

	Catalog No.	Size	
<b>Package contents</b>	14001-012	50 reactions	Kit contents
	14001-013	200 reactions	
	14001-014	1000 reactions	

<b>Storage conditions</b>	All the components of the kit can be stored at 4°C for periods up to 3 months. For longer storage, keep all components at -20°C.
---------------------------	--

<b>Required materials</b>	<ul style="list-style-type: none"> <li>Template: cDNA, genomic DNA, plasmid DNA, phage DNA</li> <li>Forward and reverse primers</li> <li>Invitrogen™ E-Gel™ EX Agarose Gels, 1% (Cat. No. G4010-01)</li> <li>Invitrogen™ E-Gel™ 1 kb Plus Express DNA Ladder (Cat. No. 10488-091)</li> <li>0.2 or 0.5-mL nuclease-free microcentrifuge tubes</li> </ul>
---------------------------	---

<b>Timing</b>	Varies depending on amplicon length.
---------------	--------------------------------------

<b>Product description</b>	<ul style="list-style-type: none"> <li>Invitrogen™ Platinum™ II Hot-Start Green PCR Master Mix (2X) contains Platinum™ II <i>Taq</i> Hot-Start DNA Polymerase premixed in an optimized Platinum™ II PCR buffer with dNTPs. The master mix is supplemented with tracking dyes for direct loading of PCR products on gels.</li> <li>Platinum™ II <i>Taq</i> Hot-Start DNA Polymerase is an engineered <i>Taq</i> DNA polymerase that shows increased resistance to reaction inhibitors originating from sample material or DNA purification steps.</li> <li>The polymerase activity is blocked at ambient temperatures and restored after the initial denaturation step at 94°C. This automatic “hot start” provides increased sensitivity, specificity, and yield, while allowing reaction assembly at room temperature.</li> <li>Due to unique composition of the Platinum™ II PCR buffer, the annealing temperature is 60°C for most primer pairs designed following general primer design rules.</li> <li>Platinum™ II <i>Taq</i> DNA polymerase extends 1 kb in 15 seconds. The extension step can be prolonged without a negative effect on specificity.</li> <li>The enzyme has a template independent terminal transferase activity that adds a single deoxyadenosine (A) to the 3' ends of PCR products. Like standard <i>Taq</i>, it has both 5' to 3' polymerase and 5' to 3' exonuclease activities, but lacks 3' to 5' exonuclease activity.</li> <li>The tracking dyes (a blue dye and a yellow dye) in the master mix do not interfere with PCR performance and are compatible with downstream applications such as ligation or restriction digestion.</li> </ul>
----------------------------	--

<b>Online resources</b>	Visit our <a href="#">product page</a> for additional information. Find out more at <a href="http://thermofisher.com/platinumiiataq">thermofisher.com/platinumiiataq</a> . For support, visit <a href="http://thermofisher.com/support">thermofisher.com/support</a> .
-------------------------	--

## Enzyme characteristics

<b>Hot-start:</b>	Antibody
<b>Length:</b>	Up to 5 kb
<b>Fidelity vs. <i>Taq</i>:</b>	1X
<b>Format:</b>	Master mix

## PCR setup

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

Component	20-µL rxn	50-µL rxn	Custom	Final conc.
Water, nuclease-free	to 20 µL	to 50 µL	to µL	—
Platinum™ II Hot-Start Green PCR Master Mix (2X) <sup>1</sup>	10 µL	25 µL	µL	1X
10 µM forward primer	0.4 µL	1 µL	µL	0.2 µM
10 µM reverse primer	0.4 µL	1 µL	µL	0.2 µM
Template DNA <sup>2</sup>	varies	varies	µL	<500 ng/rxn
Platinum™ GC Enhancer (optional) <sup>3</sup>	4 µL	10 µL	µL	1X

<sup>1</sup> Provides 1.5 mM MgCl<sub>2</sub> in final reaction concentration.

<sup>2</sup> 0.5–500 ng genomic DNA, 1 pg–50 ng plasmid or viral DNA, or 1–5 µL of cDNA synthesis reaction per 50-µL PCR reaction.

<sup>3</sup> Recommended for targets with >65% GC sequences.

## PCR protocol

Go to page 2 for instructions to prepare and run your PCR experiment.

## Important guidelines

Click here for important PCR guidelines.

## Optimization strategies

Click here for guidelines to optimize your PCR experiment.

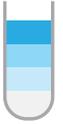
## Troubleshooting

Click here to troubleshoot your PCR experiment.

**Limited warranty, disclaimer, and licensing information**

## PCR protocol

The example procedure below shows the appropriate volumes for a single **50- $\mu$ L** reaction. For multiple reactions, prepare a master mix of components common to all reactions to minimize pipetting error, then dispense the appropriate volumes into each 0.2-mL or 0.5-mL PCR tube before adding template DNA and primers.

