Glowing products for science

## Mix-n-Stain™ Maxi HRP Antibody Labeling Kit, 1 mg Labeling

This item is discontinued.



## **Product attributes**

Conjugation	HRP
Labeling scale	1 mg
Purification method	Purification not required after labeling

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

## **Product Description**

This item is discontinued.

This datasheet was generated on November 2, 2025 at 07:41:36 PM. Visit product page to check for updated information before use. Product link:  $\frac{\text{https://biotium.com/product/mix-n-stain-maxi-hrp-antibody-labeling-kit-1-mg-labeling/}{\text{https://biotium.com/product/mix-n-stain-maxi-hrp-antibody-labeling-kit-1-mg-labeling/}}$