



Glowing products for science

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

Mix-n-Stain™ Fluorescent Protein & Tandem Dye Antibody Labeling Kits

This item is discontinued.



Product attributes

Product Description

This item is discontinued.

References

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

This datasheet was generated on June 9, 2026 at 12:04:51 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-fluorescent-protein-tandem-dyes-antibody-labeling-kit/>