

AmpliTaq Gold® DNA Polymerase

	Package Contents	<p>Catalog Number N808-0240 AmpliTaq Gold® with Buffer I 250 Units N808-0241 AmpliTaq Gold® with Buffer II 250 Units 4311806 AmpliTaq Gold® with Gold Buffer 250 Units</p> <p> Kit Contents</p>
	Storage Conditions	<ul style="list-style-type: none"> Store all contents at -20°C.
	Required Materials	<ul style="list-style-type: none"> Template: cDNA, gDNA, λDNA 10 mM dNTP mix (Cat. no. 18427-088) Forward and reverse gene-specific primers Autoclaved, distilled water E-Gel® General Purpose Gels, 1.2% (Cat. no. G5018-01) TrackIt™ 1 Kb Plus DNA Ladder (Cat. no. 10488-085) 0.2 or 0.5-mL nuclease-free microcentrifuge tubes
	Timing	Varies depending on amplicon length
	Selection Guide	<p>PCR Enzymes and Master Mixes Go online to view related products.</p>
	Product Description	<ul style="list-style-type: none"> AmpliTaq Gold® is derived from recombinant, thermostable, 94 kDa DNA polymerase, encoded by a modified form of the <i>Thermus aquaticus</i> DNA polymerase gene. It is available with 3 different buffers, which provide preferred pH and ionic strength for PCR amplification reactions. With AmpliTaq Gold® DNA Polymerase, Hot Start and Time Release PCR can be introduced into existing amplification systems by modifying cycling parameters or reaction conditions for increased specificity, sensitivity, and yield. AmpliTaq Gold® is provided in an inactive state and can be completely or partially activated in a pre-PCR heat step, allowing flexibility and assembly of reactions at room temperature.
	Important Guidelines	<ul style="list-style-type: none"> Take precautions to avoid cross-contamination by using aerosol-resistant barrier tips and analyzing PCR products in a separate area from PCR assembly. If the samples contain EDTA or other chelators, raise the MgCl₂ concentration in the reaction mix proportionately. If proteases are present in the sample DNA (e.g. impure genomic DNA), inactivate the proteases by heating samples to 95°C for 5 minutes before adding AmpliTaq Gold®.



Online Resources

Visit our [product page](#) for additional information and protocols. For support, visit www.lifetechnologies.com/support.



Enzyme Characteristics

Hot-start:	Chemical
Length:	Up to 5 kb
Fidelity vs. Taq:	1X
Format:	Separate components

PCR Reaction Setup

Use the measurements below to prepare your PCR experiment, or enter your own parameters in the column provided.

Component	25-µL rxn	50-µL rxn	Custom	Final Conc.
Autoclaved, distilled water	to 25 µL	to 50 µL	to µL	–
10X PCR Buffer I, II, or Gold Buffer	2.5 µL	5.0 µL	µL	1X
10 mM dNTP Mix*	0.5–5.0 µL	1.0 µL	µL	0.2 mM
25 mM MgCl ₂ **	1.5 µL	3.0 µL	µL	1.5 mM
10 µM forward primer	0.5 µL	1.0 µL	µL	0.2 µM
10 µM reverse primer	0.5 µL	1.0 µL	µL	0.2 µM
Template DNA	varies	varies		< 1 µg/rxn
AmpliTaq Gold® DNA Polymerase (5 U/µL)	0.125 µL	0.25 µL	µL	1.25 U/ 50-µL rxn

* Substituting dUTP for dTTP to control PCR product carryover may require higher concentrations of dUTP (typically twice that of any other dNTP) for optimal amplification.

** Use MgCl₂ with Buffer II or Gold only. Buffer I already contains Mg.

PCR Protocol

See page 2 to view a procedure for preparing and running your PCR experiment.






Optimization Strategies

Refer to the pop-up for guidelines to optimize your PCR reactions.

Limited Warranty, Disclaimer, and Licensing Information

AmpliQa Gold® DNA Polymerase Protocol

The example PCR procedure below shows appropriate volumes for a single 50- μL reaction. For multiple reactions, prepare a master mix of components common to all reactions to minimize pipetting error, and then dispense appropriate volumes into each 0.2–0.5 mL PCR reaction tube prior to adding template DNA and primers.

