

## Goat Anti-Alpaca IgG, VHH Antibody

Goat anti-alpaca IgG antibody that recognizes the VHH domain of heavy chain alpaca IgG. Labeled with CF® Dyes and suitable for western or IF applications.



### Product Description

Goat Anti-Alpaca IgG, VHH antibody is an affinity-purified polyclonal goat IgG whole antibody that recognizes the VHH domain of heavy chain alpaca IgG, subclasses 2 and 3, and the VHH domain of llama IgG, subclasses 2 and 3. The antibody is cross-adsorbed against bovine, human, mouse, rabbit, and rat serum proteins. It is not highly cross-adsorbed against immunoglobulins from other species. The antibody is available conjugated to Biotium's bright and photostable CF® Dyes.

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

### Product attributes

<b>Clonality</b>	Polyclonal
<b>Antibody type</b>	Secondary
<b>Host species</b>	Goat
<b>Antibody reactivity (target)</b>	Alpaca IgG, VHH domain, Llama IgG, VHH domain
<b>Species reactivity</b>	Alpaca, Llama
<b>Cross adsorption</b>	Bovine, Human, Mouse, Rabbit, Rat
<b>Antibody/conjugate formulation</b>	Liquid: PBS/50% glycerol/2 mg/mL BSA/0.05% azide
<b>Concentration</b>	1 mg/mL
<b>Secondary/tag antibody applications</b>	IHC, IF (cells or tissue sections), Western blot
<b>Product origin</b>	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.

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/516 nm	100 uL (100 ug)	<a href="#">20882-100uL</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	100 uL (100 ug)	<a href="#">20883-100uL</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/615 nm	100 uL (100 ug)	<a href="#">20884-100uL</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/663 nm	100 uL (100 ug)	<a href="#">20885-100uL</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	652/668 nm	100 uL (100 ug)	<a href="#">20886-100uL</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@680</a>	681/698 nm	100 uL (100 ug)	<a href="#">20887-100uL</a>	<a href="#">CF@680 Features</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.

### 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.

This product has been made available through a collaborative agreement between Jackson ImmunoResearch and Biotium.

CF is a registered trademark of Biotium, Inc.

### References

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

This datasheet was generated on November 21, 2024 at 03:07:36 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/goat-anti-alpaca-igg-vhh/>