 \times 10^{-6}$	16.0

<sup>a</sup> Fidelity is measured using a published PCR forward mutation assay that is based on the *lacI* target gene.<sup>2</sup>

<sup>b</sup> The error rate equals mutation frequency per base pair per duplication.

<sup>c</sup> The percentage of mutated PCR products after amplification of a 1-kb target sequence for 20 effective cycles ( $2^{20}$ - or  $10^6$ -fold amplification).

# CRITICAL OPTIMIZATION PARAMETERS FOR *PfuTurbo* HOTSTART DNA POLYMERASE-BASED PCR

All PCR amplification reactions require optimization to achieve the highest product yield and specificity. Critical optimization parameters for successful PCR using *PfuTurbo* hotstart DNA polymerase are outlined in Table III and are discussed in the following section. These parameters include the use of an extension time that is adequate for full-length DNA synthesis, sufficient enzyme concentration, adequate primer–template purity and concentration, optimal primer design, and appropriate nucleotide concentration.

**TABLE III**  
**Optimization Parameters and Suggested Reaction Conditions**  
**(50  $\mu$ l reaction volume)**

<b>Parameter</b>	<b>Targets: &lt; 10 kb vector DNA or &lt; 6 kb genomic DNA</b>	<b>Targets: &gt; 10 kb vector DNA or &gt; 6 kb genomic DNA</b>
Extension time	1 min per kb	2 min per kb
<i>PfuTurbo</i> hotstart DNA polymerase	2.5 U	5.0 U
Input template	50–100 ng genomic DNA <sup>o</sup>	200–250 ng genomic DNA <sup>o</sup>
Primers (each)	100–200 ng (0.2–0.5 $\mu$ M)	200 ng (0.5 $\mu$ M)
dNTP concentration	100–250 $\mu$ M each dNTP (0.4–1.0 mM total)	500 $\mu$ M each dNTP (2 mM total)
Final reaction buffer concentration	1.0 $\times$	1.5 $\times$ (genomic DNA targets) 1.0 $\times$ (vector DNA targets)
Denaturing temperature	95°C	92°C
Extension temperature	72°C	68°C

<sup>o</sup> See *Primer-Template Purity and Concentration* for recommended amounts of other forms of template DNA.

## Extension Time

Extension time is one of the most critical parameters affecting the yield of PCR product obtained using *PfuTurbo* hotstart DNA polymerase. For optimal yield, use an extension time of 1.0 minute per kb for vector targets up to 10 kb and genomic targets up to 6 kb. When amplifying vector targets greater than 10 kb or genomic targets greater than 6 kb in length, use an extension time of 2.0 minutes per kb.

## Enzyme Concentration

The concentration of *PfuTurbo* hotstart DNA polymerase required for optimal PCR product yield and specificity depends on the individual target system to be amplified. Most amplifications are successful with 2.5 U of enzyme per 50  $\mu$ l reaction for vector targets up to 19 kb and for genomic targets up to 6 kb.

## Primer-Template Purity and Concentration

The most successful PCR results are achieved when the amplification reaction is performed using purified primers and templates that are essentially free of extraneous salts. Gel-purified primers, generally >18 nucleotides in length, are strongly recommended for use in *PfuTurbo* hotstart DNA polymerase-based PCR.

Additionally, an adequate concentration of primers and template should be used to ensure a good yield of the desired PCR products. When DNA of known concentration is available, amounts of 50–100 ng of DNA template per 50- $\mu$ l reaction are typically used for amplifications of single-copy chromosomal targets. When amplifying genomic targets greater than 6 kb, increase the template amount to 200–250 ng. The amplification of a single-copy target from complex genomic DNA is generally more difficult than amplification of a fragment from a plasmid or phage. Less DNA template can be used for amplification of lambda (1–30 ng) or plasmid (100 pg–10 ng) PCR targets or for amplification of multicopy chromosomal genes (10–100 ng).<sup>5</sup>

Use primers at a final concentration of 0.2–0.5  $\mu$ M, which is equivalent to ~100–200 ng of an 18- to 25-mer oligonucleotide primer in a 50- $\mu$ l reaction volume.

