



**Contents and storage**

Catalog No.	Amount	Storage
15966005	500 U	-20°C
15966025	5 × 500 U	-20°C

Kit contents



**Product description**

- The Invitrogen™ Platinum™ Taq DNA Polymerase, DNA-free, is an ideal choice for PCR-based applications requiring the highest sensitivity without false-positive results from reagent-borne contamination.
- Platinum™ Taq DNA Polymerase, DNA-free, is manufactured using closed and single-use system technology to minimize DNA contamination risk. It contains ≤0.01 genome equivalent of bacterial DNA and ≤0.001 copy of human genomic DNA per enzyme unit. Go to [thermofisher.com/dna-free](http://thermofisher.com/dna-free) for additional information.
- The enzyme is a recombinant Taq polymerase complexed with a proprietary antibody that blocks polymerase activity at ambient temperatures.
- Activity is restored after the initial 94°C denaturation step during PCR, providing an automatic “hot start” and offering increased sensitivity, specificity, and yield, while allowing reaction assembly at room temperature.



**Required materials**

- Template: cDNA, gDNA
  - Forward and reverse gene-specific primers
  - 10 mM dNTP mix (Cat. No. 18427-088)
  - Water, nuclease-free
  - 0.2-mL or 0.5-mL nuclease-free microcentrifuge tubes
- for PCR:*
- E-Gel™ General Purpose Gels, 1.2% (Cat. No. G5018-01)
  - TrackIt™ 1 kb Plus DNA Ladder (Cat. No. 10488-085)
- for qPCR:*
- 100 μM qPCR probe
  - ROX Reference Dye



**Online resources**

- Visit our [product page](#) for protocols, safety, and additional product information.
- For support, visit [thermofisher.com/support](http://thermofisher.com/support).



**Important guidelines**

**IMPORTANT:** Always maintain a DNA-free environment when handling vials with polymerase and other PCR reagents to prevent DNA contamination.

[Click here for important PCR guidelines.](#)

**Enzyme characteristics**

<b>Hot-start:</b>	Antibody
<b>Amplicon size:</b>	Up to 4 kb
<b>Fidelity vs. Taq:</b>	1X
<b>Product overhang:</b>	3' A
<b>Exonuclease activity:</b>	5' → 3'

**PCR setup**

Use the following volumes to prepare your PCR reaction, or enter your own parameters in the column provided. For qPCR set-up, see page 2.

Component	25-μL rxn	50-μL rxn	Custom	Final conc. in 25-μL rxn
Water, nuclease-free	to 25 μL	to 50 μL	to.....μL	—
10X PCR Buffer ( - MgCl <sub>2</sub> ), DNA-free	2.5 μL	5 μL	μL	1X
50 mM MgCl <sub>2</sub> , DNA-free	0.75 μL	1.5 μL	μL	1.5 mM
10 mM dNTP mix	0.5 μL	1 μL	μL	0.2 mM each
10 μM forward primer	0.5 μL	1 μL	μL	0.2 μM
10 μM reverse primer	0.5 μL	1 μL	μL	0.2 μM
Template DNA	varies	varies	μL	≤500 ng/rxn
Platinum™ Taq DNA Polymerase, DNA-free (5 U/μL)	0.25 μL	0.5 μL	μL	1.25 U/rxn




**Optimization strategies**



[Click here for guidelines to optimize your PCR experiment.](#)

**Purchaser notification**

[Click here for Limited Warranty, Disclaimer, and Licensing information.](#)

The example procedure below shows appropriate volumes for a single **25- $\mu$ L** 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 tube or well of a MicroAmp™ EnduraPlate™ Optical 96- or 384-well plate prior to adding template DNA and primers. For 384-well plates, we recommend a maximum reaction volume of 10  $\mu$ L per well.

