# **Certificate of Analysis**

# GoTaq<sup>®</sup> MDx DNA Polymerase:

For Laboratory Use. Supplied with:

Cat.#	GoTaq® MDx DNA Polymerase	5X Green GoTaq® Flexi Buffer	5X Colorless GoTaq® Flexi Buffer	Magnesium Chloride Solution, 25mM
D4001	100 units (D400A)	1ml (M891A)	1ml (M890A)	0.75ml (A351B)
D4005	500 units (D400B)	4 × 1ml (M891A)	4 × 1ml (M890A)	3 × 0.75ml (A351B)
D4006	5 × 500 units (D400B)	20 × 1ml (M891A)	20 × 1ml (M890A)	15 × 0.75ml (A351B)

**Description:** GoTaq<sup>®</sup> MDx DNA Polymerase<sup>(a)</sup> contains GoTaq<sup>®</sup> MDx DNA Polymerase, 5X Green GoTaq<sup>®</sup> Flexi Buffer, 5X Colorless GoTaq<sup>®</sup> Flexi Buffer and 25mM MgCl<sub>2</sub>. The enzyme is supplied in a proprietary formulation containing 50% glycerol with buffers designed for enhanced amplification. The enzyme is a full-length form of *Taq* DNA polymerase that exhibits  $5' \rightarrow 3'$  exonuclease activity. The 5X Green GoTaq<sup>®</sup> Flexi Buffer contains two dyes (a blue dye and a yellow dye) that separate during electrophoresis to indicate migration progress. Use the colorless buffer when direct fluorescence or absorbance readings are required without prior purification of the amplified DNA from the PCR. The Flexi Buffers do not contain magnesium, allowing easy optimization of magnesium concentration in amplification reactions. GoTaq<sup>®</sup> MDx DNA Polymerase is manufactured under cGMP.

Biological Source: The enzyme is derived from bacteria.

Concentration: See product label for measured unit activity.

Storage Conditions: Store at -30°C to -10°C.

Expiration Date: See product label for expiration date.

**5X Green GoTaq® Flexi Buffer (Part# M891A):** Proprietary formulation supplied at pH 8.5 containing blue dye and yellow dye. The blue dye migrates at the same rate as a 3–5kb DNA fragment in a 1% agarose gel. The yellow dye migrates at a rate faster than primers (<50bp) in a 1% agarose gel. Green GoTaq® Flexi Buffer also increases the density of the sample, so it will sink into the well of the agarose gel, allowing reactions to be loaded directly onto gels without loading dye. This buffer does not contain magnesium.

5X Colorless GoTaq<sup>®</sup> Flexi Buffer (Part# M890A): Proprietary formulation supplied at pH 8.5. This buffer does not contain magnesium.

Magnesium Chloride Solution, 25mM (Part# A351B): Provided to allow you to optimize MgCl<sub>2</sub> concentration according to your individual requirements. Vortex the MgCl<sub>2</sub> thoroughly after thawing and prior to use.

# **Quality Control Assays**

Test		Specification		Result
Concentration		5.2–7.6u/µl One unit is defined as the amount of enzyme required to catalyze the incorporation of		Pass See attached
		2 nanomoles of dNTPs in 30 minutes at 55°C in a fluorescent extension assay.		
Purity	DNA Contamination	Bacterial DNA	One unit of enzyme contains less than 10 copies of bacterial genomic DNA determined by quantitative amplification of a 16S rRNA gene.	Pass
		Fungal DNA	One unit of enzyme contains less than 1 genome equivalent of fungal genomic DNA by quantitative amplification of a 18S rRNA gene.	Pass
		Mammalian DNA	One unit of enzyme contains less than 1 genome equivalent of mammalian gDNA by quantitative amplification of mitochondrial genomic DNA.	Pass
	Nuclease Contamination	Endonuclease/ Nicking	No observable nicking of 0.5µg of supercoiled DNA after incubation for 8 hours at 22°C, followed by 8 hours at 45°C in the presence of 15 units of enzyme.	Pass
		Exonuclease	No observable degradation of 1.0µg of Lambda DNA/HindIII markers after incubation for 8 hours at 22°C, followed by 8 hours at 45°C in the presence of 15 units of enzyme.	Pass
		Ribonuclease	No observable degradation of RNA target after incubation for 1 hour at 37°C in the presence of 10 units of enzyme.	Pass
Function		1 unit of enzyme quantitatively amplifies mitochondrial genomic DNA.		Pass



