



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Cheetah™ FLEX HotStart Taq DNA Polymerase



Product attributes

Storage Conditions Store at -10 to -35 °C

A fast-activating, antibody-free hot-start Taq polymerase engineered for fast-cycling qPCR and advanced applications like proximity extension assays (PEA).

Product Description

Cheetah™ FLEX HotStart Taq is Biotium's proprietary chemically modified Taq polymerase designed for reducing non-specific DNA amplification in PCR. Cheetah™ Taq features a novel chemical modification that completely suppresses its activity prior to activation, and offers a much faster activation time than AmpliTaq Gold® or HotStartTaq®. Cheetah™ FLEX HotStart Taq is validated for fast cycling PCR and proximity extension assays (PEA).

Note: This product is available for licensing and bulk purchase; contact us at techsupport@biotium.com for evaluation samples and to discuss custom size, packaging, or formulation options.

A superior HotStart Taq for FLEXible applications

Antibody-based hot-start Taq polymerases, such as AmpliTaq Gold®, can suffer from inconsistent activation, limited pH tolerance, reduced shelf life, and contamination risk from animal DNA. Cheetah™ FLEX HotStart Taq solves these problems with a proprietary chemical modification that ensures complete inactivity until heat activation—no antibodies, no animal DNA. It offers a rapid 2-minute activation at 95 °C—up to 10x faster than AmpliTaq Gold® or HotStarTaq®. Cheetah™ FLEX is also more stable due to reduced risk of intramolecular cross-linking that can occur in AmpliTaq® Gold. It delivers the same fidelity and efficiency as native Taq polymerase.

Cheetah™ FLEX HotStart Taq is validated for fast-cycling PCR and also third-party validated for qPCR-based proximity extension assay (PEA) applications for protein detection. Cheetah™ FLEX HotStart Taq is available as a ready-to-use kit or standalone enzyme.

Also explore our [Forget-Me-Not™ qPCR Master Mixes](#) featuring Cheetah™ Taq, and [Cheetah™ PEA-PLEX](#), specially formulated for NGS-based, high-throughput PEA.

Product Name	Format	Size	Cat. No.	Price
Cheetah™ FLEX HotStart Taq	500 U Taq with Buffer Pack	Kit (100 uL of 5 U/uL)	29079-KIT	\$250
Enzyme Only	1 mL (5 U/uL)	29079-1ML	\$2,250	
10 mL (5 U/uL)	29079-10ML	\$20,000		
100 mL (5 U/uL)	29079-100ML	\$125,000		
Cheetah™ PEA-PLEX HotStart Taq	Enzyme Only	100 uL (5 U/uL)	29138-100UL	\$275
1 mL (5 U/uL)	29138-1ML	\$2,500		
10 mL (5 U/uL)	29138-10ML	\$22,500		
100 mL (5 U/uL)	29138-100ML	\$150,000		

Please contact us at techsupport@biotium.com for purchasing of bulk 10 mL and 100 mL sizes, or to discuss sampling, custom packaging, or custom formulation options.

AmpliTaq Gold

This datasheet was generated on June 19, 2026 at 08:32:21 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cheetah-flex-hotstart-taq-dna-polymerase/>