

# 2X PCR Solution

## PrimeSTAR® HS (Premix)

Code No. R040A  
Size: 500  $\mu$ l x 5  
(for 100 PCR reactions)

Shipping at  $-20^{\circ}\text{C}$   
Store at  $-20^{\circ}\text{C}$

**Storage: Excessive freezing and thawing may decrease enzyme activity.**

Lot No.

Expiration Date:

For a detailed protocol, refer to the product manual supplied with this product.

### Description:

PrimeSTAR HS (Premix) is a 2X PCR enzyme mixture composed of PrimeSTAR HS DNA Polymerase, a high-fidelity DNA polymerase, reaction buffer, and dNTP mixture. The premix formulation allows for quick preparation of reactions and reduces the risk of contamination. This product is also useful for high-throughput applications.

PrimeSTAR HS DNA Polymerase is a high-fidelity enzyme that has a very strong 3'-5' exonuclease activity. It also has a higher amplification efficiency than standard *Taq* DNA Polymerase.

### Composition:

PrimeSTAR HS DNA Polymerase\* : 1.25 units/25  $\mu$ l  
dNTP Mixture : 2X conc. ; ea. 0.4 mM  
PrimeSTAR Buffer : 2X conc. ; including 2 mM  $\text{Mg}^{2+}$

### \* : Specification of PrimeSTAR HS DNA Polymerase (Cat. #R010A)

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

### Reaction mixture for unit definition:

100 mM	Tris-HCl (pH8.3 at 37°C)
10 mM	KCl
2 mM	$\text{MgCl}_2$
6 mM	$(\text{NH}_4)_2\text{SO}_4$
0.1%	Triton X-100
200 $\mu$ M	each dATP, dGTP, dCTP
100 $\mu$ M	$[\text{^3H}]$ TTP
0.001%	BSA
0.4 mg/ml	activated salmon sperm DNA

**Purity:** Nicking, endonuclease, and exonuclease activity were not detected after incubation of 0.6  $\mu$ g of supercoiled pBR322 DNA, or 0.6  $\mu$ g of  $\lambda$ -*Hind* III digest with 10 units of this enzyme for 1 hour at 74°C.

### Applications:

High fidelity DNA amplification by hot start PCR

### PCR product:

A significant percentage of PCR products obtained using PrimeSTAR HS DNA polymerase will have blunt-ends. Thus, the obtained PCR products can be directly cloned into blunt-end vectors. If necessary, phosphorylate PCR products before cloning.

### PCR test:

Good enzyme performance was confirmed by PCR amplification using both  $\lambda$  DNA (amplified fragments: 8, 10, 12 kb) and human genomic DNA (amplified fragment: 0.5, 1, 2, 4, 6, 8 kb) as templates.

### General reaction mixture for PCR (total 50 $\mu$ l):

PrimeSTAR HS (Premix)	25 $\mu$ l
Primer 1	0.2 - 0.3 $\mu$ M (final conc.)
Primer 2	0.2 - 0.3 $\mu$ M (final conc.)
Template	< 200 ng
Sterile distilled water	to 50 $\mu$ l

### Recommended template amount

Human genomic DNA	5 ng - 200 ng (< 200 ng)
<i>E. coli</i> genomic DNA	100 pg - 100 ng
cDNA library	1 ng - 200 ng
$\lambda$ DNA	10 pg - 10 ng
plasmid DNA	10 pg - 1 ng

### PCR conditions:

98°C	10 sec.	} 30 cycles
55°C	5 sec. or 15 sec.	
72°C	1 min./kb	
	or	
98°C	10 sec.	} 30 cycles
68°C	1 min./kb	

**Note:** Because this enzyme has an extremely high priming efficiency, a short annealing time is strongly recommended. Please refer to the user manual supplied with this product for detailed cycling recommendations.

### NOTICE TO PURCHASER: LIMITED LICENSE

#### [P1] PCR Notice

Use of this product is covered by one or more of the following US patents and corresponding patent claims outside the US: 5,789,224, 5,618,711, 6,127,155 and claims outside the US corresponding to expired US Patent No. 5,079,352. 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, no right to perform any patented method, 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.

#### [M54] PrimeSTAR HS DNA Polymerase

This product is covered by the claims of U.S. Patent No. 7,704,713 and its foreign counterparts.

PrimeSTAR is a registered 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 77 543 7247 or from our website at [www.takara-bio.com](http://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.

