



Glowing products for science

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

Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits

Our revolutionary technology allows antibody labeling with Cyanine 555 or Cyanine 647 in 15 minutes without a purification step. The labeling procedure tolerates many common buffer components and antibody stabilizers.



Product attributes

Product Description

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

Unrivaled 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. | <ul style="list-style-type: none"> • Optimized for Nanobodies® with single exposed cysteine residue |
| Mix-n-Stain™ Nanobody Thiol Labeling Kits | CF® Dyes | ~ 2 hours | | |
| 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 June 20, 2026 at 03:22:19 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cyanine-dye-antibody-labeling-kits/>