



Glowing products for science

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

## VivoBrite™ Rapid Antibody Labeling Kits for Small Animal In Vivo Imaging

Kits for labeling & purification of near-IR antibody conjugates for small animal *in vivo* imaging, featuring Biotium's industry-leading near-IR CF® Dyes.



### Product attributes

## Product Description

This labeling kit contains reagents for labeling antibodies with our superior near-infrared amine-reactive CF® dyes for small animal *in vivo* imaging. Learn more about Biotium's superior [Near-IR CF® dyes](#).

See our full line of [Antibody & Protein Labeling Kits](#).

## VivoBrite™ Rapid Antibody Labeling Kits for Small Animal In Vivo Imaging

| Product  | Ex/Em      | Size               | Catalog No.           |
|--|------------|--------------------|-----------------------|
| CF®680 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging | 681/698 nm | 3 x 1 mg labelings | <a href="#">92160</a> |
| CF®750 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging | 755/777 nm | 3 x 1 mg labelings | <a href="#">92161</a> |
| CF®770 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging | 770/797 nm | 3 x 1 mg labelings | <a href="#">92162</a> |
| CF®790 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging | 784/806 nm | 3 x 1 mg labelings | <a href="#">92163</a> |

## 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<br>20-50 ug IgG<br>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<br>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<br>Cyanine Dyes   | 1 mg IgG                                      | ~ 30 min.     |  |
| <a href="#">Mix-n-Stain™ CF® Dye IgM Antibody Labeling Kits</a> | CF® Dyes<br>FITC           | 25 ug IgM<br>100 ug IgM                       | 15-30 min.    | <ul style="list-style-type: none"><li>• Rapid, simple labeling</li><li>• No purification</li></ul>   |

| Product Name   | Label Options      | Labeling Scale                          | Labeling Time  | Features   |
|--|--------------------|---|--|--|
| <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> </ul>   |
| <a href="#">Mix-n-Stain™ Nanobody Labeling Kits</a>                                      | CF® Dyes           | 5-20 ug Nanobody®<br>20-50 ug Nanobody® | ~ 30 min.  | <ul style="list-style-type: none"> <li>• No purification</li> <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<br>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>  |

## References

1. J Immunother Cancer (2019) 7, 55. <https://doi.org/10.1186/s40425-019-0537-9>
2. Biomolecules (2020) 10, 51. [doi:10.3390/biom10010051](https://doi.org/10.3390/biom10010051)

Download a list of [CF® dye references](#).

This datasheet was generated on January 6, 2026 at 05:15:03 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/vivobrite-rapid-antibody-labeling-kits-for-small-animal-in-vivo-imaging/>