Alpaca Anti-Mouse (Fab) Kappa, rVHH (N2507.MFAB) - MiniMab™

VHH single-domain antibody that binds to the mouse kappa light chain, part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.

Biotum

Product Description

Alpaca Anti-Mouse (Fab) Kappa, rVHH (N2507.MFAB) is a single domain antibody (SdAb, VHH, also known as Nanobody®) that recognizes mouse antigen-binding fragment (Fab) kappa light chain. This SdAb is part of our MiniMab™ series of highly optimized conjugated probes. It has been validated for immunofluorescence microscopy, flow cytometry, and western blot, and is available conjugated to CF® Dves.

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
- 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>CytoLinerTM Fixed Cell Membrane Stains</u> 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

Call us: 800-304-5357

Fiduuci attributes				
Antibody number	N005			
Antibody type	MiniMab™ SdAb (VHH)			
Clonality	Recombinant single-domain antibody			
Host species	Alpaca			
Clone	N2507.MFAB			
Isotype	VHH			
Antibody reactivity (target)	Mouse Fab (Kappa)			
Synonyms	Mouse immunoglobulin kappa light chain			
Species reactivity	Mouse			
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, 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 $^{\circ}\text{C}$ to 8 $^{\circ}\text{C},$ Protect fluorescent conjugates from light			
Regulatory status	For research use only (RUO)			
Product origin	Animal, Recombinant alpaca VHH produced in E.coli, Recombinant BSA produced in Chinese hamster ovary cells			
Shipping condition	Room temperature			

Email: btinfo@biotium.com

Alpaca Anti-Mouse (Fab) