Mix-n-Stain™ Nanobody Thiol Labeling

Product attributes

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

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 Description

Kits

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 lgG 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 <u>Product Protocol</u> to see our Kit Compatibility guidelines.

For labeling IgG antibodies, please use one of our Mix-n-StainTM Antibody Labeling Kits. We also offer Mix-n-StainTM Small Ligand Labeling Kits, for labeling SNAP-tag®, CLIP-tagTM, and HaloTag® ligands, or other small molecules. See Mix-n-StainTM 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
CF®488A	490/516 nm	5-20 ug	<u>92585</u>	CF®488A Features
		20-50 ug	<u>92586</u>	
<u>CF®568</u>	562/584 nm	5-20 ug	<u>92587</u>	CF®568 Features
		20-50 ug	<u>92588</u>	
CF®583R	585/609 nm	5-20 ug	<u>92589</u>	CF®583R Features
		20-50 ug	<u>92590</u>	
CF®594	593/615 nm	5-20 ug	<u>92591</u>	CF®594 Features
		20-50 ug	<u>92592</u>	
<u>CF®647</u>	652/668 nm	5-20 ug	<u>92593</u>	CF®647 Features
		20-50 ug	<u>92594</u>	
<u>CF®680</u>	681/698 nm	5-20 ug	<u>92595</u>	CF®680 Features
		20-50 ug	<u>92596</u>	
<u>CF®740</u>	742/767 nm	5-20 ug	<u>92597</u>	CF®740 Features
		20-50 ug	<u>92598</u>	

Note: Before selecting a kit, please download the Product Protocol to see Kit Compatibility guidelines.

Antibody & Protein Labeling Kits

Antibody Labeling Kits

· · · · · · · · · · · · · · · · · · ·						
Product Name	Label Options	Labeling Scale	Labeling Time	Features		
Mix-n-Stain™ CF® Dye Antibody Labeling Kits	CF® Dyes	≤5-20 ug lgG 20-50 ug lgG 50-100 ug lgG	~ 15 min.	Rapid, simple labeling		
				 No purification 		
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			Tolerates BSA & other additives		
Mix-n-Stain™ Cyanine	Cyanine 555			additives		
Dye Antibody Labeling Kits	Cyanine 647			Dyes for super-resolution, spectral flow, & NIR detection		
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin					
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxygenin (DIG)					
Mix-n-Stain™ DNP	DNP					

Product Name	Label Options	Labeling Scale	Labeling Time	Features
Mix-n-Stain™ Maxi Antibody Labeling Kits	CF® Dyes Cyanine Dyes	1 mg lgG	~ 30 min.	
Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits	CF® Dyes FITC	25 ug IgM 100 ug IgM	15-30 min.	Rapid, simple labelingNo purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	Optimized to provide low 1-2.5 DOLNo purification
Mix-n-Stain™ Nanobody Labeling Kits	CF® Dyes	5-20 ug Nanobody® 20-50 ug Nanobody®	~ 30 min.	Optimized for Nanobodies® Tolerates BSA & other additives
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.	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 lgG	~ 2 hours	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	Dyes, buffers, & spin vials for labeling + purification

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 $\underline{\mathsf{CF} \$}$ Dye and Mix-n-Stain $\!\!\!\!\mathsf{TM}$ references.

This datasheet was generated on November 9, 2025 at 10:49:45 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mix-n-stain-nanobody-thiol-labeling-kits/