

## Alpaca Anti-Mouse IgG1 (Fc), rVHH (N2504.M1FC) - MiniMab™

VHH single-domain antibody targeting the heavy chain (Fc region) of mouse IgG1, part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.



### Product Description

Alpaca Anti-Mouse IgG1 (Fc), rVHH (N2504.M1FC) is a secondary antibody that binds to the heavy chain (Fc region) of mouse IgG1. 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, flow cytometry, and western blot, 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.

### Product attributes

Antibody number	N004
Antibody type	MiniMab™ SdAb (VHH)
Clonality	Recombinant single-domain antibody
Host species	Alpaca
Clone	N2504.M1FC
Isotype	VHH
Antibody reactivity (target)	Mouse IgG1, Fc
Synonyms	IgG1 Fc Fragment, immunoglobulin G1 fragment; immunoglobulin G1 Fc region
Species reactivity	Human (weak), Mouse, Rat
Verified antibody applications	Flow (surface) (verified), IF (verified), Western Blot (verified)
Antibody application notes	Immunofluorescence: 0.1 ug/mL; Flow cytometry: 0.1 ug per million cells/0.1 mL; Western blot: 250 ng/mL; Optimal concentration to be determined by end-user.
Antibody/conjugate formulation	Conjugates: PBS/0.1% rBSA/0.05% azide
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot, STORM
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates from light
Regulatory status	For research use only (RUO)
Product origin	Recombinant alpaca VHH produced in E.coli, Recombinant BSA produced in Chinese hamster ovary cells
Shipping condition	Room temperature

## Alpaca Anti-Mouse IgG1 (Fc), rVHH (N2504.M1FC) – MiniMab™

Conjugation	Ex/Em	Conc.	STORM <sup>1</sup> compatibility	Catalog No.	Dye Features
<a href="#">CF@405S</a>	404/431 nm	100 ug/mL	No	<a href="#">N004-405S-500UL</a>	<a href="#">CF@405S Features</a>
<a href="#">CF@488A</a>	490/516 nm	100 ug/mL	Yes	<a href="#">N004-488A-500UL</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	100 ug/mL	Yes	<a href="#">N004-568-500UL</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@583R</a>	585/609 nm	100 ug/mL	Yes	<a href="#">N004-583R-500UL</a>	<a href="#">CF@583R Features</a>
<a href="#">CF@594</a>	593/615 nm	100 ug/mL	No	<a href="#">N004-594-500UL</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/663 nm	100 ug/mL	No	<a href="#">N004-640R-500UL</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	652/668 nm	100 ug/mL	Yes	<a href="#">N004-647-500UL</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@660R</a>	662/682 nm	100 ug/mL	No	<a href="#">N004-660R-500UL</a>	<a href="#">CF@660R Features</a>
<a href="#">CF@680</a>	681/698 nm	100 ug/mL	Yes	<a href="#">N004-680-500UL</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@680R</a>	680/701 nm	100 ug/mL	Yes	<a href="#">N004-680R-500UL</a>	<a href="#">CF@680R Features</a>
<a href="#">CF@740</a>	742/767 nm	100 ug/mL	No	<a href="#">N004-740-500UL</a>	<a href="#">CF@740 Features</a>
<a href="#">CF@770</a>	770/797 nm	100 ug/mL	No	<a href="#">N004-770-500UL</a>	<a href="#">CF@770 Features</a>

<sup>1</sup> 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 August 20, 2025 at 02:02:08 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/alpaca-anti-mouse-igg1-fc-rvhh-n2504-m1fc-minimab/>