

Cat. # RR070A

For Research Use

TAKARA

**SpeedSTAR™ HS
DNA Polymerase**

Product Manual

v201704Da

Table of Contents

I.	Description.....	3
II.	Components	3
III.	Storage	3
IV.	Supplied buffers	3
V.	General reaction mixture	4
VI.	Thermal cycling conditions.....	4
VII.	Optimization of parameters.....	4
VIII.	Experimental examples	5
IX.	Amplified products.....	7
X.	Troubleshooting.....	8

I. Description

SpeedSTAR HS DNA Polymerase is designed for high speed PCR and includes two optimized buffers, Fast Buffer I and II. SpeedSTAR HS DNA Polymerase is compatible with an extension time as fast as 10 sec/kb, compared to general PCR enzymes that typically require 1 min/kb. SpeedSTAR HS DNA polymerase utilizes a monoclonal antibody-mediated hot start formulation, thereby preventing non-specific amplification due to mispriming and/or formation of primer dimers during reaction assembly. This enzyme can be used with routine PCR reaction conditions since the monoclonal antibody is denatured in the initial DNA-denaturation step.

II. Components (for 200 reactions ; 50 μ l reactions)

SpeedSTAR HS DNA Polymerase (5 units/ μ l)*1	50 μ l
10X Fast Buffer I (Mg ²⁺ plus)*2	1 ml
10X Fast Buffer II (Mg ²⁺ plus)*2	1 ml
dNTP Mixture (2.5 mM each)	800 μ l

*1 【 Storage Buffer 】

20 mM	Tris-HCl (pH 8.0)
100 mM	KCl
0.1 mM	EDTA
1 mM	DTT
0.5%	Tween 20
0.5%	Nonidet P-40
50%	Glycerol

【 Unit definition 】

One unit is the amount of enzyme that will incorporate 10 nmol of dNTP into acid-insoluble products in 30 minutes at 74°C with activated salmon sperm DNA as the template-primer.

*2 Mg²⁺ concentration: 10X Fast Buffer I, 30 mM ; 10X Fast Buffer II, 20 mM

III. Storage — 20°C

IV. Supplied buffers

Two Fast Buffers are supplied with this product. Select the appropriate buffer depending on the target amplification size. For amplicons up to 2 kb, Fast Buffer I should be used. Either Fast Buffer I or II can be used for amplicons between 2 - 4 kb. For amplification of targets over 4 kb, Fast Buffer II is preferred. The use of Fast Buffer I for longer amplification or Fast Buffer II for shorter amplification can result in lower amplification efficiency.

V. General reaction mixture (for 50 μ l reactions)

Reagent	Volume/Amount	Final Conc.
SpeedSTAR HS DNA Polymerase (5 units/ μ l)	0.25 μ l	1.25 units/50 μ l
dNTP Mixture (2.5 mM each)	4 μ l	200 μ M
Primer 1	10 - 50 pmol	0.2 - 1 μ M
Primer 2	10 - 50 pmol	0.2 - 1 μ M
Template	< 500 ng	
10X Fast Buffer I or II	5 μ l	1X
Sterile purified water	up to 50 μ l	

NOTE: The reaction mixture can be prepared at room temperature. Keep reaction components on ice while assembling the reaction mixture.

VI. Thermal cycling conditions

SpeedSTAR polymerase can be used for both 2 step and 3 step PCR. For fast PCR, try the 2 step PCR method first. When using short primers, the 3 step PCR method is recommended. For long amplification, we recommend using a longer extension time. In some cases, using SpeedSTAR polymerase for long amplifications may result in smearing of the amplified products.

(A) 2 step PCR

- Amplification up to 4 kb or 6 kb (with Fast Buffer I or II)

95°C	5 sec] 30 cycles
65°C	10 sec (or up to 20 sec)/kb	

- Amplification of 6 kb or longer (with Fast Buffer II)

98°C	5 sec] 30 cycles
68°C	10 sec (or up to 20 sec)/kb	

NOTE: Efficient amplification can be achieved by optimizing the temperature of each step, depending on the amplicon size.

