

**Package contents**

Catalog No.	Size
12594025	25 reaction
12594100	100 reaction
12595025	25 reaction (with ezDNase™ Enzyme)
12595100	100 reaction (with ezDNase™ Enzyme)

 Kit contents

**Storage conditions**

Store all contents at -20°C (non-frost-free)

**Required materials**

- Template: RNA
- Gene specific primers (forward and reverse)

**Timing**

- Preparation time: 5 minutes
- Total cycling time: varies with target size (1–3.5 hours)

**Product description**

- The Invitrogen™ SuperScript™ IV One-Step RT-PCR System is designed for sensitive end-point detection and analysis of RNA by RT-PCR. The convenient formulation enables both cDNA synthesis and PCR amplification to be performed in a single reaction tube using gene-specific primers.
- The system can amplify RNA targets up to 14 kb in length, and is compatible with multiplex RT-PCR.
- SuperScript™ IV RT Mix allows high efficiency cDNA synthesis due to high thermostability, processivity, and ability to synthesize cDNA efficiently from a variety of RNA samples.
- 2X Platinum™ SuperFi™ RT-PCR Master Mix contains Platinum™ SuperFi™ DNA Polymerase which has high specificity, offers high yield, and is ideally suited for PCR applications that require sequence accuracy. The master mix uses a buffer system optimized for both RT and PCR reactions in the same tube.
- ezDNase™ enzyme (available separately as [Cat. No. 11766051](#)) is included with the SuperScript™ IV One-Step RT PCR System with ezDNase™ Kits for protocols that require removal of residual gDNA contamination.

**Online resources**

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

**Guidelines for RNA samples**

- This kit is optimized for use with 0.01 pg to 1 µg of total RNA, and is compatible with total RNA, mRNA, viral RNA, or *in vitro* transcribed RNA.
- High-quality, intact RNA is essential for RT-PCR, particularly for long targets. RNA must be devoid of RNase contamination and handled using aseptic conditions.
- Isolate total RNA with [TRIZOL™ Reagent](#), the [PureLink™ RNA Mini Kit](#), or the [MagMAX™-96 Total RNA Isolation Kit](#).
- Determine RNA quality using a bioanalyzer or by agarose gel electrophoresis.

**Guidelines for primers**

- Use gene-specific primers (GSPs) with the SuperScript™ IV One-Step RT-PCR System. Oligo(dT) or random primers are not recommended, because nonspecific products can be generated, thereby reducing the amount of target RT-PCR product.
- A final concentration of 0.5 µM for each primer is recommended, but further optimization may be necessary.
- Design primers that anneal to the mRNA sequence in exons on both sides of an intron or exon/exon boundary, to allow differentiation between the amplified cDNA and potential contaminating genomic DNA. If this approach is not feasible, use the protocol that involves residual gDNA removal with ezDNase™ enzyme.
- Ensure that primers are not self-complementary or complementary to each other at the 3' ends.
- To calculate primer T<sub>m</sub> and estimate appropriate annealing temperatures for PCR, use the T<sub>m</sub> calculator at [thermofisher.com/tmcalculator](http://thermofisher.com/tmcalculator)

** Limited product warranty****Important licensing information**

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**Disclaimer:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

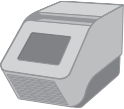


Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.



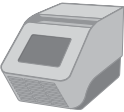


## Guidelines for cDNA synthesis and PCR

- Keep all components, reaction mixes, and samples on ice.
- After preparing reaction mixes, transfer them to the preheated thermal cycler and start the RT-PCR program.
- The optimal conditions for RT-PCR depend on primer and target sequences.
- Efficient cDNA synthesis can be accomplished by a 10-minute incubation at 45–60°C. A 50°C incubation is recommended as a general starting point.
- For GC-rich or structurally complex RNA templates, increasing the cDNA synthesis incubation temperatures up to 55–60 °C may improve RT-PCR results.
- Use 40 cycles of amplification for RT-PCR products ≤3 kb, and 35 cycles for RT-PCR products >3 kb.
- For very low RNA input or rare targets, increasing the number of PCR cycles to 40 may improve results.
- The PCR extension time varies with the size of the amplicon (recommended extension time is approximately 30 seconds per 1 kb of amplicon).

