

2X Taq Master Mix
Technical Manual No. 0228Cat. No. E00019
Version 10112010

I	Description.....	1
II	Key Features.....	1
III	Contents.....	1
IV	Stability.....	2
V	Storage.....	2
VI	Storage.....	2
VII	General PCR Protocol Using 2X Taq DNA Master Mix.....	2
VII	Order Information.....	3

I. DESCRIPTION

2X Taq Master Mix is a premixed 2X concentrated solution of Taq DNA Polymerase (GenScript, Cat. No. E00007), reaction buffer, MgCl₂ and dNTPs. 2X Taq Master Mix contains all components for PCR*, except DNA template and primers. The mixture is optimized for consistent and efficient routine PCR amplifications. It can amplify up to 8 kb fragment from lambda DNA. For a 50 µl reaction, simply add 25 µl of 2X Taq Master Mix to primers, DNA template and PCR-Qualified H₂O.

II. KEY FEATURES

- Taq DNA Polymerase in ready-to-use mix
- Low contamination risk
- Low risk of pipetting errors

III. CONTENTS

- 0.1 U/µl Taq DNA Polymerase (GenScript, Cat. No. E00007)
- Reaction buffer
- 3 mM MgCl₂
- 0.4 mM dNTPs

IV. STABILITY

1. Freeze-thaw stability of 2X *Taq* master mix: Following 25 freeze-thaw cycles, no effect on performance is observed.

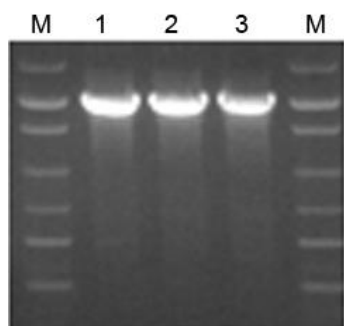


Fig. 1 Stability after 25 freeze-thaw cycles.

2 kb fragment amplification.

Lane 1 25 µl of 2X *Taq* Master Mix

Lane 2 25 µl of 2X *Taq* Master Mix after 25 freeze-thaw cycles

Lane 3 2.5 U *Taq* DNA Polymerase (GenScript, Cat. No. E00007)

2. Stability at 4°C: No effect on performance is observed after storage at 4°C for 2 months.

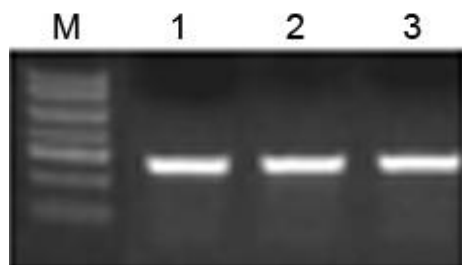


Fig. 2 Stability at 4°C

0.5 kb fragment amplification.

Lane 1 25 µl of 2X *Taq* Master Mix

Lane 2 25 µl of 2X *Taq* Master Mix storage at 4°C for 2 months

Lane 3 2.5 U *Taq* DNA Polymerase (GenScript, Cat. No. E00007)

V. STORAGE

This product is shipped on blue ice. Store the product at –20°C.

VI. GENERAL PCR PROTOCOL USING 2X TAQ MASTER MIX

This is a general PCR amplification protocol, optimization may be needed to get satisfactory results.

1. Thaw the 2X *Taq* Master Mix at room temperature. Vortex the 2X *Taq* Master Mix and then spin it briefly in a microcentrifuge to collect the material in the bottom of the tube.
2. Prepare one of the following reaction mixes on ice:

- **For a 25 µl reaction volume:**

Component	Volume	Final Concentration
2X <i>Taq</i> Master Mix	12.5 µl	1X
Upstream Primer, 10 µM	0.5 µl	0.1–1.0 µM
Downstream Primer, 10 µM	0.5 µl	0.1–1.0 µM
DNA Template	1-5 µl	<500 ng
Nuclease-Free Water to	25 µl	

- **For a 50 µl reaction volume:**

Component	Volume	Final Concentration
2X <i>Taq</i> Master Mix	25 µl	1X
Upstream Primer, 10 µM	1 µl	0.1–1.0 µM
Downstream Primer, 10 µM	1 µl	0.1–1.0 µM
DNA Template	1-5 µl	<500 ng
Nuclease-Free Water to	50 µl	

3. Gently mix the reaction and spin down in microcentrifuge.

4. Set up cycling conditions for a routine PCR reactions:

- Initial Denaturation 94-95°C for 1-5 minutes
- 25-40 cycles 94-96°C for 30 sec.
 45-70°C for 10-30 seconds
 72°C for X minutes (1 min/kb)
- Final extension: 72°C for 7 minutes
- Final soak 4-10°C

VII. ORDER INFORMATION

2X *Taq* Master Mix Cat. No. E00019

GenScript USA Inc.

860 Centennial Ave., Piscataway, NJ 08854

Tel: 732-885-9188, 732-885-9688

Fax: 732-210-0262, 732-885-5878

Email: product@genscript.com

Web: <http://www.genscript.com>

For Research Use Only

* The PCR process is covered by US. Patent numbers 4683195 and 4683202 issued to Cetus and owned by Hoffman-La Roche Inc. GenScript does not encourage or support the unauthorized use of the PCR process. Use of this product is recommended for persons that either have a license to perform PCR or are not required to obtain a license.