

## Thermo Scientific DreamTaq Hot Start PCR Master Mix (2X)

Pub. No. MAN0015976 Rev. Date 29 July 2016 (Rev. A.00)

## **Ordering Information**

Component	<b>#K9011</b> 200 rxns of 50 μL	<b>#K9012</b> 1000 rxns of 50 μL
DreamTaq Hot Start PCR Master Mix (2X)	4 × 1.25 mL	20 × 1.25 mL
Water, nuclease-free	$4 \times 1.25 \text{ mL}$	$20\times1.25~\text{mL}$

## Store at -20°C

www.thermofisher.com

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

## DESCRIPTION

Thermo Scientific™ DreamTaq™ Hot Start PCR Master Mix (2X) is a ready-to-use solution containing DreamTaq Hot Start DNA polymerase, optimized DreamTaq buffer, MgCl₂, and dNTPs. This pre-mixed formulation saves time and reduces contamination due to a reduced number of pipetting steps required for PCR set up.

Thermo Scientific™ DreamTaq™ Hot Start DNA Polymerase ensures higher sensitivity, specificity, and yields compared to conventional hot start *Taq* DNA polymerase. It is capable of amplifying long amplicons such as 6 kb genomic DNA and 20 kb λ DNA.

DreamTaq Hot Start DNA Polymerase combines *Taq* DNA polymerase and a specific antibody that inhibits the DNA polymerase activity at ambient temperatures, thus preventing the amplification of non-specific products. At polymerization temperatures, the antibody molecule is released, rendering the polymerase fully active. Due to the antibody-based hot start, the reactions can be set up at room temperature.

## **FEATURES**

- · High specificity due to antibody based hot start.
- Robust amplification with minimal optimization.
- High yields of PCR products.
- Higher sensitivity compared to conventional hot start Taq DNA polymerases.
- Amplification of long targets up to 6 kb from genomic DNA and up to 20 kb from viral DNA.
- Generates 3'-dA overhangs.

## **APPLICATIONS**

- Routine PCR amplification of DNA fragments up to 6 kb from genomic DNA and up to 20 kb from viral DNA.
- RT-PCR.
- Genotyping.
- · Generation of PCR products for TA cloning.

## DREAMTAQ HOT START PCR MASTER MIX (2X) COMPOSITION

DreamTaq Hot Start PCR Master Mix contains all the necessary reaction components except for template DNA and primers. It includes DreamTaq Hot Start DNA Polymerase, 2X DreamTaq buffer, dATP, dCTP, dGTP, and dTTP, 0.4 mM each, and 4 mM MgCl<sub>2</sub>. DreamTaq buffer is a proprietary formulation optimized for robust PCR performance.

## **PROTOCOL**

- Gently vortex and briefly centrifuge DreamTaq Hot Start PCR Master Mix (2X) after thawing.
- For each 50-µL reaction, add the following components into a separate thin-walled PCR tube:

DreamTaq Hot Start PCR Master Mix (2X)	25 µL	
Forward primer	0.1–1.0 µM	
Reverse primer	0.1–1.0 µM	
Template DNA	10 pg–1 μg	
Water, nuclease-free	to 50 µL	
Total volume	50 μL	

- Gently vortex the samples and briefly centrifuge.
- 4. When using a thermal cycler that does not contain a heated lid, overlay the reaction mixture with 25  $\mu L$  of mineral oil.
- Place the reactions in a thermal cycler. Perform PCR using the recommended thermal cycling conditions outlined below:

Step	Temperature, °C	Time	Number of cycles
Initial denaturation	95	1–3 min	1
Denaturation	95	30 s	
Annealing	Tm	30 s	25–40
Extension*	72	1 min	
Final Extension	72	5–15 min	1

<sup>\*</sup> The recommended extension step is 1 minute for PCR products up to 2 kb. For longer products, the extension time should be prolonged by 1 minute/kb.

## GUIDELINES FOR PREVENTING CONTAMINATION OF PCR REACTION

During PCR more than 10 million copies of template DNA are generated. Therefore, care must be taken to avoid contamination with other templates and amplicons that may be present in the laboratory environment. Follow the general recommendations below to lower the risk of contamination.

- Prepare your DNA sample, set up the PCR mixture, perform thermal cycling and analyze PCR products in separate areas.
- Set up PCR mixtures in a laminar flow cabinet equipped with an UV lamp.
- Wear fresh gloves for DNA purification and reaction set up.
- Use reagent containers dedicated for PCR. Use positive displacement pipettes, or use pipette tips with aerosol filters to prepare DNA samples and perform PCR set up.
- Always perform "no template control" (NTC) reactions to check for contamination.

## **GUIDELINES FOR PRIMER DESIGN**

Use special design software or follow the general recommendations for PCR primer design as outlined below to design optimal primers:

- Use PCR primers that are 15–30 nucleotides long.
- Optimal GC content of the primer is 40–60%. Ideally, C and G nucleotides should be distributed uniformly along the primer.
- Avoid placing more than three G or C nucleotides at the 3'-end to lower the risk of non-specific priming.
- If possible, the primer should terminate with a G or C at the 3'-end.
- Avoid self-complementary primer regions, and complementarities between the primers and direct primer repeats to prevent hairpin formation and primer dimerization.
- Check for possible sites of undesired complementarity between primers and template DNA.
- When designing degenerate primers, place at least 3 conserved nucleotides at the 3'-end.
- Differences in melting temperatures (Tm) between the two primers should not exceed 5°C.

