



***PfuUltra* High-Fidelity DNA Polymerase**

Instruction Manual

Catalog #600380, #600382, and #600384

Revision F0

Laboratory Reagent.

600380-12



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.

TECHNICAL SERVICES

United States and Canada

Email: techservices@agilent.com

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

All Other Locations

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

***PfuUltra* High-Fidelity DNA Polymerase**

CONTENTS

Materials Provided	1
Storage Conditions	1
Additional Materials Required	1
Notices to Purchaser	1
Introduction	2
Critical Optimization Parameters for <i>PfuUltra</i> High-Fidelity DNA Polymerase-Based PCR	3
Extension Time.....	3
Enzyme Concentration.....	3
Primer-Template Purity and Concentration.....	4
Primer Design.....	4
Deoxynucleoside Triphosphates.....	5
Reaction Buffer.....	5
PCR Cycling Parameters.....	5
Amplification of Genomic Targets >6 kb.....	6
Additional Optimization Parameters	6
Order of Addition of Reaction Mixture Components.....	6
Magnesium Ion Concentration.....	6
Adjuncts and Cosolvents.....	7
Application Notes	8
Thermostability.....	8
Inherent “Hot Start” Properties.....	8
Terminal Transferase Activity.....	8
Reverse Transcriptase Activity.....	8
PCR Protocol for <i>PfuUltra</i> High-Fidelity DNA Polymerase	9
Troubleshooting	11
References	12
MSDS Information	12

***PfuUltra* High-Fidelity DNA Polymerase**

MATERIALS PROVIDED

Material provided	Quantity		
	Catalog #600380	Catalog #600382	Catalog #600384
<i>PfuUltra</i> HF DNA polymerase (2.5 U/ μ l)	100 U	500 U	1000 U
10 \times <i>PfuUltra</i> HF reaction buffer	1 ml	2 \times 1 ml	4 \times 1 ml

STORAGE CONDITIONS

All components: -20°C

ADDITIONAL MATERIALS REQUIRED

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

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.

INTRODUCTION

The *PfuUltra* high-fidelity DNA polymerase formulation features a genetically engineered mutant of *Pfu* DNA polymerase and the ArchaeMaxx polymerase-enhancing factor. *PfuUltra* high-fidelity DNA polymerase exhibits an average error rate three-fold lower than *PfuTurbo* DNA polymerase and 18-fold lower than *Taq* DNA polymerase, making it the highest fidelity enzyme available (See Table I). Amplification of a 500-bp fragment using *PfuUltra* high-fidelity DNA polymerase results in errors in less than 0.5% of full-length PCR products, making *PfuUltra* high-fidelity DNA polymerase the ideal enzyme for PCR cloning.

In addition to high accuracy, *PfuUltra* high-fidelity DNA polymerase provides robust amplification of long, complex genomic targets. A key component of *PfuUltra* high-fidelity DNA polymerase is the 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.¹ Once incorporated, dU-containing DNA inhibits *Pfu* and most archaeal proofreading DNA polymerases, such as Vent and Deep Vent DNA polymerases, limiting their efficiency.¹ The ArchaeMaxx factor functions as a dUTPase, converting poisonous dUTP to harmless dUMP and inorganic pyrophosphate, resulting in improved overall PCR performance.

TABLE I

Comparison of Thermostable DNA Polymerases Using a *lacIOZ α* -Based Fidelity Assay^a

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

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

^b The error rate equals mutation frequency per base pair per duplication.

^c 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 *PfuUltra* HIGH-FIDELITY 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 *PfuUltra* high-fidelity DNA polymerase are outlined in Table II 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 II
Optimization Parameters and Suggested Reaction Conditions
(50- μ l reaction volume)

Parameter	Targets: < 10 kb vector DNA or < 6 kb genomic DNA	Targets: > 10 kb vector DNA or > 6 kb genomic DNA
Extension time	1 min per kb	2 min per kb
<i>PfuUltra</i> high-fidelity DNA polymerase	2.5 U	5.0 U
Input template	50–100 ng genomic DNA ^a	200–250 ng genomic DNA ^a
Primers (each)	100–200 ng (0.2–0.5 μ M)	200 ng (0.5 μ M)
dNTP concentration	200–250 μ M each dNTP (0.8–1.0 mM total)	500 μ 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

