

TECHNICAL BULLETIN

GoTaq[®] PCR Core System

Instructions for Use of Product
M7660



GoTaq[®] PCR Core System

All technical literature is available at: www.promega.com/protocols/
Visit the web site to verify that you are using the most current version of this Technical Bulletin.
E-mail Promega Technical Services if you have questions on use of this system: techserv@promega.com

| | |
|--|----|
| 1. Description..... | 2 |
| 2. Product Components and Storage Conditions | 2 |
| 3. PCR Protocol | 3 |
| 3.A. Protocol..... | 3 |
| 3.B. Analysis | 5 |
| 4. Optimization of PCR..... | 5 |
| 4.A. Routine PCR | 5 |
| 4.B. Magnesium Concentration | 6 |
| 4.C. Buffer Considerations | 6 |
| 4.D. Enzyme Concentration | 7 |
| 4.E. Primer Design | 7 |
| 4.F. Template Considerations..... | 7 |
| 4.G. Primer Annealing Temperature..... | 9 |
| 4.H. Extension Temperature | 9 |
| 4.I. Nucleic Acid Cross-Contamination | 9 |
| 5. Troubleshooting..... | 10 |
| 6. References..... | 12 |
| 7. Related Products..... | 13 |
| 8. Summary of Changes | 16 |



1. Description

The GoTaq[®] PCR Core System I^(a,b) provides all of the reagents necessary for the exponential amplification of specific regions of DNA using the polymerase chain reaction (PCR; 1). The GoTaq[®] PCR Core System I is supplied with GoTaq[®] DNA Polymerase, a proprietary formulation of *Taq* DNA polymerase designed for enhanced amplification, and GoTaq[®] Flexi Buffers, magnesium-free proprietary buffer formulations supplied at pH 8.5. The 5X Green GoTaq[®] Flexi Buffer increases the density of the sample and contains two dyes (a blue dye and a yellow dye) that separate during electrophoresis, allowing reactions to be loaded directly onto an agarose gel without the addition of loading dye. During electrophoresis in a 1% agarose gel, the blue dye migrates at the same rate as a 3–5kb DNA fragment, and the yellow dye migrates slightly faster than primers (<50bp). The 5X Colorless GoTaq[®] Flexi buffer allows fluorescence or absorbance measurements to be made directly on the amplified DNA without the need to purify it from the amplification reaction.

Although simple in theory, PCR can benefit from optimization of several parameters. This technical bulletin provides important and convenient information about performing PCR and includes protocols and troubleshooting tips for successful PCR amplifications.

Comprehensive information on the PCR process is available online in the PCR Chapter of the *Protocols and Applications Guide* at: www.promega.com/paguide

2. Product Components and Storage Conditions

| PRODUCT | CAT.# |
|--------------------------------------|-------|
| GoTaq [®] PCR Core System I | M7660 |

For Laboratory Use. Includes:

- 250u GoTaq[®] DNA Polymerase
- 2 × 1ml 5X Colorless GoTaq[®] Flexi Buffer, Mg-Free
- 2 × 1ml 5X Green GoTaq[®] Flexi Buffer, Mg-Free
- 1.2ml MgCl₂, 25mM
- 200µl PCR Nucleotide Mix, 10mM

Storage Conditions: Store all components at –20°C. See product label for expiration date.

3. PCR Protocol

To facilitate optimization, troubleshooting and validation of PCR, we strongly recommend performing concurrent positive and negative control reactions.

If working with multiple samples, assemble a master mix of water, $MgCl_2$, 5X buffer, primers, PCR Nucleotide Mix and GoTaq[®] DNA Polymerase. Combine the appropriate multiples of these components (except template) and dispense aliquots of the appropriate volume, minus the specific volume to be occupied by the template DNA, to each reaction tube. Initiate the reaction by adding the template. Use individual pipette tips for all additions, being careful not to cross-contaminate the samples.

If possible, start with $>10^4$ copies of the target sequence to obtain a signal in 25–30 cycles. Excess template is not beneficial to the reaction. Always ensure that the final DNA concentration is $\leq 10ng/\mu l$. Less than 10 copies of a target can be amplified (2), but more cycles may be required to detect a signal by gel electrophoresis. Additional cycles may increase nonspecific amplification, evidenced by multiple bands when analyzed by gel electrophoresis.

