

Product Information

SYBR® Green JumpStart™ Taq ReadyMix™ without MgCl₂

Catalog Number **S5193**

Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

SYBR Green JumpStart Taq ReadyMix without MgCl₂ combines JumpStart Taq Antibody for hot start^{1,2} PCR with SYBR Green I fluorescent dye, and provides the convenience of an easy-to-use ReadyMix solution. This mixture, provided as a 2× concentrate, is ideal for performing high-throughput, quantitative PCR, containing SYBR Green I dye, JumpStart Taq DNA polymerase, 99% pure deoxynucleotides, glass passivator, and reaction buffer. Simply add 25 μL of the 2× mix to DNA template, primers, the optimal amount of magnesium, and water. The JumpStart Taq antibody inactivates the DNA polymerase at room temperature. When the temperature is raised above 70 °C in the first denaturation step of the cycling process, the complex dissociates and the polymerase becomes fully active. There are no special preparations or protocol changes required to activate this hot start.

Features and Benefits

- The perfect ReadyMix for high throughput, quantitative PCR applications.
- SYBR Green I dye is ideal for quantifying any DNA sequence³. The dye binds to double-stranded DNA and detection is monitored by measuring the increase in fluorescence throughout cycling.
- The hot start mechanism, using JumpStart Taq antibody, prevents non-specific product formation and allows assembled PCR reactions to be placed at room temperature up to 2 hours without compromising performance.
- Magnesium chloride is provided separately for reaction optimization.
- An Internal Reference Dye is provided for reaction normalization. Maximum excitation of this dye is 586 nm and maximum emission is 605 nm.
- When performing large numbers of PCR reactions, the ReadyMix can save a significant amount of preparation time, reduce the risk of contamination from multiple pipetting steps, and provide consistent batch-to-batch and reaction-to-reaction performance.

Reagents

Sufficient for 100 or 400 PCR reactions (50 μL reaction volume)

- SYBR Green JumpStart Taq ReadyMix, Catalog Number S3443; 20 mM Tris-HCl, pH 8.3, 100 mM KCl, 0.4 mM each dATP, dCTP, dGTP, TTP, glass passivator, stabilizers, 0.05 unit/μL Taq DNA Polymerase, JumpStart Taq antibody, and SYBR Green I dye.
- Internal Reference Dye, Catalog Number R4526; 100x dye, provided in a 0.3 mL package size.
- Magnesium chloride solution, 25 mM, Catalog Number. M8787; 1 vial of 1.5 ml provided for 100 reactions, 1 vial of 5 ml provided for 400 reactions

Materials and reagents required but not provided

- Water, PCR Reagent, Catalog Number W1754
- Primers
- DNA template
- Thermal cycler for Quantitative PCR

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

Products can be stored at $-20\text{ }^{\circ}\text{C}$ for up to one year.

Preliminary Considerations

DNA Preparation

The single most important step in assuring success with PCR is high quality DNA preparation. Integrity and purity of DNA template is essential. Quantitative PCR involves multiple rounds of enzymatic reactions and is therefore more sensitive to impurities such as proteins, phenol/chloroform, salts, EDTA, and other chemical solvents. Contaminants can also interfere with fluorescence detection. The ratio of absorbance values at 260 nm and 280 nm gives an estimate of DNA purity. Pure DNA has an A_{260}/A_{280} ratio of 1.8-2.0. Lower ratios indicate the presence of contaminants such as proteins.

Primer Design

Specific primers for PCR should be designed with the aid of primer design software to eliminate the complications introduced with primer-dimers and secondary structures. Lower primer concentrations decrease the accumulation of primer-dimer formation and nonspecific product formation, which is critical in using SYBR Green I dye in quantitative PCR.

Magnesium Concentration

Lower magnesium chloride concentrations usually result in the formation of fewer nonspecific products. This ReadyMix solution contains no magnesium chloride. A separate tube of a 25 mM magnesium chloride solution is provided for optimization.

Internal Reference Dye

An internal reference dye is included for reaction normalization. Maximum excitation of this dye is 586 nm and maximum emission is 605 nm. Standard instrument settings for ROX reference dye are satisfactory for the measurement of the internal reference dye. This internal reference dye is necessary for ABI Sequence Detection Systems and does not interfere with the Roche LightCycler® SYBR Green quantitative PCR results.

Glass Passivator

A glass passivator is present in the solution due to the Roche LightCycler quantitative PCR reaction requirements. The glass passivator does not interfere with quantitative PCR performed on ABI Sequence Detection Systems.