## SuperScript™ IV One-Step RT-PCR System protocol

Step	Action	Procedure details																								
<b>1</b> 	<b>Program thermal cycler</b>	<table border="1"> <thead> <tr> <th>Step</th> <th>Temp</th> <th>Time</th> <th>No. cycles</th> </tr> </thead> <tbody> <tr> <td>Reverse transcription</td> <td>45–60°C</td> <td>10 min</td> <td rowspan="2">1</td> </tr> <tr> <td>RT inactivation/initial denaturation</td> <td>98°C</td> <td>2 min</td> </tr> <tr> <td rowspan="2">Amplification</td> <td>98°C</td> <td>10 s</td> <td rowspan="3">35–40 <sup>[2]</sup></td> </tr> <tr> <td>55–72°C <sup>[1]</sup></td> <td>10 s</td> </tr> <tr> <td></td> <td>72°C</td> <td>30 s/kb</td> </tr> <tr> <td>Final extension</td> <td>72°C</td> <td>5 min</td> <td>1</td> </tr> </tbody> </table>	Step	Temp	Time	No. cycles	Reverse transcription	45–60°C	10 min	1	RT inactivation/initial denaturation	98°C	2 min	Amplification	98°C	10 s	35–40 <sup>[2]</sup>	55–72°C <sup>[1]</sup>	10 s		72°C	30 s/kb	Final extension	72°C	5 min	1
		Step	Temp	Time	No. cycles																					
		Reverse transcription	45–60°C	10 min	1																					
		RT inactivation/initial denaturation	98°C	2 min																						
		Amplification	98°C	10 s	35–40 <sup>[2]</sup>																					
			55–72°C <sup>[1]</sup>	10 s																						
	72°C	30 s/kb																								
Final extension	72°C	5 min	1																							
<p>[1] <b>IMPORTANT!</b> Use the <math>T_m</math> calculator at <a href="http://thermofisher.com/tmcalculator">thermofisher.com/tmcalculator</a> to determine actual annealing temperature.</p> <p>[2] Use 40 cycles for short (≤3 kb) templates only.</p>																										
<b>2</b> 	<b>Prepare RT-PCR reaction mix</b>	<p>a. Combine the following components in a 0.2-mL, nuclease-free, thin-walled PCR tube on ice.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>Volume</th> </tr> </thead> <tbody> <tr> <td>2X Platinum™ SuperFi™ RT-PCR Master Mix</td> <td>25 µL</td> </tr> <tr> <td>Forward primer (10 µM)</td> <td>2.5 µL</td> </tr> <tr> <td>Reverse primer (10 µM)</td> <td>2.5 µL</td> </tr> <tr> <td>SuperScript™ IV RT Mix <sup>[1]</sup></td> <td>0.5 µL</td> </tr> <tr> <td>Template RNA (0.01 pg to 1 µg total RNA)</td> <td>varies</td> </tr> <tr> <td>Nuclease-free water</td> <td>to 50 µL</td> </tr> </tbody> </table>	Component	Volume	2X Platinum™ SuperFi™ RT-PCR Master Mix	25 µL	Forward primer (10 µM)	2.5 µL	Reverse primer (10 µM)	2.5 µL	SuperScript™ IV RT Mix <sup>[1]</sup>	0.5 µL	Template RNA (0.01 pg to 1 µg total RNA)	varies	Nuclease-free water	to 50 µL										
		Component	Volume																							
		2X Platinum™ SuperFi™ RT-PCR Master Mix	25 µL																							
		Forward primer (10 µM)	2.5 µL																							
		Reverse primer (10 µM)	2.5 µL																							
		SuperScript™ IV RT Mix <sup>[1]</sup>	0.5 µL																							
		Template RNA (0.01 pg to 1 µg total RNA)	varies																							
Nuclease-free water	to 50 µL																									
<p>[1] Use all components except the SuperScript™ IV RT Mix for no RT control reactions.</p>																										
<p>b. Mix gently and ensure all the components are at the bottom of the amplification tube. Centrifuge briefly if needed.</p>																										
<b>3</b> 	<b>Run thermal cycler</b>	<p>Place the reaction in the pre-heated thermal cycler and run program set up in step 1.</p>																								

