










# Platinum® *Taq* PCR<sub>x</sub> DNA Polymerase

invitrogen<sup>®</sup>  
by life technologies™

	<b>Package Contents</b>	Catalog Number 11509-015	Size 500 units	 Kit Contents
	<b>Storage Conditions</b>	<ul style="list-style-type: none"> <li>Store all contents at -20°C.</li> </ul>		
	<b>Required Materials</b>	<ul style="list-style-type: none"> <li>Template: cDNA, gDNA, λDNA</li> <li>Forward and reverse gene-specific primers</li> <li>10 mM dNTP mix (Cat. no. 18427-088)</li> <li>Autoclaved, distilled water</li> <li>E-Gel® General Purpose Gels, 1.2% (Cat. no. G5018-01)</li> <li>TrackIt™ 1 Kb Plus DNA Ladder (Cat. no. 10488-085)</li> <li>0.2 or 0.5-mL nuclease-free microcentrifuge tubes</li> </ul>		
	<b>Timing</b>	Varies depending on amplicon length		
	<b>Selection Guide</b>	<a href="#">PCR Enzymes and Master Mixes</a> Go online to view related products.		
	<b>Product Description</b>	<ul style="list-style-type: none"> <li>Platinum® <i>Taq</i> PCR<sub>x</sub> DNA Polymerase is recombinant <i>Taq</i>, complexed with a proprietary antibody that blocks polymerase activity at ambient temperatures, providing an automatic “hot-start” upon initial denaturation.</li> <li>An optimized buffer system is also supplied, including 10X PCR<sub>x</sub> Amplification Buffer, 50 mM MgSO<sub>4</sub>, and 10X PCR<sub>x</sub> Enhancer Solution, a novel PCR cosolvent that simplifies amplification of problematic and/or GC-rich templates.</li> <li>Take precautions to avoid cross-contamination by using aerosol-resistant barrier tips and analyzing PCR products in a separate area from PCR assembly.</li> <li>Two buffers are provided: The standard 10X PCR Buffer and 50 mM MgCl<sub>2</sub> are recommended for PCR of routine templates (30–50% GC content). Substitute standard components with 10X PCR<sub>x</sub> Amplification Buffer and MgSO<sub>4</sub> for more robust amplification.</li> <li>Use 10X PCR<sub>x</sub> Enhancer with 10X PCR<sub>x</sub> Amplification Buffer and MgSO<sub>4</sub> for GC-rich templates to widen reaction optima and increase PCR success rates.</li> <li>The optimal concentration of 10X PCR<sub>x</sub> Enhancer Solution will vary depending on GC content, Mg<sup>++</sup> concentration, and annealing temperature.</li> </ul>		
	<b>Important Guidelines</b>			
	<b>Online Resources</b>	Visit our <a href="#">product page</a> for additional information and protocols. For support, visit <a href="http://www.lifetechnologies.com/support">www.lifetechnologies.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>	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, Minus Mg	2.5 μL	5 μL	μL	1X
50 mM MgCl <sub>2</sub>	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		< 500 ng
Platinum® <i>Taq</i> DNA Polymerase (5 U/μL)	0.25 μL	0.5 μL	μL	2.5 U

- Incubate in a thermal cycler at 94°C for 2 minutes to denature the template and activate the enzyme.
- Perform 25 to 35 cycles of PCR amplification:
  - Denature—94°C for 15 seconds
  - Anneal—~55°C depending on Primer T<sub>m</sub> for 30 seconds
  - Extend—72°C for 1 minute/kb

## Optimization PCR Protocol for Problematic/GC-Rich

-  See page 2 to view a procedure for optimizing your PCR experiment by testing multiple concentrations of the PCR<sub>x</sub> Enhancer Solution in your reactions.

## Optimization Strategies


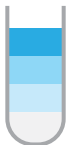



-  Refer to the pop-up for guidelines to optimize your PCR reactions.

 **Limited Warranty, Disclaimer, and Licensing Information**

life  
technologies™

## Platinum® Taq PCR<sub>x</sub> DNA Polymerase Protocol

The example PCR procedure below, for GC-rich or problematic templates, is designed to test six varying concentrations (0X to 4X) of PCR<sub>x</sub> Enhancer Solution by preparing enough master mix for seven 50-μL reactions. You can also prepare individual PCR mixtures according to the guidelines below by selecting a final concentration of PCR<sub>x</sub> Enhancer Solution to use in your reaction set-up.

Timeline	Steps
1 	Thaw reagents
2 	Prepare PCR master mix
3 	Add enhancer solution
4 	Incubate reactions in a thermal cycler
5 	Analyze with gel electrophoresis

Procedure Details							
Thaw, mix, and briefly centrifuge each component before use.							
Add the following components to each PCR tube. Mix and briefly centrifuge the contents.							
<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.							
Component		Seven-Rxn Master Mix	One 50-μL Rxn	Final Concentration			
Autoclaved, distilled water		to 210 μL	to 30 μL				
10X PCR <sub>x</sub> Amplification Buffer		35 μL	5 μL	1X			
10 mM dNTP Mixture		7 μL	1 μL	0.2 mM each			
50 mM MgSO <sub>4</sub>		10.5 μL	1.5 μL	1.5 mM			
10 μM Forward Primer		7 μL	1 μL	0.2 μM			
10 μM Reverse Primer		7 μL	1 μL	0.2 μM			
Template DNA		≥ 7 μL	≥ 1 μL	< 500 ng			
Platinum® Taq DNA Polymerase (5 U/μL)		3.5 μL	0.5 μL	2.5 U/rxn			
Add the following reagents to <b>six</b> microcentrifuge tubes:							
Component		PCR <sub>x</sub> Enhancer Concentration					
		0X	0.5X	1X	2X	3X	4X
Master Mix from step 2		30	30	30	30	30	30
10X PCR <sub>x</sub> Enhancer Solution		-	2.5	5	10	15	20
Autoclaved, distilled water		20	17.5	15	10	5	-
Cap each tube, mix, and then briefly centrifuge the contents.							
Step		Temperature [°C]			Time		
Initial Denaturation		95			2 minutes		
25–35 PCR Cycles	Denature	95			45 seconds		
	Anneal	~55 (depending on primer T <sub>m</sub> )			30 seconds		
	Extend	68			1 minute/kb		
Hold		4			indefinitely		
Analyze 10 μL using agarose gel electrophoresis.							
Use your PCR reaction immediately for down-stream applications, or store it at -20°C.							