Step	Action	Procedure details																																								
1 	<b>Thaw reagents</b>	<p>a. UV irradiate your work area prior to beginning work to ensure a DNA-free environment.</p> <p>b. Thaw, mix, and briefly centrifuge each component before use.</p>																																								
2 	<b>Prepare PCR master mix</b>	<p>a. Add the following components to each PCR 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>25-<math>\mu</math>L rxn</th> <th>Custom</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>Water, nuclease-free</td> <td>to 25 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>—</td> </tr> <tr> <td>10X PCR Buffer (– MgCl<sub>2</sub>), DNA-free</td> <td>2.5 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>1X</td> </tr> <tr> <td>50 mM MgCl<sub>2</sub>, DNA-free</td> <td>0.75 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>1.5 mM</td> </tr> <tr> <td>10 mM dNTP mix</td> <td>0.5 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.2 mM each</td> </tr> <tr> <td>Platinum™ Taq DNA Polymerase, DNA-free (5 U/<math>\mu</math>L)</td> <td>0.25 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>1.25 U/rxn</td> </tr> </tbody> </table> <p>b. Mix and then briefly centrifuge the components.</p>	Component	25- $\mu$ L rxn	Custom	Final conc.	Water, nuclease-free	to 25 $\mu$ L	$\mu$ L	—	10X PCR Buffer (– MgCl <sub>2</sub> ), DNA-free	2.5 $\mu$ L	$\mu$ L	1X	50 mM MgCl <sub>2</sub> , DNA-free	0.75 $\mu$ L	$\mu$ L	1.5 mM	10 mM dNTP mix	0.5 $\mu$ L	$\mu$ L	0.2 mM each	Platinum™ Taq DNA Polymerase, DNA-free (5 U/ $\mu$ L)	0.25 $\mu$ L	$\mu$ L	1.25 U/rxn																
Component	25- $\mu$ L rxn	Custom	Final conc.																																							
Water, nuclease-free	to 25 $\mu$ L	$\mu$ L	—																																							
10X PCR Buffer (– MgCl <sub>2</sub> ), DNA-free	2.5 $\mu$ L	$\mu$ L	1X																																							
50 mM MgCl <sub>2</sub> , DNA-free	0.75 $\mu$ L	$\mu$ L	1.5 mM																																							
10 mM dNTP mix	0.5 $\mu$ L	$\mu$ L	0.2 mM each																																							
Platinum™ Taq DNA Polymerase, DNA-free (5 U/ $\mu$ L)	0.25 $\mu$ L	$\mu$ L	1.25 U/rxn																																							
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 25 <math>\mu</math>L.</p> <p><b>For PCR<sup>1</sup>:</b></p> <table border="1"> <thead> <tr> <th>Component</th> <th>25-<math>\mu</math>L rxn</th> <th>Custom</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>10 <math>\mu</math>M forward gene-specific primer</td> <td>0.5 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>10 <math>\mu</math>M reverse gene-specific primer</td> <td>0.5 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>Template DNA</td> <td>varies</td> <td><math>\mu</math>L</td> <td><math>\leq</math>500 ng/rxn (human gDNA)</td> </tr> </tbody> </table> <p><sup>1</sup> See “Optimization strategies”, page 1.