

<b>Package contents</b>	<b>Sample Kit No.</b> EP204SMP	<b>Size</b> 1 kU	Kit contents
-------------------------	-----------------------------------	---------------------	--------------

<b>Storage conditions</b>	Store all contents at $-20^{\circ}\text{C}$ . Product is designed to withstand at least 10 freeze-thaw cycles.		
---------------------------	--	--	--

<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>▪ qPCR probe</li> <li>▪ Water, nuclease-free</li> <li>▪ 0.2-mL 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™ Lyo-ready Platinum™ II Taq Hot-Start DNA Polymerase is an engineered Taq DNA polymerase that shows increased resistance to reaction inhibitors originating from sample material or DNA purification steps. The lyo-ready enzyme formulation combines feasibility to lyophilize, while retaining all favourable standard enzyme (with glycerol) properties.</li> <li>▪ The polymerase activity is blocked at ambient temperatures and restored after the initial denaturation step at <math>95^{\circ}\text{C}</math>. This automatic “hot start” provides increased sensitivity, specificity, and yield, while allowing reaction assembly at room temperature.</li> <li>▪ Lyo-ready Platinum™ II Taq Hot-Start 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 Taq, it has both 5' to 3' polymerase and 5' to 3' exonuclease activities, but lacks 3' to 5' exonuclease activity.</li> </ul>
----------------------------	---

<b>Online resources</b>	Find out more at <a href="#">Lyo-ready PCR Enzymes</a> . For further information, contact <a href="mailto:LCSVilnius@thermofisher.com">LCSVilnius@thermofisher.com</a> .
-------------------------	---

## Enzyme characteristics

<b>Hot-start:</b>	Antibody
<b>Fidelity vs. Taq:</b>	1X
<b>Format:</b>	Separate components

## Experiment setup

Use the following amounts to prepare your qPCR experiment. Note that the table shows the appropriate volumes for a single **25- $\mu\text{L}$**  reaction.

Component	Volume	Final conc.
Water, nuclease-free	to 25 $\mu\text{L}$	—
5X Lyo-ready Platinum™ II PCR Buffer 2, w/o $\text{MgCl}_2$	5 $\mu\text{L}$	1X
10 mM dNTP mix	0.5 $\mu\text{L}$	0.2 mM each
$\text{MgCl}_2$ (50 mM)	0.75 $\mu\text{L}$	1.5 mM
10 $\mu\text{M}$ forward primer	0.75 $\mu\text{L}$	0.3 $\mu\text{M}$
10 $\mu\text{M}$ reverse primer	0.75 $\mu\text{L}$	0.3 $\mu\text{M}$
10 $\mu\text{M}$ primer probe	0.5 $\mu\text{L}$	0.2 $\mu\text{M}$
(Optional) ROX Reference Dye <sup>1</sup>	varies	varies
Template DNA <sup>1</sup>	varies	$\leq 500$ ng/rxn
Lyo-ready Platinum™ II Taq Hot-Start DNA Polymerase (5 U/ $\mu\text{L}$ )	0.2 $\mu\text{L}$	1 U/rxn

<sup>1</sup> See “Optimization strategies”, below.

## Protocol





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

## Optimization strategies

Click here for guidelines to optimize your qPCR experiment.