Materials to Be Supplied by the User

- Mineral Oil
- Nuclease-Free Water (Cat.# P1193)
- template DNA
- upstream primer
- downstream primer

3.A. Protocol

1. Combine the components, as listed in Table 1, in sterile, 0.5–0.6ml microcentrifuge tubes on ice. Amplification reactions may be scaled up or down as necessary. Use of a master mix, as described above, greatly facilitates the reaction setup and decreases tube-to-tube variability of the reaction components. Thaw the 25mM $MgCl_2$ solution and GoTaq[®] Flexi Buffer completely and vortex thoroughly before use. We strongly advise optimizing the $MgCl_2$ concentration.
2. If using a thermal cycler without a heated lid, overlay the reaction mix with 1–2 drops (approximately 50 μ l) of Mineral Oil to prevent evaporation during thermal cycling. Centrifuge the reaction mix in a microcentrifuge for 5 seconds.
3. Place the reactions in a thermal cycler that has been preheated to 95°C. We recommend heating the samples at 95°C for 2 minutes to ensure that the target DNA is completely denatured. Incubation longer than 2 minutes at 95°C is unnecessary and may reduce the yield.
4. Start the thermal cycling program. The cycling profile given in Table 2 may be used as a guideline. We recommend that you optimize the amplification profile for each primer/target combination.

3.A. Protocol (continued)

Table 1. Recommended Reaction Volumes and Final Concentrations of the GoTaq® PCR Core System Components.

| Component | Component Volume (25µl) | Component Volume (50µl) | Final Concentration |
|---|--------------------------------|--------------------------------|-----------------------------|
| MgCl ₂ , 25mM Solution (see Section 4.B) | 1.0–4.0µl | 2.0–8.0µl | 1.0–4.0mM |
| 5X Colorless GoTaq® Flexi Buffer OR 5X Green GoTaq® Flexi Buffer | 5µl | 10µl | 1.0X |
| PCR Nucleotide Mix, 10mM each | 0.5µl | 1µl | 200µM each |
| upstream primer | 2.5–25pmol | 5–50pmol | 0.1–1.0µM |
| downstream primer | 2.5–25pmol | 5–50pmol | 0.1–1.0µM |
| GoTaq® DNA Polymerase, 5u/µl | 0.125µl | 0.25µl | 0.625u/25µl 1.25u/50µl |
| template DNA | variable | variable | <0.25µg/25µl <0.5µg/50µl |
| Nuclease-Free Water to a final volume of | 25µl | 50µl | |



Thaw the 25mM MgCl₂ solution and GoTaq® Flexi Buffer completely and vortex thoroughly before use. We strongly advise optimizing the MgCl₂ concentration.

Table 2. Thermal Cycling Guidelines for PCR Amplification.

| Step | Temperature | Time | Number of Cycles |
|----------------------|--------------------|--------------|-------------------------|
| Initial Denaturation | 95°C | 2 minutes | 1 cycle |
| Denaturation | 95°C | 0.5–1 minute | 25–35 cycles |
| Annealing | 42–65°C* | 0.5–1 minute | |
| Extension | 72°C | 1 minutes/kb | |
| Final Extension | 72°C | 5 minutes | 1 cycle |
| Soak | 4°C | indefinite | 1 cycle |

These guidelines apply to target sequences between 200–2,000bp and may need to be adapted for your thermal cycler.

*The annealing temperature for a specific amplification reaction will depend upon the sequences of the two primers. See Section 4.G for discussions on how to determine optimal annealing temperatures for PCR amplification.

3.B. Analysis

1. Analyze PCR products by agarose gel electrophoresis. The products should be readily visible in an ethidium bromide-stained gel under UV light. Reactions containing Green GoTaq® Buffer do not need loading dye added before electrophoresis.
2. Store PCR products at -20°C until needed. The PCR products can be further purified using a number of procedures including the Wizard® SV Gel and PCR Clean-Up System (Cat.# A9281).