</p> <p><b>For qPCR<sup>1</sup>:</b></p> <table border="1"> <thead> <tr> <th>Component</th> <th>25-<math>\mu</math>L rxn</th> <th>Custom</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>10 <math>\mu</math>M forward gene-specific primer</td> <td>0.75 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.3 <math>\mu</math>M</td> </tr> <tr> <td>10 <math>\mu</math>M reverse gene-specific primer</td> <td>0.75 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.3 <math>\mu</math>M</td> </tr> <tr> <td>100 <math>\mu</math>M qPCR probe</td> <td>0.05 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>30 <math>\mu</math>M ROX Reference Dye</td> <td>0.025 <math>\mu</math>L</td> <td><math>\mu</math>L</td> <td>30 nM<sup>2</sup></td> </tr> <tr> <td>Template DNA</td> <td>varies</td> <td><math>\mu</math>L</td> <td><math>\leq</math>500 ng/rxn (human gDNA)</td> </tr> </tbody> </table> <p><sup>1</sup> See “Optimization strategies”, page 1.</p> <p><sup>2</sup> The recommended final ROX concentration depends on the instrument (see “Important guidelines”, page 1).</p> <p>b. Cap each tube, mix, and then briefly centrifuge the contents.</p>	Component	25- $\mu$ L rxn	Custom	Final conc.	10 $\mu$ M forward gene-specific primer	0.5 $\mu$ L	$\mu$ L	0.2 $\mu$ M	10 $\mu$ M reverse gene-specific primer	0.5 $\mu$ L	$\mu$ L	0.2 $\mu$ M	Template DNA	varies	$\mu$ L	$\leq$ 500 ng/rxn (human gDNA)	Component	25- $\mu$ L rxn	Custom	Final conc.	10 $\mu$ M forward gene-specific primer	0.75 $\mu$ L	$\mu$ L	0.3 $\mu$ M	10 $\mu$ M reverse gene-specific primer	0.75 $\mu$ L	$\mu$ L	0.3 $\mu$ M	100 $\mu$ M qPCR probe	0.05 $\mu$ L	$\mu$ L	0.2 $\mu$ M	30 $\mu$ M ROX Reference Dye	0.025 $\mu$ L	$\mu$ L	30 nM <sup>2</sup>	Template DNA	varies	$\mu$ L	$\leq$ 500 ng/rxn (human gDNA)
Component	25- $\mu$ L rxn	Custom	Final conc.																																							
10 $\mu$ M forward gene-specific primer	0.5 $\mu$ L	$\mu$ L	0.2 $\mu$ M																																							
10 $\mu$ M reverse gene-specific primer	0.5 $\mu$ L	$\mu$ L	0.2 $\mu$ M																																							
Template DNA	varies	$\mu$ L	$\leq$ 500 ng/rxn (human gDNA)																																							
Component	25- $\mu$ L rxn	Custom	Final conc.																																							
10 $\mu$ M forward gene-specific primer	0.75 $\mu$ L	$\mu$ L	0.3 $\mu$ M																																							
10 $\mu$ M reverse gene-specific primer	0.75 $\mu$ L	$\mu$ L	0.3 $\mu$ M																																							
100 $\mu$ M qPCR probe	0.05 $\mu$ L	$\mu$ L	0.2 $\mu$ M																																							
30 $\mu$ M ROX Reference Dye	0.025 $\mu$ L	$\mu$ L	30 nM <sup>2</sup>																																							
Template DNA	varies	$\mu$ L	$\leq$ 500 ng/rxn (human gDNA)																																							

