

iScript™ One-Step RT-PCR Kit With SYBR® Green

170-8892 50 x 50 µl reactions
 170-8893 200 x 50 µl reactions

For research purposes only.

Storage at –20 °C Guaranteed for 12 months at –20 °C in a constant temperature freezer (multiple freeze/thaw cycles not recommended)

Storage at +4 °C Guaranteed for 4 months at 4 °C (protected from light)

The iScript one-step RT-PCR kit with SYBR® Green is a convenient and highly sensitive solution for real-time quantitative PCR of RNA templates. 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 without compromising fluorescent signal. The proprietary reaction buffer has been specifically formulated to maximize activities of both the reverse transcriptase and the iTaq™ DNA polymerase, while minimizing the potential for primer-dimer and other nonspecific PCR artifacts.

2x SYBR® Green RT-PCR reaction mix 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. Highly specific amplification is essential for successful qRT-PCR with SYBR® Green I technology, since this dye binds to any double-stranded DNA generated during amplification.

Kit Contents

Reagents

iScript reverse transcriptase for one-step RT-PCR (yellow cap)
 2x SYBR® Green RT-PCR reaction mix (green cap)

Nuclease-free water

Description

Optimized 50x formulation of iScript MMLV for one-step RT-PCR procedures
 2x reaction buffer containing 0.4 mM of each dNTP (dATP, dCTP, dGTP, dTTP), magnesium ions, iTaq DNA polymerase, 20 nM fluorescein, SYBR® Green I dye, stabilizers

Reaction Set Up

To maximize specificity, reactions should be assembled on ice.

Component

2x SYBR® Green RT-PCR reaction mix
 Forward primer (300 nM final concentration)
 Reverse primer (300 nM final concentration)
 Nuclease-free water
 RNA template (1 pg to 100 ng total RNA)
 iScript reverse transcriptase for one-step RT-PCR

Volume per Reaction

25 µl
 Variable
 Variable
 Variable
 Variable
 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
Reverse transcriptase inactivation	5 min at 95 °C
PCR cycling and detection (30 to 45 cycles)	10 sec at 95 °C
	30 sec at 55 °C-60 °C (data collection step)
Melt curve analysis (optional)	1 min at 95 °C
	1 min at 55 °C
	10 sec at 55 °C-95 °C (80 cycles, increasing by 0.5 °C each cycle)

Recommendations for Optimal Results Using the iScript One-Step RT-PCR Kit With SYBR® Green

Primers should be designed according to standard PCR guidelines with a length of 18-25 nucleotides, and a GC content of 40%-65%. Primer design should avoid internal secondary structure and complementarity at the 3' ends within each primer and primer pair. Optimal results may require titration of primer concentration between 100 and 500 nM. A final concentration of 300 nM per primer is effective for most reactions. In general, reaction efficiency and/or specificity can be optimized using equal concentrations of each primer. For best results, amplicon size should be limited to 50–200 bp.

Suggested input quantities of template are: 1 pg to 100 ng 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 min incubation at 50°C. Incubation at temperatures higher than 50°C can delay or eliminate the detection of some nonspecific amplification artifacts. However, this may also delay the C_t for detection of specific targets.

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.

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.

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

www.bio-rad.com/amplification.

NOTICE TO PURCHASER: LIMITED LICENSE

Practice of the patented polymerase chain reaction (PCR) process requires a license. The Bio-Rad real-time 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. AuroM, iScript, iTaq, iQ, iCycler, and MyiQ are trademarks of Bio-Rad Laboratories. SYBR Green is a registered trademark of Invitrogen Corporation.

This product is provided under an agreement between Molecular Probes, Inc. (a wholly owned subsidiary of Invitrogen Corporation) and Bio-Rad Laboratories and the manufacture, use, sale or import of this product is subject to one or more of U.S. Patents and corresponding international equivalents, owned by Molecular Probes, Inc. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer, where such research does not include testing, analysis or screening services for any third party in return for compensation on a per test basis. The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing, except for quality control or quality assurance purposes; (2) use of the product or its components to provide a service, information, or data, except for quality control or quality assurance purposes; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, USA. Tel: (541) 465-8300; Fax: (541) 335-0354.

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 to 23 only), 5,733,258 (claims 1 and 6 only), 5,407,800, 5,322,770, 5,310,652, 5,994,056, 6,171,785, 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. No right under any other patent claim (such as apparatus or system claims in US Patent No. 6,814,934) 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 by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.