

AMPLIFICATION: PCR REAGENTS

iScript Reverse Transcription Kits

Get the Most Out of Your cDNA Synthesis






Choose the best cDNA synthesis kit for your application, with formulations optimized for ease, yield, flexibility, fidelity, or genomic DNA (gDNA) clearance. Bio-Rad's iScript Reverse Transcription (RT) Kits provide fast and sensitive first-strand cDNA synthesis with high reproducibility.

- Highly efficient cDNA synthesis with high-quality results
- RNase H+ for accurate representation of gene expression
- RNase inhibition prevents RNA degradation



RNase H function. RNase H works to degrade RNA that is hybridized to DNA. During cDNA synthesis, the RNase H domain of the Moloney murine leukemia virus (MMLV) reverse transcriptase degrades the template RNA as the cDNA is synthesized. This activity ensures the one-to-one conversion of RNA into cDNA molecules.

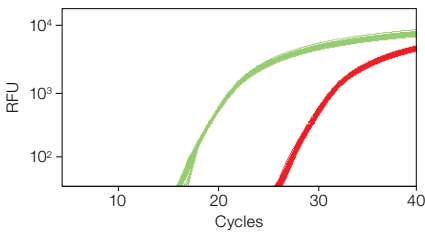
Which iScript Is My iScript?

Feature	 Minimum Setup Time MOST POPULAR	 Effective gDNA Elimination before RT	 Maximum Input RNA for High cDNA Yields	 Flexible Priming Options and High Fidelity	 Reliable Value Solution
Product	iScript Reverse Transcription Supermix for RT-qPCR	iScript gDNA Clear cDNA Synthesis Kit	iScript Advanced cDNA Synthesis Kit for RT-qPCR	iScript Select cDNA Synthesis Kit	iScript cDNA Synthesis Kit
Applications	<ul style="list-style-type: none"> ▪ Gene expression ▪ RNA quantification 	<ul style="list-style-type: none"> ▪ Gene expression ▪ RNA quantification 	<ul style="list-style-type: none"> ▪ Gene expression ▪ RNA quantification 	<ul style="list-style-type: none"> ▪ Cloning ▪ Sequencing and next-generation sequencing ▪ Gene expression ▪ RNA quantification 	<ul style="list-style-type: none"> ▪ Gene expression ▪ RNA quantification
Total input RNA	1 µg–1 pg	1 µg–1 pg	7.5 µg–100 fg	1 µg–1 pg	1 µg–100 fg
Format	1 tube	3 tubes	2 tubes	5 tubes	2 tubes
Kit contents	<ul style="list-style-type: none"> ▪ 5x iScript RT Supermix ▪ No-RT control supermix 	<ul style="list-style-type: none"> ▪ 5x iScript RT Supermix ▪ No-RT control supermix ▪ DNase ▪ DNase buffer 	<ul style="list-style-type: none"> ▪ iScript Reverse Transcriptase ▪ 5x iScript Advanced Reaction Mix 	<ul style="list-style-type: none"> ▪ iScript Reverse Transcriptase ▪ 5x iScript Reaction Mix ▪ 3 priming options 	<ul style="list-style-type: none"> ▪ iScript Reverse Transcriptase ▪ 5x iScript Reaction Mix
Time to produce cDNA	26 min	36 min	21 min	15–65 min	26 min
RNase H+ activity	✓	✓	✓	✓	✓

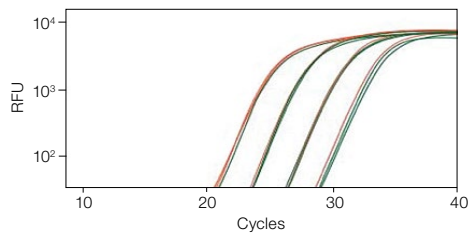
Visit bio-rad.com/myiScriptRTkit for more information.



