Glowing products for science

# Mix-n-Stain™ Enzyme Antibody Labeling Kits

This item is discontinued.

#### **Product attributes**

Product origin

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

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 9, 2025 at 05:06:47 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mix-n-stain-enzyme-antibody-labeling-kit/">https://biotium.com/product/mix-n-stain-enzyme-antibody-labeling-kit/</a>