## ESTIMATION OF PRIMER MELTING TEMPERATURE

For primers containing less than 25 nucleotides, the approximate melting temperature (Tm) can be calculated using the following equation:

$$Tm = 4 (G + C) + 2 (A + T),$$

where G, C, A, T represent the number of respective nucleotides in the primer.

If the primer contains more than 25 nucleotides, we recommend using specialized computer programs to account for interactions of adjacent bases, effect of salt concentration, etc.

## COMPONENTS OF THE REACTION MIXTURE

#### Template DNA

Optimal amount of template DNA for a 50- $\mu$ L reaction volume is 1 pg–1 ng for both plasmid and phage DNA, and 100 pg–1  $\mu$ g for genomic DNA. Higher amounts of template increase the risk of non-specific PCR products. Lower amounts of template reduce the accuracy of the amplification.

All routine DNA purification methods are suitable for template preparation; e.g., Thermo Scientific™ GeneJET™ Genomic DNA Purification Kit (#K0721) or GeneJET Plasmid Miniprep Kit (#K0502). Trace amounts of certain agents used for DNA purification, such as phenol, EDTA, and proteinase K, can inhibit DNA polymerases. Ethanol precipitation and repeated washes of the DNA pellet with 70% ethanol normally removes trace contaminants from DNA samples.

## **Primers**

The recommended concentration range of the PCR primers is  $0.1-1~\mu M$ . Excessive primer concentrations increase the probability of mispriming and generation of non-specific PCR products.

For degenerate primers and primers used for long PCR, we recommend higher primer concentrations in the range of 0.3–1  $\mu$ M.

## CYCLING PARAMETERS

## Initial DNA denaturation and enzyme activation

DreamTaq Hot Start DNA polymerase is inactive at room temperature during the reaction set up and is activated during the 1–3 minute initial denaturation/enzyme activation step.

It is essential to completely denature the template DNA at the beginning of the PCR run to ensure efficient utilization of the template during the first amplification cycle. If the GC content of the template is 60% or less, an initial 1–3 minute denaturation at 95°C is sufficient. For GC-rich templates, this step can be prolonged.

## Denaturation

A DNA denaturation time of 30 seconds per cycle at 95°C is normally sufficient. For GC-rich DNA templates, this step can be prolonged to 3–4 minutes. DNA denaturation can also be enhanced by the addition of 5–10% glycerol, 5% DMSO, 1% formamide, or 1–1.5 M betaine. The melting temperature of the primer-template complex decreases significantly in the presence of these reagents. Therefore, the annealing temperature has to be adjusted accordingly.

## Primer annealing

The annealing temperature should be equal to the melting temperature (Tm) of the primers. Annealing for 30 seconds is normally sufficient. If non-specific PCR products appear, the annealing temperature should be optimized stepwise in 1–2°C increments. When additives that change the melting temperature of the primer-template complex are used (glycerol, DMSO, formamide and betaine), the annealing temperature must also be adjusted.

#### Extension

The optimal extension temperature for DreamTaq Hot Start DNA Polymerase is 70–75°C. The recommended extension step is 1 minute at 72°C for PCR products up to 2 kb. For longer products, the extension time should be increased by 1 minute/kb. For amplification of templates >6 kb, we recommend reducing the extension temperature to 68°C.

## Number of cycles

The number of cycles may vary depending on the amount of template DNA in the PCR mixture and the expected PCR product yield.

If less than 10 copies of the template is present in the reaction, about 40 cycles are required. For higher template amounts, 25–35 cycles are sufficient.

## Final extension

After the last cycle, we recommend incubating the PCR mixture at 72°C for an additional 5–15 minutes to fill in any possible incomplete reaction products. If the PCR product will be cloned into TA vectors such as the Thermo Scientific™ InsTAclone™ PCR Cloning Kit (#K1213), the final extension step may be prolonged to 15 minutes to ensure the complete 3'-dA tailing of the PCR product. If the PCR product will be used for cloning using Thermo Scientific™ CloneJET™ PCR Cloning Kit (#K1231), the final extension step can be omitted.

## **TROUBLESHOOTING**

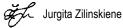
For troubleshooting, visit www.thermofisher.com.

## **CERTIFICATE OF ANALYSIS**

## **Functional Assay**

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

Quality authorized by:



## LIMITED USE LABEL LICENSE No. 593: Newcastle License for Modified DNA Polymerase

Notice to Purchaser: This product is licensed under patents owned by University of Newcastle upon Tyne

# LIMITED USE LABEL LICENSE No. 599: Internal Research and Development Use Only

Notice to Purchaser: 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., 5791 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.

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

Refer to www.thermofisher.com/support for the Safety Data Sheets.

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.