

# Product Information

## CF® Dye/HRP Conjugated Rabbit Anti-DNP Antibodies

Conjugate	Ex/Em (nm)	Catalog no.	Unit size
CF@405S	404/431	20865-50uL	50 uL
		20865-250uL	250 uL
CF@488A	490/515	20866-50uL	50 uL
		20866-250uL	250 uL
CF@568	562/583	20867-50uL	50 uL
		20867-250uL	250 uL
CF@594	593/614	20868-50uL	50 uL
		20868-250uL	250 uL
CF@640R	642/662	20869-50uL	50 uL
		20869-250uL	250 uL
CF@680R	680/701	20870-50uL	50 uL
		20870-250uL	250 uL
CF@750	755/777	20875-50uL	50 uL
		20875-250uL	250 uL
HRP	---	20871-100uL	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 conjugates: 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

Rabbit anti-DNP, whole antibody is an affinity purified rabbit antibody that reacts with dinitrophenol (DNP) hapten for detection of DNP conjugated proteins and nucleic acids. The antibodies are conjugated to Biotium's bright and photostable CF® dyes and 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](http://www.biotium.com).

### Recommended Dilution Range

Fluorescence microscopy: 1-2 ug/mL

Flow cytometry: 1 ug/10<sup>6</sup> 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 Number	Product
92325-92327	Mix-n-Stain™ DNP Antibody Labeling Kit
96019	DNP Tyramide
30015	DAB Substrate Kit
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23014	TrueBlack® Plus Lipofuscin Autofluorescence Quencher
23007	TrueBlack® Lipofuscin Autofluorescence Quencher
22030	AntiFix™ Universal Antigen Retrieval Buffer
40061	RedDot™2 Far Red Nuclear Counterstain
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses
22003	Mini-Cell Scrapers
23006	Flow Cytometry Fixation/Permeabilization Kit
22015	Fixation Buffer
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
22023	Paraformaldehyde, 4% in PBS, Ready-To-Use Fixative

Please visit [www.biotium.com](http://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; 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.