

PRODUCT INFORMATION

Thermo Scientific

Phire Hot Start II PCR Master Mix

Pub. No. MAN0013122

Rev. Date 21 November 2016 (Rev. B.00)

#_
 Lot _ Expiry Date _

Store at -20 °C

Ordering information

Component	#F-125S 200 rxns	#F-125L 1000 rxns
	200 × 50 µL rxns	1000 × 50 µL rxns
2X Phire Hot Start II PCR Master Mix	4 × 1.25 mL	20 × 1.25 mL
100% DMSO	1 × 0.5 mL	4 × 0.5 mL
Water, nuclease free	4 × 1.25 mL	20 × 1.25 mL

www.thermofisher.com

For Research Use Only. Not for use in diagnostic procedures.

1. Introduction

Thermo Scientific™ Phire™ Hot Start II DNA Polymerase is a unique DNA polymerase designed for use in all routine and high throughput PCR applications. A special DNA-binding domain linked to the Phire DNA Polymerase enhances the processivity of the polymerase, enabling short extension times and improved yields. The polymerase is also capable of amplifying long DNA fragments, such as the 7.5 kb genomic DNA used in Thermo Scientific quality control assays. Phire Hot Start II DNA Polymerase provides 2-fold fidelity compared to *Taq* DNA polymerase.

The hot start modification of the polymerase is based on the Affibody™ inactivation method.^{1,2} It inhibits DNA polymerase activity at ambient temperatures, thus preventing amplification of non-specific products. At polymerization temperatures the Affibody molecule is released, rendering the polymerase fully active. Phire Hot Start II DNA Polymerase generates blunt ends in the amplification products. It does not possess the 5'→3' exonuclease activity needed for hydrolysis experiments.

Phire Hot Start II PCR Master Mix is a convenient 2X mix containing Phire Hot Start II DNA Polymerase, nucleotides and optimized reaction buffer including MgCl₂. Only template and primers need to be added by the user.

2. Important Notes

- Use 98 °C for denaturation. (See 5.1 & 5.2)
- The annealing rules are different from many common DNA polymerases (such as *Taq* DNA polymerases). Read Section 5.3 carefully before setting annealing temperature.
- Use 10–15 s/kb for extension. (See 5.4)
- Phire Hot Start II DNA Polymerase produces blunt end DNA products.

3. Setting up PCR reactions using Phire Hot Start II PCR Master Mix

Carefully mix and centrifuge all tubes before opening to ensure homogeneity and improve recovery. When using Phire Hot Start II PCR Master Mix, it is not necessary to perform the PCR setup on ice.

Due to the unique nature of Phire Hot Start II DNA Polymerase, optimal reaction conditions may differ from standard enzyme protocols. Phire Hot Start II DNA Polymerase tends to work better at elevated denaturation and annealing temperatures due to higher salt concentrations in its buffer. Please pay special attention to the conditions listed in section 5 when running your reactions. Following the guidelines will ensure optimal enzyme performance.

Table 1. Pipetting instructions: add items in this order.

Component	20 µL rxn	50 µL rxn	Final conc.
H ₂ O	add to 20 µL	add to 50 µL	
2X Phire HS II Master Mix	10 µL	25 µL	1X
Forward primer*	X µL	X µL	0.5 µM
Reverse primer*	X µL	X µL	0.5 µM
Template DNA	X µL	X µL	
(DMSO**, optional)	(0.6 µL)	(1.5 µL)	(3%)

* The recommendation for final primer concentration is 0.5 µM, but it can be varied in a range of 0.2–1.0 µM, if needed.

** Addition of DMSO is recommended for GC-rich amplicons. DMSO is not recommended for amplicons with very low GC %.

Table 2. Cycling instruction

Cycle step	2-step protocol		3-step protocol		Cycles
	Temp.	Time	Temp.	Time	
Initial Denaturation	98 °C	30 s	98 °C	30 s	1
Denaturation	98 °C	5 s	98 °C	5 s	25–35
Annealing (see 5.3)	–	–	X °C	5 s	
Extension (see 5.4)	72 °C	10–15 s/kb	72 °C	10–15 s/kb	
Final extension	72 °C 4 °C	1 min hold	72 °C 4 °C	1 min hold	1

4. Notes about reaction components

4.1. 2X Phire Hot Start II PCR Master Mix

Phire Hot Start II Master Mix contains all the necessary reaction components except for template DNA and primers. The master mix provides 1.5 mM MgCl₂ and 200 µM of each dNTP in final reaction concentration. Phire Hot Start II DNA Polymerase concentration is optimized to give good results in most reactions.

4.2. Template

For low complexity DNA (e.g. plasmid, lambda or BAC DNA) it is recommended to use 2.5 pg–25 ng per 50 µL reaction volume. For high complexity genomic DNA, the amount of DNA template should be 25–250 ng per 50 µL reaction volume.

If cDNA synthesis reaction mixture is used as a source of template, the volume of the template should not exceed 10% of the final PCR reaction volume.