Controls

A negative control containing no template is necessary to determine the presence of contamination. A positive control is helpful in accessing functionality of the kit components.

Data Analysis

Follow the recommendations of the real time instrument manufacturer to perform quantitative PCR using SYBR Green I dye. Generally, the log of relative fluorescence is plotted against the number of cycles to determine the threshold cycle (C_t) or crossing point. The C_t value is used to determine the amount of template in each sample. Consider the following points when determining the C_t :

- C_t is the first detectable increase in fluorescence due to PCR product formation
- Cycles before the C_t are the baseline cycles
- The threshold can be adjusted manually
- Threshold should always be set using a logarithmic amplification plot
- Threshold should be set in the most exponential phase of the reaction, not after reaching the plateau.

Melting Curves

Performing a melting curve analysis at the end of the run will help analyze only the PCR product. Follow the real time instrument manufacturer's instructions for melting curve analysis. After running a melting curve, any additional runs involving the same PCR product can be done with data collected in an additional detection step to eliminate primer-dimer and other misprimed product signal.

Methods of Quantification

Standard curves are necessary for both absolute and relative quantification. When generating standard curves, at least five different concentrations of the DNA should be amplified. Each concentration should be run in duplicate.

This product may be used to quantitate target DNA using either absolute or relative quantification. Absolute quantification determines the absolute amount of target DNA whereas relative quantification determines the ratio between the amount of target DNA and a reference amplicon. The reference DNA amplicon can be a housekeeping gene.

Absolute quantification uses external standards to determine the absolute amount of target nucleic acid. These external standards contain sequences that are the same as the target sequence or which vary only slightly from the target sequence. The primer binding sites of the external standards are always identical to the target sequence. The similarity between the external standard sequence and the target sequence is necessary for amplification efficiencies between the two to be essentially equivalent. Equivalent amplification efficiencies between the target and external standard are necessary for absolute quantification. A standard curve of external standard dilutions is generated and used to determine the concentrations of unknown target samples.

Relative quantification calculates the ratio between the amount of target template and a reference template in a sample. The relative amount of gene expression is a common application for relative quantification. The reference gene, usually a housekeeping gene, must not vary in concentration in different experimental conditions or tissue states for relative quantification to be possible. Amplification of the target and reference template dilutions in the sample should be performed in separate tubes. SYBR Green PCR quantification does not allow for multiplexing. If the reference template and the target template have different amplification efficiencies, then two standard curves need to be generated. The ratio of the resulting amounts of target and reference in the sample of interest can then be determined from these two standard curves. If the reference template and the target template have very similar amplification efficiencies, then only one standard curve for the reference template needs to be generated to determine the ratio of the amounts of target and reference in the sample.

Determination of PCR Reaction Efficiencies

The PCR efficiency between a reference sample and a target sample is determined by preparing a dilution series for each target. The C_t values from either the reference or target is then subtracted from the other. The difference in C_t values is then plotted against the log of the template amount. If the resulting slope of the straight line is less than 0.1, the amplification efficiencies are similar.

References

1. Dieffenbach, C. and Dveksler, G (Eds). *PCR Primer: A Laboratory Manual*, Cold Spring Harbor Laboratory Press, (Cold Spring Harbor Press, New York, 1995). (Catalog Number Z364118).
2. Kellogg, D. E., et al., *TaqStart Antibody: "hot start" PCR facilitated by a neutralizing monoclonal antibody directed against Taq DNA polymerase. *BioTechniques* 16, 1134-1137 (1994).*
3. Morrison, T. B., et al., Quantification of Low-Copy Transcripts by Continuous SYBR[®] Green I Monitoring during Amplification. *BioTechniques* 24, 954-962 (1998).
4. Sambrook, J, et al. *Molecular Cloning: A Laboratory Manual*, Second Edition, (Cold Spring Harbor Laboratory Press, New York, 1989). (Catalog Number M8265).
5. Rees, W. *et al.*, *Biochemistry*, 32, 137-144 (1993).

Procedure

Note: Because SYBR Green I binds to all double-stranded DNA, it is important to test primers and cycling conditions to insure that the PCR product is a single band, or the results will be uninterpretable. It is best to insure PCR specificity by checking the reaction on a normal (non-quantitative) thermocycler and analyzing the result using agarose gel separation.⁴

For best results, optimal concentrations of primers, MgCl₂, KCl and PCR adjuncts need to be determined. Testing various combinations of primer concentrations (50-1000 nM) is most efficient for primer optimization. If maximum sensitivity is not required and your PCR target is abundant, satisfactory results for SYBR Green based qPCR are often obtained with final concentrations of both primers 200-400 nM.