# SuperScript™ IV One-Step RT-PCR System with ezDNase™ Enzyme protocol

Step	Action	Procedure details																							
1 	<b>Prepare gDNA digestion reaction mix</b>	<p>Prepare a 10 µL gDNA digestion reaction mix for each RT-PCR reaction. Mix the following components in a 0.2-mL, nuclease-free, thin-walled PCR tube on ice.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>Volume</th> </tr> </thead> <tbody> <tr> <td>10X ezDNase™ Buffer</td> <td>1 µL</td> </tr> <tr> <td>ezDNase™ Enzyme</td> <td>1 µL</td> </tr> <tr> <td>Template RNA (0.01 pg to 1 µg total RNA)</td> <td>varies</td> </tr> <tr> <td>Nuclease-free Water</td> <td>to 10 µL</td> </tr> </tbody> </table>	Component	Volume	10X ezDNase™ Buffer	1 µL	ezDNase™ Enzyme	1 µL	Template RNA (0.01 pg to 1 µg total RNA)	varies	Nuclease-free Water	to 10 µL													
Component	Volume																								
10X ezDNase™ Buffer	1 µL																								
ezDNase™ Enzyme	1 µL																								
Template RNA (0.01 pg to 1 µg total RNA)	varies																								
Nuclease-free Water	to 10 µL																								
2 	<b>Digest gDNA</b>	<p>a. Gently mix and incubate at 37°C for 5 minutes.            b. (Optional) For long RT-PCR products (&gt;3 kb) add 2 µL of 13.2 mM EDTA, then incubate at 65°C for 10 minutes.            c. Briefly centrifuge the reaction mix and place the tube on ice.</p>																							
3 	<b>Program thermal cycler</b>	<table border="1"> <thead> <tr> <th>Step</th> <th>Temp</th> <th>Time</th> <th>No. cycles</th> </tr> </thead> <tbody> <tr> <td>Reverse transcription</td> <td>45–60°C</td> <td>10 min</td> <td rowspan="2">1</td> </tr> <tr> <td>RT inactivation/initial denaturation</td> <td>98°C</td> <td>2 min</td> </tr> <tr> <td rowspan="3">Amplification</td> <td>98°C</td> <td>10 s</td> <td rowspan="3">35–40 [2]</td> </tr> <tr> <td>55–72°C [1]</td> <td>10 s</td> </tr> <tr> <td>72°C</td> <td>30 s/kb</td> </tr> <tr> <td>Final extension</td> <td>72°C</td> <td>5 min</td> <td>1</td> </tr> </tbody> </table> <p>[1] <b>IMPORTANT!</b> Use the <math>T_m</math> calculator at <a href="http://thermofisher.com/tmcalculator">thermofisher.com/tmcalculator</a> to determine actual annealing temperature.            [2] Use 40 cycles for short (<math>\leq 3</math> kb) templates only.</p>	Step	Temp	Time	No. cycles	Reverse transcription	45–60°C	10 min	1	RT inactivation/initial denaturation	98°C	2 min	Amplification	98°C	10 s	35–40 [2]	55–72°C [1]	10 s	72°C	30 s/kb	Final extension	72°C	5 min	1
Step	Temp	Time	No. cycles																						
Reverse transcription	45–60°C	10 min	1																						
RT inactivation/initial denaturation	98°C	2 min																							
Amplification	98°C	10 s	35–40 [2]																						
	55–72°C [1]	10 s																							
	72°C	30 s/kb																							
Final extension	72°C	5 min	1																						
4 	<b>Prepare RT-PCR reaction mix</b>	<p>a. Combine the following components in the tube on ice containing the template RNA.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>Volume</th> </tr> </thead> <tbody> <tr> <td>Template RNA (from step 2)</td> <td>10 µL</td> </tr> <tr> <td>2X Platinum™ SuperFi™ RT-PCR Master Mix</td> <td>25 µL</td> </tr> <tr> <td>Forward primer (10 µM)</td> <td>2.5 µL</td> </tr> <tr> <td>Reverse primer (10 µM)</td> <td>2.5 µL</td> </tr> <tr> <td>SuperScript™ IV RT Mix [1]</td> <td>0.5 µL</td> </tr> <tr> <td>Nuclease-free water</td> <td>to 50 µL</td> </tr> </tbody> </table> <p>[1] Use all components except the SuperScript™ IV RT Mix for no RT control reactions.</p> <p>b. Mix gently and ensure all the components are at the bottom of the amplification tube. Centrifuge briefly if needed.</p>	Component	Volume	Template RNA (from step 2)	10 µL	2X Platinum™ SuperFi™ RT-PCR Master Mix	25 µL	Forward primer (10 µM)	2.5 µL	Reverse primer (10 µM)	2.5 µL	SuperScript™ IV RT Mix [1]	0.5 µL	Nuclease-free water	to 50 µL									
Component	Volume																								
Template RNA (from step 2)	10 µL																								
2X Platinum™ SuperFi™ RT-PCR Master Mix	25 µL																								
Forward primer (10 µM)	2.5 µL																								
Reverse primer (10 µM)	2.5 µL																								
SuperScript™ IV RT Mix [1]	0.5 µL																								
Nuclease-free water	to 50 µL																								
5 	<b>Run thermal cycler</b>	Place the reaction in the pre-heated thermal cycler and run program set up in step 3.																							