# 2 × PCR Solution PrimeSTAR® HS (Premix)

Code No. R040A

Size: 500  $\mu$ l × 5

(for 100 PCR reactions)

Shipping at - 20°C

Store at - 20°C

保存：過剰に凍結融解を繰り返すと活性が低下する場合がありますのでご注意ください。

Lot No. (英文面をご覧ください。)

品質保証期限： (英文面をご覧ください。)

本製品の使用方法については、製品添付の取扱説明書をご確認ください。

## ●製品説明

PrimeSTAR HS (Premix) は、タカラバイオが独自に開発した High-Fidelity PCR 酵素、PrimeSTAR HS DNA Polymerase と反応用バッファー、dNTP Mixture をあらかじめ 2 倍濃度で混合したプレミックス PCR 酵素である。反応液調製の手間が大幅に減少し、また、コンタミネーションの危険性も軽減するため、ハイスループットな実験系にも有用である。本製品に使用している PrimeSTAR HS DNA Polymerase は非常に強力な 3'→5' exonuclease 活性を有し、DNA 合成において抜群の校正力を示す一方、Taq DNA Polymerase に優る高い増幅効率も示す。

## ●組成

PrimeSTAR HS DNA Polymerase \* : 1.25 units/25  $\mu$ l  
dNTP Mixture : 2 × conc. ; 各 0.4 mM  
PrimeSTAR Buffer : 2 × conc. ; 2 mM Mg<sup>2+</sup> を含む

## \* : PrimeSTAR HS DNA Polymerase (製品コード R010A)

### ○活性の定義

活性化サケ精子 DNA を鋳型/プライマーとして用い、下記の活性測定用反応液中にて 74°C において、30 分間に 10 nmol の全ヌクレオチドを酸不溶性沈殿物に取り込む活性を 1U とする。

### ○活性測定用反応液組成

100 mM	Tris-HCl 緩衝液 (pH8.3 at 37°C)
10 mM	KCl
2 mM	MgCl <sub>2</sub>
6 mM	(NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub>
0.1%	Triton X-100
各 200 $\mu$ M	dATP、dGTP、dCTP
100 $\mu$ M	[ <sup>3</sup> H] TTP
0.001%	BSA
0.4 mg/ml	活性化サケ精子 DNA

### ○純度

- 10 U の本酵素と 0.6  $\mu$ g の  $\lambda$ -Hind III 分解物を 74°C、1 時間反応させても DNA の電気泳動パターンに変化は起こらない。
- 10 U の本酵素と 0.6  $\mu$ g の supercoiled pBR322 DNA を 74°C、1 時間反応させても DNA の電気泳動パターンに変化は起こらない。

## ●用途

hot start PCR 法による DNA 増幅

## ●PCR 産物

PrimeSTAR HS DNA Polymerase を用いて増幅した PCR 産物のほとんどは平滑末端である。したがって、その PCR 産物をそのまま (必要に応じてリン酸化を行って) 平滑末端のベクターにクローニングすることが可能である。

## ●PCR 検定

1.  $\lambda$  DNA を鋳型とした PCR 反応 (増幅産物 8、10、12 kb) において良好な増幅が見られることを確認している。
2. ヒトゲノム DNA を鋳型とした PCR 反応 (増幅産物 0.5、1、2、4、6、8 kb) において良好な増幅が見られることを確認している。

## ●PCR 反応例 (total 50 $\mu$ l PCR)

PrimeSTAR HS (Premix)	25 $\mu$ l
Primer 1	0.2 ~ 0.3 $\mu$ M (final conc.)
Primer 2	0.2 ~ 0.3 $\mu$ M (final conc.)
Template	< 200 ng
滅菌蒸留水	up to 50 $\mu$ l

### ※ Template 量の目安

ヒトゲノム DNA	5 ng ~ 200 ng (< 200 ng)
<i>E. coli</i> ゲノム DNA	100 pg ~ 100 ng
cDNA ライブラリー	1 ng ~ 200 ng
$\lambda$ DNA	10 pg ~ 10 ng
プラスミド DNA	10 pg ~ 1 ng

## ●PCR 条件

98°C	10 sec.	} 30 cycles
55°C	5 sec. または 15 sec.	
72°C	1 min./kb	
or		
98°C	10 sec.	} 30 cycles
68°C	1 min./kb	

※ 本酵素は高いプライミング効率を有しているため、アニーリング時間を短かく設定することをお勧めします。ここに挙げた条件はあくまで一例ですので、PCR 条件の詳細な設定方法については、添付の取扱説明書をご確認ください。

## ●注意

本製品は研究用として販売しております。ヒト、動物への医療、臨床診断用には使用しないようご注意ください。また、食品、化粧品、家庭用品等として使用しないでください。タカラバイオの承認を得ずに製品の再販・譲渡、再販・譲渡のための改変、商用製品の製造に使用することは禁止されています。ライセンスに関する最新の情報は弊社ウェブカタログをご覧ください。本データシートに記載されている会社名および商品名などは、各社の商号、または登録済みもしくは未登録の商標であり、これらは各所有者に帰属します。

v201403Da