

RFP Polyclonal Rabbit Antibody

Polyclonal rabbit anti-RFP antibody labeled with our superior CF® dyes.



Product Description

This is a polyclonal rabbit anti-RFP antibody labeled with our superior CF® dyes. The antibody is designed to detect red fluorescent protein (RFP) and its variants: mCherry, tdTomato, mBanana, mOrange, mPlum, and mStrawberry. To minimize cross reactivity, the antibodies are cross-adsorbed against human, mouse, and rat proteins.

- Cross-adsorbed for specific staining with minimal background
- Available in 7 bright and photostable CF® dyes
- Suitable for western, immunofluorescence, and immunohistology

See our full selection of [anti-tag and anti-hapten antibody conjugates](#).

Product attributes

Clonality	Polyclonal
Antibody type	Tag Antibody
Concentration	1 mg/mL
Host species	Rabbit
Antibody reactivity (target)	RFP
Cross adsorption	Human, Mouse, Rat
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide, Lyophilized: PBS/15 mg/mL BSA/20 mg/mL trehalose after reconstitution
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

RFP Polyclonal Rabbit Antibody

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF@488A	490/515 nm	100 uL (100 ug)	20421	CF@488A Features
CF@543	541/560 nm	100 uL (100 ug)	20476	CF@543 Features
CF@568	562/583 nm	100 uL (100 ug)	20477	CF@568 Features
CF@594	593/614 nm	100 uL (100 ug)	20422	CF@594 Features
CF@640R	642/662 nm	100 uL (100 ug)	20423	CF@640R Features
CF@660R	663/682 nm	100 uL (100 ug)	20478	CF@660R Features
CF@680R	680/701 nm	100 uL (100 ug)	20479	CF@680R Features

View our full selection of [Secondary Antibodies](#), or search our catalog using our [Antibody Finder](#). Alternatively, you can view our [secondary antibody product listings](#) with catalog numbers.

CF® Dyes offer exceptional brightness and photostability. For more information see our [CF® Dye technology page](#).

Storage and Handling

Liquid format: 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.

Lyophilized format: Store at -20°C, protected from light. Product is stable for at least 6 months from date of receipt when stored as recommended. Reconstitute antibodies in water using the indicated volumes below:

CF® Dye and biotin conjugates: add 0.5 mL dH₂O

HRP or DNP conjugates: add 1 mL dH₂O

Add the indicated volume of water directly to the vial containing the lyophilized antibody and mix gently to dissolve. Store reconstituted antibody at -20°C and protect from light. Aliquot to avoid repeated freeze/thaw cycles. Alternatively, an equal volume of glycerol can be mixed with the reconstituted antibody so that it will remain liquid at -20°C.

Optional: A preservative such as 0.05% sodium azide (final concentration) can be added to CF® Dye and biotin conjugates. Do not add sodium azide to HRP conjugates.

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

CF is a registered trademark of Biotium, Inc.

References

Download a list of [CF® dye references](#).

This datasheet was generated on November 21, 2024 at 02:37:45 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/rabbit-anti-rfp/>