

## Mix-n-Stain™ Nanobody Thiol Labeling Kits

Label 5-50 ug of VHH (also known as Nanobodies®) containing a single free cysteine with Biotium's bright and photostable CF® Dyes, with no purification after labeling.



### Product attributes

### Product Description

Mix-n-Stain™ Nanobody Thiol Labeling Kits are designed for labeling VHH (also known as Nanobody® or single-domain antibodies) that contain a single free cysteine. The labeling is done within 2 hours in ambient conditions, and requires minimal hands-on time and no purification.

**Note:** This kit is for labeling VHH via a single cysteine (1x Cys) only. Please see our [Mix-n-Stain™ Nanobody Labeling Kits](#) for standard labeling of VHH, or [view our full selection of labeling kits for labeling IgG antibodies or other biomolecules](#).

### Superior CF® Dyes

These kits feature a choice of Biotium's next-generation CF® Dyes, which were designed for advantages in brightness and photostability compared to other commercially available fluorescent dyes. Learn more about [CF® Dyes](#).

### Choosing the Right Kit for Your Application

Before selecting a Mix-n-Stain™ Nanobody Thiol Labeling Kit, please download the [Product Protocol](#) to see our Kit Compatibility guidelines.

For labeling IgG antibodies, please use one of our Mix-n-Stain™ Antibody Labeling Kits. We also offer Mix-n-Stain™ Small Ligand Labeling Kits, for labeling SNAP-tag®, CLIP-tag™, and HaloTag® ligands, or other small molecules. See Mix-n-Stain™ Product Lines below for our full selection of antibody, protein, and ligand labeling kits.

## Mix-n-Stain™ Nanobody Thiol Labeling Kits

Conjugation	Ex/Em	Labeling size (ug Nanobody®)	Catalog No.	Dye Features
<a href="#">CF@488A</a> 20-50 ug	490/516 nm <a href="#">92586</a>	5-20 ug	<a href="#">92585</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a> 20-50 ug	562/584 nm <a href="#">92588</a>	5-20 ug	<a href="#">92587</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@583R</a> 20-50 ug	585/609 nm <a href="#">92590</a>	5-20 ug	<a href="#">92589</a>	<a href="#">CF@583R Features</a>
<a href="#">CF@594</a> 20-50 ug	593/615 nm <a href="#">92592</a>	5-20 ug	<a href="#">92591</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@647</a> 20-50 ug	652/668 nm <a href="#">92594</a>	5-20 ug	<a href="#">92593</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@680</a> 20-50 ug	681/698 nm <a href="#">92596</a>	5-20 ug	<a href="#">92595</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@740</a> 20-50 ug	742/767 nm <a href="#">92598</a>	5-20 ug	<a href="#">92597</a>	<a href="#">CF@740 Features</a>

Note: Before selecting a kit, please download the [Product Protocol](#) to see Kit Compatibility guidelines.

## Antibody & Protein Labeling Kits

Product Name	Label Options	Labeling Scale	Labeling Time	Features
<a href="#">Mix-n-Stain™ CF® Dye Antibody Labeling Kits</a>	CF® Dyes	≤5-20 ug IgG 20-50 ug IgG 50-100 ug IgG	~ 15 min.	<ul style="list-style-type: none"> <li>• Rapid, simple labeling</li> <li>• No purification</li> <li>• Tolerates BSA &amp; other additives</li> <li>• Dyes for super-resolution, spectral flow, &amp; NIR detection</li> </ul>

[Mix-n-Stain™ FITC Antibody Labeling Kits](#)

FITC

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

Nanobody is a registered trademark of Ablynx, a Sanofi company. Alexa Fluor and DyLight are registered trademarks of Thermo Fisher Scientific. Cy Dye is a registered trademark of Cytiva. SNAP-tag and CLIP-tag are trademarks or registered trademarks of New England BioLabs; HaloTag is a registered trademark of Promega Corporation.

## References

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

This datasheet was generated on June 9, 2026 at 03:34:55 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mix-n-stain-nanobody-thiol-labeling-kits/>