

## iScript™ cDNA Synthesis Kits

Highly sensitive, convenient, and flexible, the iScript cDNA synthesis kit and the iScript Select cDNA synthesis kit command respect with startling efficiency.

- Purified MMLV RNase H<sup>+</sup> reverse transcriptase for sensitive detection
- Optimized, preblended components simplify reaction setup, save time, enhance reproducibility, and minimize the risk of cross-contamination
- Consistent performance from your choice of 2 complete kits — one with a flexible selection of primers, the other with convenient preblended primers



iScript cDNA Synthesis Kit



iScript Select cDNA Synthesis Kit

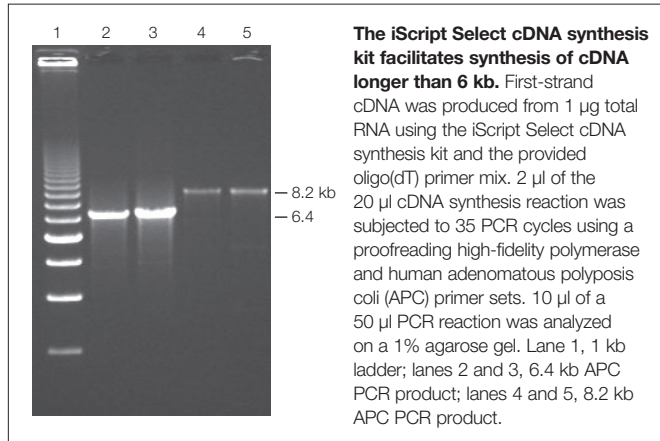
For more information, visit us on the Web at [www.bio-rad.com/iscript](http://www.bio-rad.com/iscript).

**BIO-RAD**

## High Quality and High Sensitivity

iScript cDNA synthesis kits are unique and sensitive reverse transcription systems that offer simplified handling and superior performance. Unlike other reverse transcription kits with multiple reaction components that must be combined during reaction setup, these streamlined kits feature an optimized, preblended reaction mix. Preblended reaction mixes reduce the number of steps required for reaction setup. This saves time, enhances reproducibility, and minimizes cross-contamination risk.

Choose between two optimized kits. For greatest convenience and sensitivity, choose the original iScript cDNA synthesis kit with oligo(dT) and random primers preblended into the reaction mix. For the ultimate combination of flexibility and sensitivity, choose the iScript Select cDNA synthesis kit, which allows first-strand priming with oligo(dT), random, or gene-specific primers.



## Ordering Information

Catalog #	Description
170-8890	<b>iScript cDNA Synthesis Kit</b> , 25 x 20 µl reactions, includes 5x iScript reaction mix, iScript reverse transcriptase, nuclease-free water
170-8891	<b>iScript cDNA Synthesis Kit</b> , 100 x 20 µl reactions
170-8896	<b>iScript Select cDNA Synthesis Kit</b> , 25 x 20 µl reactions, includes 5x iScript Select reaction mix, iScript reverse transcriptase, oligo(dT) mix, random primer mix, gene-specific primer (GSP) enhancer solution, nuclease-free water
170-8897	<b>iScript Select cDNA Synthesis Kit</b> , 100 x 20 µl reactions
170-8892	<b>iScript™ One-Step RT-PCR Kit With SYBR® Green</b> , 50 x 50 µl reactions
170-8893	<b>iScript™ One-Step RT-PCR Kit With SYBR® Green</b> , 200 x 50 µl reactions
170-8894	<b>iScript One-Step RT-PCR Kit for Probes</b> , 50 x 50 µl reactions
170-8895	<b>iScript One-Step RT-PCR Kit for Probes</b> , 200 x 50 µl reactions

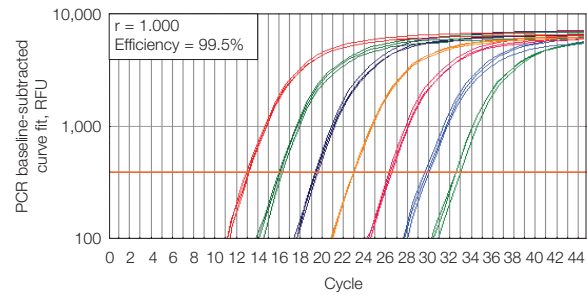
SYBR is a trademark of Molecular Probes, Inc. Bio-Rad Laboratories, Inc. is licensed by Molecular Probes, Inc. to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Practice of the patented 5' Nuclease Process requires a license from Applied Biosystems. The purchase of these products includes an immunity from suit under patents specified in the product insert to use only the amount purchased for the purchaser's own internal research when used with the separate purchase of Licensed Probe. No other patent rights are conveyed expressly, by implication, or by estoppel. Further information on purchasing licenses may be obtained from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

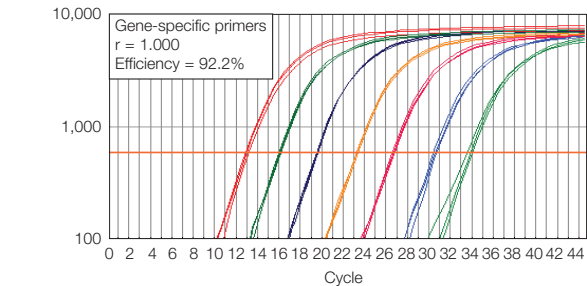
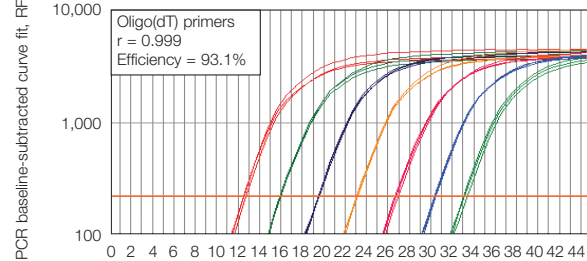
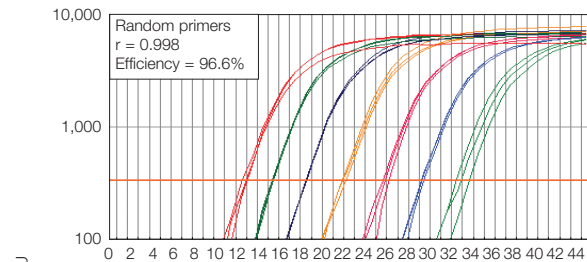
Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

## iScript cDNA Synthesis Kit



## iScript Select cDNA Synthesis Kit



**iScript cDNA synthesis kits perform reliably over 6 orders of magnitude using any first-strand priming approach.** Human total RNA from 1 pg to 1 µg was reverse transcribed using the iScript cDNA synthesis kit or the iScript Select cDNA synthesis kit and the first-strand primers indicated above. The resulting cDNA was amplified in triplicate using iQ™ SYBR® Green supermix with β-actin primers and detected in the iCycler iQ® real-time detection system. RFU, relative fluorescence units.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600  
Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000