4.4. PCR additives

The recommended reaction conditions for GC-rich templates include 3% DMSO as a PCR additive, which aids in the denaturing of templates with high GC contents. For further optimization DMSO should be varied in 2% increments. In some cases DMSO may also be required for supercoiled plasmids to relax for denaturation. If high DMSO concentration is used, the annealing temperature must be decreased, as DMSO affects the melting point of the primers. It has been reported that 10% DMSO decreases the annealing temperature by 5.5–6.0 °C.³

5. Notes about cycling conditions

5.1. Initial denaturation

Denaturation should be performed at 98 °C. Due to the high thermostability of Phire Hot Start II DNA Polymerase even higher than 98 °C denaturation temperatures can be used. We recommend 30 seconds initial denaturation at 98 °C for most templates. Some templates may require longer initial denaturation time and the length of the initial denaturation time can be extended up to 3 minutes.

5.2. Denaturation

Keep the denaturation as short as possible. Usually 5 seconds at 98 °C is enough for most templates.

Note: The denaturation time and temperature may vary depending on the ramp rate and temperature control mode of the cyclers.

5.3. Primer annealing

The optimal annealing temperature for Phire Hot Start II DNA Polymerase may be significantly different than annealing temperature with other DNA polymerases. Always use the T_m calculator and instructions on our website (www.thermofisher.com/tmcalculator) to determine the T_m values of your primers and optimal annealing temperature.

As a basic rule, for primers > 20 nt, anneal for 5 seconds at a T_m +3 °C of the lower T_m primer. For primers ≤ 20 nt, use an annealing temperature equal to the T_m of the lower T_m primer. If necessary, use a temperature gradient to find the optimal annealing temperature for each template-primer pair combination. The annealing gradient should extend up to the extension temperature (two-step PCR).

Two-step cycling without annealing step is also recommended for high T_m primer pairs.

5.4. Extension

The extension should be performed at 72 °C. Extension time of 10 seconds per 1 kb is recommended for most templates. However, higher yields may be obtained using extension time of 15 s/kb with challenging primer-template pairs.

6. Cloning recommendations

Blunt end cloning is recommended when cloning DNA fragments amplified with Phire Hot Start II PCR Master Mix. If TA cloning is required, it is necessary to add A overhangs to the PCR product (with Thermo Scientific *Taq* DNA Polymerase, for example). A detailed protocol for TA cloning of fragments amplified with Phire Hot Start II PCR Master Mix can be found on our website www.thermofisher.com.

(continued on reverse page)

7. Troubleshooting

No product at all or low yield

- | | |
|---|---|
| <ul style="list-style-type: none">• Repeat and make sure that there are no pipetting errors.• Titrate template amount.• Template DNA may be damaged. Use carefully purified template.• Increase extension time.• Increase the number of cycles.• Optimize annealing temperature. | <ul style="list-style-type: none">• Titrate DMSO (2–8%) in the reaction.• Denaturation temperature may be too low. Optimal denaturation temperature for most templates is 98 °C or higher.• Optimize denaturation time.• Check the purity and concentration of the primers.• Check primer design. |
|---|---|

Non-specific products - High molecular weight smears

- | | |
|---|--|
| <ul style="list-style-type: none">• Shorten extension time.• Reduce the total number of cycles.• Increase annealing temperature or try 2-step protocol. | <ul style="list-style-type: none">• Vary denaturation temperature• Decrease primer concentration. |
|---|--|

Non-specific products - Low molecular weight discrete bands

- | | |
|---|--|
| <ul style="list-style-type: none">• Increase annealing temperature• Shorten extension time.• Titrate template amount. | <ul style="list-style-type: none">• Decrease primer concentration.• Design new primers. |
|---|--|

8. References

1. Nord K. et al. (1997) *Nature Biotechnol.* 15: 772–777.
2. Wikman M. et al. (2004) *Protein Eng., Des. Sel.* 17: 455–462.
3. Chester N. & Marshak D.R. (1993) *Analytical Biochemistry* 209: 284–290.

CERTIFICATE OF ANALYSIS

DNA amplification assay

Performance in PCR is tested by the amplification of a 7.5 kb fragment of genomic DNA.

Quality authorized by:  Jurgita Zilinskiene

NOTICE TO PURCHASE:

The purchase price of this product includes a limited, non-transferable license under U.S. and foreign patents owned by BIO-RAD Laboratories, Inc., to use this product. No other license under these patents is conveyed expressly or by implication to the purchaser by the purchase of this product.

LIMITED USE LABEL LICENSE: Internal Research and Development Use Only.

The purchase of this product conveys to the buyer the limited, non-exclusive, non-transferable right (without the right to resell, repackage, or further sublicense) to use this product for internal research and development purposes. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of the product does not include or carry any right or license to use, develop, or otherwise exploit this product commercially and no rights are conveyed to the buyer to use the product or components of the product for purposes including but not limited to provision of services to a third party, generation of commercial databases or clinical diagnostics. This product is sold pursuant to authorization from Thermo Fisher Scientific and Thermo Fisher Scientific reserves all other rights. For information on purchasing a license for uses other than internal research and development purposes, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Inc., 5781 Van Allen Way, Carlsbad, California 92008.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively for research purposes and in vitro use only. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to www.thermofisher.com for Material Safety Data Sheet of the product.

© 2016 Thermo Fisher Scientific, Inc. All rights reserved. Affibody is a registered trademark of Affibody AB, Sweden. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.