

Cheetah™ Taq

for hot-starting PCR

September 27, 2010

Speed, Power, and More! **Cheetah™ Taq, a Chemically-modified Hot-start DNA Polymerase for PCR**

- **Fast activation:** *Takes less than 2 minutes to activate, much faster than AmpliTaq® Gold and HotStar™ Taq.*
- **Excellent activity recovery:** *Better recoveries of both 5'-3'-polymerase and 5'-exonuclease activities following activation than AmpliTaq® Gold.*
- **Improved stability:** *Better stability than AmpliTaq® Gold during storage.*
- **pH compatibility:** *Can be activated at alkaline pH (8.5-9) for optimal Taq activity and more specific target amplification.*

Cheetah™ Taq is a chemically modified hot-start Taq DNA polymerase useful for preventing or minimizing nonspecific DNA amplification in PCR. Modified using a novel modifying reagent, Cheetah™ Taq represents a major improvement over AmpliTaq® Gold and other similar chemically-modified hot-start enzymes by having a faster activation time, better shelf-life and alkaline activation condition optimal for both Taq activity and specific amplification. Under standard hot-start conditions (*i.e.*, 94 °C in pH 8-9 Tris), Cheetah™ Taq regains both the 5'-3' polymerase and 5'-exonuclease activities within two minutes. Cheetah™ Taq is also superior to antibody-based hot-start DNA polymerases due to its complete suppression of enzyme activity at room temperature, near full recovery of enzyme activity following activation and lack of animal DNA contaminants. In addition, Cheetah™ Taq is also significantly more economical to manufacture than antibody-based hot-start enzymes.

Cheetah™ Taq is a patent-pending invention. We welcome you to explore licensing opportunities from us.



Cheetah™ Taq Is Ideal for Fast PCR

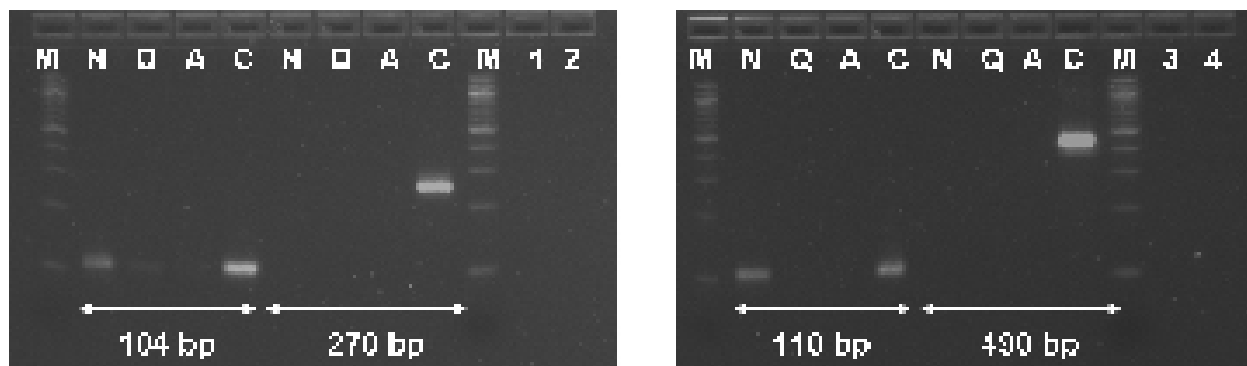


Figure 1. Comparison of PCR products in agarose gel among amplifications using different hot-start Taq DNA polymerases. PCR experiments were performed on four separate fragments of human genomic DNA (104 bp, 270 bp, 110 bp and 490 bp, respectively) using a fast PCR protocol with a 2-minute enzyme activation at 95 °C. Lanes N, Q and A are for enzymes from vendor N, vendor Q and vendor A, respectively. Lanes C are for Cheetah Taq. Lanes 1, 2, 3 and 4 are nontemplate controls (NTC) for Cheetah Taq in the amplifications of fragments 104 bp, 270 bp, 110 bp and 490 bp, respectively. The data show that under the activation condition only Cheetah Taq is capable of producing specific product with high efficiency for all four templates. The enzyme from vendor N produced weak bands for the small fragments (104 bp and 110 bp) but no band at all for the large fragments (270 bp and 490 bp). The enzymes from vendor Q and vendor A exhibited no or insignificant activity as shown by the absence of any band.

Cheetah™ Taq Can Be Activated at pH 8-9

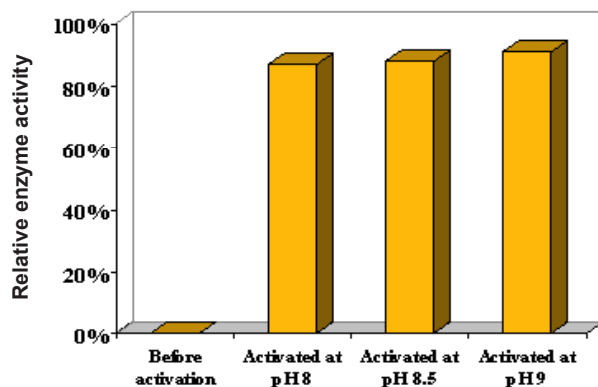


Figure 2. Relative enzyme activity of Cheetah™ Taq following activation at 94 °C in Tris buffer of pH 8.0, pH 8.5 and pH 9.0, respectively. Enzyme activity assays were performed in pH 8.0 Tris buffer at 60 °C. For comparison, the modified Taq without being activated was assayed under identical conditions. The data demonstrates that Cheetah™ Taq can be reliably activated within the pH range from 8 to 9.

Cheetah™ Taq Is Faster to Activate Than Ampli-Taq Gold

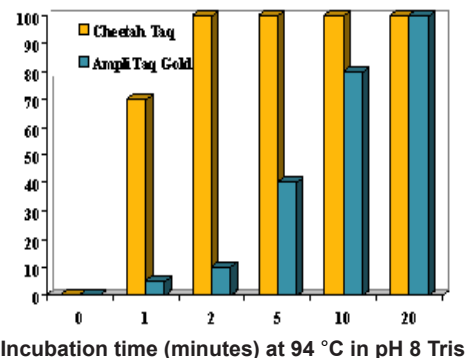


Figure 3. Relative activity of Cheetah™ Taq and Ampli-Taq Gold following different incubation times. All incubations took place with 50 nM of the enzyme in 50 mM pH 8.0 Tris at 94 °C.

Table 1. Cheetah™ Taq Product

Cat.#	Product Name	Unit Size	Unit Price (\$)
29050	Cheetah™ Hot-start Taq DNA Polymerase (Low DNA), 5 U/uL	500 units (100 uL)	210.00
		5 x 500 units (5 x 100 uL)	945.00
		25 x 500 units (25 x 100 uL)	4,200.00

Biotium products are high-quality reagents and materials intended for research purposes only. Some products are potentially hazardous chemicals - please read the Material Safety Data Sheet for additional information regarding handling potentially hazardous chemicals. Several of Biotium products and product applications are covered by U.S. and international patents and pending patents. EvaGreen® dye and applications are covered under patent nos. 7,601,498 and 7,776,567. Our products are not available for resale or other commercial uses without a specific agreement from Biotium, Inc. We welcome inquiries about licensing the use of our dyes, trademarks or technologies. Please submit inquiries by e-mail to btinfo@biotium.com. Names containing the designation ™ are trademarks of Biotium, Inc. ™ Practicing real-time PCR may require additional licensing from Roche or Applied Biosystems, Inc.

Biotium
Glowing Products for Science™

Biotium, Inc.
Tel: (510)265-1027
web: www.biotium.com