

# HA (Hemagglutinin Tag) Polyclonal Rabbit Antibody

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



## Product attributes

<b>Clonality</b>	Polyclonal
<b>Antibody type</b>	Tag Antibody
<b>Concentration</b>	1 mg/mL
<b>Host species</b>	Rabbit
<b>Antibody reactivity (target)</b>	HA Tag
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
<b>Secondary/tag antibody applications</b>	Flow cytometry, IHC, IF (cells or tissue sections), Western blot

## 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](#).

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](#).

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