

iScript™ RT-qPCR Sample Preparation Reagent

Bio-Rad's iScript RT-qPCR sample preparation reagent delivers efficient cell lysis, RNA stabilization, and removal of genomic DNA for sensitive qPCR without RNA purification. This novel buffer accelerates and streamlines RT-qPCR analysis of cultured cells by eliminating the need for RNA purification and enabling PCR or real-time PCR directly from cell lysates. This system is ideal for rapid, high-throughput gene expression analysis, such as validation of siRNA-mediated gene knockdown.

- Rapid protocol efficiently removes genomic DNA and stabilizes RNA in 5–10 min
- Sensitive detection of high-, medium-, and low-copy gene targets directly from cell lysates
- Reagent enables multiplex real-time detection of up to 4 targets from as few as 10 cells
- Isolated RNA is stable for up to 6 months when frozen

For more information, visit us on the web at www.bio-rad.com/iscript/.

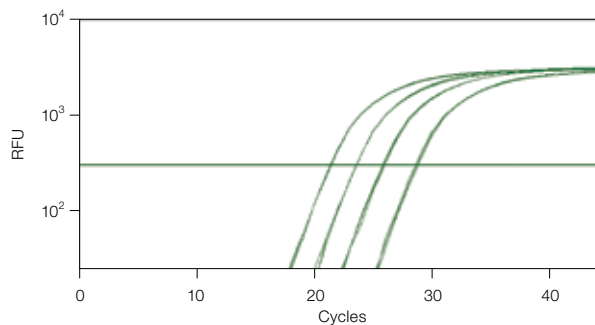
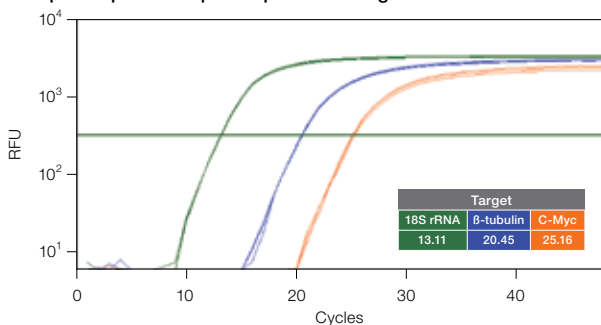


BIO-RAD

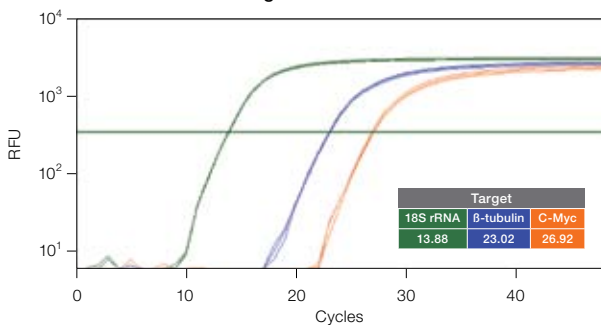
iScript RT-qPCR Sample Preparation Reagent

iScript RT-qPCR sample preparation reagent provides a simple and rapid procedure for isolating total RNA from cultured cells. It represents an entirely new approach to rapid, high-throughput gene expression analysis by eliminating the numerous, time-consuming steps required to purify RNA that is free of contaminating genomic DNA. This one buffer formulation contains all of the necessary components to perform cell lysis, RNA stabilization, and genomic DNA removal for sensitive, quantitative gene expression analysis.

iScript RT-qPCR Sample Preparation Reagent



PureZOL™ RNA Isolation Reagent



iScript RT-qPCR sample preparation reagent generates earlier threshold cycles (C_T s) compared to PureZOL RNA isolation reagent.

HeLa cells were treated with each reagent (125 cells/ μ l) according to the recommended guidelines and analyzed for 18S rRNA (■), β -tubulin (■), and c-Myc (■) expression levels using iScript cDNA synthesis kit and iQ™ SYBR® Green supermix on the CFX96™ real-time PCR detection system. iScript RT-qPCR sample preparation reagent delivers earlier C_T s as compared to PureZOL treatment. Average C_T values for triplicate reactions are shown for each gene target (inset tables). RFU, relative fluorescence units.

iScript RT-qPCR sample preparation reagent generates linear results over varying input cell amounts.

HeLa cells were treated with each reagent (125, 25, 5, and 1 cells/ μ l) according to the recommended guidelines and analyzed for *GAPDH* expression levels using iScript cDNA synthesis kit and iQ™ SYBR® Green supermix on the CFX96 real-time PCR detection system. iScript RT-qPCR sample preparation reagent delivers excellent efficiency and reproducibility for the detection of *GAPDH* from cell extracts. *GAPDH* efficiency = 94.4%, R^2 = 0.996. RFU, relative fluorescence units.

Ordering Information

Catalog #	Description
170-8898	iScript RT-qPCR Sample Preparation Reagent , 100 reactions, 10 ml, contains RNase inhibitors and RNA stabilizers
170-8899	iScript RT-qPCR Sample Preparation Reagent , 500 reactions, 5 x 10 ml, contains RNase inhibitors and RNA stabilizers

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.

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 Applied's United States Patent 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 Nos. 6,767,512 and 7,074,367.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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