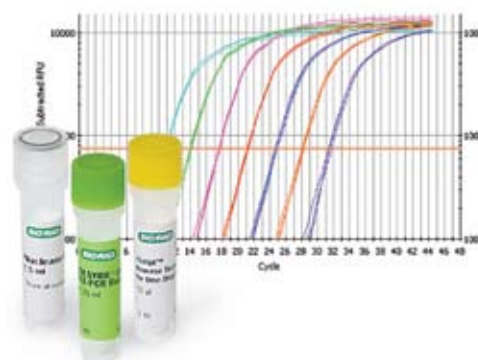


## iScript™ One-Step RT-PCR Kits

iScript one-step quantitative RT-PCR kits from Bio-Rad: two kits completely in sync with your research — one step does it all.

- Versatile — kits for use with either SYBR® Green or probe-based detection chemistries
- Sensitive — powerful combination of enzymes that accurately detect 100 fg of input RNA
- Convenient — 1-tube reaction setup minimizes handling and contamination risk

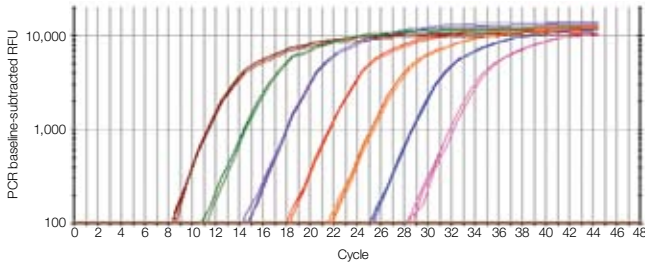
For more information, visit us on the Web at [www.bio-rad.com/iscript/](http://www.bio-rad.com/iscript/).



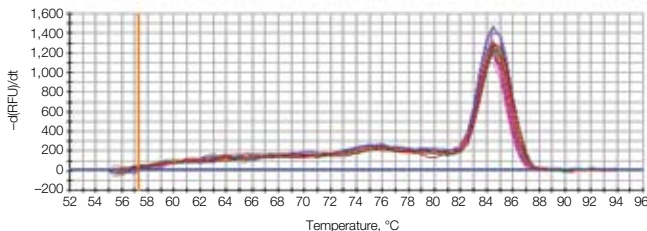
**BIO-RAD**

## iScript One-Step Quantitative RT-PCR Kits

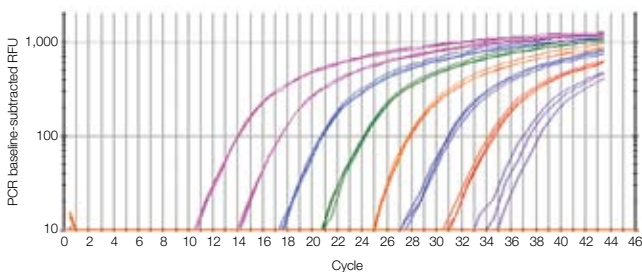
The iScript™ one-step RT-PCR kit with SYBR® Green and the iScript one-step RT-PCR kit for probes have been designed for optimal speed and convenience. Built on the powerful combination of iScript reverse transcriptase and iTaq™ DNA polymerase, these kits also allow sensitive detection down to 100 fg of input RNA. Discover quantitative RT-PCR kits that are completely in sync with your research. Just one step does it all.



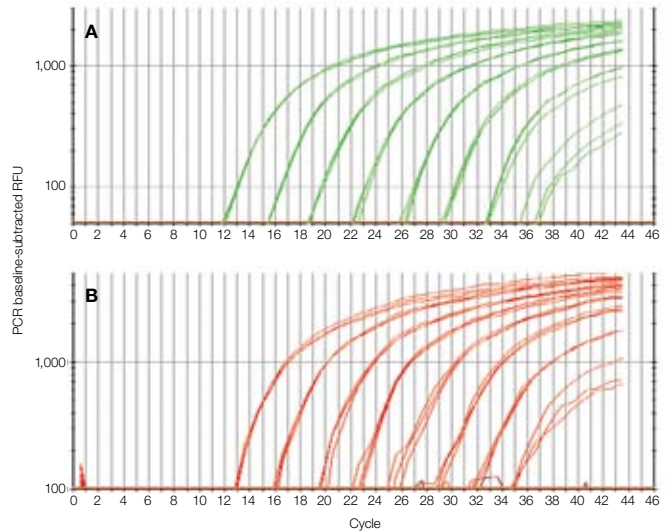
The iScript™ one-step RT-PCR kit with SYBR® Green provides high reproducibility and sensitivity across a broad range of concentrations. Reactions were performed in triplicate, along with no-template controls, using GAPDH primers and 100 ng to 100 fg of total HeLa RNA. Reactions were carried out on the iCycler iQ® real-time detection system. Standard curve had  $r = 1.000$ , slope =  $-3.466$ , and efficiency = 95%.



Melt-curve analysis of RT-PCR reactions performed using the iScript™ one-step RT-PCR kit with SYBR® Green shows absence of nonspecific PCR artifacts. Reactions were performed as for figure above. The melt-curve protocol was performed after RT-PCR, revealing a single clean melting peak for all samples tested.



The iScript one-step RT-PCR kit for probes delivers unparalleled results over an extremely wide dynamic range. RNA (1 µg to 100 fg) isolated from HeLa cells using the Aurum™ total RNA kit was reverse transcribed and amplified using primers to  $\beta$ -actin and a FAM-labeled detection probe. Each dilution was performed in triplicate, and RT-PCR was carried out on the iCycler iQ detection system. Standard curve had  $r = 1.000$ , slope =  $-3.39$ , and efficiency = 97.2%.



The iScript one-step RT-PCR kit for probes yields compelling results in a duplex real-time reaction. Tubulin and  $\beta$ -actin gene targets from HeLa total RNA were reverse transcribed and amplified in duplex over a 10-fold serial dilution. Each dilution was performed in triplicate, and RT-PCR was carried out on the iCycler iQ system. A dual-labeled FAM probe was used to detect the  $\beta$ -actin gene product (panel A) in the first dye layer. This reaction had a standard curve with  $r = 1.000$ , slope =  $-3.53$ , and efficiency = 92%. In the second dye layer, a dual-labeled Cy5 probe was used to detect the tubulin gene product (panel B). That reaction had a standard curve with  $r = 1.000$ , slope =  $-3.27$ , and efficiency = 102%.

### Ordering Information

Catalog #	Description
170-8892	iScript™ One-Step RT-PCR Kit With SYBR® Green, 50 x 50 µl reactions
170-8893	iScript™ One-Step RT-PCR Kit With SYBR® Green, 200 x 50 µl reactions
170-8894	iScript One-Step RT-PCR Kit for Probes, 50 x 50 µl reactions
170-8895	iScript One-Step RT-PCR Kit for Probes, 200 x 50 µl reactions

Cy is a trademark of GE Healthcare group companies. SYBR is a trademark of Invitrogen Corporation. Bio-Rad Laboratories, Inc. is licensed by Invitrogen Corporation to sell reagents containing SYBR Green I for use in real-time PCR, for research purposes only.

Purchase of this product 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. No other patent rights (such as 5' Nuclease Process patent rights) are conveyed expressly, by implication, or by estoppel. 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.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applied's United States Patent No. 6,814,934 B1 for use in research and for all other fields except the fields of human diagnostics and 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 Nos. 6,767,512 and 7,074,367.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 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