Glowing products for science

Cheetah™ HotStart Taq DNA Polymerase

This item is discontinued.



Product attributes

Storage Conditions Store at -10 to -35 °C

Concentration 5U/uL

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

This item is discontinued, please see our replacement products <u>Cheetah™ FLEX HotStart Taq</u> and <u>Cheetah™ PEA-PLEX HotStart Taq</u>.

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 November 6, 2025 at 01:33:33 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cheetahtm-hotstart-taq-dna-polymerase/