

Product Information

Isotype-specific goat anti-mouse antibodies, CF® dye conjugates

Highly cross-adsorbed for multiple labeling (min X Bv, Hu, Rb)

Antibody	CF® dye/catalog number													
	CF350	CF405S	CF488A	CF543	CF555	CF568	CF594	CF633	CF640R	CF647	CF680	CF680R	CF750	CF770
Goat Anti-Mouse IgG1 (γ1)	20245	20380	20246	20325	20247	20248	20249	20250	20251	20252	20253			20254
Goat Anti-Mouse IgG2a (γ2a)	20255	20381	20256	20326	20257	20258	20259	20260	20261	20262	20263	20842		20264
Goat Anti-Mouse IgG2b (γ2b)	20265	20382	20266	20327	20267	20268	20269	20270	20271	20272	20273		20430	20274
Goat Anti-Mouse IgG3 (γ3)													20431	

Unit Size:

50 uL or 0.25 mL

Concentration:

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.

Spectral Properties

($\lambda_{abs}/\lambda_{em}$ maxima in pH 7.4 PBS buffer)

Conjugate	Abs (nm)	Em (nm)
CF@350	347	448
CF@405S	404	431
CF@488A	490	515
CF@543	541	560
CF@555	555	565
CF@568	562	583
CF@594	593	614
CF@633	630	650
CF@640R	642	662
CF@647	650	665
CF@680	681	698
CF@680R	680	701
CF@770	770	797

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. Antibodies contain 50% glycerol and will not freeze at -20°C.

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

Product Description

Staining of mouse tissues with anti-mouse secondary antibodies results in background staining due to reactivity with endogenous mouse immunoglobulins. Isotype-specific anti-mouse secondary antibodies are useful for staining mouse tissues because IgG is present at relatively low levels in mouse tissues. Biotium offers goat anti-mouse isotype specific antibodies conjugated to a wide selection of

our outstanding series of CF® dyes developed by Biotium. CF® dyes are superior to leading commercial dyes for antibody labeling, with combined advantages in brightness, photostability, specificity and novel features ideal for in vivo imaging (Please visit our website for details). A concentration of 1 - 10 ug/mL is typical for most immunofluorescence staining applications, however, optimal conditions should be determined empirically.

Related Products

Cat.#	Product Name	Unit Size
40061-T	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO, Trial Size (15-20 tests)	25 uL
23001	EverBrite™ Mounting Medium	10 mL
23002	EverBrite™ Mounting Medium with DAPI	10 mL
23003	EverBrite™ Hardset Mounting Medium	10 mL
23004	EverBrite™ Hardset Mounting Medium with DAPI	10 mL
23005	CoverGrip™ Coverslip Sealant	15 mL
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses	1 pen
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses	1 pen
22015	Fixation Buffer	100 mL
22016	Permeabilization Buffer	100 mL
22017	Permeabilization and Blocking Buffer	100 mL
22010	10X Fish Gelatin Blocking Agent	100 mL
22011	Fish Gelatin Powder	2 x 50 g
22014	30% Bovine Serum Albumin Solution	100 mL
22002	Tween®-20	50 mL

Please visit www.biotium.com to view our full selection of products featuring bright and photostable fluorescent CF® dyes, including secondary antibodies and Mix-n-Stain™ antibody labeling kits, and R-PE dye conjugates. Biotium also offers a variety of apoptosis and cell viability assays for flow cytometry analysis, including mitochondrial membrane potential dyes, fluorescent Annexin V conjugates, and NucView® 488 Caspase-3 Substrate for live cells.

CF® dye technology is covered by pending US and international patents. TWEEN is a registered trademark of Uniqema Americas LLC.

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