



Contents and storage

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

SuperScript™ IV UniPrime™ One-Step RT-PCR System			
Component	Cat. No. 12596025 (25 Rx ^[1])	Cat. No. 12596100 (100 Rx ^[1])	Cat. No. 12596500 (500 Rx ^[1])
SuperScript™ IV RT Mix	50 µL	200 µL	5 × Cat. No. 12596100
2X UniPrime™ RT-PCR Master Mix	625 µL	2 × 1.25 mL	
Water, Nuclease-free	1.25 mL	2 × 1.25 mL	
SuperScript™ IV UniPrime™ One-Step RT-PCR System, Colored			
Component	Cat. No. 12597025 (25 Rx ^[1])	Cat. No. 12597100 (100 Rx ^[1])	Cat. No. 12597500 (500 Rx ^[1])
SuperScript™ IV RT Mix, Red	50 µL	200 µL	5 × Cat. No. 12597100
2X UniPrime™ RT-PCR Master Mix, Blue	625 µL	2 × 1.25 mL	
Water, Nuclease-free	1.25 mL	2 × 1.25 mL	

[1] Number of reactions is based on 50 µL RT-PCR reaction volumes.



Product description

The Invitrogen™ SuperScript™ UniPrime™ 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 provides sensitivity down to 0.01 pg, can amplify RNA targets up to 13 kb in length, and is compatible with multiplex RT-PCR.
- A two-phase hot start activation mechanism allows easy room-temperature setup and benchtop stability of assembled reactions.
- The SuperScript™ IV RT Mix allows highly efficient reverse transcription step due to high thermostability, processivity, and ability to synthesize cDNA from variety of RNA samples.



Online resources

- Visit the product page at thermofisher.com for additional information and protocols.
- For support, visit thermofisher.com/support.



Product description, continued

- The 2X UniPrime™ RT-PCR Master Mix contains modified Platinum™ SuperFi™ II DNA Polymerase, which provides high specificity, high yields, and is ideally suited for PCR applications that require sequence accuracy. Novel master mix formulation enables universal primer annealing at 60°C for most primer pairs designed following the general design rules.
- A colored version of the kit makes it easy to track pipetting steps and enables direct sample loading on gels.



Required materials

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



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, and *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 primer design

- Use gene-specific primers (GSPs) with the SuperScript™ IV UniPrime™ 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 for single-plex reactions 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, or perform gDNA removal on template RNA (See “(Optional) Perform gDNA removal”).
- Ensure that primers are not self-complementary or complementary to each other at the 3' ends.

Guidelines for cDNA synthesis and PCR

- Efficient cDNA synthesis can be accomplished by a 10-minute incubation at 50°C. This incubation is recommended as a general starting point.
- For GC-rich or structurally complex RNA templates, increasing the cDNA synthesis incubation temperatures up to 65°C may improve RT-PCR results.
- Use 30–40 cycles of amplification for RT-PCR products ≤3 kb, and up to 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. The recommended extension time is 30 seconds per 1 kb of amplicon (for a faster protocol, optimization of up to 20 seconds per 1 kb is possible).
- Prepared reactions can be stored at room temperature for up to 24 hours, when the target is ≤3 kb in length. For amplification of longer targets, preassembled reactions can be stored for up to 4 hours.
- In the colored format (Cat. No. 12597025, 12597100, 12597500), when the red SuperScript™ IV RT mix is added to the blue UniPrime™ RT-PCR Master Mix, the final reaction mix turns purple. Colored buffers allow the user to track pipetting steps, and also enable direct loading of RT-PCR products onto the gel.
- RT-PCR products obtained with both colored and colorless formats can be successfully used in further downstream applications such as sequencing, cloning, and genotyping.
- Certain components of the SuperScript™ IV UniPrime™ One-Step RT-PCR System may form precipitates after performing the RT-PCR reaction. These precipitates do not influence RT-PCR results or usability of the RT-PCR product for downstream purposes.