## Primer Design

Primer pairs that exhibit similar melting temperatures and are completely complementary to the template are recommended. Depending on the primer design and the desired specificity of the PCR amplification reaction, melting temperatures between 55° and 80°C generally yield the best results.<sup>5</sup> The following formula<sup>6</sup> is commonly used for estimating the melting temperature ( $T_m$ ) of the primers:

$$T_m(^{\circ}\text{C}) \cong 2(N_A + N_T) + 4(N_G + N_C)$$

where  $N$  equals the number of primer adenine (A), thymidine (T), guanine (G), or cytosine (C) bases. Several other articles present additional equations for estimating the melting temperature of primers.<sup>7, 8</sup> Finally, care must be taken when using degenerate primers. Degenerate primers should be designed with the least degeneracy at the 3' end. Optimization of degenerate primer concentration is necessary.

**Note** *Because of the unique composition of the Pfu buffer, the actual primer  $T_m$  may be 3°–5°C lower than that estimated by this formula.*



## Deoxynucleoside Triphosphates

For *PfuTurbo* hotstart DNA polymerase-based PCR, use a dNTP concentration range of 100–250  $\mu\text{M}$  each dNTP (0.4–1.0 mM total) for optimal product yield. Supplying dNTPs in this concentration range generally results in the optimal balance of product yield (greatest at high dNTP concentrations) versus specificity and fidelity (highest at low dNTP concentration).<sup>2, 5</sup> The yield produced from genomic targets >6 kb in length can be improved by increasing nucleotide concentration to 500  $\mu\text{M}$  (each dNTP) and the reaction buffer to 1.5 $\times$  final concentration.<sup>1</sup> The use of a balanced pool of dNTPs (equimolar amounts of each dNTP) ensures the lowest rate of misincorporation errors.

## Reaction Buffer

The reaction buffer provided with this enzyme has been formulated for optimal PCR yield and fidelity when performing PCR amplification using *PfuTurbo* hotstart DNA polymerase. Use this reaction buffer in standard PCR amplification reactions. If alterations are made to the buffer, significant increases in the error rate of *Pfu* hotstart DNA polymerase can be avoided by maintaining the  $\text{Mg}^{2+}$  concentration above 1.5 mM, the total dNTP concentration at 0.4–1.0 mM, and the pH of Tris-based buffers above pH 8.0 when measured at 25°C.<sup>2</sup>

To improve yields of genomic targets >6 kb, increase the final concentration of reaction buffer from 1 $\times$  to 1.5 $\times$ .<sup>1</sup>

## Magnesium Ion Concentration

Magnesium ion concentration affects primer annealing and template denaturation, as well as enzyme activity and fidelity. Generally, excess  $\text{Mg}^{2+}$  results in accumulation of nonspecific amplification products, whereas insufficient  $\text{Mg}^{2+}$  results in reduced yield of the desired PCR product.<sup>9</sup> PCR amplification reactions should contain *free*  $\text{Mg}^{2+}$  in excess of the total dNTP concentration. For *PfuTurbo* hotstart DNA polymerase-based PCR, yield is optimal when the *total*  $\text{Mg}^{2+}$  concentration is  $\sim 2$  mM in a standard reaction mixture, and  $\sim 3$  mM for amplification of cDNA. A 2 mM total  $\text{Mg}^{2+}$  concentration is present in the final 1 $\times$  dilution of the 10 $\times$  cloned *Pfu* DNA polymerase reaction buffer. For the amplification of cDNA,  $\text{Mg}^{2+}$  should be added to the PCR reaction to a final concentration of 3 mM.

## PCR Cycling Parameters

Standard PCR amplification reactions typically require 25–30 cycles to obtain a high yield of PCR product. Thermal cycling parameters should be chosen carefully to ensure (1) the shortest denaturation times to avoid template damage, (2) adequate extension times to achieve full-length target synthesis, and (3) the use of annealing temperatures near the primer melting temperature to improve yield of the desired PCR product.