^a 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 *PfuUltra* high-fidelity 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 *PfuUltra* high-fidelity 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- μ l reaction for vector targets up to 17 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 *PfuUltra* high-fidelity 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- μ 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).³

We suggest using primers at a final concentration of 0.2–0.5 μ M, which is equivalent to ~100–200 ng of an 18- to 25-mer oligo-nucleotide primer in a 50- μ 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.³ The following formula⁴ 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.^{5, 6} 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 *PfuUltra* high-fidelity DNA polymerase-based PCR, We recommend a dNTP concentration range of 200–250 μM each dNTP (0.8–1.0 mM total) for optimal product yield. Supplying dNTPs at 100–175 μM each dNTP generally results in good product yield; however, at these dNTP concentrations, yields achieved in amplifications performed using *PfuUltra* high-fidelity DNA polymerase may be lower compared to identical amplifications performed using *PfuTurbo* DNA polymerase. The yield of genomic targets >6 kb in length can be improved by increasing nucleotide concentration to 500 μM (each dNTP) and the reaction buffer to 1.5 \times final concentration.⁷ 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 *PfuUltra* high-fidelity DNA polymerase. We strongly recommend use of the reaction buffer provided with *PfuUltra* high-fidelity DNA polymerase in standard PCR amplification reactions. If alterations in these buffers are made, significant increases in the error rate of *Pfu* DNA polymerase can be avoided by maintaining the 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.²

To improve yields of genomic targets >6 kb, we recommend increasing the final concentration of reaction buffer from 1 \times to 1.5 \times .⁷

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 suggest 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 target 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, we recommend increasing the amount of *PfuUltra* high-fidelity DNA polymerase from 2.5 U to 5.0 U, and increasing the final concentration of reaction buffer from 1× to 1.5×.⁷ 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.

ADDITIONAL OPTIMIZATION PARAMETERS

Order of Addition of Reaction Mixture Components

Because *PfuUltra* high-fidelity DNA polymerase exhibits 3'- to 5'-exonuclease activity that enables the polymerase to proofread nucleotide misincorporation errors, it is critical that the enzyme is the last component added to the PCR mixture (i.e., **after** the dNTPs). In the absence of dNTPs, the 3'- to 5'-exonuclease activity of proofreading DNA polymerases may degrade primers. When primers and nucleotides are present in the reaction mixture at recommended levels (i.e., primer concentrations of 0.2–0.5 μM and nucleotide concentrations of 200–250 μM each dNTP), primer degradation is not a concern.

Magnesium Ion Concentration

Magnesium ion concentration affects primer annealing and template denaturation, as well as enzyme activity and fidelity. Generally, excess Mg²⁺ results in accumulation of nonspecific amplification products, whereas insufficient Mg²⁺ results in reduced yield of the desired PCR product.⁸ PCR amplification reactions should contain free Mg²⁺ in excess of the total dNTP concentration. For *PfuUltra* DNA polymerase-based PCR, yield is optimal when the total Mg²⁺ concentration is ~2 mM in a standard reaction mixture, and ~3 mM for amplification of cDNA. A 2 mM total Mg²⁺ concentration is present in the final 1× dilution of the 10× *PfuUltra* HF reaction buffer. For the amplification of cDNA, Mg²⁺ should be added to the PCR reaction to a final concentration of 3 mM.³

Adjuncts and Cosolvents

The adjuncts or cosolvents listed in the following table may be advantageous with respect to yield when used in the PCR buffer. Fidelity may or may not be affected by the presence of these adjuncts or cosolvents.

Adjunct or cosolvent	Optimal PCR final concentration
Formamide	1.25–10%
Dimethylsulfoxide (DMSO)	1–10%
Glycerol	5–20%
Perfect Match PCR enhancer	1 U per 50- μ l reaction (genomic DNA template) 0.01–1 U per 50- μ l reaction (plasmid DNA template)

Formamide

Formamide facilitates certain primer–template annealing reactions and also lowers the denaturing temperature of melt-resistant DNA.⁹

Dimethylsulfoxide and Glycerol

Cosolvents, such as DMSO and glycerol, improve the denaturation of GC-rich DNA and help overcome the difficulties of polymerase extension through secondary structures. Studies indicate that the presence of 1–10% DMSO in PCR may be essential for the amplification of the retinoblastoma gene¹⁰ and may also enhance amplification of *Herpes simplex* virus (HSV) sequences.¹¹ Glycerol is known to improve the yield of amplification products and also serves as an enzyme stabilizer.¹¹

