



Glowing products for science

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

Mix-n-Stain™ DNP Antibody Labeling Kit

Label your antibody with DNP in just 15 minutes without a purification step. Labeling tolerates many common buffer components including BSA and ascites.



Product attributes

Product Description

Mix-n-Stain™ DNP Antibody Labeling Kits allow you to label ≤ 5 ug up to 100 ug of your antibody with dinitrophenol (DNP) in just 15 minutes, with minimal hands-on time and no purification.

Unrivaed Convenience

Mix-n-Stain™ labeling tolerates common antibody buffer formulations. An ultrafiltration centrifuge vial is provided to quickly remove interfering substances like glycerol, or to concentrate your antibody. With a slight modification in protocol, antibodies can be labeled in the presence of BSA, gelatin, or ascites fluid. The presence of other proteins like BSA or gelatin in the labeling reaction has minimal effect on background fluorescence, because any labeled non-antibody proteins readily wash away during immunofluorescence staining.

Choose the Right Labeling Kit for Your Antibody

Mix-n-Stain™ Antibody Labeling Kits are very simple to use (see the workflow overview above). But before you begin, you must check that your antibody meets the compatibility requirements for labeling, and choose the right labeling protocol. Download the updated and expanded [Product Protocol](#) for a labeling check list to help you select the right kit size and labeling protocol for your antibody. Also see our [Kit Compatibility and Protocol Selection Flowchart](#).

We offer Mix-n-Stain™ Kits for labeling antibodies, nanobodies, and small ligands with a wide selection of fluorescent dyes, enzymes, and other labels. See our full selection of Mix-n-Stain™ Antibody Labeling Kits below.

Antibody & Protein Labeling Kits

Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ CF® Dye Antibody Labeling Kits	CF® Dyes	≤5-20 ug IgG 20-50 ug IgG 50-100 ug IgG	~ 15 min.	<ul style="list-style-type: none">• Rapid, simple labeling• No purification• Tolerates BSA & other additives• Dyes for super-resolution, spectral flow, & NIR detection
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			
Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits	Cyanine 555 Cyanine 647			
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxigenin (DIG)			
Mix-n-Stain™ DNP Antibody Labeling Kits	DNP			
Mix-n-Stain™ Maxi Antibody Labeling Kits	CF® Dyes Cyanine Dyes	1 mg IgG	~ 30 min.	
Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits	CF® Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	<ul style="list-style-type: none">• Rapid, simple labeling• No purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	<ul style="list-style-type: none">• Optimized to provide low 1-2.5 DOL• No purification• Optimized for Nanobodies®• Tolerates BSA & other additives
Mix-n-Stain™ Nanobody Labeling Kits	CF® Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	
Mix-n-Stain™ Nanobody Thiol Labeling Kits	CF® Dyes	~ 2 hours	• Optimized for Nanobodies® with single exposed cysteine residue	
Mix-n-Stain™ Small Ligand Labeling Kits	CF® Dyes	0.1 umol small ligand	~ 30 min.	<ul style="list-style-type: none">• Label SNAP®, CLIP™, HaloTag® & other small ligands• Dye options for surface or intracellular targets
VivoBrite™ Rapid Antibody Labeling Kits for Small Animal In Vivo Imaging	Near-IR CF® Dyes	3 x 1 mg IgG	~ 2 hours	<ul style="list-style-type: none">• Dyes, buffers, & spin vials for labeling + purification• Sterile syringes, filters and storage vials for in vivo use
CF® Dye & Biotin SE Protein Labeling Kits	CF® Dyes Biotin	3 x 1 mg protein	~ 2 hours	<ul style="list-style-type: none">• Dyes, buffers, & spin vials for labeling + purification

References

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

This datasheet was generated on January 27, 2026 at 06:28:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mix-n-stain-dnp-antibody-labeling-kit/>