



PRODUCT INFORMATION

Thermo Scientific

Phire Green Hot Start II PCR Master Mix

Pub. No. MAN0013357

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

#\_
Lot \_ Expiry Date \_

Store at -20 °C

Ordering information

Table with 3 columns: Component, #F-126S 200 rxns, #F-126L 1000 rxns. Rows include 2X Phire Green Hot Start II PCR Master Mix, 100% DMSO, and Water, nuclease free.

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.

The hot start modification of the polymerase is based on the Affibody™ inactivation method. It inhibits DNA polymerase activity at ambient temperatures, thus preventing amplification of non-specific products.

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.

The master mix also includes a density reagent and two tracking dyes for direct loading of PCR products on a gel. The dyes do not interfere with PCR performance and are compatible with downstream applications such as DNA sequencing, ligation and restriction digestion.

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 Green 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.

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

Table with 4 columns: Component, 20 µL rxn, 50 µL rxn, Final conc. Rows include H2O, 2X Phire Green HS II Master Mix, Forward primer, Reverse primer, and Template DNA (DMSO optional).

\* 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

Table with 6 columns: Cycle step, 2-step protocol (Temp, Time), 3-step protocol (Temp, Time), Cycles. Rows include Initial Denaturation, Denaturation, Annealing, Extension, and Final extension.

4. Notes about reaction components

4.1. 2X Phire Green Hot Start II PCR Master Mix

Phire Green Hot Start II PCR 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.

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.

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.

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.

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 cyler.

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 Tm calculator and instructions on our website (www.thermofisher.com/tmcalculator) to determine the Tm values of your primers and optimal annealing temperature.

As a basic rule, for primers > 20 nt, anneal for 5 seconds at a Tm +3 °C of the lower Tm primer. For primers ≤ 20 nt, use an annealing temperature equal to the Tm of the lower Tm primer. If necessary, use a temperature gradient to find the optimal annealing temperature for each template-primer pair combination.

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 Green 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 Green Hot Start II PCR Master Mix can be found on our website www.thermofisher.com.

## 7. Troubleshooting

<b>No product at all or low yield</b>	
<ul style="list-style-type: none"><li>• Repeat and make sure that there are no pipetting errors.</li><li>• Titrate template amount.</li><li>• Template DNA may be damaged. Use carefully purified template.</li><li>• Increase extension time.</li><li>• Increase the number of cycles.</li><li>• Optimize annealing temperature.</li></ul>	<ul style="list-style-type: none"><li>• Titrate DMSO (2–8%) in the reaction.</li><li>• Denaturation temperature may be too low. Optimal denaturation temperature for most templates is 98 °C or higher.</li><li>• Optimize denaturation time.</li><li>• Check the purity and concentration of the primers.</li><li>• Check primer design.</li></ul>
<b>Non-specific products - High molecular weight smears</b>	
<ul style="list-style-type: none"><li>• Shorten extension time.</li><li>• Reduce the total number of cycles.</li><li>• Increase annealing temperature or try 2-step protocol.</li></ul>	<ul style="list-style-type: none"><li>• Vary denaturation temperature</li><li>• Decrease primer concentration.</li></ul>
<b>Non-specific products - Low molecular weight discrete bands</b>	
<ul style="list-style-type: none"><li>• Increase annealing temperature</li><li>• Shorten extension time.</li><li>• Titrate template amount.</li></ul>	<ul style="list-style-type: none"><li>• Decrease primer concentration.</li><li>• Design new primers.</li></ul>

## 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: LIMITED LICENSE

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](mailto: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](http://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.