

## User Guide

# JumpStart™ REDAccuTaq® LA DNA Polymerase

Long and accurate hot-start *Taq* with inert dye, 10X buffer included

**D1313**

## Product Description

JumpStart™ REDAccuTaq® LA DNA Polymerase is a combination of AccuTaq™ LA DNA Polymerase, the *Taq* neutralizing JumpStart™ *Taq* antibody, and an inert dye.

The AccuTaq™ LA DNA Polymerase can perform 3'→5' exonuclease activity necessary for proofreading and thus has the ability to amplify genomic targets ≥ 20 kb. The JumpStart™ *Taq* antibody binds to *Taq* DNA polymerase, inactivating it until the complex is dissociated at ≥70 °C, providing a simple and efficient hot start PCR. Because JumpStart™ REDAccuTaq® DNA polymerase is inactive at room temperature, reaction mixtures can be prepared on the bench and transported directly to the thermocycler as the enzyme will become active in the first denaturation step of the cycling process. The red tracer provides quick recognition of reactions to which the enzyme has been added as well as visual confirmation of complete mixing. It also serves as a tracking dye, co-migrating at the same rate as a 125 bp fragment in a 1% agarose gel.

The enzyme is supplied at 1 unit/μL and comes with an optimized 10X reaction buffer.

## Features

- Allows for room-temperature PCR set up
- Amplifies 0.25 to 22 kb for complex genomic DNA and up to 40 kb for less complex templates
- 3'-5' exonuclease proofreading activity provides higher fidelity (up to 6.5x) than standard *Taq* DNA polymerase
- Ready for direct gel loading after PCR
- Inert dye does not interfere with DNA polymerase, ligase, restriction enzymes, or transformation

## Applications

JumpStart™ REDAccuTaq® can be used for high-fidelity amplification of:

- Genomic DNA
- cDNA
- Multiple targets (multiplex PCR)
- Long amplicons

## Unit Definition

One unit incorporates 10 nmol of total deoxyribonucleoside triphosphates into acid precipitable DNA in 30 minutes at 74 °C.

## Reagents Provided

- JumpStart™ REDAccuTaq® LA DNA Polymerase - 1 unit/μL AccuTaq™ LA DNA polymerase with JumpStart™ *Taq* antibody in 20 mM Tris-HCl, pH 8.0, 37 mM KCl, 0.1 mM EDTA, 7 mM DTT, 0.1% TWEEN® 20, stabilizers, 50% glycerol, inert dye (D1938)
- AccuTaq™ LA 10× Buffer, 500 mM Tris-HCl, 150 mM ammonium sulfate (pH 9.3, adjusted with NH<sub>4</sub>OH), 25 mM MgCl<sub>2</sub>, 1% TWEEN® 20 (B0174)

## Materials and Reagents Required

(Not included)

- Deoxynucleotide (dNTP) Mix, containing 10 mM each of dATP, dCTP, dGTP, and dTTP sodium salts
- Nuclease-free water
- Custom ordered primers specific to gene target
- PCR tubes or plates
- Sample containing template DNA
- Thermal cycler

## Precautions and Disclaimer

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

## Storage/Stability

Store JumpStart™ REDAccuTaq® LA DNA Polymerase and 10X Buffer at -20 °C.

## Directions for Use

For best reproducibility, assemble a master mix of PCR reagents by multiplying the number of reactions needed (plus 10% to account for pipetting error) by the suggested volumes in the table below. Aliquot reaction mixture into PCR tubes.



### Assemble Reaction Mix

Reagent	Final Concentration	Amount per 20 µL reaction
AccuTaq™ LA 10X buffer	1X	2 µL
JumpStart™ REDAccuTaq® LA DNA Polymerase (1 U/µL)	0.05 U/µL	1 µL
dNTP Mix, 10 mM	200 µM	0.4 µL
Primers	0.1–0.5 µM	Variable
Template	Variable	Variable
Nuclease-free Water	-	To 20 µL total

**Note:** JumpStart™ REDAccuTaq® LA is a magnesium ion-dependent enzyme, optimal concentrations of template DNA, primers, and MgCl<sub>2</sub> will be target-specific. The supplied AccuTaq™ LA 10X Buffer contains 25 mM MgCl<sub>2</sub>, for a final concentration of 2.5 mM. The recommended MgCl<sub>2</sub> concentration range for endpoint PCR is 1.5–3.5 mM, thus the reaction can be supplemented accordingly if necessary (see [Product Ordering](#)).



### Add Template

Recommended input template is 80–200 ng gDNA. Less DNA is needed for less complex templates, such as lambda DNA.

A suggested thermocycling protocol using JumpStart™ REDAccuTaq® is provided below:



### Amplify

Initial denaturation		94 °C	2 min
40 cycles	Denaturation	94 °C	15 sec
	Annealing	60 °C or 5 °C below lowest primer T <sub>m</sub>	30 sec
	Extension	72 °C	1 min/kb
Final extension		72 °C	1 min
Hold		4 °C	∞

Amplification parameters will vary depending on primers, template, and instrument used. For tips on optimizing PCR conditions as well as a 2-step cycling protocol please see [Technical Guide](#).



