

## 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

### 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 [Mix-n-Stain™ Small Ligand Labeling Kit Flyer](#) and the [Product Information Sheet](#). We also offer [TMP-PEG3-Amine](#) for making fluorescent TMP ligands. Also see our [Mix-n-Stain™ Antibody Labeling Kits](#) and other [Reactive Dyes and Biotin Reagents](#).

### Mix-n-Stain™ Small Ligand Labeling Kits

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

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 January 7, 2026 at 04:08:24 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/>