

# Alpaca Anti-Rabbit IgG (Fc), rVHH (N2507.RFC) - MiniMab™

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

# Bietum

## **Product Description**

Alpaca Anti-Rabbit IgG (Fc), rVHH (N2507.RFC) is a secondary antibody that binds to the heavy chain (Fc region) of rabbit 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, 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 <u>CF® Dyes for super-resolution</u> as well as our innovative reagents for <u>immunofluorescence microscopy</u>; this includes our <u>NucSpot® Nuclear Stains</u> for bright and nuclear-specific staining in a wide color selection, and <u>CytoLiner™ Fixed Cell Membrane Stains</u> for robust membrane staining in formaldehyde-fixed cells.

<u>View our full selection of primary and secondary antibodies</u> available with bright CF® Dyes and other labels.

#### Product attributes

Call us: 800-304-5357

Antibody number	N006			
Antibody reactivity (target)	Rabbit IgG, Fc			
Antibody type	MiniMab™ SdAb (VHH)			
Host species	Alpaca			
Clonality	Recombinant single-domain antibody			
Clone	N2507.RFC			
Isotype	VHH			
Synonyms	Rabbit IgG FC Fragment; fragment crystallizable region; Fc region			
Verified antibody applications	Flow (surface) (verified), IF (verified), Western Blot (verified)			
Species reactivity	Rabbit IgG, Fc			
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.			
Shipping condition	Room temperature			
Storage Conditions	Store conjugates at 2 $^{\circ}\text{C}$ to 8 $^{\circ}\text{C},$ Protect fluorescent conjugate 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% rBSA/0.05% azide			
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot, STORM			
Product origin	Recombinant alpaca VHH produced in E.coli, Recombinant BSA produced in Chinese hamster ovary cells			

## Alpaca Anti-Rabbit IgG (Fc), rVHH (N2507.RFC) - MiniMab™

Conjugation	Ex/Em	Conc.	STORM¹ compatibility	Catalog No.	Dye Features
<u>CF®405S</u>	404/431 nm	100 ug/mL	No	N006-405S-500UL	CF®405S Features
CF®488A	490/516 nm	100 ug/mL	Yes	N006-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	Yes	N006-568-500UL	CF®568 Features
CF®583R	585/609 nm	100 ug/mL	Yes	N006-583R-500UL	CF®583R Features
CF®594	593/615 nm	100 ug/mL	No	N006-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	No	N006-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	Yes	N006-647-500UL	CF®647 Features
CF®660R	662/682 nm	100 ug/mL	No	N006-660R-500UL	CF®660R Features
CF®680	681/698 nm	100 ug/mL	Yes	N006-680-500UL	CF®680 Features
CF®680R	680/701 nm	100 ug/mL	Yes	N006-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	No	N006-740-500UL	CF®740 Features
CF®770	770/797 nm	100 ug/mL	No	N006-770-500UL	CF®770 Features

<sup>1</sup> STORM: Stochastical optical reconstruction microscopy. Learn more about CF® Dyes for super-resolution. NANOBODY is a registered trademark of ABLYNX.

This datasheet was generated on November 25, 2025 at 03:32:53 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/alpaca-anti-rabbit-igg-fc-rvhh-n2507-rfc-minimab/">https://biotium.com/product/alpaca-anti-rabbit-igg-fc-rvhh-n2507-rfc-minimab/</a>