Timeline		Steps	Procedure Details																																																	
1		Thaw reagents	<p>Thaw, mix, and briefly centrifuge each component before use. Keep components on ice.</p> <p>Add the following components into appropriate wells or tubes. Note: Consider the volumes for all components listed in steps 2 and 3 to determine the correct amount of water required to reach your final reaction volume.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>50-μL rxn</th> <th>Final Concentration</th> </tr> </thead> <tbody> <tr> <td>Autoclaved, distilled water</td> <td>to 50 μL</td> <td></td> </tr> <tr> <td>10X PCR Buffer I, II, or Gold Buffer</td> <td>5.0 μL</td> <td>1X</td> </tr> <tr> <td>10 mM dNTP mix</td> <td>1.0 μL</td> <td>0.2 mM each</td> </tr> <tr> <td>25 mM MgCl_2 (with Buffer II or Gold only)</td> <td>3.0 μL</td> <td>1.5 mM</td> </tr> <tr> <td>AmpliQa Gold® DNA Polymerase (5 U/μL)</td> <td>0.25 μL</td> <td>1.25 U/rxn*</td> </tr> </tbody> </table> <p>* The amount of AmpliQa Gold® needed for the typical PCR amplification depends on cycling parameters. Start with 1.25 U/reaction.</p> <p>Mix and briefly centrifuge the components.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>50-μL rxn</th> <th>Final Concentration</th> </tr> </thead> <tbody> <tr> <td>10 μM forward primer</td> <td>1.0 μL</td> <td>0.2 μM</td> </tr> <tr> <td>10 μM reverse primer</td> <td>1.0 μL</td> <td>0.2 μM</td> </tr> <tr> <td>Template DNA</td> <td>varies</td> <td>< 1 μg/rxn*</td> </tr> </tbody> </table> <p>* Preferably > 10^4 copies of template but < 1 μg DNA/reaction.</p> <p>Cap each tube, mix, and then briefly centrifuge the contents.</p> <p>Note: The pre-PCR heat step can completely activate AmpliQa Gold®. If using the Time Release method, skip or reduce the duration of the initial denaturation step.</p> <table border="1"> <thead> <tr> <th>Step</th> <th>Temperature ($^{\circ}\text{C}$)</th> <th>Time</th> </tr> </thead> <tbody> <tr> <td>Initial Denaturation</td> <td>95</td> <td>10 minutes</td> </tr> <tr> <td rowspan="3">25–35 PCR Cycles</td> <td>Denature</td> <td>95</td> </tr> <tr> <td>Anneal</td> <td>~55 (depending on primer T_m)</td> </tr> <tr> <td>Extend</td> <td>72</td> </tr> <tr> <td>Final Extension</td> <td>72</td> <td>5 minutes</td> </tr> <tr> <td>Hold</td> <td>4</td> <td>indefinitely</td> </tr> </tbody> </table>	Component	50- μL rxn	Final Concentration	Autoclaved, distilled water	to 50 μL		10X PCR Buffer I, II, or Gold Buffer	5.0 μL	1X	10 mM dNTP mix	1.0 μL	0.2 mM each	25 mM MgCl_2 (with Buffer II or Gold only)	3.0 μL	1.5 mM	AmpliQa Gold® DNA Polymerase (5 U/ μL)	0.25 μL	1.25 U/rxn*	Component	50- μL rxn	Final Concentration	10 μM forward primer	1.0 μL	0.2 μM	10 μM reverse primer	1.0 μL	0.2 μM	Template DNA	varies	< 1 μg /rxn*	Step	Temperature ($^{\circ}\text{C}$)	Time	Initial Denaturation	95	10 minutes	25–35 PCR Cycles	Denature	95	Anneal	~55 (depending on primer T_m)	Extend	72	Final Extension	72	5 minutes	Hold	4	indefinitely
Component	50- μL rxn	Final Concentration																																																		
Autoclaved, distilled water	to 50 μL																																																			
10X PCR Buffer I, II, or Gold Buffer	5.0 μL	1X																																																		
10 mM dNTP mix	1.0 μL	0.2 mM each																																																		
25 mM MgCl_2 (with Buffer II or Gold only)	3.0 μL	1.5 mM																																																		
AmpliQa Gold® DNA Polymerase (5 U/ μL)	0.25 μL	1.25 U/rxn*																																																		
Component	50- μL rxn	Final Concentration																																																		
10 μM forward primer	1.0 μL	0.2 μM																																																		
10 μM reverse primer	1.0 μL	0.2 μM																																																		
Template DNA	varies	< 1 μg /rxn*																																																		
Step	Temperature ($^{\circ}\text{C}$)	Time																																																		
Initial Denaturation	95	10 minutes																																																		
25–35 PCR Cycles	Denature	95																																																		
	Anneal	~55 (depending on primer T_m)																																																		
	Extend	72																																																		
Final Extension	72	5 minutes																																																		
Hold	4	indefinitely																																																		
2		Prepare PCR master mix																																																		
3		Add template DNA and primers																																																		
4		Incubate reactions in a thermal cycler																																																		
5		Analyze with gel electrophoresis	<p>Analyze 10 μL using agarose gel electrophoresis. Use your PCR reaction immediately for down-stream applications, or store it at -20°C.</p>																																																	