

iScript™ One-Step RT-PCR Kit for Probes

50 x 50 µl reactions	170-8894
200 x 50 µl reactions	170-8895

For research purposes only.
Store at –20°C.

The iScript one-step RT-PCR kit for probes is a convenient and highly sensitive solution for real-time quantitative PCR of RNA templates using hybridization-based probe detection chemistries such as 5'-hydrolysis probes or molecular beacons. cDNA synthesis and PCR amplification are carried out in the same tube. This kit is optimized to deliver maximum RT-PCR efficiency, sensitivity and specificity. The reaction buffer has been specifically formulated to maximize activities of both the iScript reverse transcriptase and the iTaq™ DNA polymerase, while minimizing the potential for primer-dimer and other non-specific PCR artifacts.

The 2x RT-PCR reaction mix for probes includes the iTaq antibody-mediated hot-start DNA polymerase that sequesters activity prior to the initial PCR denaturation step. Upon heat activation, the antibody denatures irreversibly, releasing fully active and unmodified iTaq DNA polymerase.

Storage and Stability

Store the iScript one-step RT-PCR kit for probes at –20°C in a constant temperature freezer. When stored under these conditions, the kit components are stable for a minimum of one year after ship date. For extended stability, the iScript reverse transcriptase for one-step RT-PCR may be stored at –70°C. For convenience, the 2x RT-PCR reaction mix for probes may be stored unfrozen at 2°C to 8°C for up to 6 months. Repeated freeze-thaw cycles should be avoided.

Kit Contents

Reagent

iScript reverse transcriptase
for one-step RT-PCR
(yellow cap)

Description

Optimized 50x formulation of iScript MMLV reverse
transcriptase for one-step RT-PCR procedures

2x RT-PCR reaction
mix for probes
(blue cap)

2x reaction buffer containing 0.5 mM of each dNTP
(dATP, dCTP, dGTP, dTTP), magnesium ions,
iTaq DNA polymerase, stabilizers

Nuclease-free water

Reaction Set Up

To avoid primer-dimers, reactions should be assembled on ice.

Component	Volume per reaction
2x RT-PCR reaction mix for probes	25 µl
Forward primer, 100 to 500 nM final concentration	variable
Reverse primer, 100 to 500 nM final concentration	variable
Probe, 50 to 200 nM final concentration	variable
Nuclease-free water	variable
RNA template	variable
iScript reverse transcriptase for one-step RT-PCR	1 µl
Total Volume	50 µl

Reaction Protocol

Incubate complete reaction mix in a real-time thermal detection system as follows:

cDNA synthesis:	10 min at 50°C
iScript reverse transcriptase inactivation:	5 min at 95°C
PCR cycling and detection (30 to 45 cycles):	10 to 15 sec at 95°C
	30 sec at 55°C to 60°C (data collection step)

Recommendations for optimal results using the iScript One-Step RT-PCR Kit for Probes

Probe and primers should be designed according to standard qPCR guidelines.

Suggested input quantities of template are: 1 pg to 1 µg total RNA; 10 fg to 100 ng polyA(+) RNA.

First strand synthesis can be performed between 40°C and 52°C. Optimal results are generally obtained with a 10-minute incubation at 50°C. Incubation at temperatures higher than 50°C can delay or eliminate the detection of some non-specific amplification artifacts. However, this may also delay the C_t for detection of specific targets. We also recommend a 5-minute incubation at 95°C to fully inactivate the reverse transcriptase prior to PCR cycling.

Thaw all components, except the iScript reverse transcriptase, at room temperature. Mix gently, but thoroughly, and then centrifuge at 4°C to collect contents to the bottom of the tube. Chill on ice before using. Centrifuge again briefly at 4°C if needed.

Preparation of a reaction cocktail is crucial in quantitative PCR applications to reduce pipetting errors and maximize assay precision and accuracy. Assemble the reaction cocktail with all required components except sample template (total RNA) and dispense equal aliquots into each reaction tube. Add target sample to each reaction as the final step. Addition of sample as 5–10 µl volumes will improve assay precision. Replicate samples should be assembled as a master mix with a single addition of sample template.

Reagents and Materials Not Supplied

Gene-specific primers and probe

Pipet tips, aerosol barrier tips, such as:

Xcluda® Style B, 211-2006

Nuclease-free tubes or plates, such as:

0.2 ml thin-wall tubes, 223-9473 or plates, 223-9441

RNA purification kit, such as:

Aurum™ total RNA mini kit, 732-6820, or

Aurum total RNA kit, 2 x 96 well, 732-6800

To learn more about Bio-Rad's complete solution for Amplification, visit our website:

www.bio-rad.com/genomics

NOTICE TO PURCHASER: LIMITED LICENSE

Practice of the patented polymerase chain reaction (PCR) process requires a license. The Bio-Rad real-time detection systems include a licensed thermal cycler and may be used with PCR licenses available from Applied Biosystems. Its use with authorized reagents also provides a limited PCR license in accordance with the label rights accompanying such reagents. Some applications may require licenses from other parties. Aurum, iScript, iTaq, iQ, iCycler, Xcluda, and MyiQ are trademarks of Bio-Rad Laboratories. SYBR Green is a registered trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

A license to perform the patented 5' Nuclease Process for research is obtained by the purchase of (i) both Authorized 5' Nuclease Core Kit and Licensed Probe, (ii) a Licensed 5' Nuclease Kit, or (iii) license rights from Applied Biosystems.

This product is an Authorized 5' Nuclease Core Kit. Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,079,352, 5,789,224, 5,618,711, 6,127,155, 5,677,152, (Claims 1-23 only), 5,773,258, (claims 1 and 6 only), 5,407,800, 5,322,770, 5,310,652, 5,210,015, 5,487,972, and claims outside the US corresponding to US Patent No. 4,889,818. The purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. Separate purchase of a Licensed Probe would convey rights under the applicable claims on US Patents Nos. 5,538,848, 5,723,591, 5,876,930, 6,030,787, 6,258,569, 5,804,375 (claims 1-12 only), and 6,214,979, and corresponding claims outside the United States. No right under any other patent claim and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under Roche patents require a separate license from Roche. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.