4. Optimization of PCR

4.A. Routine PCR

As originally developed, the PCR process amplifies short (approximately 100–500bp) segments of a longer DNA molecule (1). A typical amplification reaction includes the target DNA, a thermostable DNA polymerase, two oligonucleotide primers, deoxynucleotide triphosphates (dNTPs), reaction buffer, magnesium and optional additives. The components of the reaction are mixed and placed in a thermal cycler, an automated instrument that ‘cycles’ the reaction through a predetermined series of specific temperatures and times. One cycle of amplification is defined by the series of temperature and time adjustments. Each cycle of PCR after the first cycle theoretically doubles the amount of targeted template sequence (amplimer). Therefore, ten cycles theoretically multiplies the amplimer by a factor of about one thousand; 20 cycles, by a factor of more than one million. PCR amplification can be completed in as little as 2 hours.

Each cycle of PCR amplification consists of a defined number of reaction steps. The steps are designed using temperature and duration time to denature the template, anneal the two oligonucleotide primers and extend the new complementary DNA strands by polymerization. These steps can be optimized for each template and primer pair combination. The target DNA is denatured by heating to 95°C or higher for 15–120 seconds. During denaturation, the two intertwined strands of DNA separate from one another, producing the necessary single-stranded DNA (ssDNA) template for primer annealing and polymerization (extension) by a thermostable polymerase. To anneal the oligonucleotide primers, the temperature of the next step in the cycle is reduced to approximately $42\text{--}65^{\circ}\text{C}$. At this temperature, the oligonucleotide primers can anneal to the ssDNA strands and serve as primers for DNA synthesis by the polymerase. This step requires approximately 30–60 seconds. Finally, to extend from the primer-bound template DNA, the reaction temperature is raised to the optimum for most thermostable DNA polymerases, which is approximately 72°C . Extension of the primer by the thermostable polymerase requires approximately 1 minute per kilobase to be amplified. Extension completes one cycle, and the next cycle begins by returning the reaction to 95°C for denaturation. After 20–40 cycles, the amplified nucleic acid may then be analyzed (e.g., for size, quantity or sequence) or it may be used in further experimental procedures (e.g., cloning or mutagenesis).

4.B. Magnesium Concentration

Magnesium concentration is a crucial factor affecting the performance of thermostable enzymes. Reaction components, including template DNA, chelating agents present in the sample (e.g., EDTA or citrate), dNTPs and proteins, can affect the amount of free magnesium. In the absence of adequate free magnesium, GoTaq® DNA Polymerase is inactive. Conversely, excess free magnesium reduces enzyme fidelity (3) and may increase the level of nonspecific amplification (4,5). For these reasons, it is important to empirically determine the optimal MgCl₂ concentration for each reaction. This is accomplished by preparing a series of reactions containing 1.0–4.0mM Mg²⁺, in increments of 0.5mM, by adding 2, 3, 4, 5, 6, 7 and 8µl of a 25mM MgCl₂ stock to 50µl reactions.

The GoTaq® PCR Core System includes 5X Green GoTaq® Flexi Buffer, Mg-Free; 5X Colorless GoTaq® Flexi Buffer, Mg-Free; and a solution of 25mM MgCl₂. The buffers allow you to adjust the Mg²⁺ concentration to the level that is optimal for each reaction using the 25mM MgCl₂.

Two important steps will ensure the reaction contains the appropriate amount of Mg²⁺: Magnesium solutions must be thawed completely and vortexed for several seconds prior to use because magnesium chloride solutions can form concentration gradients when frozen. Thawing and vortexing is required to render the solution uniform with respect to magnesium salts. These two simple steps can eliminate a major source of many failed experiments.

4.C. Buffer Considerations

We recommend using the 5X Green GoTaq® Flexi Buffer in any amplification reaction that will be visualized by agarose gel electrophoresis followed by ethidium bromide staining. The 5X Green GoTaq® Flexi Buffer is not recommended for any downstream applications using absorbance or fluorescence excitation, as the yellow and blue dyes in the reaction buffer may interfere with these applications. The dyes absorb between 225–300nm, making standard A₂₆₀ readings to determine DNA concentration unreliable. Also, the dyes have excitation peaks at 488nm and between 600–700nm that correspond to the excitation wavelengths commonly used in fluorescence detection instrumentation. However, for some instrumentation, such as a fluorescent gel scanner that uses a 488nm excitation wavelength, there will be minimal interference, since it is the yellow dye that absorbs this wavelength. Gels scanned by this method will have a light gray dye front below the primers corresponding to the yellow dye front. The Green and Colorless GoTaq® Flexi Buffers give approximately equivalent amplification yields. Balanced amplifications between the two buffers may require further optimization.