(B) 3 step PCR (with Fast Buffer I or II)

98°C	5 sec] 30 cycles
55°C	10 - 15 sec	
72°C	5 - 10 sec/kb	

NOTE: Denaturation conditions vary depending on the thermal cyler and tubes used for PCR. We recommend 5 - 10 sec at 98°C or 20 - 30 sec at 95°C.

VII. Optimization of parameters

The following parameters should be optimized for best performance of SpeedSTAR HS DNA polymerase.

1) Enzyme amount:

Use 1.25 units for a 50 μ l PCR. This amount may be adjusted slightly depending on the template amount or amplicon size. Excess enzyme can result in non-specific amplification causing smeared bands on a gel; too little enzyme can lower the amplification efficiency.

2) Template DNA:
Excess template can result in non-specific amplification or smeared bands. Refer to the following recommendations for the amount of template for a 50 μl reaction.

Human genomic DNA	5 ng - 500 ng
<i>E. coli</i> genomic DNA	50 pg - 100 ng
Plasmid DNA	10 pg - 1 ng

3) Concentration of dNTPs and Mg²⁺:
The supplied Fast Buffer I includes Mg²⁺ at a final concentration of 3 mM, and Fast Buffer II includes Mg²⁺ at a final concentration of 2 mM. The best results will be obtained when dNTPs are used at a final concentration of 200 μM each.

4) Primers:
Design primers using a commercially available primer design software (e.g. OLIGO Primer Analysis Software, Molecular Biology Insights)

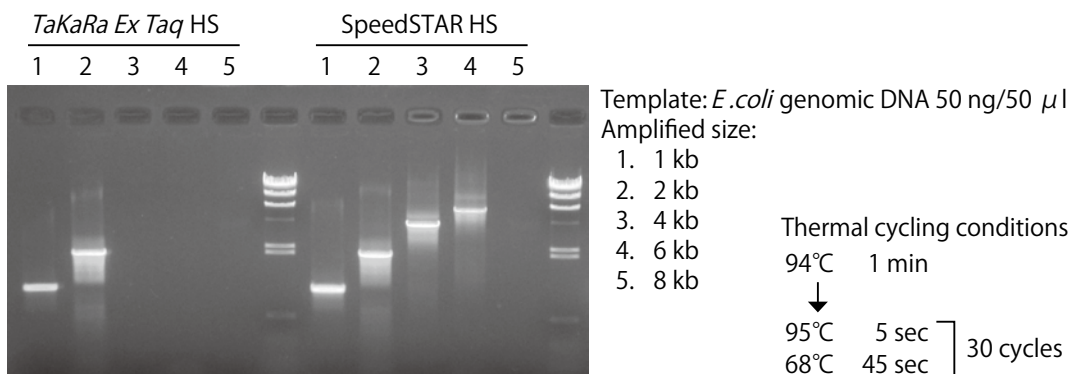
- Primer length
In general, 20 - 25 -mer primers provide good results. For longer amplifications, 25 - 30 -mer primers may be better.
- GC content
Primer GC content should be 40 - 60%. GC residues should be evenly distributed throughout each primer. The 3' termini should not be GC-rich.
- Tm value
Tm values of the upstream and downstream primers should be the same.
- Primer concentration
The concentration should be empirically determined within the range of 0.2 μM - 1.0 μM. Low primer concentration can result in lower yield of amplified products. In contrast, high primer concentration can cause non-specific amplification which can inhibit specific amplification.