Guidelines for multiplex RT-PCR

Due to its high processivity and specificity, as well as the specially formulated reaction buffer, SuperScript™ IV UniPrime™ One-Step RT-PCR System is suitable for multiplex RT-PCR applications without the need for significant optimization.

Template

- Recommended RNA range for detection is from 0.1 ng up to 1 µg.

Primers

- For multiplex RT-PCR, reduce primer concentration to 0.1 µM each. If required, primer concentration may be optimized in the range of 0.1–0.5 µM.
- Special attention to primer design parameters is critical for the success of multiplex RT-PCR.
- Optimal primers for multiplexing should be 21–34 nt in length and have 40–60% GC content. Ideally, G and C nucleotides should be distributed uniformly along the primer.
- Avoid significant homology between the primers and self-complementary primer regions, as well as three or more G or C nucleotides at the 3' end.
- If possible, the primer should terminate with a G or C at the 3' end.

Amplicons



- For product analysis on gel, up to 12 RNA targets can be amplified in a single reaction. Maximal amplicon length is up to 13 kb.
- For effective separation in agarose gel electrophoresis, ensure that the amplicon sizes differ sufficiently.

Cycling



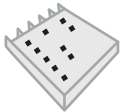
- For amplification of multiple targets, an annealing time of 1 minute and an extension time of 30 seconds/kb at 72°C is recommended.
- In order to achieve more output, the extension time can be extended to 60 seconds/kb, for for targets up to 1 kb. For longer targets prolonged extension time is not recommended.
- 35 cycles for multiplex reaction is recommended.

Prepare RT-PCR reaction mix

Note: If necessary, perform a gDNA digestion reaction with the template RNA before use in the RT-PCR reaction mix (See “(Optional) Perform gDNA removal” on page 5).


Step	Action	Procedure details																												
1 	Thaw reagents	Thaw reagents on ice. Mix and briefly centrifuge all reagents to ensure homogeneity before use.																												
2 	Prepare RT-PCR reaction mix	<p>a. Add the following components to each PCR tube.</p> <table border="1"> <thead> <tr> <th>Component</th> <th>20-μL rxn</th> <th>50-μL rxn</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>2X UniPrime™ RT-PCR Master Mix or 2X UniPrime™ RT-PCR Master Mix, Blue</td> <td>10 μL</td> <td>25 μL</td> <td>1 X</td> </tr> <tr> <td>Forward primer (10 μM)</td> <td>1 μL</td> <td>2.5 μL</td> <td>0.5 μM</td> </tr> <tr> <td>Reverse primer (10 μM)</td> <td>1 μL</td> <td>2.5 μL</td> <td>0.5 μM</td> </tr> <tr> <td>25X SuperScript™ IV RT Mix 25X SuperScript™ IV RT Mix, Red ^[1]</td> <td>0.8 μL</td> <td>2.0 μL</td> <td>1 X</td> </tr> <tr> <td>Template RNA</td> <td>0.01 μg to 0.4 μg</td> <td>0.01 μg to 1 μg</td> <td>0.01 μg to 1 μg total RNA</td> </tr> <tr> <td>Water, nuclease-free</td> <td>Fill to 20 μL</td> <td>Fill to 50 μL</td> <td>—</td> </tr> </tbody> </table> <p>[1] For no RT control reactions, add all components except the SuperScript™ IV RT Mix.</p> <p>b. Mix gently and ensure all the components are at the bottom of the amplification tube. Centrifuge briefly if needed. Note: Assembled reactions can be stored for up to 24 h at room temperature prior amplification of targets up to 3 kb. For targets \geq3 kb, do not keep assembled reactions at room temperature for more than 4 h. With RNA samples of suboptimal purity, proceed immediately to thermal cycling.</p> <p>c. Proceed to “Perform RT-PCR and analyze results” on page 4.</p>	Component	20- μ L rxn	50- μ L rxn	Final concentration	2X UniPrime™ RT-PCR Master Mix or 2X UniPrime™ RT-PCR Master Mix, Blue	10 μ L	25 μ L	1 X	Forward primer (10 μ M)	1 μ L	2.5 μ L	0.5 μ M	Reverse primer (10 μ M)	1 μ L	2.5 μ L	0.5 μ M	25X SuperScript™ IV RT Mix 25X SuperScript™ IV RT Mix, Red ^[1]	0.8 μ L	2.0 μ L	1 X	Template RNA	0.01 μ g to 0.4 μ g	0.01 μ g to 1 μ g	0.01 μ g to 1 μ g total RNA	Water, nuclease-free	Fill to 20 μ L	Fill to 50 μ L	—
Component	20- μ L rxn	50- μ L rxn	Final concentration																											
2X UniPrime™ RT-PCR Master Mix or 2X UniPrime™ RT-PCR Master Mix, Blue	10 μ L	25 μ L	1 X																											
Forward primer (10 μ M)	1 μ L	2.5 μ L	0.5 μ M																											
Reverse primer (10 μ M)	1 μ L	2.5 μ L	0.5 μ M																											
25X SuperScript™ IV RT Mix 25X SuperScript™ IV RT Mix, Red ^[1]	0.8 μ L	2.0 μ L	1 X																											
Template RNA	0.01 μ g to 0.4 μ g	0.01 μ g to 1 μ g	0.01 μ g to 1 μ g total RNA																											
Water, nuclease-free	Fill to 20 μ L	Fill to 50 μ L	—																											

