

Product Information

CF® Dye/HRP Conjugated Goat Anti-Digoxigenin (DIG) Antibodies

Conjugation	Ex/Em (nm)	Catalog No.	Unit Size
CF®405S	404/431	20858-50uL	50 uL
		20858-250uL	250 uL
CF®488A	490/515	20859-50uL	50 uL
		20859-250uL	250 uL
CF®568	562/583	20860-50uL	50 uL
		20860-250uL	250 uL
CF®594	593/614	20861-50uL	50 uL
		20861-250uL	250 uL
CF®640R	642/662	20862-50uL	50 uL
		20862-250uL	250 uL
CF®680R	680/701	20863-50uL	50 uL
		20863-250uL	250 uL
HRP	--	20864-100uL	100 uL

Concentration:

CF® dye conjugates: 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.

HRP conjugate: 1 mg/mL in pH ~7.4 PBS containing 50% glycerol and 15 mg/ml bovine serum albumin (IgG-free and protease-free).

Storage and Handling

Product is stable for at least 6 months at -20°C as an undiluted liquid. Storage of the antibody for more than a day at final working dilution is not recommended. Protect from light.

Product Description

Goat anti-DIG, whole antibody, is an affinity purified goat antibody that reacts with Digoxigenin and is conjugated to Biotium's bright and photostable CF® dyes or HRP.

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

Catalog No.	Product
40078	Digoxigenin-dUTP, Alkali stable
92328-92330	Mix-n-Stain Digoxigenin labeling kit
40061	RedDot™2 Far Red Nuclear Counterstain
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23014	TrueBlack® Plus Lipofuscin Autofluorescence Quencher
23007	TrueBlack® Lipofuscin Autofluorescence Quencher
23005	CoverGrip™ Coverslip Sealant
22005	Mini Super ^{HT} Pap Pen 2.5 mm tip, ~400 uses
22006	Super ^{HT} Pap Pen 4 mm tip, ~800 uses
22015	Fixation Buffer
22016	Permeabilization Buffer
22017	Permeabilization and Blocking Buffer
22010	10% Fish Gelatin Blocking Buffer
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 streptavidin and other bioconjugates, secondary antibodies, Mix-n-Stain™ antibody labeling kits, and R-PE 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 U.S. and international patents.

Alexa Fluor is a registered trademark of Molecular Probes; Cy dye is a registered trademark of GE Healthcare; DyLight is a registered trademark of Pierce Biotechnology; 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.