PCR Satisfaction Guarantee

Promega's PCR Systems, enzymes and reagents are proven in PCR to ensure reliable, high performance results. Your success is important to us. Our products are backed by a worldwide team of Technical Support scientists. Please contact them for applications or technical assistance. If you are not completely satisfied with any Promega PCR product we will send a replacement or refund your account. That's Our PCR Guarantee! Product must be within expiration date and have been stored and used in accordance with product literature. See Promega Product Insert for specific tests performed.

For Wheeler

R. Wheeler, Quality Assurance





**O** Promega

#### **Promega Corporation**

2800 Woods Hollow Road	ł
Madison, WI 53711-5399	) USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

#### PRODUCT USE LIMITATIONS, WARRANTY DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DIS-CLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PAR-TICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifica-

(a)U.S. Pat. No. 6,242,235, Australian Pat. No. 761757, Canadian Pat. No. 2,335,153, Chinese Pat. No. ZL99808861.7, Hong Kong Pat. No. HK 1040262, Japanese Pat. No. 3673175, European Pat. No. 1088060 and other patents pending.

© 2015, 2016, 2018 Promega Corporation. All Rights Reserved.

GoTaq is a registered trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PID400

Printed in USA Revised 5/18.

Signed by:



## 1. Standard Application

## Reagents to Be Supplied by the User

- dNTP Mix (Cat.# U1511)
- upstream primer
  template DNA
- downstream primer
  mineral oil (optional)
- 1. In a sterile, nuclease-free microcentrifuge tube, combine the following components at room temperature:

Nuclease-Free Water (Cat.# P1193)

Component	Final Volume	Final Conc.			
5X GoTaq <sup>®</sup> Flexi Buffer <sup>1</sup>					
(Green or Colorless)	10µI	1X			
MgCl <sub>2</sub> Solution, 25mM <sup>1</sup>	2–8µI	1.0-4.0mM			
dNTP Mix, 10mM each	1µI	0.2mM each dNTP			
upstream primer	XμI	0.1–1.0µM			
downstream primer	YμI	0.1–1.0µM			
GoTaq® MDx DNA Polymerase	ZμI	1.25u <sup>2</sup>			
template DNA	<u> </u>	<0.5µg/50µl			
Nuclease-Free Water to	50µl				
<sup>1</sup> Thaw completely and vortex thoroughly prior to use.					
2Recommended optimization of enzyme quantity specific to assay.					

- If using a thermal cycler without a heated lid, overlay the reaction with 1–2 drops (approximately 50µl) of mineral oil to prevent evaporation during thermal cycling.
- 3. Place reactions in a thermal cycler that has been heated to 94–95°C, and begin PCR.

## 2. General Guidelines for Amplification by PCR

### A. Denaturation

 Following the initial 2-minute 94–95°C denaturation step, denaturation steps should be between 15 seconds and 1 minute.

#### B. Annealing

- Optimize the annealing conditions by performing the reaction starting approximately 5°C below the calculated melting temperature of the primers and increasing the temperature in increments of 1°C to the annealing temperature.
- The annealing step is typically 30 seconds to 1 minute.

## C. Extension

- The extension reaction is typically performed at the optimal temperature for Taq DNA polymerase, which is 72–74°C.
- Allow approximately 1 minute for every 1kb of DNA to be amplified.
- A final extension of 5 minutes at 72-74°C is recommended.