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

The following example procedure shows the appropriate volumes for a single **25- $\mu$ L** qPCR 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 or well of a MicroAmp™ EnduraPlate™ Optical 96-well Fast Plate 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 25-<math>\mu</math>L rxn</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>Water, nuclease-free</td> <td>to 25 <math>\mu</math>L</td> <td>—</td> </tr> <tr> <td>5X Lyo-ready Platinum™ II PCR Buffer 2, w/o MgCl<sub>2</sub></td> <td>5 <math>\mu</math>L</td> <td>1X</td> </tr> <tr> <td>10 mM dNTP mix (10 mM each)</td> <td>0.5 <math>\mu</math>L</td> <td>0.2 mM each</td> </tr> <tr> <td>MgCl<sub>2</sub> (50 mM)</td> <td>0.75 <math>\mu</math>L</td> <td>1.5 mM</td> </tr> <tr> <td>Lyo-ready Platinum™ II Taq Hot-Start DNA Polymerase</td> <td>0.2 <math>\mu</math>L</td> <td>0.04 U/<math>\mu</math>L</td> </tr> </tbody> </table> <p>b. Mix, then briefly centrifuge the components.</p>	Component	Volume for 25- $\mu$ L rxn	Final concentration	Water, nuclease-free	to 25 $\mu$ L	—	5X Lyo-ready Platinum™ II PCR Buffer 2, w/o MgCl <sub>2</sub>	5 $\mu$ L	1X	10 mM dNTP mix (10 mM each)	0.5 $\mu$ L	0.2 mM each	MgCl <sub>2</sub> (50 mM)	0.75 $\mu$ L	1.5 mM	Lyo-ready Platinum™ II Taq Hot-Start DNA Polymerase	0.2 $\mu$ L	0.04 U/ $\mu$ L
Component	Volume for 25- $\mu$ L rxn	Final concentration																		
Water, nuclease-free	to 25 $\mu$ L	—																		
5X Lyo-ready Platinum™ II PCR Buffer 2, w/o MgCl <sub>2</sub>	5 $\mu$ L	1X																		
10 mM dNTP mix (10 mM each)	0.5 $\mu$ L	0.2 mM each																		
MgCl <sub>2</sub> (50 mM)	0.75 $\mu$ L	1.5 mM																		
Lyo-ready Platinum™ II Taq Hot-Start DNA Polymerase	0.2 $\mu$ L	0.04 U/ $\mu$ L																		
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> <table border="1"> <thead> <tr> <th>Component</th> <th>Volume for 25-<math>\mu</math>L rxn</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>10 <math>\mu</math>M forward primer</td> <td>0.75 <math>\mu</math>L</td> <td>0.3 <math>\mu</math>M</td> </tr> <tr> <td>10 <math>\mu</math>M reverse primer</td> <td>0.75 <math>\mu</math>L</td> <td>0.3 <math>\mu</math>M</td> </tr> <tr> <td>10 <math>\mu</math>M primer probe</td> <td>0.5 <math>\mu</math>L</td> <td>0.2 <math>\mu</math>M</td> </tr> <tr> <td>(Optional) ROX Reference Dye</td> <td>varies</td> <td>varies</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 25- $\mu$ L rxn	Final concentration	10 $\mu$ M forward primer	0.75 $\mu$ L	0.3 $\mu$ M	10 $\mu$ M reverse primer	0.75 $\mu$ L	0.3 $\mu$ M	10 $\mu$ M primer probe	0.5 $\mu$ L	0.2 $\mu$ M	(Optional) ROX Reference Dye	varies	varies	Template DNA	varies	<500 ng/rxn
Component	Volume for 25- $\mu$ L rxn	Final concentration																		
10 $\mu$ M forward primer	0.75 $\mu$ L	0.3 $\mu$ M																		
10 $\mu$ M reverse primer	0.75 $\mu$ L	0.3 $\mu$ M																		
10 $\mu$ M primer probe	0.5 $\mu$ L	0.2 $\mu$ M																		
(Optional) ROX Reference Dye	varies	varies																		
Template DNA	varies	<500 ng/rxn																		
4 	<b>Incubate reactions in a thermal cycler</b>	<table border="1"> <thead> <tr> <th>Step</th> <th>Temperature</th> <th>Time</th> </tr> </thead> <tbody> <tr> <td>Initial denaturation</td> <td>95°C</td> <td>2 minutes</td> </tr> <tr> <td rowspan="2">40 cycles</td> <td>Denature</td> <td>95°C</td> </tr> <tr> <td>Anneal/Extend<sup>1</sup></td> <td>60°C</td> </tr> <tr> <td></td> <td></td> <td>15 seconds</td> </tr> </tbody> </table> <p><sup>1</sup> Data acquisition should be performed during the annealing/extension step for probe-based assays.</p> <p><b>Note:</b> Refer to “Optimization strategies”, page 1, for guidelines to optimize cycling conditions.</p>	Step	Temperature	Time	Initial denaturation	95°C	2 minutes	40 cycles	Denature	95°C	Anneal/Extend <sup>1</sup>	60°C			15 seconds				
Step	Temperature	Time																		
Initial denaturation	95°C	2 minutes																		
40 cycles	Denature	95°C																		
	Anneal/Extend <sup>1</sup>	60°C																		
		15 seconds																		