For reactions going directly from a thermal cycler to an application using absorbance or fluorescence, the 5X Colorless GoTaq® Flexi Buffer is recommended. If both agarose gel analysis and further downstream applications involving absorbance or fluorescence will be used, the two dyes can be removed from the Green GoTaq® Flexi reactions using standard PCR cleanup systems like the Wizard® SV Gel and PCR Clean-Up System (Cat.# A9281) or the Wizard® SV 96 PCR Clean-Up System (Cat.# A9341).

Both reaction buffers are compatible with common PCR additives such as DMSO and betaine. These additives do not change the color of the Green GoTaq® Flexi Buffer or affect dye migration.

4.D. Enzyme Concentration

We recommend that 1.25 units of GoTaq® DNA Polymerase be used per 50µl amplification reaction. For most applications, enzyme will be in excess; the inclusion of more enzyme will not significantly increase product yield. Increased amounts of enzyme and excessively long extension times increase the likelihood of generating artifacts. Artifacts generally can be seen as smearing of bands in ethidium bromide-stained agarose gels (6–8).

The most frequent cause of excessive enzyme levels is pipetting error. Accurate dispensing of submicroliter (<1µl) volumes of enzyme solutions in 50% glycerol is nearly impossible. We strongly recommend the use of reaction master mixes sufficient for the number of reactions being performed to obviate this problem. A master mix increases the volumes of pipetted reagents and reduces pipetting errors.

4.E. Primer Design

PCR primers generally range in length from 15–30 bases and are designed to flank the region of interest. Primers should contain 40–60% (G + C), and care should be taken to avoid sequences that might produce internal secondary structure. To avoid the production of primer-dimers, the 3' ends of the primers should not be complementary. Primer dimers unnecessarily deplete primers from the reaction and result in an unwanted polymerase reaction that competes with the desired reaction. Avoid three G or C nucleotides in a row near the 3' end of the primer, as this may result in nonspecific primer annealing, increasing the synthesis of undesirable reaction products. Ideally, both primers should have nearly identical melting temperatures (T_m); in this manner, the two primers should anneal roughly at the same temperature. In any case, the annealing temperature of the reaction is dependent upon the primer with the lowest T_m .

The sequence of the primers can also include regions at the 5' ends that may prove useful for downstream applications. For example, restriction enzyme sites can be designed into the primer pair for ease in downstream manipulations such as cloning. Regardless of primer choice, the final concentration of the primer in the reaction must be optimized. We recommend adding 50pmol of each primer (1µM final concentration in a 50µl reaction) as a starting point for optimization. Generally, nanograms of primer DNA equivalent to 50pmol is: $16.3\text{ng} \times b$, where b is the number of bases in the primer.

4.F. Template Considerations

Successful amplification of the region of interest depends upon the amount and quality of the template DNA. Reagents commonly used to purify nucleic acids (e.g., salts, guanidine, proteases, organic solvents and SDS) are potent inhibitors of DNA polymerases. A final ethanol precipitation of the nucleic acid sample will eliminate most of the inhibitory agents. Spiking a control DNA fragment and the appropriate primer pair into the DNA preparation may be useful in verifying the purity of the DNA sample.

The amount of template required for successful amplification depends upon the complexity of the DNA sample. For example, whereas a 4kb plasmid containing a 1kb insert equates to 25% of the DNA being the target of interest, a 1kb gene in human genomic DNA (genome of 3.3×10^9 bp) represents approximately 0.00003% of the input DNA. Therefore, approximately 1,000,000-fold more human genomic DNA is required to maintain the same number of target copies per reaction.

4.F. Template Considerations (continued)

Two common mistakes encountered when trying to amplify target DNA are using too much plasmid DNA and too little genomic DNA. Table 3 lists the correlation of molecules per microgram of nucleic acids from some common RNA and DNA targets. Table 4 shows typical yields of genomic DNA from a variety of source materials. As a general guide for how much template DNA to use, start with a minimum of 10^4 copies of the target sequence to obtain a signal in 25–30 cycles, but keep the final DNA concentration of the reaction $\leq 10\text{ng}/\mu\text{l}$.