When performing PCR on a new target system, we suggests using an annealing temperature 5°C below the lowest primer melting temperature.

For best results, PCR primers should be designed with similar melting temperatures ranging from 55° to 80°C. The use of primers with melting temperatures within this range reduces false priming and ensures complete denaturation of unextended primers at 94–95°C (see also *Primer-Template Purity and Concentration* and *Primer Design*).

See Table IV for suggested PCR cycling parameters, depending upon template size and complexity. Optimized cycling parameters are not necessarily transferable between thermal cyclers designed by different manufacturers. Therefore, each manufacturer's recommendations for optimal cycling parameters should be consulted.

### **Amplification of Genomic Targets >6 kb**

To improve yields of genomic targets >6 kb, increase the amount of *PfuTurbo* hotstart DNA polymerase from 2.5 U to 5.0 U, and increasing the final concentration of reaction buffer from 1× to 1.5×.<sup>1</sup> Use 200–250 ng of genomic template DNA, 200 ng of each primer, and 500 μM each dNTP. Use a denaturing temperature of 92°C, an extension temperature of 68°C, and an extension time of 2.0 minutes per kilobase. Finally, overlay each reaction with ~50 μl of RNase/DNase-free mineral oil prior to thermal cycling.

## **APPLICATION NOTES**

---

### **Thermostability**

*PfuTurbo* hotstart DNA polymerase is a highly thermostable enzyme, retaining 94–99% of its polymerase activity after 1 hour at 95°C. The half life of *Pfu* DNA polymerase is approximately 19 hours at 95°C. Unlike *Taq* DNA polymerase, denaturing temperatures up to 98°C can be used successfully with *PfuTurbo* hotstart DNA polymerase to amplify GC-rich regions.<sup>10,11</sup>

### **Terminal Transferase Activity**

Studies demonstrate that thermostable DNA polymerases, with the exception of *Pfu* DNA polymerase, exhibit terminal deoxynucleotidyl-transferase (TdT) activity, which is characterized by the addition of nontemplate-directed nucleotide(s) at the 3' end of PCR-generated fragments.<sup>12,13</sup> *Pfu* DNA polymerase is devoid of TdT activity and *PfuTurbo* hotstart DNA polymerase generates blunt-ended PCR products exclusively. Therefore, *PfuTurbo* hotstart DNA polymerase is optimal for use with the PCR-Script Amp SK(+) cloning kit<sup>14</sup> and the PCR-Script Cam SK(+) cloning kit.<sup>15</sup> Alternatively, *PfuTurbo* hotstart DNA polymerase can be used to remove 3' overhangs (polishing) or to fill-in 5' overhangs with greater efficiencies than either Klenow polymerase or T4 DNA polymerase.<sup>16,17</sup>

### **Reverse Transcriptase Activity**

*PfuTurbo* hotstart DNA polymerase lacks detectable reverse transcriptase activity.

## PCR PROTOCOL USING *PfuTurbo* HOTSTART DNA POLYMERASE

---

1. Prepare a reaction mixture for the appropriate number of samples to be amplified. Add the components *in order* while mixing gently. The table shown below provides an example of a reaction mixture for the amplification of a typical single-copy chromosomal target. The recipe listed in the table below is for one reaction and must be adjusted for multiple samples. The final volume of each sample reaction is 50  $\mu$ l.