Perform RT-PCR and analyze results

Step	Action	Procedure details																								
1 	Program thermal cycler	<p>a. Add the following components to each PCR tube. If preparing a master mix, multiply the volumes of components to all.</p> <table border="1"> <thead> <tr> <th>Step</th> <th>Temperature</th> <th>Time</th> <th>No. cycles</th> </tr> </thead> <tbody> <tr> <td>Reverse transcription</td> <td>50°C</td> <td>10 minutes</td> <td>1</td> </tr> <tr> <td>RT inactivation/initial denaturation</td> <td>98°C</td> <td>2 minutes</td> <td>1</td> </tr> <tr> <td rowspan="3">Amplification</td> <td>98°C</td> <td>10 seconds</td> <td rowspan="3">30–40 ^[1]</td> </tr> <tr> <td>60°C</td> <td>10 seconds</td> </tr> <tr> <td>72°C</td> <td>30 seconds/kb</td> </tr> <tr> <td>Final extension</td> <td>72°C</td> <td>5 minutes</td> <td>1</td> </tr> </tbody> </table> <p>[1] Use up to 35 cycles for targets >3 kb.</p>	Step	Temperature	Time	No. cycles	Reverse transcription	50°C	10 minutes	1	RT inactivation/initial denaturation	98°C	2 minutes	1	Amplification	98°C	10 seconds	30–40 ^[1]	60°C	10 seconds	72°C	30 seconds/kb	Final extension	72°C	5 minutes	1
Step	Temperature	Time	No. cycles																							
Reverse transcription	50°C	10 minutes	1																							
RT inactivation/initial denaturation	98°C	2 minutes	1																							
Amplification	98°C	10 seconds	30–40 ^[1]																							
	60°C	10 seconds																								
	72°C	30 seconds/kb																								
Final extension	72°C	5 minutes	1																							
2 	Run thermal cycler	<p>a. Dispense the prepared reaction mix to separate PCR tubes, 8-well strips, or individual wells of a 96-well PCR plate.</p> <p>b. Place the reaction in the pre-heated thermal cycler and run program.</p>																								
3 	Analyze with gel electrophoresis	<p>a. Add gel loading buffer to PCR product, mix, and briefly centrifuge the contents.</p> <p>b. Reactions in the colored format do not need to be stained, load the purple sample directly into the gel well. Note: The blue and red tracking dyes that resolve on the gel correspond to PCR products of about 4 and 2 kb in length.</p> <p>c. For optimal separation using E-Gel™ precast agarose gels, dilute the sample 2- to 30-fold as directed in the recommendations for use of E-Gel™ agarose gels.</p> <p>d. After electrophoresis use your PCR product immediately in downstream applications, or store it at –20°C.</p>																								

(Optional) Perform gDNA removal

If not using primers that allow differentiation between the amplified cDNA and potential contaminating gDNA, perform a gDNA digestion reaction with the template RNA. Use Invitrogen™ ezDNase™ Enzyme for fast, easy, and RNA-friendly gDNA removal. Prepare a 10 µL gDNA digestion reaction mix for each RT-PCR reaction following the ezDNase™ Enzyme protocol.