Step	Action	Procedure details																																								
<p data-bbox="73 544 94 576">4</p> 	<p data-bbox="304 503 556 609"><b>Incubate reactions in a thermal cycler</b></p>	<p data-bbox="598 105 903 138"><b>PCR cycling parameters:</b></p> <table border="1" data-bbox="598 138 1764 414"> <thead> <tr> <th data-bbox="598 138 840 186">Step</th> <th data-bbox="840 138 1501 186">Temperature</th> <th data-bbox="1501 138 1764 186">Time</th> </tr> </thead> <tbody> <tr> <td data-bbox="598 186 840 235">Initial denaturation</td> <td data-bbox="840 186 1501 235">94°C</td> <td data-bbox="1501 186 1764 235">2 minutes</td> </tr> <tr> <td data-bbox="598 235 840 365" rowspan="3">25–40 PCR cycles</td> <td data-bbox="840 235 1501 284">Denature</td> <td data-bbox="1501 235 1764 284">94°C</td> </tr> <tr> <td data-bbox="840 284 1501 332">Anneal</td> <td data-bbox="1501 284 1764 332">~60°C (depending on primer T<sub>m</sub>)</td> </tr> <tr> <td data-bbox="840 332 1501 365">Extend</td> <td data-bbox="1501 332 1764 365">72°C</td> </tr> <tr> <td data-bbox="598 365 840 414">Hold</td> <td data-bbox="840 365 1501 414">4°C</td> <td data-bbox="1501 365 1764 414">hold</td> </tr> </tbody> </table> <p data-bbox="598 430 1239 462"><b>qPCR cycling parameters (Two-step cycling protocol):</b></p> <table border="1" data-bbox="598 462 1764 641"> <thead> <tr> <th data-bbox="598 462 840 511">Step</th> <th data-bbox="840 462 1501 511">Temperature</th> <th data-bbox="1501 462 1764 511">Time</th> </tr> </thead> <tbody> <tr> <td data-bbox="598 511 840 560">Initial denaturation</td> <td data-bbox="840 511 1501 560">94°C</td> <td data-bbox="1501 511 1764 560">2 minutes</td> </tr> <tr> <td data-bbox="598 560 840 641" rowspan="2">25–40 PCR cycles</td> <td data-bbox="840 560 1501 609">Denature</td> <td data-bbox="1501 560 1764 609">94°C</td> </tr> <tr> <td data-bbox="840 609 1501 641">Anneal/Extend<sup>1</sup></td> <td data-bbox="1501 609 1764 641">~60°C (depending on primer T<sub>m</sub>)</td> </tr> </tbody> </table> <p data-bbox="598 657 1659 690"><sup>1</sup> Data acquisition should be performed during the annealing/extension step for probe-based assays.</p> <p data-bbox="598 706 1260 738"><b>qPCR cycling parameters (Three-step cycling protocol):</b></p> <table border="1" data-bbox="598 738 1764 966"> <thead> <tr> <th data-bbox="598 738 840 787">Step</th> <th data-bbox="840 738 1501 787">Temperature</th> <th data-bbox="1501 738 1764 787">Time</th> </tr> </thead> <tbody> <tr> <td data-bbox="598 787 840 836">Initial denaturation</td> <td data-bbox="840 787 1501 836">94°C</td> <td data-bbox="1501 787 1764 836">2 minutes</td> </tr> <tr> <td data-bbox="598 836 840 966" rowspan="3">25–40 PCR cycles</td> <td data-bbox="840 836 1501 885">Denature</td> <td data-bbox="1501 836 1764 885">94°C</td> </tr> <tr> <td data-bbox="840 885 1501 933">Anneal</td> <td data-bbox="1501 885 1764 933">~60°C (depending on primer T<sub>m</sub>)</td> </tr> <tr> <td data-bbox="840 933 1501 966">Extend<sup>1</sup></td> <td data-bbox="1501 933 1764 966">72°C</td> </tr> </tbody> </table> <p data-bbox="598 982 1543 1015"><sup>1</sup> Data acquisition should be performed during the extension step for probe-based assays.</p>	Step	Temperature	Time	Initial denaturation	94°C	2 minutes	25–40 PCR cycles	Denature	94°C	Anneal	~60°C (depending on primer T <sub>m</sub> )	Extend	72°C	Hold	4°C	hold	Step	Temperature	Time	Initial denaturation	94°C	2 minutes	25–40 PCR cycles	Denature	94°C	Anneal/Extend <sup>1</sup>	~60°C (depending on primer T <sub>m</sub> )	Step	Temperature	Time	Initial denaturation	94°C	2 minutes	25–40 PCR cycles	Denature	94°C	Anneal	~60°C (depending on primer T <sub>m</sub> )	Extend <sup>1</sup>	72°C
Step	Temperature	Time																																								
Initial denaturation	94°C	2 minutes																																								
25–40 PCR cycles	Denature	94°C																																								
	Anneal	~60°C (depending on primer T <sub>m</sub> )																																								
	Extend	72°C																																								
Hold	4°C	hold																																								
Step	Temperature	Time																																								
Initial denaturation	94°C	2 minutes																																								
25–40 PCR cycles	Denature	94°C																																								
	Anneal/Extend <sup>1</sup>	~60°C (depending on primer T <sub>m</sub> )																																								
Step	Temperature	Time																																								
Initial denaturation	94°C	2 minutes																																								
25–40 PCR cycles	Denature	94°C																																								
	Anneal	~60°C (depending on primer T <sub>m</sub> )																																								
	Extend <sup>1</sup>	72°C																																								
<p data-bbox="73 1104 94 1136">5</p> 	<p data-bbox="315 1104 546 1136"><b>Analyze results</b></p>	<ol data-bbox="598 1047 2026 1193" style="list-style-type: none"> <li>Check the specificity of the PCR/qPCR products by agarose gel electrophoresis. Before loading, add gel loading buffer to 10 µL of the PCR/qPCR sample, mix, and briefly centrifuge the contents.</li> <li>Analyze qPCR results following your real-time instrument manufacturer’s guidelines.</li> <li>You can store your samples overnight at 2–8°C, or at –20°C for longer period.</li> </ol>																																								