### Evaluate

Amplified DNA can be evaluated by directly loading 8–10 µL of PCR product to a 0.8–1% agarose gel for electrophoresis and subsequent stain with nucleic acid dye.

**Note:** A minimum of 0.6 units of JumpStart™ REDAccuTaq® LA DNA polymerase must be added per 20 µL reaction to ensure enough glycerol is present for direct gel loading.

## Technical Guide

### Considerations for Primer Design

Thoughtful primer design is essential for PCR efficiency and specificity. For successful amplification consider the following:<sup>1</sup>

- Select an 18-30 nucleotide-long sequence with 40-60% G/C content and even distribution of all 4 bases.
- Avoid inverted repeat or self-complementary sequences >3 bp.
- Primer pairs should not differ in length by >3 bp and should not contain complementarity to one another.
- Maintain calculated primer  $T_m$  between 55-60 °C, permitting only 2-3 °C variation between primer pairs.
- Priming efficiency can be increased by including a terminal G at the 3' end; however, the number of Gs or Cs in the last 5 bases of the primer sequence should be no more than 3.
- Ensure each primer sequence is unique to the gene of interest and is absent in other genes in the gDNA sample or within the vector.

### Optimization of PCR Conditions

PCR involves the cycling of denaturing, annealing, and extension steps for DNA synthesis by a polymerase enzyme. To obtain the best product yield and accuracy, each step must be optimized.

- The **denaturing** step (94-96 °C) activates the JumpStart™ REDAccuTaq® DNA polymerase and separates double-stranded DNA strands, making it accessible to primers.<sup>1</sup> The duration of this step should be long enough to denature DNA but not so long that it compromises Taq DNA polymerase integrity.<sup>1</sup> High salt conditions, GC-rich (>55%) templates, and gDNA templates may require longer denaturation times and/or higher temperatures. For maximum retention of JumpStart™ REDAccuTaq® activity during thermocycling, use 94 °C for denaturation.
- The **annealing** temperature can be calculated by subtracting 5 °C from the lowest reaction primer  $T_m$ . The annealing time should be long enough for the primer to anneal to the template but not too long for non-specific annealing to occur.<sup>1</sup>
- The optimal **extension** temperature for REDAccuTaq® DNA polymerase is 72 °C; however, lower temperatures may be used for some reactions. Extension time depends on length and complexity of the target sequence. For complex templates, use 1 minute/kb, with 15 seconds added if the PCR product is >2 kb. Short or non-complex templates may be amplified with extension times of 30 seconds/kb.

- The **number of cycles** needed for amplification depends on the amount of template input, with higher amount of input requiring less cycling.<sup>1</sup> Generally, 25-30 cycles are sufficient to produce detectable product; however, low concentration templates may require up to 45 cycles.
- To maintain **enzyme fidelity**, or accuracy of nucleotide incorporation, limit the number of PCR cycles and use an equimolar concentration of each dNTP.<sup>1</sup> Magnesium ion concentration also affects enzyme fidelity at a concentration that varies by reaction conditions and should thus be optimized.<sup>1</sup>

### Handling gDNA Templates

To prevent genomic DNA (gDNA) shearing, add template last and mix gently using a wide pore pipet tip. DO NOT VORTEX!

### Multiplex PCR

When performing multiplex PCR, competition between products for reagents may occur. Consider adjusting the following for optimization:<sup>2</sup>

- Proportion of primer pair concentration: if a target sequence produces a relatively “weaker” signal, the amount of primer used may be increased to compensate. For sequences with low copy numbers, or high-complexity, primer concentration can be used at 0.3-0.5 µM.
- Primer concentration can also be decreased for target sequences producing “stronger” signal to achieve balance. For high copy number or low-complexity sequences, primer concentration can be used at 0.04-0.4 µM.
- dNTP: Perform a stepwise increase of dNTP to a concentration ≤ 400 µM. Keep MgCl<sub>2</sub> concentration constant for this optimization.
- PCR Buffer: Use 2X buffer concentration for the reaction instead of 1X.
- Supplement the reaction with additional MgCl<sub>2</sub>, or PCR-enhancing additives.

## PCR-Enhancing Additives

When optimizing PCR conditions for a new experiment, the following can be added to the reaction mix individually. After performing PCR amplification, samples with and without additive can be compared using agarose gel electrophoresis or other standard methods to look for improved product specificity and yield.

Additive	Purpose
Bovine Serum Albumin (10-100 µg/mL)	<i>Taq</i> DNA polymerase stabilization <sup>3</sup>
Formamide (1.25-10%)	Increases specificity in G/C rich regions <sup>4</sup>
DMSO (Up to 5%)	Accelerates strand renaturation <sup>5</sup>
	Nucleic acid thermal stability against depurination <sup>5</sup>
Glycerol (Up to 10%)	Increases thermal stability of the polymerase and lowers the temperature necessary for strand separation <sup>5</sup>
Ammonium sulfate (15-30 mM)	Affects the denaturing and annealing temperatures of the DNA <sup>6</sup>
Single strand binding protein (0.7-1.5 µg)	Inhibits formation of secondary structures, improving fidelity and <i>Taq</i> processivity <sup>7</sup>
Betaine (0.8-1.6 M)	Reduces base pair composition dependence of DNA melting <sup>8</sup>

## Two-Step PCR Amplification

Application of a two-step PCR process is possible when the annealing and extension temperatures are similar.