VIII. Experimental examples

< Fast PCR >

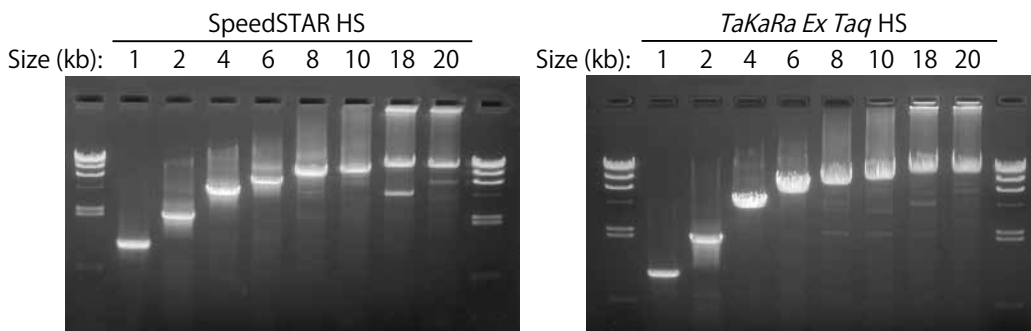
Comparison of SpeedSTAR HS DNA Polymerase and *TaKaRa Ex Taq*® Hot Start Version: Amplification of products of various sizes was compared for the enzymes using 2 step PCR with extension time of 45 sec. Under these conditions, SpeedSTAR HS was able to amplify fragments up to 6 kb, while *TaKaRa Ex Taq* HS was only able to amplify products up to 2 kb.*

* *TaKaRa Ex Taq* HS can amplify 6 kb products when optimized conditions are used.



< **High Performance** >

Comparison of reaction time between SpeedSTAR polymerase and *TaKaRa Ex Taq* HS with varying amplification target sizes:



Thermal cycling conditions

**Amplification of 1 kb, 2 kb
(with Fast Buffer I)**

94°C 1 min
↓
95°C 5 sec
65°C 20 sec] 30 cycles

Total reaction time: approx. 33 min

**Amplification of 4 kb, 6 kb
(with Fast Buffer II)**

94°C 1 min
↓
95°C 5 sec
65°C 60 sec] 30 cycles

Total reaction time: approx. 53 min

**Amplification of 8 kb, 10 kb
(with Fast Buffer II)**

94°C 1 min
↓
98°C 5 sec
68°C 2 min] 30 cycles

Total reaction time: approx. 83 min

**Amplification of 18 kb, 20 kb
(with Fast Buffer II)**

94°C 1 min
↓
98°C 5 sec
68°C 5 min] 35 cycles
↓
72°C 5 min

Total reaction time: approx. 3 hr 29 min

Amplification of 1 kb, 2 kb

94°C 1 min
↓
98°C 10 sec
68°C 2 min] 30 cycles

Total reaction time: approx. 96 min

Amplification of 4 kb, 6 kb

94°C 1 min
↓
98°C 10 sec
68°C 6 min] 30 cycles
↓
72°C 10 min

Total reaction time: approx. 3 hr 46 min

Amplification of 8 kb, 10 kb

94°C 1 min
↓
98°C 10 sec
68°C 10 min] 30 cycles
↓
72°C 10 min

Total reaction time: approx. 5 hr 46 min

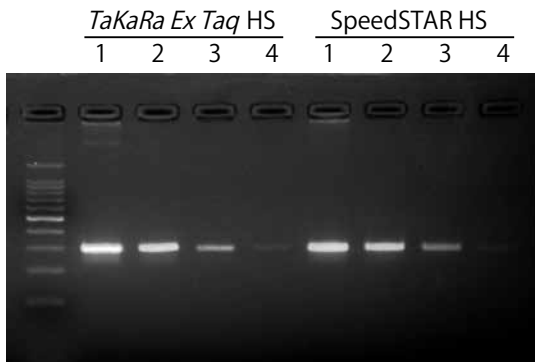
Amplification of 18 kb, 20 kb

94°C 1 min
↓
98°C 10 sec
68°C 15 min] 30 cycles
↓
72°C 10 min

Total reaction time: approx. 8 hr 16 min

< **Fast PCR with high sensitivity and performance** >

Comparison of detection sensitivity between SpeedSTAR polymerase and *TaKaRa Ex Taq* HS with varying amplification product sizes:



