



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## DNP Polyclonal Rabbit Antibody

Polyclonal rabbit anti-DNP antibody labeled with our superior CF® dyes and HRP.



### Product Description

This is a polyclonal rabbit anti-DNP antibody labeled with our superior CF® dyes and HRP. This antibody recognizes the dinitrophenol (DNP) hapten for detection of DNP conjugated proteins and nucleic acids.

- Available in 7 bright and photostable CF® dyes and HRP
- Suitable for western, immunofluorescence, and immunohistology in FFPE tissues

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

### Rabbit Anti-DNP Tag

Conjugation	Ex/Em	Size	Catalog No.
CF®405S	404/431 nm	50 uL (100 ug)	<a href="#">20865-50uL</a>
250 uL (500 ug)	<a href="#">20865-250uL</a>		
CF®488A	490/515 nm	50 uL (100 ug)	<a href="#">20866-50uL</a>
250 uL (500 ug)	<a href="#">20866-250uL</a>		
CF®568	562/583 nm	50 uL (100 ug)	<a href="#">20867-50uL</a>
250 uL (500 ug)	<a href="#">20867-250uL</a>		
CF®594	593/614 nm	50 uL (100 ug)	<a href="#">20868-50uL</a>
250 uL (500 ug)	<a href="#">20868-250uL</a>		
CF®640R	642/662 nm	50 uL (100 ug)	<a href="#">20869-50uL</a>
250 uL (500 ug)	<a href="#">20869-250uL</a>		
CF®680R	680/701 nm	50 uL (100 ug)	<a href="#">20870-50uL</a>
250 uL (500 ug)	<a href="#">20870-250uL</a>		
CF®750	755/777 nm	50 uL (100 ug)	<a href="#">20875-50uL</a>
250 uL (500 ug)	<a href="#">20875-250uL</a>		
HRP	N/A	100 uL (100 ug)	<a href="#">20871-100uL</a>

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.

### References

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

This datasheet was generated on December 14, 2025 at 12:22:08 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-hrp-rabbit-anti-dnp-antibody/>

#### Product attributes

Clonality	Polyclonal
Antibody type	Tag Antibody
Concentration	1 mg/mL (HRP conjugate), 2 mg/mL
Host species	Rabbit
Antibody reactivity (target)	DNP
Antibody/conjugate formulation	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
Cross adsorption	Not cross-adsorbed
Secondary/tag antibody applications	ELISA, 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.