



Glowing products for science

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 [MiniMab™ single domain antibodies](#) paired with Biotium's industry-leading [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-Mouse IgG1 (Fc), rVHH (N2504.M1FC) – MiniMab™

Conjugation	Ex/Em	Conc.	STORM ¹ compatibility	Catalog No.	Dye Features
CF®405S	404/431 nm	100 ug/mL	No	N004-405S-500UL	CF®405S Features
CF®488A	490/516 nm	100 ug/mL	Yes	N004-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	Yes	N004-568-500UL	CF®568 Features
CF®583R	585/609 nm	100 ug/mL	Yes	N004-583R-500UL	CF®583R Features
CF®594	593/615 nm	100 ug/mL	No	N004-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	No	N004-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	Yes	N004-647-500UL	CF®647 Features
CF®660R	662/682 nm	100 ug/mL	No	N004-660R-500UL	CF®660R Features
CF®680	681/698 nm	100 ug/mL	Yes	N004-680-500UL	CF®680 Features
CF®680R	680/701 nm	100 ug/mL	Yes	N004-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	No	N004-740-500UL	CF®740 Features
CF®770	770/797 nm	100 ug/mL	No	N004-770-500UL	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 January 29, 2026 at 12:07:17 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/>

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

Product attributes

Antibody number	N004
Antibody type	MiniMab™ SdAb (VHH), Secondary
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), WB (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