

Product Information

Llama Anti-Mouse IgG (H+L), Whole Antibody, CF® Dye Conjugates

| CF® Dye | Ex/Em | Catalog No. | Size |
|---------|------------|-------------|--------------------|
| CF®488A | 490/515 nm | 20454-50uL | 50 uL |
| | | 20454-500uL | 500 uL |
| | | 20454-1mg | 1 mg (lyophilized) |
| CF®568 | 562/583 nm | 20455-50uL | 50 uL |
| | | 20455-500uL | 500 uL |
| | | 20455-1mg | 1 mg (lyophilized) |
| CF®594 | 593/614 nm | 20456-50uL | 50 uL |
| | | 20456-500uL | 500 uL |
| | | 20456-1mg | 1 mg (lyophilized) |
| CF®640R | 630/650 nm | 20457-50uL | 50 uL |
| | | 20457-500uL | 500 uL |
| | | 20457-1mg | 1 mg (lyophilized) |
| CF®647 | 650/665 nm | 20458-50uL | 50 uL |
| | | 20458-500uL | 500 uL |
| | | 20458-1mg | 1 mg (lyophilized) |

Concentration

Liquid format: 2 mg/mL in pH ~7.4 PBS containing 50% glycerol, 2 mg/ml bovine serum albumin (IgG-free and protease-free) and 0.05% sodium azide.

Lyophilized format (after reconstitution): 2 mg/mL in pH ~7.4 PBS containing 15 mg/mL bovine serum albumin (IgG-free and protease-free) and 20 mg/mL trehalose.

Storage and Handling

Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Liquid format antibodies contain 50% glycerol and will not freeze at -20°C.

Reconstitution (lyophilized format only): add 0.5 mL dH₂O and mix gently to dissolve. Store at -20°C, protected from light. Aliquot to avoid freeze-thaw cycles. Alternatively, glycerol can be added to the antibody so that it will not freeze at -20°C: add 0.25 mL dH₂O to the lyophilized antibody and mix gently to dissolve, then add 0.25 mL glycerol and mix well. Optional: a preservative may be added, such as 0.05% (final concentration) sodium azide.

Note: storage of the antibody for more than a day at final working dilution is not recommended.

Product Description

Llama anti-mouse IgG (H+L), whole antibody is an affinity purified llama antibody that recognizes mouse IgG. This antibody is not highly-cross adsorbed against other species. Visit www.biotium.com to see our highly cross-adsorbed secondary antibodies for multiple labeling.

CF® dyes are superior to Alexa Fluor®, DyLight®, and Cy® dyes for antibody labeling by having combined advantages in brightness, photostability, specificity and novel features ideal for in vivo imaging. For more information about CF® dyes, download our CF® Dye Selection Guide at www.biotium.com.

Recommended Dilution Range

Fluorescence microscopy: 1-2 ug/mL

Flow cytometry: 1 ug/10⁶ cells

Near-infrared western detection: 50-100 ng/mL

These concentrations are provided as a starting point for optimization; appropriate dilutions should be determined empirically. Generally IgG conjugates are used in the range of 1-10 ug/mL.

For more detailed protocols for immunofluorescence staining for fluorescence microscopy, flow cytometry, or fluorescence-based western detection, please visit our website.

Related Products

| Cat. No. | Product Name | Features |
|----------|--|---|
| 23001 | EverBrite™ Mounting Medium | <ul style="list-style-type: none"> • Excellent protection from photobleaching • Compatible with a wide variety of dyes including cyanine dyes • Available in wet-set or hardset formulations • With or without DAPI |
| 23002 | EverBrite™ Mounting Medium with DAPI | |
| 23003 | EverBrite™ Hardset Mounting Medium | |
| 23004 | EverBrite™ Hardset Mounting Medium with DAPI | |
| 23005 | CoverGrip™ Coverslip Sealant | <ul style="list-style-type: none"> • Replaces nail polish for coverslip sealing • Won't mix with aqueous mounting media |
| 23007 | TrueBlack™ Lipofuscin Autofluorescence Quencher | <ul style="list-style-type: none"> • Eliminates lipofuscin autofluorescence with less background than Sudan Black B • Reduces background from other sources • Can be used before or after IF staining |
| 40061 | RedDot™2 Far Red Nuclear Counterstain | <ul style="list-style-type: none"> • Far-red nuclear dye for the Cy®5 channel • More specific than Draq®7 |
| 22005 | Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses | <ul style="list-style-type: none"> • Create hydrophobic barriers around tissue sections • Heat-stable to 120°C • Insoluble in aqueous buffers, detergents, alcohol and acetone; can be removed with xylene |
| 22006 | Super ^{HT} Pap Pen 4 mm tip, ~800 uses | |
| 22003 | Mini-Cell Scrapers | <ul style="list-style-type: none"> • Collect cells from 96-well to 24-well plates • Polypropylene, sterile, disposable |
| 23006 | Flow Cytometry Fixation/Permeabilization Kit | <ul style="list-style-type: none"> • Ready-to-use fix/perm buffers for intracellular staining |
| 22015 | Fixation Buffer | <ul style="list-style-type: none"> • Convenient buffers and blocking agents for immunofluorescence or western |
| 22016 | Permeabilization Buffer | |
| 22017 | Permeabilization and Blocking Buffer | |
| 22010 | 10X Fish Gelatin Blocking Agent | |
| 22011 | Fish Gelatin Powder | |
| 22014 | 30% Bovine Serum Albumin Solution | |
| 22002 | Tween®-20 | |

Please visit www.biotium.com to view our full selection of products featuring bright and photostable fluorescent CF® dyes, including Mix-n-Stain™ antibody labeling kits, primary antibody conjugates, streptavidin, phalloidin, and other bioconjugates, as well as conjugates of biotin, HRP, AP, R-PE, APC and PerCP.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

CF dye technology is covered by pending US and international patents. Alexa Fluor is a registered trademark of Molecular Probes; CYDYE is a registered trademark of GE Healthcare; Draq7 is a registered trademark of Biostatus Ltd; DyLight is a registered trademark of Pierce Biotechnology; TWEEN is a registered trademark of Uniqema Americas LLC.