Protocol Pub. No. MAN0000948 Rev. A.0

Platinum[®] *Taq* DNA Polymerase High Fidelity

S	Package Contents	Catalog Number 11304-011 11304-029 11304-102	Size 100 rxns 500 rxns 5,000 rxns	i Kit Contents			
	Storage Conditions	 Store all contents at -20°C. 					
	Required Materials	 Template: cDNA, gDNA, λDNA Forward and reverse gene-specific primers 10 mM dNTP mix (Cat. no. 18427-088) Autoclaved, distilled water E-Gel[®] General Purpose Gels, 1.2% (Cat. no. G5018-01) TrackIt[™] 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					
R	Selection Guide	PCR Enzymes and Master Mixes Go online to view related products.					
¢	Product Description	 Platinum[®] <i>Taq</i> DNA Polymerase High Fidelity contains recombinant <i>Taq</i> DNA polymerase, <i>Pyrococcus species</i> GB-D polymerase, and Platinum[®] <i>Taq</i> Antibody. This enzyme allows amplification of simple and complex DNA templates over a large range of target sizes and provides 6X higher fidelity over <i>Taq</i>. 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. 					
	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. For gDNA or cDNA, use a primer concentration of 0.2 μM. For Plasmid or λDNA, increase to 0.4 μM. Do not perform the initial denaturation for more than 30 seconds if the target is greater than 12 kb. 					
	Online Resources	Visit our product pa information and provisit www.lifetechn	age for additional otocols. For suppor ologies.com/supp	rt, P			

Enzyme Characteristics

Hot-start:	Antibody
Length:	Up to 20 kb
Fidelity vs. <i>Taq</i> :	6X
Format:	Separate components
	1 1

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	5-μL rxn 50-μL rxn		Final Conc.	
Autoclaved, distilled water	to 25 µL	to 50 µL	to µL	-	
10X High Fidelity PCR Buffer	2.5 µL	5 µL	μL	1X	
50 mM MgSO_4	1 µL	2 µL	μL	2.0 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 High Fidelity (5 U/µL)	0.1 µL	0.2 μL	μL	1 U/rxn	

PCR Protocol

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



invitrogen

by *life* technologies"

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

Platinum[®] Taq DNA Polymerase High Fidelity Protocol

The example PCR procedure below shows appropriate volumes for a single **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	Procedure Details					
1		Thaw reagents	Thaw, mix, and briefly centrifuge each component before use					
		Prepare PCR master mix	Add the following components to each PCR reaction tube.					
			correct amount of water required to reach your final reaction volume.					
			Component 50-µL rxn Final Concentration					
			Autoclaved,	Autoclaved, distilled water		to 50 µL		
0			10X High Fic	10X High Fidelity PCR Buffer		5 µL	1X	
2			10 mM dNTI	10 mM dNTP mix		1 µL	0.2 mM each	
			50 mM MgS0	50 mM MgSO ₄		2 µL	2 mM	
			Platinum [®] Taq	DNA High Fide	lity Polymerase (5 U/µL)	0.2 µL	1 U*	
			* 1 U is sufficient for amplifying most targets. In some cases, more enzyme may be required (up to 2.5 units). Use 1 U of enzyme for targets > 12 kb.					
			Mix and briefly centrifuge the components.					
		Add template DNA and primers	Add your template DNA and primers to each tube for a final reaction volume of 50 μ L.					
			Component 50-µL rxn Final Concentration				Concentration	
3	200		10 µM forwa	10 μM forward primer 1–2 μL		0.2 μM (Ge 0.4 μM (Pl	0.2 μ M (Genomic and cDNA)– 0.4 μ M (Plasmid and λ DNA)	
			10 µM revers	10 µM reverse primer 1–2		0.2 μ M (Genomic and cDNA)– 0.4 μ M (Plasmid and λ DNA)		
			Template DNA varies		< 500 ng			
			Cap each tube, mix, and then briefly centrifuge the contents.					
		Incubate reactions in a thermal cycler	St	ep	Temperature (°C)	Time	
			Initial Denaturation		94	30	seconds-2 minutes*	
			DE DE	Denature	94		15 seconds	
4			25–35 PCR	Anneal	~55 (depending on primer T _m)		30 seconds	
	Y		Cycles	Extend	68	1 minute/kb		
			H	Hold 4			indefinitely	
			* For targets > 12 kb, do not exceed 30 seconds for initial denaturation.					
	KARAA .	Analyze with gel	Analyze 10 µL using agarose gel electrophoresis.					
5		electrophoresis	Use your PCR reaction immediately for down-stream applications, or store it at -20°C.					
			For support, visit www.lifetechnologies.com/support					