Step	Action	Procedure details										
<p>1</p> 	<p>gDNA removal</p>	<p>a. Mix the following components in a 0.2-mL, nuclease-free, thin-walled PCR tube on ice.</p> <table border="1" data-bbox="579 297 1398 537"> <thead> <tr> <th>Component</th> <th>10-µL rxn</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</td> <td>0.01 µg to 1 µg total RNA</td> </tr> <tr> <td>Water, nuclease-free</td> <td>Fill to 10 µL</td> </tr> </tbody> </table> <p>b. Gently mix and incubate at 37°C for 2 minutes.</p> <p>c. <i>(Optional)</i> If the RNA sample is to be used for RT-PCR of fragments ≥3 kb, incubate the sample at 55°C for 5 minutes in the presence of 10 mM DTT to inactivate the enzyme.</p> <p>d. Chill the tube on ice to bring the sample to room temperature, then briefly centrifuge and place the tube on ice.</p> <p>e. Proceed to “Prepare RT-PCR reaction mix” on page 3.</p> <p>Note: Depending on the RT-PCR reaction volume, use all or part of the prepared RNA volume after gDNA removal. Do not exceed 20% of final reaction volume (up to 10 µL per 50 µL RT-PCR reaction or up to 4 µL per 20 µL RT-PCR reaction).</p>	Component	10-µL rxn	10X ezDNase™ Buffer	1 µL	ezDNase™ Enzyme	1 µL	Template RNA	0.01 µg to 1 µg total RNA	Water, nuclease-free	Fill to 10 µL
Component	10-µL rxn											
10X ezDNase™ Buffer	1 µL											
ezDNase™ Enzyme	1 µL											
Template RNA	0.01 µg to 1 µg total RNA											
Water, nuclease-free	Fill to 10 µL											

Related products and equipment

Product	Cat. No.
TRIzol™ Reagent	15596026
PureLink™ RNA Mini Kit	12183020
MagMAX™-96 Total RNA Isolation Kit	AM1830
ezDNase™ Enzyme	11766051
RNaseZap™ RNase Decontamination Solution	AM9784
RNaseZap™ RNase Decontamination Wipes	AM9786
Water, nuclease-free	R0581
E-Gel™ EX Agarose Gels, 1%	G401001
E-Gel™ EX Agarose Gels, 2%	G401002
E-Gel™ EX Double Comb Agarose Gels, 1%	A42345
E-Gel™ EX Double Comb Agarose Gels, 2%	A42346
E-Gel™ Power Snap Plus Electrophoresis Systems with Extended Warranty	G9301
E-Gel™ Power Snap Plus Electrophoresis Systems, Starter Kit with Extended Warranty (E-Gel 48 1% with SYBR™ Safe)	G9341
E-Gel™ Power Snap Plus Electrophoresis Systems, Starter Kit with Extended Warranty (E-Gel 48 2% with SYBR™ Safe)	G9342
MicroAmp™ Optical 8-Tube Strip with Attached Optical Caps, 0.2 mL	A30588
EasyStrip™ Plus Tube Strip with Attached Ultra Clear Caps	AB2005
ProFlex™ 96-well PCR System	4484075
ProFlex™ 3 × 32-well PCR System	4484073
PCR plastics	thermofisher.com/us/en/home/life-science/pcr/pcr-plastics/plastics-selection-guide.html

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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.

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.

 **Manufacturer:** Thermo Fisher Scientific Baltics UAB | V. A. Graiciuno 8 | LT-02241 Vilnius, Lithuania

© 2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

For support, visit thermofisher.com/support.