

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

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 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.

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](#).

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

Product attributes

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