Initial denaturation	94 °C	2 min
40 Cycles	Denaturation	94 °C 3 sec
	Annealing/extension	60 °C* 15-30 sec
Final extension	72 °C	1 min
Hold	4 °C	∞

\*Consult primer  $T_m$  regarding temperature selection. Extension time is target-dependent, with larger targets requiring more than the recommended time.

## Product Ordering

Description	Catalog Number
1 kb DNA Ladder (0.5-10 Kb)	D0428
Betaine solution	B0300
Bovine Serum Albumin solution	B8667
DMSO	D8418
Single strand binding protein	S3917
Mineral Oil	M5904
Magnesium chloride solution	M8787
REDAccu <i>Taq</i> ® LA DNA Polymerase	D4812
GenElute™ Bacterial Genomic DNA Kit	NA2120
GenElute™ Gel Extraction Kit	NA1111
GenElute™ Mammalian Genomic DNA Miniprep Kit	G1N10
GenElute™ PCR Clean-Up Kit	NA1020
GenElute™ Plant Genomic DNA Miniprep Kit	G2N70
GenElute™-E Single Spin DNA Cleanup Kit	EC600
	P5722
	P5972
Precast Agarose Gels	P6097
	P5472
	P6222
Water, Microbial DNA-free	MBD0025
Nuclease-Free Water, for Molecular Biology	W4502

## Troubleshooting Guide

Problem	Suggestions
No or low product amplification	<ul style="list-style-type: none"> <li>• Titrate <math>MgCl_2</math> concentration in 0.5 mM increments using molecular grade <math>MgCl_2</math> (see <a href="#">Product Ordering</a>). Each amplicon target must be optimized individually.</li> <li>• Adjust the annealing temperature in 2-3 °C increments or use a gradient PCR to find the optimal annealing temperature.</li> <li>• Increase the number of amplification cycles. If currently using 25-30 cycles, increase the cycle number to 35-40.</li> <li>• For complex templates like human genomic DNA, increase the initial denaturation time by 1-2 minutes and/or increase the denaturation temperature to 95 °C to overcome denaturation difficulties.</li> <li>• Check concentration of input template. For complex templates like intact eukaryotic genomic DNA, 1000 genome copies may be required for amplification of difficult targets. For highly concentrated templates, such as purified plasmid, consider diluting 1:1000 to improve amplification.</li> <li>• Assess DNA quality to ensure absence of PCR inhibitors in sample. If presence of inhibitors is suspected, DNA can be diluted 1:10-1:100. Alternatively, lysis and DNA purification can be performed using the GenElute™ genomic DNA miniprep kits.</li> <li>• Refer to "<a href="#">PCR-Enhancing Additives</a>" section of the <a href="#">Technical Guide</a> to improve amplification.</li> <li>• If yield is too low for downstream applications, increase the reaction volume to 50-75 µL.</li> </ul>
Amplification of nonspecific product(s)	<ul style="list-style-type: none"> <li>• Raise the annealing temperature in 2-3 °C increments or use a gradient PCR to find the optimal annealing temperature. Raising the temperature improves the specificity of binding by the primers; however, it may also result in reduced binding and extension of the primers.<sup>1</sup> If raising the annealing temperature causes reduced yield of the specific product without eliminating side reaction products, it may be necessary to redesign the primers to improve specificity.</li> <li>• Take precautions to avoid crossover contamination of PCR with both specific and nonspecific PCR products, including primer-dimer artifacts.<sup>9</sup></li> <li>• The use of more than 5% v/v DMSO with JumpStart™ REDAccuTaq® is not recommended as it may interfere with the enzyme-antibody complex. Other co-solvents, salts, and extremes in pH can also reduce the affinity of the JumpStart™ Taq antibody for the AccuTaq™ LA DNA Polymerase and compromise its effectiveness for hot start PCR.</li> </ul>

## References

1. Green MR, Sambrook J., Cold Spring Harb Protoc. 6, 436-456(2019).
2. Markoulatos, P., et al., J. Clin. Lab. Anal. 16, 47-51 (2002).
3. Paabo, S., et al., Nucleic Acids Res., 16, 9775-9787 (1988).
4. Sarkar, G., et al., Nucleic Acids Res., 18, 7465 (1990).
5. Cheng, S., et al., Proc. Natl. Acad. Sci. USA, 91, 5695-5699 (1994).
6. Grunenwald H. PCR Protocols. Methods in Molecular Biology™. 226, 89-99 (2003).
7. Chou, Q., Nucleic Acids Res., 20, 4371 (1992).
8. Rees, W. A., et al., Biochemistry, 32, 137-144 (1993).
9. Kwok, S., and Higuchi, R., Nature 339, 237-238 (1989).

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

### Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

### Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

### Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck, JumpStart, REDAccuTaq, AccuTaq, GenElute and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2022 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

D1313ug Rev 11/22