### Reaction Mixture for a Typical Single-Copy Chromosomal Locus PCR Amplification

Component	Amount per reaction
Distilled water (dH <sub>2</sub> O)	40.6 $\mu$ l
10 $\times$ Cloned <i>Pfu</i> reaction buffer <sup>a</sup>	5.0 $\mu$ l
dNTPs (25 mM each dNTP)	0.4 $\mu$ l
DNA template (100 ng/ $\mu$ l)	1.0 $\mu$ l <sup>b</sup>
Primer #1 (100 ng/ $\mu$ l)	1.0 $\mu$ l <sup>c</sup>
Primer #2 (100 ng/ $\mu$ l)	1.0 $\mu$ l <sup>c</sup>
<i>PfuTurbo</i> hotstart DNA polymerase (2.5 U/ $\mu$ l)	1.0 $\mu$ l (2.5 U) <sup>d</sup>
Total reaction volume	50 $\mu$ l

- <sup>a</sup> The 10 $\times$  buffer provides a final 1 $\times$  Mg<sup>2+</sup> concentration of 2 mM. To amplify cDNA, Mg<sup>2+</sup> may need to be added to a final 1 $\times$  concentration of 3 mM.
  - <sup>b</sup> The amount of DNA template required varies depending on the type of DNA being amplified. Generally 50–100 ng of genomic DNA template is recommended. Less DNA template (typically 0.1–30 ng) can be used for amplification of lambda or plasmid PCR targets or 10–100 ng for amplification of multicopy chromosomal genes.
  - <sup>c</sup> Primer concentrations between 0.2 and 0.5  $\mu$ M are recommended (this corresponds to 100–250 ng for typical 18- to 25-mer oligonucleotide primers in a 50- $\mu$ l reaction volume).
  - <sup>d</sup> The amount of *PfuTurbo* hotstart DNA polymerase varies depending on the length of the template to be amplified. The standard amount for vector targets up to 19 kb and genomic targets up to 6 kb in length is 1  $\mu$ l (2.5 U).
2. Aliquot 50  $\mu$ l of the reaction mixture into the appropriate number of sterile thin-wall PCR tubes or standard 0.5-ml microcentrifuge tubes.
  3. If the extension times are >15 minutes, overlay each reaction with DNase-, RNase-, and protease-free mineral oil (Sigma, St. Louis, Missouri) even if the temperature-cycler is equipped with a heated cover.
  4. Perform PCR using optimized cycling conditions (see also *PCR Cycling Parameters*). Suggested cycling parameters are indicated in Table IV.
  5. Analyze the PCR amplification products on a 0.7–1.0% (w/v) agarose gel.

**TABLE IV****PCR Cycling Parameters for *PfuTurbo* DNA Polymerase with Single-Block Temperature Cyclers <sup>a,b</sup>****A. Targets <10 kb vector DNA or <6 kb genomic DNA**

Segment	Number of cycles	Temperature	Duration
1	1	95°C <sup>c</sup>	2 minutes
2	30	95°C	30 seconds
		Primer $T_m - 5^\circ\text{C}^d$	30 seconds
		72°C	1 minute for targets ≤1 kb 1 minute per kb for targets >1 kb <sup>e</sup>
3	1	72°C	10 minutes

**B. Targets >10 kb vector DNA or >6 kb genomic DNA**

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

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

<sup>b</sup> Optimized cycling parameters are not necessarily transferable between thermal cyclers designed by different manufacturers; therefore, each manufacturer's recommendations for optimal cycling parameters should be consulted.

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

<sup>d</sup> The annealing temperature may require optimization. Typically annealing temperatures will range between 55° and 72°C.<sup>5</sup>

<sup>e</sup> The annealing temperature may require optimization. Typical annealing temperatures will range between 60 and 65°C.

