



Glowing products for science

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

HotStart Polymerase Modification Kit

A simple kit for reversibly modifying thermostable DNA polymerases, rendering the enzyme inactive and preventing primer dimers during reaction set-up.



Product attributes

Storage Conditions

Store at -10 to -35 °C

Product Description

The HotStart Polymerase Modification Kit provides an easy way to reversibly modify lysine residues of thermostable DNA polymerase, rendering the enzyme inactive. The modification is reversed after heating to >90 °C. HotStart modification of DNA polymerase for PCR prevents amplification of non-specific PCR products due to low stringency annealing of primers at low temperature during reaction assembly.

This kit is sold to U.S customers only.

Biotium also offers the WarmStart™ Enzyme Modification Kit ([29053](#)) for temperature control of non-thermostable enzymes.

This datasheet was generated on June 5, 2026 at 02:23:04 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/hotstart-polymerase-modification-kit/>