The following procedure serves as a guideline to establish optimal primer concentrations. Further optimization may be necessary due to primer specificity. For more optimization information, please read the qPCR user guide available online at www.sigmaaldrich.com.

Note: The use of up to 5% (v/v) dimethyl sulfoxide (DMSO) will not disturb the enzyme-antibody complex. Other co-solvents, solutes (salts) and extremes in pH or other reaction conditions may reduce the affinity of the JumpStart *Taq* antibody for the *Taq* polymerase and thereby compromise its effectiveness.

A. Optimizing Primer Concentrations

1. Prepare and dispense diluted primers (Fig 1).
 - a. Prepare 60 μ L of 8 μ M working solutions of both forward (fwd) and reverse (rev) primers in the first tubes of 2 separate 8-tube strips.
 - b. Dispense 30 μ L of water into tubes 2-5.
 - c. Transfer 30 μ L of the 8 μ M primer solution from tube 1 into tube 2. Mix thoroughly by pipetting up and down at least 5 times.
 - d. Repeat transfer and mixing from tube 2 to 3, 3 to 4, and 4 to 5.
 - e. Using a multichannel pipettor, transfer 5 μ L from the strip-tubes containing diluted fwd primer into the first 5 wells down columns 1-5 of a 96-well PCR plate. After adding fwd primer, PCR mix and template, final concentrations of fwd primer will be 1000, 500, 250, 125, 62.5 nM.

- f. Similarly transfer 5 μ L from the strip-tubes containing diluted rev primer into the first 5 wells across rows A-E. After adding PCR mix and template, final concentrations of rev primer will be 1000, 500, 250, 125 and 62.5 nM

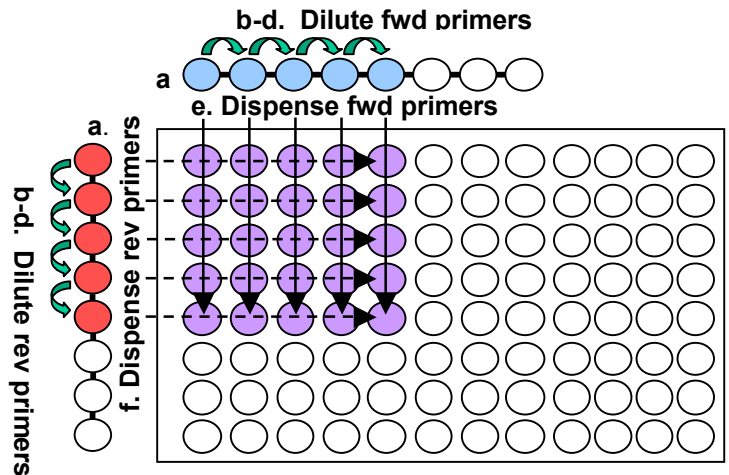


Fig 1: Follow steps 1a – 1f using diagram above

2. Prepare qPCR master mix:

Add reagents below in an appropriate sized DNase-free tube. Mix gently by vortexing and briefly centrifuge to collect all components at the bottom of the tube.

Volume	Reagent	Final Concentration
520 μ L	2 \times SYBR Green JumpStart <i>Taq</i> ReadyMix	1.25 units <i>Taq</i> DNA polymerase, 10 mM Tris-HCl, 50 mM KCl, 0.2 mM dNTP, stabilizers
(7 μ L)	Reference Dye* (optional)	1 \times
--- μ L	25 mM MgCl ₂ (optional)	
q.s. to 676 μ L	Water	
676 μ L	Total Volume	

*Use 0.1 \times for ABI 7500 and Stratagene instruments; replace with FITC for BioRad iCycler.

3. Aliquot 26 μL master mix into all wells in the PCR plate that contain primers (A1-E5)
4. Mix Thoroughly and transfer 18 μL from each of wells A1 through E5 to wells A8 through E12.
5. Add 2 μL template DNA (10-50 ng genomic DNA or 0.1-1 ng plasmid) to one set of reactions (columns 1-5) and 2 μL of water to the other columns (8-12).
6. Mix gently by vortexing and briefly centrifuge to collect all components at the bottom of the tube.
7. Perform Thermal cycling:

Optimal cycling parameters vary with primer design and thermal cycler. Consult your thermal cycler manual. It may be necessary to optimize the cycling parameters to achieve maximum product yield and/or quality.