Table 3. Conversion of Nucleic Acids from Microgram Amount to Number of Molecules.

| Nucleic Acid | Amount | # of Molecules |
|------------------------------|-----------------|-----------------------|
| 1kb RNA | 1 μg | 1.8×10^{12} |
| 1kb dsDNA | 1 μg | 9.18×10^{11} |
| pGEM [®] Vector DNA | 1 μg | 2.9×10^{11} |
| lambda (λ) DNA | 1 μg | 1.9×10^{10} |
| <i>E. coli</i> genomic DNA | 1 μg | 2×10^8 |
| human genomic DNA | 1 μg | 3.0×10^5 |

Table 4. DNA Yields from Different Human Tissue Sources.

| Amount of Material | Source of DNA Typically Used | Typical Yield |
|---------------------------|-------------------------------------|----------------------|
| Whole blood | 30 μl | 0.5–1 μg |
| Blood spot | 1/2 of a 5mm spot | 1–3 μg |
| Cell suspension | 5×10^5 cells | 2–5 μg |
| Buccal cells | Single mouth rinse | 0.1–1 μg |
| Chorionic villus biopsy | Small frond | 1–3 μg |
| Semen | 30 μl | 5–10 μg |
| Hair root | Single root | 10–200ng |
| Tissue block | 50mg | 0.1–10 μg |

4.G. Primer Annealing Temperature

The sequences of the primers are a major consideration in determining the optimal temperature of the PCR amplification cycles. For primers with a high T_m , it may be advantageous to increase the annealing temperature. Higher temperatures minimize nonspecific primer annealing, increase the amount of specific product produced and reduce the amount of primer-dimer formation.

Numerous formulas exist to determine the theoretical T_m of nucleic acids (9,10), and these may serve as a starting point for annealing conditions. However, it is best to optimize the annealing conditions by performing the reaction at several temperatures starting approximately 5°C below the calculated T_m .

For a T_m calculator that can calculate melting temperatures for primers in a GoTaq® reaction, visit our web site at: www.promega.com/biomath/

The formula below can be used to estimate the melting temperature for any oligonucleotide:

$$T_m = 81.5 + 16.6 \times (\log_{10}[\text{Na}^+]) + 0.41 \times (\%G+C) - 675/n,$$

where $[\text{Na}^+]$ is the molar concentration of monovalent cations and
 n = number of bases in the oligonucleotide.

Example: To calculate the melting temperature of a 22mer oligonucleotide with 60% G+C in 50mM KCl:

$$\begin{aligned} T_m &= 81.5 + 16.6 \times (\log_{10}[0.05]) + 0.41 \times (60) - 675/22 \\ &= 81.5 + 16.6 \times (-1.30) + 24.60 - 30.68 \\ &= 53.84^\circ\text{C} \end{aligned}$$

4.H. Extension Temperature

During the extension step, allow approximately 1 minute for every 1kb to be amplified (minimum extension time of 1 minute). Generally, 25–40 cycles are sufficient for most reactions.

4.I. Nucleic Acid Cross-Contamination

It is important to minimize cross-contamination between samples and prevent carryover of RNA and DNA from one experiment to the next. Use separate work areas and pipettors for pre- and post-amplification steps. Use positive displacement pipettes or aerosol-resistant tips to reduce cross-contamination during pipetting. Wear gloves and change them often. Consider using a contamination control technique (11) to prevent DNA carryover to subsequent reactions.



5. Troubleshooting

For questions not addressed here, please contact your local Promega Branch Office or Distributor. Contact information available at: www.promega.com. E-mail: techserv@promega.com

Symptoms

Low yield or no amplification

Causes and Comments

Insufficient number of cycles. Return reactions to thermal cycler for 5 or more cycles.

Template degraded. Verify the integrity of the DNA by electrophoresis after incubation in the presence of Mg^{2+} .

Thermal cycler programmed incorrectly. Verify that times and temperatures are correct. Use step cycles, not hold segments.

Temperature too low in some positions of thermal cycler. Perform a set of control reactions to determine whether certain positions in the thermal cycler give low yields.

