



Glowing products for science

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

Cheetah™ HotStart Taq DNA Polymerase

This item is discontinued.



Product attributes

Storage Conditions	Store at -10 to -35 °C
Concentration	5U/μL

Product Description

This item is discontinued, please see our replacement products [Cheetah™ FLEX HotStart Taq](#) and [Cheetah™ PEA-PLEX HotStart Taq](#).

Note: Existing stock of Cheetah™ Hot Start Taq DNA Polymerase (Cat. No. 29050) may be available for special order for existing users. Contact us at techsupport@biotium.com to inquire.

This datasheet was generated on January 7, 2026 at 04:08:22 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cheetahm-hotstart-taq-dna-polymerase/>