



Glowing products for science

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

## AntiFix™ Universal Antigen Retrieval Buffer, 10X

AntiFix™ Universal Antigen Retrieval Buffer provides excellent heat unmasking for a wide variety of targets at neutral pH, reducing the need for buffer optimization.



### Product attributes

### Product Description

AntiFix™ Universal Antigen Retrieval Buffer is designed for heat-induced epitope retrieval (HIER) of formalin-fixed, paraffin embedded (FFPE) tissues. The buffer uses a proprietary reagent that catalyzes the reversal of formaldehyde cross-links in FFPE tissue to restore antigen reactivity.

#### Features:

- Excellent antigen retrieval for a wide variety of targets at neutral pH
- Simplifies HIER optimization
- Validated with fluorescence staining using tyramide amplification
- Non-toxic and non-flammable

Fixation of tissue with formalin (formaldehyde) preserves tissue morphology, but the cross-linking of proteins by formaldehyde can modify antibody binding sites or epitopes so they are no longer detectable by immunohistochemistry. Traditionally, epitope reactivity is recovered by using protease treatment, or by heating slides in either low pH citrate buffer or high pH Tris buffer. The optimal method for unmasking varies for different epitopes, requiring multiple buffers to be used for different targets.

AntiFix™ Universal Antigen Retrieval Buffer has been formulated to provide excellent heat unmasking for a wide variety of targets at neutral pH, reducing the need for buffer optimization. It has been validated in immunofluorescence staining of FFPE tissue with our next-generation fluorescent CF® dyes using HRP-mediated tyramide signal amplification.

This datasheet was generated on December 25, 2025 at 09:01:45 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/antifix-universal-antigen-retrieval-buffer-10x/>