Top of thermal cycler is open. The top must be closed for correct heating and cooling.

Inhibitor present in DNA template. Reduce the volume of sample in the reaction. Ethanol precipitate template DNA to remove inhibitor.

Improper reaction conditions. Reduce the annealing temperature and/or allow longer extension times for longer amplifiers. Also, optimize Mg^{2+} concentration.

Missing reaction component. Check the reaction components and repeat the reaction.

Mineral oil problem. The reaction must be overlaid with high-quality, nuclease-free light mineral oil. **Do not use autoclaved mineral oil.**

Poor primer design. Make sure primers are not self-complementary or complementary to each other.

Incorrect primer specificity. Verify that the primers are complementary to the appropriate strands.

Primer concentration too low. Verify primer concentration in the reaction and increase the concentration as needed.

Suboptimal reaction conditions. Optimize Mg^{2+} concentration, annealing temperature and extension time. Always vortex the Mg^{2+} . Verify that primers are present in equal concentrations.

Symptoms
Causes and Comments

Low yield or no amplification (continued)

Degraded nucleotides or primers. Keep nucleotides and primers frozen in aliquots, thaw quickly and keep on ice once thawed. Avoid multiple freeze-thaw cycles.

Target sequence not present in target DNA. Redesign experiment or try other sources of target DNA.

Low yield or no amplification in mouse tail genotyping applications

Adjust annealing temperature. Reaction buffer composition affects the melting properties of DNA. See the BioMath Calculator to calculate the melting temperature for primers in the GoTaq[®] reaction (www.promega.com/biomath). The salt-adjusted and base-stacking T_m s are lower in GoTaq[®] reaction buffer than in *Taq* DNA polymerase reaction buffer.

Reduce volume of template DNA in reaction or dilute template DNA prior to adding to reaction. Many DNA preparations for genotyping applications result in the copurification of amplification inhibitors. Reducing the amount of inhibitors present in the reaction may improve results. Diluting samples even 1:10,000 has been shown to be effective, depending on initial DNA concentration.

Ethanol precipitate template DNA. Include an ethanol precipitation and wash step prior to amplification to remove inhibitors that copurified with the DNA.

Add PCR enhancers. Adding PCR-enhancing agents may improve yields. General stabilizing agents such as BSA (Sigma Cat.# A7030; final concentration 0.16mg/ml) may also help to overcome amplification failure.

Multiple, nonspecific amplification products

Suboptimal reaction conditions. Optimize MgCl₂ concentration, annealing temperature, size, extension time and cycle number to minimize nonspecific priming.

Poor primer design. Make sure primers are not self-complementary or complementary to each other, especially near the 3' ends. Try a longer primer. Avoid use of three consecutive G or C nucleotides at the 3' end of the primers.

Primer concentration too high. Verify primer concentration in the reaction. Try a lower concentration in the reaction.

5. Troubleshooting (continued)

| <u>Symptoms</u> | <u>Causes and Comments</u> |
|--|--|
| Multiple, nonspecific amplification products (continued) | <p>Contamination by another target DNA.</p> <ul style="list-style-type: none"> • Use positive displacement pipettes or aerosol-resistant tips to reduce cross-contamination during pipetting. • Use a separate work area and pipettor for pre- and post-amplification. • Wear gloves and change them often. • Use UNG (11) or another contamination control technique to prevent DNA carryover to subsequent reactions. <hr/> <p>Multiple target sequences exist in target DNA. Design new primers with higher specificity to target sequence.</p> |

