

Alpaca Anti-Human IgG (Fc), rVHH (N2508.HFC) - MiniMab™

VHH single-domain antibody targeting the heavy chain (Fc region) of human IgG, part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.



Product Description

Alpaca Anti-Human IgG (Fc), rVHH (N2508.HFC) is a secondary antibody that binds to the heavy chain (Fc region) of human IgG. This high-affinity single-domain antibody (SdAb), also known as camelid VHH or Nanobody®, is part of our MiniMab™ series of highly optimized conjugated probes. The SdAb has been validated for immunofluorescence microscopy and flow cytometry, and is available conjugated to CF® Dyes.

Features of MiniMab™ single-domain antibodies

- Superior to conventional antibodies: deeper tissue penetration, higher solubility and stability, and faster staining
- Minimal epitope-dye displacement—perfect for super-resolution imaging
- Specifically developed and optimized for immunofluorescence
- Labeled with bright, photostable CF® Dyes, including near-infrared CF®740 option
- Available as conjugates with Biotium's best-in-class dyes for STORM

Learn more about [CF® Dyes for super-resolution](#) as well as our innovative reagents for [immunofluorescence microscopy](#); this includes our [NucSpot® Nuclear Stains](#) for bright and nuclear-specific staining in a wide color selection, and [CytoLiner™ Fixed Cell Membrane Stains](#) for robust membrane staining in formaldehyde-fixed cells.

[View our full selection of primary and secondary antibodies](#) available with bright CF® Dyes and other labels.

Alpaca Anti-Human IgG (Fc), rVHH (N2508.HFC) - MiniMab™

Conjugation	Ex/Em	Conc.	STORM ¹ compatibility	Catalog No.	Dye Features
CF®405S	404/431 nm	100 ug/mL	No	N007-405S-500UL	CF®405S Features
CF®488A	490/516 nm	100 ug/mL	Yes	N007-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	Yes	N007-568-500UL	CF®568 Features
CF®583R	585/609 nm	100 ug/mL	Yes	N007-583R-500UL	CF®583R Features
CF®594	593/615 nm	100 ug/mL	No	N007-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	No	N007-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	Yes	N007-647-500UL	CF®647 Features
CF®660R	662/682 nm	100 ug/mL	No	N007-660R-500UL	CF®660R Features
CF®680	681/698 nm	100 ug/mL	Yes	N007-680-500UL	CF®680 Features
CF®680R	680/701 nm	100 ug/mL	Yes	N007-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	No	N007-740-500UL	CF®740 Features
CF®770	770/797 nm	100 ug/mL	No	N007-770-500UL	CF®770 Features

¹ STORM: Stochastic optical reconstruction microscopy. [Learn more about CF® Dyes for super-resolution](#). NANOBODY is a registered trademark of ABLYNX.

This datasheet was generated on December 12, 2025 at 07:31:46 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/alpaca-anti-human-igg-fc-rvhh-n2508-hfc-minimab/>

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

Product attributes

Antibody number	N007
Antibody reactivity (target)	Human IgG, Fc
Antibody type	MiniMab™ SdAb (VHH)
Host species	Alpaca
Clonality	Recombinant single-domain antibody
Clone	N2508.HFC
Isotype	VHH
Synonyms	IgG1 Fc Fragment, immunoglobulin G1 fragment; immunoglobulin G1 Fc region
Verified antibody applications	Flow (surface) (verified), IF (verified)
Species reactivity	Human
Antibody application notes	Immunofluorescence: 0.1 ug/mL; Flow cytometry: 0.1 ug per million cells/0.1 mL; Optimal concentration to be determined by end-user.
Shipping condition	Room temperature
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/0.1% BSA/0.05% azide
Product origin	Recombinant alpaca VHH produced in E.coli, Recombinant BSA produced in Chinese hamster ovary cells
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot, STORM