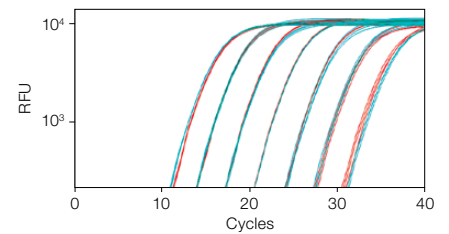
Unmatched Performance



Excellent data reproducibility. Forty-eight replicate cDNA synthesis reactions were prepared using the iScript Reverse Transcription Supermix for RT-qPCR and 100 ng (■) and 100 pg (■) of total RNA. A 160 bp fragment of the *PGK-1* gene was amplified from one-tenth of each of the resulting cDNA syntheses using SsoFast Probes Supermix on a CFX96 Real-Time PCR Detection System. Low standard deviations across the replicate cDNA synthesis reactions demonstrate exceptional data reproducibility. 100 ng RNA (■): mean Cq = 21.35 and SD = 0.123; 100 pg RNA (■): mean Cq = 31.56 and SD = 0.147. Cq, quantification cycle; RFU, relative fluorescence units; SD, standard deviation (in Cq).



Unbiased representation of 3' and 5' transcript regions. Reverse transcription of 100, 10, 1, and 0.1 ng input RNA was performed with iScript Reverse Transcription Supermix for RT-qPCR. Primer pairs were designed to the 5' (■, ~60 bp) and 3' (■, ~70 bp) ends of the *MAP* gene, and quantitative PCR (qPCR) was performed with one-tenth of the cDNA as input for the iTaq Universal SYBR® Green Supermix. There were no significant differences (<0.5 Cq difference) observed between the 5' and 3' assays. RFU, relative fluorescence units.



Broad linear dynamic range. Reverse transcription was performed from a tenfold dilution series of a human universal reference RNA (1 µg–1 pg) using the iScript cDNA Synthesis Kit (■) and the iScript Reverse Transcription Supermix for RT-qPCR (■). Amplification was performed in triplicate from one-tenth of the resulting cDNA using SsoAdvanced Universal SYBR® Green Supermix and the *GAPDH* PrimePCR Gene Expression Assay on a CFX96 Real-Time PCR Detection System. iScript cDNA Synthesis Kit: R² = 0.999, efficiency = 100.4%, and slope = -3.31; iScript Reverse Transcription Supermix for RT-qPCR: R² = 0.999, efficiency = 98.2%, and slope = -3.37. RFU, relative fluorescence units.

Need to Reliably Detect Low-Copy Targets in Limited Sample Material?

The new iScript Explore One-Step RT and PreAmp Kit generates preamplified cDNA directly from RNA in a single step, enabling unbiased target-specific preamplification of up to 100 targets.

Single Step

Go directly from RNA to preamplified cDNA in one step.



gDNA Clearance

Amplify only your target of interest.



Reaction Booster

Benefit from enhanced RT performance.



Visit bio-rad.com/iScriptExplore for more information.

Ordering Information

Kit Size	iScript Reverse Transcription Supermix for RT-qPCR	iScript gDNA Clear cDNA Synthesis Kit	iScript Advanced cDNA Synthesis Kit for RT-qPCR	iScript Select cDNA Synthesis Kit	iScript cDNA Synthesis Kit	iScript Explore One-Step RT and PreAmp Kit
25 x 20 µl reactions	1708840	1725034	1725037	1708896	1708890	—
100 x 20 µl reactions	1708841	1725035	1725038	1708897	1708891	—
500 x 20 µl reactions	1708841BUN	1725035BUN	1725038BUN	1708897BUN	1708891BUN	—
50 x 50 µl reactions	—	—	—	—	—	12004856
250 x 50 µl reactions	—	—	—	—	—	17002826

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

Bio-Rad's thermal cyclers and 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.

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

All trademarks used herein are the property of their respective owner.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00
Germany 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600
Japan 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30
Poland 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580
Sweden 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300
United Kingdom 44 01923 47 1301

