Certificate of Analysis

GoTaq® MDx Hot Start Polymerase, Glycerol-Free:

For Laboratory Use. Supplied with:

GoTaq® MDx Hot Cat.# Start Polymerase D6201 500 units (D620A) 5X Colorless GoTaq® Flexi Buffer 4 × 1ml (M890A) Magnesium Chloride Solution, 25mM 3 × 0.75ml (A351B)

Description: GoTaq[®] MDx Hot Start Polymerase, Glycerol-Free,^(a) contains GoTaq[®] MDx DNA polymerase bound to a proprietary antibody that blocks polymerase activity. The polymerase activity is restored during the initial denaturation step when amplification reactions are heated at 94–95°C for two minutes. This allows hot-start PCR in which polymerase activity is inhibited at temperatures below 70°C, allowing convenient, room-temperature reaction setup. Hot-start PCR is advantageous for some amplification targets because it may eliminate or minimize primer-dimer and secondary products. In some cases, hot-start PCR may improve yields. This product formulation is further purified to remove glycerol from the product, making is suitable for further manufacturing and lyophilization. GoTaq[®] MDx DNA polymerase is manufactured under cGMP.

Biological Source: The enzyme is derived from bacteria. The antibody is derived from murine cell culture.

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 Colorless GoTaq® 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 MgCl₂ concentration according to your individual requirements. Vortex the MgCl₂ thoroughly after thawing and prior to use.

Quality Control Assays

Test		Specification		Result
		6.5–9.5u/µl		Pass
Concentration		One unit is defined as the amount of enzyme required to catalyze the incorporation of		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.	
		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	se 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.	
		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
		Amplify a hot-start model template to produce a single 1.5kb product, eliminating extraneous amplification products. In a non-hot start PCR, this template produces an additional 410bp amplification product.		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.

Per Wheele

R. Wheeler, Quality Assurance

Part# 9PID620 Revised 5/18



O Promega

Promega Corporation

2800 Woods Hollow Road	t
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 specifications.

(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# 9PID620

Printed in USA Revised 5/18.

Signed by:



1. Standard Application

Reagents to Be Supplied by the User

- dNTP Mix (Cat.# U1511) Nuclease-Free Water (Cat.# P1193)
- upstream primer
 - downstream primer template DNA . mineral oil (optional)
- 1. In a sterile, nuclease-free microcentrifuge tube, combine the following components at room temperature:

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

- 2. If using a thermal cycler without a heated lid, overlay the reaction with 1-2 drops (approximately 50µI) of mineral oil to prevent evaporation during thermal cycling.
- 3. Place reactions in a room-temperature thermal cycler. A 2-minute initial denaturation step at 94-95°C is required to inactivate the antibody and initiate hot-start 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 30 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. Enzyme Concentration

We have found that 1.25 units of GoTag® 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.

B. 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. Primerdimers 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/

C. 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 GoTag® 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.