## TROUBLESHOOTING

Observation	Solution(s)
No product or low yield	Increase extension time to 2 minutes/kb of PCR target
	Lower the annealing temperature in 5°C increments
	Ensure that 10× cloned <i>Pfu</i> DNA polymerase reaction buffer is used
	Use higher denaturing temperatures (94–98°C) if the template DNA contains a high GC content or secondary structures (see also Reference 12)
	Use the recommended primer concentrations between 0.2 and 0.5 μM (generally 100–200 ng for typical 18- to 25-mer oligonucleotide primers in a 50-μl reaction volume)
	Use only high quality, gel purified primers (see <i>Primer–Template Purity and Concentration</i> and <i>Primer Design</i> )
	Check the melting temperature, purity, GC content, and length of the primers
	Remove extraneous salts from the PCR primers and DNA preparations. High ionic strength can inhibit reactions
	Denaturation times of 30–60 seconds at 94–95°C are usually sufficient while longer denaturation times may damage the DNA template; use the shortest denaturation time compatible with successful PCR on the thermal cycler
	Increase the amount of <i>PfuTurbo</i> hotstart DNA polymerase
	Use intact and highly purified templates at an adequate concentration (see <i>Primer–Template Purity and Concentration</i> and <i>Primer Design</i> )
	Excessive template DNA can be inhibitory. Follow the recommendations given in the manual for template amount
	Insufficient heat exchange between the reaction tubes and the thermal cycler can cause inadequate temperature ramping. Use thin-wall PCR tubes. These PCR tubes are optimized to ensure ideal contact with the multiblock design to permit more efficient heat transfer and to maximize thermal-cycling performance
Multiple bands	Increase the annealing temperature in 5°C increments to ensure sufficient primer annealing
	Use Perfect Match PCR enhancer to improve PCR product specificity
Artifactual smears	Decrease the amount of <i>PfuTurbo</i> hotstart DNA polymerase
	Reduce the extension time utilized

## PREPARATION OF MEDIA AND REAGENTS

### 10× Cloned *Pfu* DNA Polymerase Reaction Buffer

- 200 mM Tris-HCl (pH 8.8)
- 20 mM MgSO<sub>4</sub>
- 100 mM KCl
- 100 mM (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>
- 1% Triton X-100
- 1 mg/ml nuclease-free BSA

## REFERENCES

---

1. Cline, J., Braman, J. and Kretz, K. (1995) *Strategies* 8(1):24–25.
2. Cline, J., Braman, J. C. and Hogrefe, H. H. (1996) *Nucleic Acids Res* 24(18):3546-51.
3. Nielson, K. B., Cline, J., Bai, F., McMullan, D., McGowan, B. et al. (1997) *Strategies* 10(1):29–32.
4. Hogrefe, H. H., Hansen, C. J., Scott, B. R. and Nielson, K. B. (2002) *Proc Natl Acad Sci U S A* 99(2):596–601.
5. Innis, M. A., Gelfand, D. H., Sninsky, J. J. and White, T. J. (1990). *PCR Protocols: A Guide to Methods and Applications*. Academic Press, New York.
6. Thein, S. L. and Wallace, R. B. (1986). *Human Genetic Diseases: A Practical Approach*. IRL Press, Herndon, Virginia.
7. Rychlik, W., Spencer, W. J. and Rhoads, R. E. (1990) *Nucleic Acids Res* 18(21):6409-12.
8. Wu, D. Y., Ugozzoli, L., Pal, B. K., Qian, J. and Wallace, R. B. (1991) *DNA Cell Biol* 10(3):233-8.
9. Saiki, R. (1989). Chapter 1. In *PCR Technology: Principles and Applications for DNA Amplification*, H. A. Erlich (Ed.). Stockton Press, New York.
10. Chong, S. S., Eichler, E. E., Nelson, D. L. and Hughes, M. R. (1994) *Am J Med Genet* 51(4):522-6.
11. Nielson, K. B., Braman, J. and Kretz, K. (1995) *Strategies* 8(1):26.
12. Costa, G. L., Grafsky, A. and Weiner, M. P. (1994) *PCR Methods Appl* 3(6):338-45.
13. Hu, G. (1993) *DNA Cell Biol* 12(8):763-770.
14. Bauer, J. C., Deely, D., Braman, J., Viola, J. and Weiner, M. P. (1992) *Strategies* 5(3):62-64.
15. Costa, G. L., Sanchez, T. and Weiner, M. P. (1994) *Strategies* 7(2):52.
16. Costa, G. L. and Weiner, M. P. (1994) *Nucleic Acids Res* 22(12):2423.
17. Eliason, E. and Detrick, J. (1995) *Strategies* 8(1):27–28.

## MSDS INFORMATION

---

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