Perfect Match PCR Enhancer

Perfect Match PCR enhancer (Catalog #600129) improves the specificity of PCR products. This adjunct performs this function by destabilizing mismatched primer–template complexes and by helping to remove secondary structures that could impede normal extension.¹²

APPLICATION NOTES

Thermostability

Pfu DNA polymerase is a highly thermostable enzyme, retaining 94–99% of its polymerase activity after 1 hour at 95°C. Unlike *Taq* DNA polymerase, denaturing temperatures up to 98°C can be used successfully with *PfuUltra* high-fidelity DNA polymerase to amplify GC-rich regions.^{13,14}

Inherent “Hot Start” Properties

PfuUltra high-fidelity DNA polymerase exhibits optimal polymerase activity at $\geq 75^\circ\text{C}$ and only 2–8% activity between 40–50°C. *Taq* DNA polymerase, however, exhibits optimal polymerase activity between 60–70°C and 27–70% activity between 40–50°C. The minimal activity of *PfuUltra* high-fidelity DNA polymerase at lower temperatures should result in fewer mispaired primer-extension reactions than occur with *Taq* DNA polymerase during the lower temperatures encountered during PCR cycling (e.g., primer annealing). Consequently, hot start techniques, which are commonly used with *Taq* DNA polymerase to improve product yield and specificity, generally are not required for PCR amplifications with *PfuUltra* high-fidelity DNA polymerase.¹⁵

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.^{16, 17} *PfuUltra* high-fidelity DNA polymerase, like *Pfu* and *PfuTurbo* DNA polymerases, is devoid of TdT activity and generates blunt-ended PCR products exclusively. Therefore, *PfuUltra* high-fidelity, *PfuTurbo* and *Pfu* DNA polymerases are the enzymes of choice for use with the PCR-Script Amp SK(+) cloning kit¹⁸, and the PCR-Script Cam SK(+) cloning kit.¹⁹

In addition, *PfuUltra*, *PfuTurbo* or *Pfu* 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.^{20,21}

Reverse Transcriptase Activity

PfuUltra high-fidelity DNA polymerase lacks detectable reverse transcriptase activity.

PCR PROTOCOL FOR *PfuUltra* HIGH-FIDELITY DNA POLYMERASE

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

TABLE III

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

Component	Amount per reaction
Distilled water (dH ₂ O)	40.6 μ l
10 \times <i>PfuUltra</i> HF reaction buffer ^a	5.0 μ l
dNTPs (25 mM each dNTP)	0.4 μ l
DNA template (100 ng/ μ l)	1.0 μ l ^b
Primer #1 (100 ng/ μ l)	1.0 μ l ^c
Primer #2 (100 ng/ μ l)	1.0 μ l ^c
<i>PfuUltra</i> HF DNA polymerase (2.5 U/ μ l)	1.0 μ l (2.5 U) ^d
Total reaction volume	50 μ l

^a The 10 \times buffer provides a final 1 \times Mg²⁺ concentration of 2 mM. To amplify cDNA, Mg²⁺ may need to be added to a final 1 \times concentration of 3 mM.

^b 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.

^c Primer concentrations between 0.2 and 0.5 μ M are recommended (this corresponds to 100–250 ng for typical 18- to 25-mer oligonucleotide primers in a 50- μ l reaction volume).

^d The amount of *PfuUltra* HF DNA polymerase varies depending on the length of the template to be amplified. The standard amount for vector targets up to 17 kb and genomic targets up to 6 kb in length is 1 μ l (2.5 U).

2. Immediately before thermal cycling, aliquot 50 μ l of the reaction mixture into the appropriate number of sterile thin-wall PCR tubes or standard 0.5-ml microcentrifuge tubes.
3. Perform PCR using optimized cycling conditions (see also *PCR Cycling Parameters*). Suggested cycling parameters are indicated in Table IV.
4. Analyze the PCR amplification products on a 0.7–1.0% (w/v) agarose gel.

TABLE IV**PCR Cycling Parameters for *PfuUltra* High-Fidelity DNA Polymerase ^{a,b}****A. Targets <10 kb vector DNA or <6 kb genomic DNA**

Segment	Number of cycles	Temperature	Duration
1	1	95°C ^c	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 ^e
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

^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 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.