Typical cycling parameters for 100 bp – 600 bp fragments:

This protocol has been successfully tested on the following thermal cyclers: Stratagene MX 3000P, BioRad iCycler, MJ Opticon and ABI 7700.

Initial denaturation	94 °C	2 min
40 cycles:		
Denaturation	94 °C	15 sec
Annealing, extension, and read fluorescence	60 °C or 5 °C below lowest primer T_M	1 min
(Optional) Hold	4 °C - only if products will be run out on a gel	

8. Evaluate fluorescence plots (ΔR_n) for reactions containing target nucleic acid (columns 1-5). Primer combinations with the lowest C_i and the highest fluorescence will give the most sensitive and reproducible assays.

B. Procedure for Routine Analysis

1. Preparation of a reaction master mix is highly recommended to give best reproducibility. Mix all reagents but template in a common mix, using ~10% more than needed. Once template is diluted into the reaction vessel, master mix is aliquoted into the proper tube or plate for thermocycling.

Volume*	Reagent	Final Concentration
25 μL	2x JumpStart <i>Taq</i> ReadyMix	1.25 units <i>Taq</i> DNA polymerase, 10 mM Tris-HCl, 50 mM KCl, 0.2 mM dNTP, stabilizers
(0.5 μL)	Reference Dye** (optional)	1x
--- μL	Forward Primer	Optimal Conc. from Sec. A
--- μL	Reverse Primer	Optimal Conc. from Sec. A
--- μL	25 mM MgCl_2 (optional)	
--- μL	Template DNA	10-100 ng
q.s. to 50 μL	Water	
50 μL	Total Volume	

* Volume for 50 μL reaction, however component volumes may be scaled to give the desired reaction volumes.

** Use 0.1x for ABI 7500 and Stratagene instruments; replace with FITC for BioRad iCycler.

2. Mix gently by vortexing and briefly centrifuge to collect all components at the bottom of the tube.
3. Perform Thermal cycling

Typical cycling parameters for 100 bp – 600 bp fragments:

Initial denaturation	94 °C	2 min
40 cycles:		
Denaturation	94 °C	15 sec
Annealing, extension, and read fluorescence	60 °C or 5 °C below lowest primer T_M	1 min
(Optional) Hold	4 °C - only if products will be run out on a gel	

Troubleshooting Guide

Symptom	Possible Cause	Solution
No PCR product (signal) is observed	A PCR primer is missing or degraded.	A positive control should always be run to insure components are functioning. A checklist is also recommended when assembling reactions.
	Too few cycles were performed.	Increase the number of cycles.
	The annealing temperature is too high.	Decrease the annealing temperature in increments of 2-4 °C.
	The primers are not designed optimally.	Confirm the accuracy of the sequence information. If the primers are less than 27 nucleotides long, try to lengthen the primer to 27-33 nucleotides. If the primer has a GC content of less than 45%, try to redesign the primer with a GC content of 45-60%.
	There is not enough template.	After increasing the number of cycles has shown no success, repeat the reaction with a 10-fold higher concentration of template.
	The template is of poor quality.	Evaluate the template integrity by agarose gel electrophoresis. It may be necessary to repurify the template using methods that minimize shearing and nicking.
	Target template is difficult	In most cases, inherently difficult targets are due to unusually high GC content and/or secondary structure. Betaine has been reported to help amplification of high GC content templates at a concentration of 0.8-1.3 M. ⁵
Signal is independent of template dilution (Multiple products or smeared products)	The annealing temperature is too low.	Increase the annealing temperature in increments of 2-3 °C.
	The primers are not designed optimally.	Confirm the accuracy of the sequence information. If the primers are less than 27 nucleotides long, try to lengthen the primers to 27-33 nucleotides. If the primer has a GC content of less than 45%, try to redesign the primers with a GC content of 45-60%.
	The template concentration is too high.	Reduce the concentration of the template in the PCR reaction.
	The primer concentration is too high.	Reduce primer concentration in a series of two-fold dilutions (i.e. 0.1 μM, 0.05 μM, 0.025 μM and 0.0125 μM) and subject these trial reactions to PCR.

NOTICE TO PURCHASER: LIMITED LICENSE

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,994,056 and 6,171,785.. 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 (such as apparatus or system claims in US Patent No. 6,814,934) 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 by contacting the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

SYBR is a registered trademark of Molecular Probes, Inc. LightCycler is a trademark of a member of the Roche Group. JumpStart and ReadyMix are trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co/

AH,RS,PHC 09/10-1