Mix-n-Stain™ CF® Dye Small Ligand Labeling Kits

Easily label SNAP-tag®, CLIP-tag™, HaloTag®, or other ligands with fluorescent dyes, no purification needed. Cell membrane-permeant and impermeant dye options for intracellular or surface targets.



Product attributes

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

Product Description

Mix-n-Stain™ Small Ligand Labeling Kits are designed for rapid labeling of low molecular weight (MW ~150 – 5000), high affinity biological ligands or substrates.

- For SNAP-tag®, CLIP-tag™ and HaloTag®, or other ligands with an aliphatic amine
- Label 100 nmol ligand
- Reaction takes ~30 minutes with minimal hands-on time and no purification
- Choose from 15 bright & photostable CF® Dyes
- Cell membrane-permeant or impermeant options for intracellular or surface targets

For more information, please download the <u>Mix-n-Stain™ Small Ligand Labeling Kit Flyer</u> and the <u>Product Information Sheet</u>. We also offer <u>TMP-PEG3-Amine</u> for making fluorescent TMP ligands. Also see our <u>Mix-n-Stain™ Antibody Labeling Kits</u> and other <u>Reactive Dyes and Biotin Reagents</u>.

Mix-n-Stain™ Small Ligand Labeling Kits

Dye	Catalog no.	Spectrally similar dyes	Ex/Em (nm)	Staining
CF®405M	92362	Pacific Blue™, BD Horizon™ V450	408/452	Surface
<u>CF®408</u>	92356	Pacific Blue™, BD Horizon™ V450	408/450	Intracellular
CF®488A	92350	FITC, Alexa Fluor® 488	490/515	Surface
<u>CF®500</u>	92357	FITC, Alexa Fluor® 488	500/510	Intracellular
CF®540	<u>92358</u>	TAMRA, TRITC, Alexa Fluor® 546, Cy®3	540/570	Intracellular
<u>CF®555</u>	92364	Alexa Fluor® 555, Cy®3, TRITC	555/585	Intracellular
CF®568	<u>92351</u>	Rhodamine Red, Alexa Fluor® 568	562/583	Surface
<u>CF®594</u>	92352	Texas Red®, Alexa Fluor® 594	593/614	Surface
CF®633	92353	Cy®5, Alexa Fluor® 633	630/650	Surface
<u>CF®640R</u>	<u>92354</u>	Cy®5, Alexa Fluor® 647	642/662	Surface
CF®647	92359	Cy®5, Alexa Fluor® 647	650/665	Surface
<u>CF®650</u>	<u>92363</u>	Cy®5, Alexa Fluor® 647	650/670	Intracellular
CF®660C	<u>92360</u>	Alexa Fluor® 660	667/685	Surface
<u>CF®680</u>	<u>92361</u>	Cy®5.5, Alexa Fluor® 660, IRDye®680LT	681/698	Surface
<u>CF®680R</u>	<u>92355</u>	Cy®5.5, Alexa Fluor® 660, IRDye®680LT	680/701	Surface

Pacific Blue, Texas Red, and Alexa Fluor are trademarks or registered trademarks of Thermo Fisher Scientific; BD Horizon is a trademark of BD Biosciences; Cy Dye is a registered trademark of Cytiva; IRDye is a registered trademark of LI-COR Biosciences.

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 lgG 20-50 ug lgG 50-100 ug lgG	~ 15 min.	Rapid, simple labeling
	FITO			 No purification
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			 Tolerates BSA & other additives
Mix-n-Stain™ Cyanine Dye Antibody Labeling Kits	Cyanine 555 Cyanine 647			Dyes for super-resolution, spectral
Mix-n-Stain™ Biotin Antibody Labeling Kits	Biotin			flow, & NIR detection
Mix-n-Stain™ Digoxigenin Antibody Labeling Kits	Digoxygenin (DIG)			
Mix-n-Stain™ DNP Antibody Labeling Kits	DNP			
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 lgM 100 ug lgM	15-30 min.	Rapid, simple labeling
				No purification
Mix-n-Stain™ STORM CF® Dye Antibody Labeling Kits	CF® Dyes	50 ug IgG	~ 30 min.	Optimized to provide low 1-2.5 DOL
	05e D	500 N L L	00 '	No 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

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 curated CF® Dye references.

This datasheet was generated on November 4, 2025 at 08:13:34 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mix-n-stain-cf-dye-small-ligand-labeling-kit/