



iTaq™ Universal SYBR® Green One-Step Kit

Catalog #	Reverse Transcriptase Volume	Reaction Mix Volume	Kit Size
172-5150	50 µl (1 x 1 ml vials)	1 ml (1 x 1 ml vials)	100 x 20 µl reactions
172-5151	125 µl (1 x 1 ml vials)	5 ml (5 x 1 ml vials)	500 x 20 µl reactions

For research purposes only.

Storage and Stability

Guaranteed for 12 months in a constant temperature freezer at -20°C protected from light. For convenience, the reaction mix tube can be stored at 4°C for up to three months. The reverse transcriptase tube must be stored at -20°C .

Kit Contents

iTaq universal SYBR® Green one-step kit contains iScript™ reverse transcriptase, which is an RNase H+ MMLV enzyme engineered to deliver uncompromised sensitivity and true representation of target RNA level, plus a potent blend of RNase inhibitors and our patented RT inhibition reducer that prevents RNA degradation and mispriming during reaction setup and reverse transcription to ensure optimal RT efficiency.

The reaction mix is 2x concentrated and optimized for SYBR® Green-based, one-step real-time PCR on any real-time PCR instrument (ROX-independent and ROX-dependent). It contains antibody-mediated hot-start iTaq DNA polymerase, dNTPs, Mg^{++} , enhancers, stabilizers, and a blend of passive reference dyes (including ROX).

Instrument Compatibility

This supermix is compatible with all Bio-Rad and ROX-dependent Applied Biosystems real-time PCR instruments, and with the Roche LightCycler LC480, Qiagen Rotor-Gene Q, Eppendorf Mastercycler ep realplex, and Stratagene Mx real-time PCR systems.

Reaction Mix Preparation and Thermal Cycling Protocol

1. Thaw iTaq universal SYBR® Green reaction mix and other frozen reaction components to 4°C . Mix thoroughly, centrifuge briefly to collect solutions at the bottom of tubes, and then store on ice protected from light.
2. Prepare on ice enough reaction setup for all reactions by adding all required components *except* RNA according to the following recommendations (Table 1).

Component	Volume per 20 µl Reaction	Volume per 10 µl Reaction	Final Concentration
iTaq universal SYBR® Green reaction mix (2x)	10 µl	5 µl	1x
iScript reverse transcriptase	0.25 µl	0.125 µl	1x
Forward and reverse primers	Variable	Variable	300 nM** each
RNA (add at step 4)	Variable	Variable	RNA: Up to 500 ng
Nuclease-free H ₂ O	Variable	Variable	—
<i>Total reaction mix volume</i>	<i>20 µl</i>	<i>10 µl</i>	—

* Scale all components proportionally according to sample number and reaction volumes.

** To validate the optimal primer concentration, perform a primer matrix to determine final primer concentration.

3. Mix the reaction setup thoroughly to ensure homogeneity and dispense equal aliquots into each PCR tube or into the wells of a PCR plate. Use good pipetting practice to ensure assay precision and accuracy.
4. Add RNA (and nuclease-free H₂O, if needed) to the PCR tubes or wells containing the reaction setup (Table 1), seal tubes or wells with flat caps or optically transparent film, and gently vortex to ensure thorough mixing of the reaction components. Spin the tubes or plate to remove any air bubbles and collect the reaction mixture in the vessel bottom.
5. Program thermal cycling protocol on the real-time PCR instrument according to Table 2.

Table 2. Thermal Cycling Protocol							
Real-Time PCR System	Setting/ Scan Mode	Reverse Transcription Reaction	Polymerase Activation and DNA Denaturation	Amplification			Melt-Curve Analysis
				Denaturation at 95°C	Annealing/ Extension + Plate Read ** at 60°C	Cycles	
Bio-Rad® CFX96™, CFX384™, CFX96 Touch™, CFX384 Touch™, CFX Connect™ systems	SYBR®/FAM	10 min at 50°C	1 min at 95°C	10 sec	10–30 sec	35–40	65–95°C 0.5°C increment 2–5 sec/step (or use instrument default setting)
Bio-Rad® iQ™ 5, MiniOpticon™, Chromo4™, MyiQ™	Standard				15–30 sec		
AB 7500, StepOne, StepOnePlus, 7900HT, ViiA7	Standard			60 sec			
Roche LightCycler 480	Fast			15 Sec	10–30 sec		
	Standard				60 sec		
Qiagen Rotor-Gene and Stratagene Mx series	Fast			30 sec			

** Shorter annealing/extension times (1–10 sec) can be used for amplicons <100 bp. Longer annealing/extension times (30–60 sec or more) can be used for amplicons >250 bp, GC- or AT- rich targets, and crude samples, or for higher input amounts (for example, >100 ng of RNA).

- Load the PCR tubes or plate onto the real-time PCR instrument and start the RT-qPCR run.
- Perform data analysis according to the instrument-specific instructions.

Recommendations for Assay Design and Optimization

- For best qPCR efficiency, design assays targeting an amplicon size of 70–150 bp
- The iTaq universal SYBR® Green one-step kit cycling protocols have been optimized for assays with a primer melting temperature (T_m) of 60°C that were designed using the open source Primer3, Primer3Plus, or Primer-BLAST programs under default settings. If primers are designed using other programs, adjust the temperature accordingly

Quality Control

iTaq universal SYBR® Green one-step kit demonstrates high RT-qPCR efficiency and linear resolution over a wide linear dynamic range. Stringent specifications are maintained to ensure lot-to-lot consistency. This product is free of detectable DNase and RNase activities.

Related Products

- Reverse transcription reagents for two-step real-time PCR:
 - iScript advanced cDNA synthesis kit for RT-qPCR (170-8842)
 - iScript reverse transcription supermix for RT-qPCR (170-8840)
- Real-time PCR supermixes for SYBR® Green-based qPCR:
 - SsoAdvanced™ universal SYBR® Green supermix (172-5270)
 - iTaq universal SYBR® Green supermix (172-5120)

To learn more about Bio-Rad's complete solution for amplification, visit www.bio-rad.com/amplification.

NOTICE TO PURCHASER: LIMITED LICENSE: Use of this product is covered by one or more of the following U.S. patents and corresponding patent claims outside the U.S.: 5,804,375; 5,538,848; 5,723,591; 5,876,930; 6,030,787; and 6,258,569. 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 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 from the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA

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

Bio-Rad 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.

StepOne and StepOnePlus are trademarks of Applied Biosystems. Mastercycler is a trademark of Eppendorf AG. Rotor-Gene is a trademark of the Qiagen Group. LightCycler is a trademark of Roche Diagnostics. Mx is a trademark of Stratagene Corporation. ViiA is a trademark of Life Technologies Corp.