



Glowing products for science

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

This item is discontinued.

Product Description

This item is discontinued.

This datasheet was generated on January 1, 2026 at 10:30:39 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-maxi-hrp-antibody-labeling-kit-1-mg-labeling/>



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

Product attributes

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