

	Package contents	Catalog number 12351-010 12351-050 12351-250	Size 100 Units 500 Units 5 × 500 Units	Kit contents
	Storage conditions	<ul style="list-style-type: none"> Store all contents at -20°C. Template: gDNA, plasmid DNA, phage DNA, cDNA Forward and reverse gene-specific primers Invitrogen™ 10 mM dNTP Mix (Cat. No. 18427-088) Invitrogen™ E-Gel™ General Purpose Gels, 1.2% (Cat. No. G5018-01) Invitrogen™ TrackIt™ 1 kb Plus DNA Ladder (Cat. No. 10488-085) 0.2 or 0.5-mL nuclease-free microcentrifuge tubes Gel loading buffer Water, nuclease-free 		
	Required materials	<ul style="list-style-type: none"> 0.2 or 0.5-mL nuclease-free microcentrifuge tubes Gel loading buffer Water, nuclease-free 		
	Timing	Varies depending on amplicon length.		
	Selection guide	PCR Enzymes and Master Mixes Go online to view related products.		
	Product description	<ul style="list-style-type: none"> Platinum™ SuperFi™ DNA Polymerase is a proofreading DNA polymerase that combines fidelity with Platinum™ hot-start technology, and is ideally suited for cloning, mutagenesis, and other applications. Platinum™ hot-start technology inhibits DNA polymerase activity at ambient temperatures, allowing room temperature reaction setup and storage of pre-assembled PCR reactions for up to 24 hours prior to the PCR. Enzyme activity is restored after the initial denaturation step. Platinum™ SuperFi™ DNA Polymerase has 5' to 3' polymerase and 3' to 5' exonuclease activities, but lacks 5' to 3' exonuclease activity. It produces blunt end DNA products. Platinum™ SuperFi™ DNA Polymerase is supplied with a separate vial of SuperFi™ GC Enhancer designed for GC-rich templates (>65% GC). 		
	Important guidelines	Click here for important PCR guidelines.		
	Online resources	Visit our product page for additional information and protocols. For support, visit thermofisher.com/support .		

Enzyme characteristics

Hot-start:	Antibody
Length:	Up to 20 kb
Fidelity vs. Taq:	>100X
Format:	Separate components

PCR setup

Component	25- μL rxn	50- μL rxn	Custom	Final conc.
Water, nuclease-free	to 25 μL	to 50 μL	to μL	—
5X SuperFi™ Buffer ¹	5 μL	10 μL	μL	1X
10 mM dNTP mix	0.5 μL	1 μL	μL	0.2 mM each
10 μM forward primer	1.25 μL	2.5 μL	μL	0.5 μM
10 μM reverse primer	1.25 μL	2.5 μL	μL	0.5 μM
Template DNA ²	varies	varies		varies
5X SuperFi™ GC Enhancer (optional) ³	5 μL	10 μL	μL	1X
Platinum™ SuperFi™ DNA Polymerase (2 U/ μL)	0.25 μL	0.5 μL	μL	0.02 U/ μL

¹ Includes 7.5 mM MgCl_2 .

² 5–50 ng gDNA or 1 pg–10 ng plasmid DNA (see “**Optimization strategies**”, below, for more information).

³ Recommended for targets with >65% GC sequences.

PCR protocol

See page 2 and page 3 to prepare and run your PCR experiment.

Optimization strategies and troubleshooting




Click here for guidelines to optimize your PCR experiment.


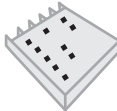
Click here for guidelines to troubleshoot your PCR experiment.

Purchaser notification

Click here for Limited Warranty, Disclaimer, and Licensing information.