Steps	Action	Procedure details												
1 	<b>Thaw reagents</b>	Thaw, mix, and briefly centrifuge each component before use.												
2 	<b>Prepare PCR master mix</b>	<p>a. Add the following components to each reaction tube.</p> <p><b>Note:</b> 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>Volume for 50-<math>\mu</math>L rxn</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>Water, nuclease-free</td> <td>to 50 <math>\mu</math>L</td> <td>—</td> </tr> <tr> <td>Platinum™ II Hot-Start Green PCR Master Mix (2X)</td> <td>25 <math>\mu</math>L</td> <td>1X</td> </tr> <tr> <td>Platinum™ GC Enhancer (<i>optional</i>)<sup>1</sup></td> <td>10 <math>\mu</math>L</td> <td>1X</td> </tr> </tbody> </table> <p><sup>1</sup> Recommended for targets with &gt;65% GC sequences.</p> <p>b. Mix, then briefly centrifuge the components.</p>	Component	Volume for 50- $\mu$ L rxn	Final concentration	Water, nuclease-free	to 50 $\mu$ L	—	Platinum™ II Hot-Start Green PCR Master Mix (2X)	25 $\mu$ L	1X	Platinum™ GC Enhancer ( <i>optional</i> ) <sup>1</sup>	10 $\mu$ L	1X
Component	Volume for 50- $\mu$ L rxn	Final concentration												
Water, nuclease-free	to 50 $\mu$ L	—												
Platinum™ II Hot-Start Green PCR Master Mix (2X)	25 $\mu$ L	1X												
Platinum™ GC Enhancer ( <i>optional</i> ) <sup>1</sup>	10 $\mu$ L	1X												
3 	<b>Add template DNA and primers</b>	<p>a. Add your template DNA and primers to each tube for a final reaction volume of 50 <math>\mu</math>L.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>Volume for 50-<math>\mu</math>L rxn</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>10 <math>\mu</math>M forward primer</td> <td>1 <math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>10 <math>\mu</math>M reverse primer</td> <td>1 <math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>Template DNA</td> <td>varies</td> <td>&lt;500 ng/rxn</td> </tr> </tbody> </table> <p>b. Cap each tube, mix, then briefly centrifuge the contents.</p>	Component	Volume for 50- $\mu$ L rxn	Final concentration	10 $\mu$ M forward primer	1 $\mu$ L	0.2 $\mu$ M	10 $\mu$ M reverse primer	1 $\mu$ L	0.2 $\mu$ M	Template DNA	varies	<500 ng/rxn
Component	Volume for 50- $\mu$ L rxn	Final concentration												
10 $\mu$ M forward primer	1 $\mu$ L	0.2 $\mu$ M												
10 $\mu$ M reverse primer	1 $\mu$ L	0.2 $\mu$ M												
Template DNA	varies	<500 ng/rxn												

Steps	Action	Procedure details																															
<p>4 </p>	<p><b>Incubate reactions in a thermal cycler</b></p>	<table border="1" data-bbox="695 126 1780 435"> <thead> <tr> <th rowspan="2">Step</th> <th colspan="2">3-step protocol</th> <th colspan="2">2-step protocol<sup>1</sup></th> </tr> <tr> <th>Temperature</th> <th>Time</th> <th>Temperature</th> <th>Time</th> </tr> </thead> <tbody> <tr> <td>Initial denaturation</td> <td>94°C</td> <td>2 minutes</td> <td>94°C</td> <td>2 minutes</td> </tr> <tr> <td rowspan="3">25–35 PCR cycles</td> <td>Denature</td> <td>94°C</td> <td>98°C</td> <td>5 seconds</td> </tr> <tr> <td>Anneal<sup>2</sup></td> <td>60°C</td> <td rowspan="2">60°C</td> <td rowspan="2">15 seconds</td> </tr> <tr> <td>Extend</td> <td>68°C</td> <td>15 seconds/kb</td> </tr> <tr> <td>Hold</td> <td>4°C</td> <td>hold</td> <td>4°C</td> <td>hold</td> </tr> </tbody> </table> <p><sup>1</sup> Recommended for simple amplicons up to 1 kb with 45–65% GC sequences. For longer, GC-rich, and complex amplicons, or cDNA targets, use the 3-step cycling protocol.</p> <p><sup>2</sup> 60°C annealing temperature works for most primers. In cases when annealing temperature requires additional optimization, we recommend performing gradient PCR or redesigning the primers.</p> <p><b>Note:</b> Refer to “Optimization strategies”, page 1, for guidelines to optimize cycling conditions.</p>	Step	3-step protocol		2-step protocol <sup>1</sup>		Temperature	Time	Temperature	Time	Initial denaturation	94°C	2 minutes	94°C	2 minutes	25–35 PCR cycles	Denature	94°C	98°C	5 seconds	Anneal <sup>2</sup>	60°C	60°C	15 seconds	Extend	68°C	15 seconds/kb	Hold	4°C	hold	4°C	hold
Step	3-step protocol			2-step protocol <sup>1</sup>																													
	Temperature	Time	Temperature	Time																													
Initial denaturation	94°C	2 minutes	94°C	2 minutes																													
25–35 PCR cycles	Denature	94°C	98°C	5 seconds																													
	Anneal <sup>2</sup>	60°C	60°C	15 seconds																													
	Extend	68°C			15 seconds/kb																												
Hold	4°C	hold	4°C	hold																													
<p>5 </p>	<p><b>Analyze with gel electrophoresis</b></p>	<p>a. Analyze the sample using agarose gel electrophoresis.</p> <p><b>Note:</b> Dilute the PCR sample 2- to 20-fold for optimal separation on E-Gel™ agarose gels.</p> <p>b. Use your PCR product immediately in down-stream applications, or store it at –20°C.</p>																															