

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)	20432	CF@488A Features
CF@543	543/563 nm	100 uL (100 ug)	20433	CF@543 Features
CF@594	593/615 nm	100 uL (100 ug)	20434	CF@594 Features
CF@640R	642/663 nm	100 uL (100 ug)	20435	CF@640R Features
CF@680	681/698 nm	50 uL (50 ug)	20973	CF@680 Features
CF@770	770/797 nm	50 uL (50 ug)	20974	CF@770 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 February 14, 2025 at 03:08:11 AM. 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 (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Host species	Rabbit
Antibody reactivity (target)	FLAG tag