FLAG Tag Polyclonal Rabbit Antibody

Polyclonal rabbit anti-FLAG antibody labeled with our superior CF® Dyes.



Product Description

This is a polyclonal rabbit anti-FLAG antibody labeled with our superior CF® Dyes. This antibody recognizes the FLAG epitope tag (DYKDDDDK) fused to either the amino- or carboxy- termini of targeted proteins.

- Available in 6 bright and photostable CF® Dyes
- Suitable for western, immunofluorescence, and immunohistology

See our full selection of anti-tag and anti-hapten antibody conjugates.

FLAG Tag Polyclonal Rabbit Antibody

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 uL (100 ug)	<u>20432</u>	CF®488A Features
CF®543	543/563 nm	100 uL (100 ug)	<u>20433</u>	CF®543 Features
CF®594	593/615 nm	100 uL (100 ug)	<u>20434</u>	CF®594 Features
<u>CF®640R</u>	642/663 nm	100 uL (100 ug)	<u>20435</u>	CF®640R Features
CF®680	681/698 nm	50 uL (50 ug)	<u>20973</u>	CF®680 Features
<u>CF®770</u>	770/797 nm	50 uL (50 ug)	<u>20974</u>	CF®770 Features

View our full selection of <u>Secondary Antibodies</u>, or search our catalog using our <u>Antibody Finder</u>. Alternatively, you can view our <u>secondary antibody product listings</u> with catalog numbers.

CF® Dyes offer exceptional brightness and photostability. For more information see our CF® Dye technology page.

Storage and Handling

<u>Liquid format</u>: 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 7, 2025 at 02:44:25 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rabbit-anti-flag/

Product attributes

Clonality	Polyclonal		
Antibody type	Tag Antibody		
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide		
Concentration	1 mg/mL		
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 (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.		
Host species	Rabbit		
Antibody reactivity (target)	FLAG tag		

Call us: 800-304-5357 Email: btinfo@biotium.com