

## 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](#). 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 dH<sub>2</sub>O HRP or DNP conjugates: add 1 mL dH<sub>2</sub>O. 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.

### Product attributes

Clonality	Polyclonal
Antibody type	Tag Antibody
Concentration	1 mg/mL
Host species	Rabbit
Antibody reactivity (parent)	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

### References

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

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