#### D. Refrigeration

• If the thermal cycler has a refrigeration or "soak" cycle, the cycling reaction can be programmed to end by holding the tubes at 4°C for several hours.

#### E. Cycle Number

 Generally, 25–30 cycles result in optimal amplification of desired products. Occasionally, up to 40 cycles may be performed, especially for detection of low-copy targets.

# 3. General Considerations

## A. Buffer Choice

We recommend using the 5X Green GoTaq<sup>®</sup> Flexi Buffer in any amplification reaction that will be visualized by agarose gel electrophoresis followed by ethidium-bromide staining. The 5X Green GoTaq<sup>®</sup> Flexi Buffer is not recommended for any downstream applications using absorbance or fluorescence excitation because the yellow and blue dyes in the reaction buffer may interfere with these applications. The dyes absorb between 225–300nm, making standard A<sub>260</sub> readings to determine DNA concentration unreliable. Also, the dyes have excitation peaks at 488nm and between 600–700nm that correspond to the excitation wavelengths commonly used in fluorescence detection instrumentation. However, for some

instrumentation, such as a fluorescent gel scanner that uses a 488nm excitation wavelength, there will be minimal interference, since it is the yellow dye that absorbs this wavelength. Gels scanned by this method will have a light gray dye front below the primers corresponding to the yellow dye front. The Green and Colorless GoTaq<sup>®</sup> Flexi Buffers give approximately equivalent amplification yields. Balanced amplifications between the two buffers may require further optimization. For reactions going directly from thermal cycler to an application using absorbance or fluorescence, the 5X Colorless GoTag<sup>®</sup> Flexi Buffer is recommended.

## B. Enzyme Concentration

We have found that 1.25 units of GoTaq<sup>®</sup> MDx Hot Start Polymerase per 50µl amplification reaction is adequate for most amplifications. However, optimization of enzyme concentration specific to the amplification reaction may be required to achieve optimal assay performance.

## C. Primer Design

PCR primers generally range in length from 15–30 bases and are designed to flank the region of interest. Primers should contain 40–60% (G + C), and care should be taken to avoid sequences that might produce internal secondary structure. The 3´-ends of the primers should not be complementary to avoid the production of primer-dimers. Primer-dimers unnecessarily deplete primers from the reaction and result in an unwanted polymerase reaction that competes with the desired reaction. Avoid three G or C nucleotides in a row near the 3´-end of the primer because this may result in nonspecific primer annealing, increasing the synthesis of undesirable reaction products. Ideally, both primers should have nearly identical melting temperatures ( $T_m$ ); in this manner, the two primers should anneal roughly at the same temperature. The annealing temperature of the reaction depends on the primer with the lowest  $T_m$ . For assistance with calculating the  $T_m$  of any primer, a  $T_m$  Calculator is provided on the BioMath page of the Promega web site at: www.promega.com/biomath/

### D. Amplification Troubleshooting

To overcome low yield or no yield in amplifications, we recommend the following suggestions:

- Adjust annealing temperature. The reaction buffer composition affects the melting properties of DNA. See BioMath Calculator to calculate the melting temperature for primers in the GoTaq<sup>®</sup> reaction (www.promega.com/biomath/).
- Minimize the effect of amplification inhibitors. Some DNA isolation procedures, particularly genomic DNA isolation, can result in the copurification of amplification inhibitors. Reduce the volume of template DNA in reaction, or dilute template DNA prior to adding to reaction. Diluting samples up to 1:10,000 has been shown to be effective in improving results, depending on initial DNA concentration.
- Increase template DNA purity. Include an ethanol precipitation and wash step prior to amplification to remove inhibitors that copurify with the DNA.
- Add PCR additives. Adding PCR-enhancing agents (e.g., DMSO or betaine) may improve yields. General stabilizing agents such as BSA (Sigma Cat.# A7030; final concentration 0.16mg/ml) also may help to overcome amplification failure.