



Glowing products for science

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

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
CF@488A 20-50 ug	490/516 nm 92586	5-20 ug	92585	CF@488A Features
CF@568 20-50 ug	562/584 nm 92588	5-20 ug	92587	CF@568 Features
CF@583R 20-50 ug	585/609 nm 92590	5-20 ug	92589	CF@583R Features
CF@594 20-50 ug	593/615 nm 92592	5-20 ug	92591	CF@594 Features
CF@647 20-50 ug	652/668 nm 92594	5-20 ug	92593	CF@647 Features
CF@680 20-50 ug	681/698 nm 92596	5-20 ug	92595	CF@680 Features
CF@740 20-50 ug	742/767 nm 92598	5-20 ug	92597	CF@740 Features

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
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

Product Name	Label Options	Labeling Scale	Labeling Time	Features
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
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® • 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.	<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

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 January 11, 2026 at 02:44:07 PM. Visit product page to check for updated information before use.

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