



Glowing products for science

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

WarmStart™ Enzyme Modification Kit

A simple kit for reversibly modifying non-thermostable enzymes like reverse transcriptase, rendering the enzyme inactive during reaction set-up. Activity is restored by heating to 45° C or higher.



Product attributes

Product Description

The WarmStart™ Enzyme Modification Kit provides an easy way to introduce temperature control of non-thermostable enzymes like reverse transcriptase. The WarmStart™ Enzyme Modification Kit reversibly modifies lysine residues, rendering the enzyme inactive at room temperature. The protein activity is restored by heating to 45° C or higher.

Potential enzymes to be modified:

- Reverse transcriptase
- Bst DNA polymerase
- E. coli DNA polymerase I
- Restriction enzymes
- Nucleases
- Proteases

The WarmStart™ Enzyme Modification Kit is similar to our HotStart Polymerase Modification Kit ([29054](#)).

This datasheet was generated on June 15, 2026 at 07:47:26 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/warmstart-enzyme-modification-kit/>