



Glowing products for science

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

Cheetah™ PEA-PLEX HotStart Taq DNA Polymerase

A fast-activating, antibody-free hot-start Taq polymerase formulated for NGS-based high-throughput proximity extension assays (PEA).



Product attributes

Storage Conditions

Store at -10 to -35 °C

Product Description

Cheetah™ PEA-PLEX HotStart Taq is a unique formulation of Biotium's chemically modified Cheetah™ Taq polymerase validated by third parties specifically for NGS-based high-throughput proximity extension assays (PEA) for protein detection. Cheetah™ PEA-PLEX is suitable for fast cycling qPCR, as well.

Features

- Unique formulation for NGS-based high-throughput PEA
- Also works for fast cycling qPCR
- Rapid hot-start activation via novel chemical modification
- Antibody-free, low-DNA Taq for cleaner results
- Superior long-term stability compared to other chemically modified Taq
- Bulk formats available

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.

Uniquely formulated for NGS-based high-throughput PEA

Cheetah™ PEA-PLEX HotStart Taq is a highly optimized formulation that was developed specifically for NGS-based high-throughput proximity extension assays (PEA). Similar to [Cheetah™ FLEX HotStart Taq](#), Cheetah™ PEA-PLEX HotStart 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®.

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™ PEA-PLEX 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 HotStartTaq®. Cheetah™ PEA-PLEX 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](#) offers a flexible formulation suitable for both fast cycling PCR and non-NGS PEA assays. Also explore our [Forget-Me-Not™ qPCR Master Mixes](#) featuring Cheetah™ Taq.

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 January 9, 2026 at 08:10:36 AM. Visit product page to check for updated information before use.

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