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

357

Product attributes

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

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 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	92160
CF®750 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging	755/777 nm	3 x 1 mg labelings	<u>92161</u>
CF®770 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging	770/797 nm	3 x 1 mg labelings	92162
CF®790 VivoBrite™ Rapid Antibody Labeling Kit for Small Animal In Vivo Imaging	784/806 nm	3 x 1 mg labelings	92163

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
, ,				 No purification
Mix-n-Stain™ FITC Antibody Labeling Kits	FITC			Tolerates BSA & other additives
Mix-n-Stain™ Cyanine Dye	Cyanine 555 Cyanine 647			
Antibody Labeling Kits	Gyariine 047			 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 TM CF® Dye	CF® Dyes FITC	25 ug lgM 100 ug lgM	15-30 min.	 Rapid, simple labeling
IgM Antibody Labeling Kits	FIIC			No purification
Mix-n-Stain™ STORM CF® Dye Antibody		50 ug IgG	~ 30 min.	Optimized to provide low 1-2.5 DOL
Labeling Kits				No purification

Product Name	Label Options	Labeling Scale	Labeling Time	Features
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

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

Download a list of CF® dye references.

This datasheet was generated on November 4, 2025 at 03:34:56 AM. 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/