# AmpliTaq Gold® 360 Master Mix

	Package
	Content

Catalog Number 4398876

Size 40 rxns **M** Kit Contents

4398881 200 rxns 4398886 2,000 rxns 4398901 2,000 rxns



Storage Conditions

Store all contents at -20°C.



Template: cDNA, gDNA, λDNA

Forward and reverse gene-specific primers

Autoclaved, distilled water

E-Gel® General Purpose Gels, 1.2% (Cat. no. G5018-01)

TrackIt<sup>™</sup> 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

PCR Enzymes and Master Mixes

Go online to view related products.



### Product **Description**

- AmpliTaq Gold<sup>®</sup> 360 Master Mix contains AmpliTaq Gold<sup>®</sup> 360 DNA Polymerase and optional 360 GC Enhancer.
- Activity is restored after the denaturation step in PCR cycling at 94°C, providing an automatic "hot start" and offering increased sensitivity, specificity, and yield, while allowing assembly of reactions at room temperature.
- AmpliTaq Gold® 360 Master Mix comes in a 2X format and was designed for 360° coverage of a full range of targets.



#### **Important Guidelines**

- Select the correct polymerase, PCR instrument, and cycling conditions for your application.
- Take precautions to avoid cross-contamination by using aerosol-resistant barrier tips and analyzing PCR products in a separate area from PCR assembly.
- Avoid creating bubbles when mixing the enzyme.
- If your application requires increased specificity, add 1–2 μL of 360 GC Enhancer per 50-µL reaction—2 to 5% (v/v) of the reaction.



#### **Online** Resources

Visit our product page for additional information and protocols. For support, visit www.lifetechnologies.com/support.





# **Enzyme Characteristics**

Chemical Chemical Up to 5 kb Length:

Fidelity vs. *Taq*: 1X

**Format:** Master mix

# **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	_
AmpliTaq Gold® 360 Master Mix	12.5 µL	25 μL	μL	1X
360 GC Enhancer (optional)*	0.5–5.0 μL	1.0–10 μL	μL	0–20%
10 μM forward primer	0.5 μL	1 μL	μL	0.2 μΜ
10 μM reverse primer	0.5 μL	1 μL	μL	0.2 μΜ
Template DNA	varies	varies	varies	< 1 µg

<sup>\*</sup> For targets with 65–75% GC, start with 5  $\mu$ L in a 50- $\mu$ L reaction (10% (v/v) of the reaction). For targets with > 75% GC, start with 10 µL in a 50-µL reaction (20% (v/v) of the reaction). For increased specificity, add 1-2 µL of 360 GC Enhancer per 50-µL reaction (2 to 5% (v/v) of the reaction).

#### PCR Protocol

1 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



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

# AmpliTaq Gold® 360 Master Mix Protocol

The example PCR procedure below shows appropriate volumes for a single, 70% GC-rich, 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	
1		Thaw reagents	
2		Prepare PCR master mix	
3	38	Add template DNA and primers	
4		Incubate reactions in a thermal cycler	
5	Wind.	Analyze with gel electrophoresis	

#### **Procedure Details**

Thaw, mix, and briefly centrifuge each component before use.

Avoid generating bubbles when mixing the MasterMix.

Add the following components to each PCR reaction tube.

**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.

Component	50-μL rxn	Final Concentration
Autoclaved, distilled water	to 50 µL	_
AmpliTaq Gold® 360 Master Mix	25 μL	1X
360 GC Enhancer (optional)	5.0 μL*	10%

<sup>\*</sup> Targets 65–75% GC, start with 5.0  $\mu$ L/rxn. Targets > 75% GC, start with 10  $\mu$ L/rxn. For increased specificity, add 1–2  $\mu$ L.

Cap each tube, mix, and then briefly centrifuge the contents.

Add your template DNA and primers to each tube for a final reaction volume of 50 µL.

Component	50-μL rxn	Final Concentration
10 μM forward primer	1.0 µL	0.2 μΜ
10 μM reverse primer	1.0 µL	0.2 μΜ
Template DNA	varies	< 1 µg/reaction

Cap each tube, mix, and then briefly centrifuge the contents.

S	tep	Temperature (°C)	Time
Initial D	enaturation	95°C	10 minutes 🚺
25–40	Denature	95°C	Amplicons > 2 kb: 15 seconds Amplicons ≤ 2 kb: 30 seconds
PCR Cycles	Anneal	~55°C (depending on primer $T_m$ )	30 seconds
	Extend	72°C	1 minute/kb
Final l	Extension	72°C	7 minutes
F	Hold	4°C	indefinitely

Analyze 10 µL using agarose gel electrophoresis.

Use your PCR reaction immediately for down-stream applications, or store it at -20°C.