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

^d The annealing temperature may require optimization. Typically annealing temperatures will range between 55° and 72°C.³

^e 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 per kb of PCR target
	Use the recommended amount of DNA template amounts. Use of excess template can reduce PCR product yield
	Lower the annealing temperature in 5°C increments
	Use a high-quality dNTP mix to supply a final concentration of $\geq 200 \mu\text{M}$ each dNTP
	Ensure that 10× <i>PfuUltra</i> HF reaction buffer is used
	Remove extraneous salts from the PCR primers and DNA preparations
	Add <i>PfuUltra</i> HF DNA polymerase last to the reaction mixture to minimize any potential primer degradation
	Use higher denaturing temperatures (94–98°C)
	Use cosolvents such as DMSO in a 1–10% (v/v) final concentration or glycerol in a 5–20% (v/v) final concentration (see <i>Dimethylsulfoxide and Glycerol</i>)
	Use the recommended primer concentrations between 0.2 and 0.5 μM (corresponding to 100–200 ng for typical 18- to 25-mer oligonucleotide primers in a 50- μl reaction volume)
	Use high-quality primers
	Check the melting temperature, purity, GC content, and length of the primers
	Consider using the adjuncts or cosolvents such as Perfect Match PCR Enhancer or formamide (see <i>Adjuncts and Cosolvents</i>)
	Perform further buffer optimization if necessary
	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>PfuUltra</i> HF DNA polymerase	
Use intact and highly purified templates at an adequate concentration (see <i>Primer-Template Purity and Concentration</i>)	
Multiple bands	Increase the annealing temperature in 5°C increments
	Use Perfect Match PCR enhancer to improve PCR product specificity
Artifactual smears	Decrease the amount of <i>PfuUltra</i> HF DNA polymerase
	Reduce the extension time utilized

REFERENCES

1. 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.
2. Cline, J., Braman, J. C. and Hogrefe, H. H. (1996) *Nucleic Acids Res* 24(18):3546-51.
3. 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.
4. Thein, S. L. and Wallace, R. B. (1986). *Human Genetic Diseases: A Practical Approach*. IRL Press, Herndon, Virginia.
5. Rychlik, W., Spencer, W. J. and Rhoads, R. E. (1990) *Nucleic Acids Res* 18(21):6409-12.
6. Wu, D. Y., Ugozzoli, L., Pal, B. K., Qian, J. and Wallace, R. B. (1991) *DNA Cell Biol* 10(3):233-8.
7. Cline, J., Braman, J. and Kretz, K. (1995) *Strategies* 8(1):24–25.
8. Saiki, R. (1989). Chapter 1. In *PCR Technology: Principles and Applications for DNA Amplification*, H. A. Erlich (Ed.). Stockton Press, New York.
9. Sarkar, G., Kapelner, S. and Sommer, S. S. (1990) *Nucleic Acids Res* 18(24):7465.
10. Hung, T., Mak, K. and Fong, K. (1990) *Nucleic Acids Res* 18(16):4953.
11. Smith, K. T., Long, C. M., Bowman, B. and Manos, M. M. (1990) *Amplifications* 5:16-17.
12. Nielsen, K. and Mather, E. (1990) *Strategies* 3(2):17-22.
13. Chong, S. S., Eichler, E. E., Nelson, D. L. and Hughes, M. R. (1994) *Am J Med Genet* 51(4):522-6.
14. Nielson, K. B., Braman, J. and Kretz, K. (1995) *Strategies* 8(1):26.
15. Nielson, K., Cline, J. and Hogrefe, H. H. (1997) *Strategies* 10(2):40-43.
16. Costa, G. L., Grafsky, A. and Weiner, M. P. (1994) *PCR Methods Appl* 3(6):338-45.
17. Hu, G. (1993) *DNA Cell Biol* 12(8):763-770.
18. Bauer, J. C., Deely, D., Braman, J., Viola, J. and Weiner, M. P. (1992) *Strategies* 5(3):62-64.
19. Costa, G. L., Sanchez, T. and Weiner, M. P. (1994) *Strategies* 7(2):52.
20. Costa, G. L. and Weiner, M. P. (1994) *Nucleic Acids Res* 22(12):2423.
21. 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. MSDS documents are not included with product shipments.