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

Steps	Action	Procedure details																		
1 	Thaw reagents	Thaw, mix, and briefly centrifuge each component before use.																		
2 	Prepare PCR master mix	<p>Add the following components to each PCR tube.</p> <p>Note: 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>50-μL rxn</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>Water, nuclease-free</td> <td>to 50 μL</td> <td></td> </tr> <tr> <td>5X SuperFi™ Buffer¹</td> <td>10 μL</td> <td>1X</td> </tr> <tr> <td>10 mM dNTP mix</td> <td>1 μL</td> <td>0.2 mM each</td> </tr> <tr> <td>5X SuperFi™ GC Enhancer (<i>optional</i>)²</td> <td>10 μL</td> <td>1X</td> </tr> <tr> <td>Platinum™ SuperFi™ DNA Polymerase</td> <td>0.5 μL</td> <td>0.02 U/μL</td> </tr> </tbody> </table> <p>¹ Includes 7.5 mM MgCl₂. ² Recommended for targets with >65% GC sequences.</p> <p>Mix and then briefly centrifuge the components.</p>	Component	50- μ L rxn	Final conc.	Water, nuclease-free	to 50 μ L		5X SuperFi™ Buffer ¹	10 μ L	1X	10 mM dNTP mix	1 μ L	0.2 mM each	5X SuperFi™ GC Enhancer (<i>optional</i>) ²	10 μ L	1X	Platinum™ SuperFi™ DNA Polymerase	0.5 μ L	0.02 U/ μ L
Component	50- μ L rxn	Final conc.																		
Water, nuclease-free	to 50 μ L																			
5X SuperFi™ Buffer ¹	10 μ L	1X																		
10 mM dNTP mix	1 μ L	0.2 mM each																		
5X SuperFi™ GC Enhancer (<i>optional</i>) ²	10 μ L	1X																		
Platinum™ SuperFi™ DNA Polymerase	0.5 μ L	0.02 U/ μ L																		
3 	Add template DNA and primers	<p>Add your template DNA and primers to each tube for a final reaction volume of 50 μL.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>50-μL rxn</th> <th>Final conc.</th> </tr> </thead> <tbody> <tr> <td>10 μM forward primer</td> <td>2.5 μL</td> <td>0.5 μM</td> </tr> <tr> <td>10 μM reverse primer</td> <td>2.5 μL</td> <td>0.5 μM</td> </tr> <tr> <td>Template DNA¹</td> <td>varies</td> <td>varies</td> </tr> </tbody> </table> <p>¹ Optimal amount of low complexity DNA (plasmid, phage, BAC DNA) is 1 pg–10 ng per 50 μL reaction, but it can be varied from 0.1 pg to 50 ng per 50 μL reaction. Optimal amount of genomic DNA is 5–50 ng per 50 μL reaction, but it can be varied from 0.1 ng to 250 ng per 50 μL reaction.</p> <p>Cap each tube, mix, and then briefly centrifuge the contents.</p>	Component	50- μ L rxn	Final conc.	10 μ M forward primer	2.5 μ L	0.5 μ M	10 μ M reverse primer	2.5 μ L	0.5 μ M	Template DNA ¹	varies	varies						
Component	50- μ L rxn	Final conc.																		
10 μ M forward primer	2.5 μ L	0.5 μ M																		
10 μ M reverse primer	2.5 μ L	0.5 μ M																		
Template DNA ¹	varies	varies																		

Steps	Action	Procedure details							
4 	Incubate reactions in a thermal cycler			2-step protocol (<10kb)		3-step protocol (<10kb)		Long PCR (>10kb)	
		Step		Temp.	Time	Temp.	Time	Temp.	Time
		Initial denaturation		98°C	30 sec	98°C	30 sec	95°C	2 min
		25–35 PCR cycles	Denature	98°C	5–10 sec	98°C	5–10 sec	95°C	10 sec
			Anneal ¹	—	—	varies	10 sec	varies	10 sec
			Extend	72°C	15–30 sec/kb	72°C	15–30 sec/kb	68°C	30 sec/kb
		Final extension		72°C	5 min	72°C	5 min	68°C	5 min
				4°C	hold	4°C	hold	4°C	hold
		¹ IMPORTANT! Always use the T_m calculator on our website at www.thermofisher.com/tmcalculator to calculate the T_m of your primers and the recommended annealing temperature.							
		Note: Refer to “ Optimization strategies ”, page 1, for guidelines to optimize cycling conditions.							
5 	Add gel loading buffer and analyze with gel electrophoresis	Add gel loading buffer to 10 µL of PCR product, mix, and briefly centrifuge the contents. Analyze the sample using agarose gel electrophoresis. Use your PCR product immediately in down-stream applications, or store it at –20°C.							