Amplified size: 300 bp
 Template: human genomic DNA
 1: 100 ng/50 µl PCR
 2: 10 ng/50 µl PCR
 3: 1 ng/50 µl PCR
 4: 0.1 ng/50 µl PCR

Thermal cycling conditions

***TaKaRa Ex Taq* HS**

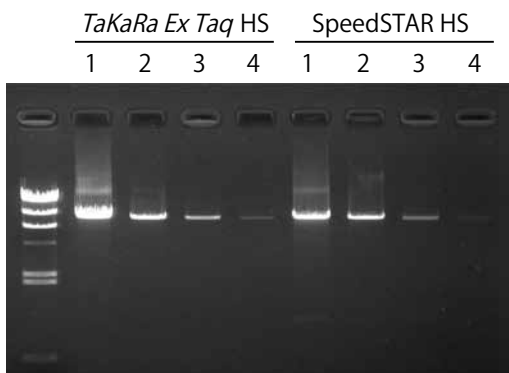
98°C 5 sec
 55°C 30 sec
 72°C 30 sec } 30 cycles

Total reaction time: approx. 67 min

SpeedSTAR HS (with Fast Buffer I)

98°C 5 sec
 55°C 10 sec
 72°C 5 sec } 30 cycles

Total reaction time: approx. 39 min



Amplified size: 8.5 kb
 Template: human genomic DNA
 1: 100 ng/50 µl PCR
 2: 10 ng/50 µl PCR
 3: 1 ng/50 µl PCR
 4: 0.1 ng/50 µl PCR

Thermal cycling conditions

***TaKaRa Ex Taq* HS**

94°C 1 min
 ↓
 98°C 5 sec
 68°C 8.5 min } 30 cycles
 ↓
 72°C 10 min

Total reaction time: approx. 4 hr 59 min

SpeedSTAR HS (with Fast Buffer II)

94°C 1 min
 ↓
 98°C 5 sec
 68°C 2 min } 35 cycles
 ↓
 72°C 3 min

Total reaction time: approx. 1 hr 40 min

IX. Amplified products

Most PCR products amplified with SpeedSTAR HS DNA Polymerase have a single A overhang at the 3'-termini. Therefore, PCR products can be directly used for cloning into a T-vector. It is also possible to clone the product using blunt-end vectors after blunting and phosphorylation of the ends.

X. Troubleshooting

Observation	Possible Cause	Solution
No amplification or low yield	Extension time	Set the extension time to 20 sec/kb.
	Annealing temperature	Lower the temperature in decrements of 2°C. Perform 3 step PCR.
	Template DNA	Repurify template DNA. For long amplification, use intact, undegraded DNA
	Primer	Re-design primers. Increase the primer amount.
Extra bands Smear bands	Extension time	Excess extension time affects the reaction. Use the following guidelines: 2 step PCR: 10 - 20 sec/kb 3 step PCR: 5 - 10 sec/kb
	Annealing temperature	Raise the temperature in increments of 2°C. Perform 2 step PCR.
	Template DNA	Use an appropriate amount of DNA. Excess template DNA affects the reaction.
	Primer	Reduce the primer amount.

TaKaRa Ex Taq is a registered trademark of TAKARA BIO INC.
SpeedSTAR is a trademark of TAKARA BIO INC.

NOTE: This product is for research use only. It is not intended for use in therapeutic or diagnostic procedures for humans or animals. Also, do not use this product as food, cosmetic, or household item, etc.

Takara products may not be resold or transferred, modified for resale or transfer, or used to manufacture commercial products without written approval from TAKARA BIO INC.

If you require licenses for other use, please contact us by phone at +81 77 565 6973 or from our website at www.takara-bio.com.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product web page. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

All trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions.