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

This item is discontinued.



## **Product attributes**

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

This item is discontinued.

## References

Download a list of Mix-n-Stain™ references.

This datasheet was generated on November 4, 2025 at 05:29:52 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mix-n-stain-fluorescent-protein-tandem-dyes-antibody-labeling-kit/">https://biotium.com/product/mix-n-stain-fluorescent-protein-tandem-dyes-antibody-labeling-kit/</a>