## HA (Hemagglutinin Tag) Polyclonal Rabbit Antibody

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

# Diotium ...

### **Product Description**

This is a polyclonal rabbit anti-HA (Hemagglutinin) Tag antibody labeled with our superior CF® Dyes. The antibody recognizes the hemagglutinin (HA) epitope tag (YPYDVPDYA).

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

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

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	50 uL (50 ug)	<u>20238</u>	CF®488A Features
<u>CF®594</u>	593/615 nm	50 uL (50 ug)	<u>20239</u>	CF®594 Features
CF®640R	642/663 nm	50 uL (50 ug)	<u>20240</u>	CF®640R Features
<u>CF®647</u>	652/668 nm	50 uL (50 ug)	<u>20486</u>	CF®647 Features
CF®680	681/698 nm	50 uL (50 ug)	<u>20975</u>	CF®680 Features
CF®770	770/797 nm	50 uL (50 ug)	<u>20976</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  $dH_2O$ 

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 November 4, 2025 at 02:15:57 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/rabbit-anti-ha-tag-antibody/">https://biotium.com/product/rabbit-anti-ha-tag-antibody/</a>

#### Product attributes

Call us: 800-304-5357

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

Email: btinfo@biotium.com