6. References

1. Saiki, R. *et al.* (1985) Enzymatic amplification of β -globin genomic sequences and restriction site analysis for diagnosis of sickle cell anemia. *Science* **230**, 1350–4.
2. Saiki, R.K. *et al.* (1988) Primer-directed enzymatic amplification of DNA with a thermostable DNA polymerase. *Science* **239**, 487–91.
3. Eckert, K.A. and Kunkel, T.A. (1990) High fidelity DNA synthesis by the *Thermus aquaticus* DNA polymerase. *Nucleic Acids Res.* **18**, 3739–44.
4. Williams, J.F. (1989) Optimization strategies for the polymerase chain reaction. *BioTechniques* **7**, 762–9.
5. Ellsworth, D.L., Rittenhouse, K.D. and Honeycutt, R.L. (1993) Artfactual variation in randomly amplified polymorphic DNA banding patterns. *BioTechniques* **14**, 214–7.
6. Longley, M.J., Bennett, S.E. and Mosbaugh, D.W. (1990) Characterization of the 5' to 3' exonuclease associated with *Thermus aquaticus* DNA polymerase. *Nucleic Acids Res.* **18**, 7317–22.
7. Sardelli, A.D. (1993) Plateau effect—understanding PCR limitations. *Amplifications* **9**, 1–6.
8. Bell, D.A. and DeMarini, D.M. (1991) Excessive cycling converts PCR products to random-length higher molecular weight fragments. *Nucleic Acids Res.* **19**, 5079.
9. Baldino, F., Chesselet, M.F. and Lewis, M.E. (1989) High-resolution in situ hybridization histochemistry. *Methods Enzymol.* **168**, 761–77.
10. Rychlik, W., Spencer, W.J. and Rhoads, R.E. (1990) Optimization of the annealing temperature for DNA amplification in vitro. *Nucleic Acids Res.* **18**, 6409–12.
11. Longo, M.C., Berninger, M.S. and Hartley, J.L. (1990) Use of uracil DNA glycosylase to control carry-over contamination in polymerase chain reactions. *Gene* **93**, 125–8.

7. Related Products

Real-Time PCR Systems

| Product | Size | Cat.# |
|--|-------------|--------------|
| GoTaq [®] Probe qPCR Master Mix | 2ml | A6101 |
| | 10ml | A6102 |
| GoTaq [®] Probe 1-Step RT-qPCR System | 2ml | A6120 |
| | 12.5ml | A6121 |
| GoTaq [®] Probe 2-Step RT-qPCR System | 2ml | A6110 |
| GoTaq [®] qPCR Master Mix (BRYT Green [®] Dye) | 5ml | A6001 |
| | 25ml | A6002 |
| GoTaq [®] 1-Step RT-qPCR System (BRYT Green [®] Dye) | 5ml | A6020 |
| GoTaq [®] 2-Step RT-qPCR System (BRYT Green [®] Dye) | 5ml | A6010 |

Endpoint Hot-Start PCR

| Product | Size | Cat.# |
|--|-----------------|--------------|
| GoTaq [®] G2 Hot Start Polymerase | 100u | M7401 |
| | 500u | M7405 |
| | 2,500u | M7406 |
| | 10,000u | M7408 |
| GoTaq [®] G2 Hot Start Green Master Mix | 100 reactions | M7422 |
| | 1,000 reactions | M7423 |
| GoTaq [®] G2 Hot Start Colorless Master Mix | 100 reactions | M7432 |
| | 1,000 reactions | M7433 |

Endpoint Standard PCR

| Product | Size | Cat.# |
|---|-------------|--------------|
| GoTaq [®] G2 DNAPolymerase | 100u | M7841 |
| | 500u | M7845 |
| | 2,500u | M7848 |
| GoTaq [®] G2 Flexi DNAPolymerase | 100u | M7801 |
| | 500u | M7805 |
| | 2,500u | M7806 |
| | 10,000u | M7808 |



7. Related Products (continued)

Endpoint Standard PCR (continued)

| Product | Size | Cat.# |
|--------------------------------|-----------------|--------------|
| GoTaq® G2 Green Master Mix | 100 reactions | M7822 |
| | 1,000 reactions | M7823 |
| GoTaq® G2 Colorless Master Mix | 100 reactions | M7832 |
| | 1,000 reactions | M7833 |
| <i>Pfu</i> DNA Polymerase | 100u | M7741 |
| | 500u | M7745 |

