V5 Tag Polyclonal Rabbit Antibody

Polyclonal rabbit anti-V5 antibody labeled with our superior CF® Dyes.



Product Description

This is a polyclonal rabbit anti-V5 antibody labeled with our superior CF® Dyes. The V5 epitope tag is derived from a small epitope (Pk) present on the P and V proteins of the paramyxovirus simian virus 5 (SV5). This antibody recognizes the V5 epitope tag (GKPIPNPLLGLDST) 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.

Product attributes

| Clonality | Polyclonal | | |
|-------------------------------------|---|--|--|
| Antibody type | Tag Antibody | | |
| Concentration | 1 mg/mL | | |
| Host species | Rabbit | | |
| Antibody reactivity (target) | V5 tag | | |
| Antibody/conjugate formulation | Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide | | |
| 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. | | |

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

V5 Tag Polyclonal Rabbit Antibody

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|----------------|------------|-----------------|--------------|------------------|
| CF®488A | 490/516 nm | 100 uL (100 ug) | 20440 | CF®488A Features |
| <u>CF®543</u> | 543/563 nm | 100 uL (100 ug) | <u>20441</u> | CF®543 Features |
| CF®594 | 593/615 nm | 100 uL (100 ug) | 20442 | CF®594 Features |
| <u>CF®640R</u> | 642/663 nm | 100 uL (100 ug) | <u>20443</u> | CF®640R Features |
| CF®680 | 681/698 nm | 50 uL (50 ug) | <u>20979</u> | CF®680 Features |
| CF®770 | 770/797 nm | 50 uL (50 ug) | <u>20980</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

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 curated CF® Dye references.

This datasheet was generated on November 3, 2025 at 04:49:17 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rabbit-anti-v5/