



**G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)**

*A Geno Technology, Inc. (USA) brand name*

## High fidelity DNA Polymerase 2X Master Mix (Cat. # 786-1968, 786-1969)



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)



|                                  |          |
|----------------------------------|----------|
| <b>INTRODUCTION .....</b>        | <b>3</b> |
| <b>FUNCTIONAL FEATURES .....</b> | <b>3</b> |
| <b>ITEM(S) SUPPLIED.....</b>     | <b>3</b> |
| <b>STORAGE CONDITIONS .....</b>  | <b>3</b> |
| <b>PROTOCOL .....</b>            | <b>4</b> |
| <b>TYPICAL QC DATA .....</b>     | <b>4</b> |





## Introduction

G-Biosciences high-fidelity DNA Polymerase 2x master mix is ready to use premix which contains high fidelity DNA Polymerase, dNTPs and optimized buffer with stabilizers. The high-fidelity DNA Polymerase possesses 5' → 3' polymerase and 3' → 5' exonuclease (proofreading) activities which has the fidelity, sensitivity and processivity with an error rate  $\sim 2.8 \times 10^{-2}$ -fold lower than Taq DNA polymerase, and significantly lower than other proofreading DNA polymerases.

## Functional features

- High fidelity PCR
- Cloning
- Long and high GC content template DNA amplification

## Items Supplied

| Cat #                                      | 786-1968   | 786-1969    |
|--------------------------------------------|------------|-------------|
| High-fidelity DNA Polymerase 2X master mix | 2 x 1.25ml | 10 x 1.25ml |

## Storage temperature

- 20°C.



## Protocol

1. Set up PCR reaction as follows:

| PCR reaction set up:                       |              |
|--------------------------------------------|--------------|
| Template                                   | ~1-20 ng     |
| Forward primer (3.2 $\mu$ M)               | 1.0 $\mu$ l  |
| Reverse primer (3.2 $\mu$ M)               | 1.0 $\mu$ l  |
| High fidelity DNA Polymerase 2X master mix | 10.0 $\mu$ l |
| H <sub>2</sub> O up to                     | 20.0 $\mu$ l |

2. Run PCR with the following program:

| PCR cycling conditions: |             |           |        |
|-------------------------|-------------|-----------|--------|
| Steps                   | Temperature | Time      | Cycles |
| Initial denaturation    | 98°C        | 2 min     | 1      |
| Denaturation            | 98°C        | 20 sec    | 30-35  |
| Annealing               | 56°C-60 °C  | 20 sec    |        |
| Extension               | 72°C        | 30 sec/kb |        |
| Final extension         | 72°C        | 5 min     | 1      |
| Hold                    | 4°C         |           |        |

- Take 7.5  $\mu$ l samples and add loading dye
- Run 1% agarose gel electrophoresis
- Scan the gel and analyze the data.

## Typical QC data

The enzyme activity was tested with lambda DNA as a template using 1 to 5kb PCR primers.