Reverse Transcription Enzymes and Systems

| Product | Size | Cat.# |
|---|-------------------------|--------------|
| GoScript™ Reverse Transcription System | 50 reactions (of 20µl) | A5000 |
| | 100 reactions (of 20µl) | A5001 |
| GoScript™ Reverse Transcriptase | 100 reactions (of 20µl) | A5003 |
| | 500 reactions (of 20µl) | A5004 |
| GoScript™ Reverse Transcriptase Mix, Oligo(dT) | 50 reactions (of 20µl) | A2790 |
| | 100 reactions (of 20µl) | A2791 |
| GoScript™ Reverse Transcriptase Mix, Random Primers | 50 reactions (of 20µl) | A2800 |
| | 100 reactions (of 20µl) | A2801 |
| AccessQuick™ RT-PCR System | 20 reactions | A1701 |
| | 100 reactions | A1702 |
| | 500 reactions | A1703 |
| Access RT-PCR System | 100 reactions | A1250 |
| | 500 reactions | A1280 |
| Access RT-PCR Introductory System | 20 reactions | A1260 |

BenchTop DNA Markers

| Product | Size | Cat.# |
|-----------------------------------|-------------|--------------|
| BenchTop ΦX174 DNA/HaeIII Markers | 250µl | G7511 |
| BenchTop pGEM® DNA Markers | 250µl | G7521 |
| BenchTop PCR Markers | 300µl | G7531 |
| BenchTop 1kb DNA Ladder | 600µl | G7541 |
| BenchTop 100bp DNA Ladder | 300µl | G8291 |

Molecular Weight Markers

| Product | Size | Cat.# |
|-----------------------|-------------|--------------|
| 1kb DNA Step Ladder | 90µg | G6941 |
| 200bp DNA Step Ladder | 100µg | G6961 |
| 100bp DNA Step Ladder | 100µg | G6951 |
| 50bp DNA Step Ladder | 90µg | G4521 |
| 25bp DNA Step Ladder | 100µg | G4511 |
| 10bp DNA Step Ladder | 32.5µg | G4471 |
| 1kb DNA Ladder | 500µl | G5711 |

PCR Nucleotide Mix

| Product | Size | Cat.# |
|---------------------------------------|--------------|--------------|
| PCR Nucleotide Mix, 10mM | 200µl | C1141 |
| | 1,000µl | C1145 |
| PCR Nucleotide Mix, 25mM | 200µl | U1431 |
| | 1,000µl | U1432 |
| dNTP Mix | 200µl | U1511 |
| | 1,000µl | U1515 |
| dNTP Sets (dATP, dCTP, dGTP and dTTP) | 10µmol each | U1330 |
| | 25µmol each | U1420 |
| | 40µmol each | U1240 |
| | 200µmol each | U1410 |
| dNTP Sets (dATP, dCTP, dGTP and dUTP) | 10µmol each | U1335 |
| | 40µmol each | U1245 |

PCR Clean-Up System

| Product | Size | Cat.# |
|--|-------------|--------------|
| Wizard® SV Gel and PCR Clean-Up System | 50 preps | A9281 |
| | 250 preps | A9282 |

Additional size available



7. Related Products (continued)

PCR Cloning

| Product | Size | Cat.# |
|---|--------------|-------|
| pGEM [®] -T Easy Vector System I | 20 reactions | A1360 |
| pGEM [®] -T Easy Vector System II (includes JM109 Competent Cells) | 20 reactions | A1380 |
| pGEM [®] -T Vector System I | 20 reactions | A3600 |
| pGEM [®] -T Vector System II (includes JM109 Competent Cells) | 20 reactions | A3610 |

Staining Dye

| Product | Size | Cat.# |
|---------------------------------------|-------|-------|
| Diamond [™] Nucleic Acid Dye | 500µl | H1181 |

8. Summary of Changes

The following changes were made to the 5/19 revision of this document:

1. The document design was updated.
2. Product and related product information were updated.

^(a)Use of this product for basic PCR is outside of any valid US patents assigned to Hoffman La-Roche or Applera. This product can be used for basic PCR in research, commercial or diagnostic applications without any license or royalty fees.

^(b)U.S. Pat. No. 6,242,235, Australian Pat. No. 761757, Canadian Pat. No. 2,335,153, Chinese Pat. No. ZL99808861.7, Hong Kong Pat. No. HK 1040262, Japanese Pat. No. 3673175, European Pat. No. 1088060 and other patents pending.

© 1997–2019 Promega Corporation. All Rights Reserved.

BRYT Green, GoTaq, pGEM and Wizard are registered trademarks of Promega Corporation. AccessQuick, Diamond and GoScript are trademarks of Promega Corporation.

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.