



Glowing products for science

Mix-n-Stain™ Enzyme Antibody Labeling Kits

This item is discontinued.

Product Description

This item is discontinued.

References

Download a list of [Mix-n-Stain™ references](#).

This datasheet was generated on June 13, 2026 at 03:02:43 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-enzyme-antibody-labeling-kit/>

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



Product attributes

Product origin	Alkaline Phosphatase: Recombinant, E. coli, Glucose Oxidase: Aspergillus sp., HRP: Horseradish (Armoracia rusticana)
----------------	--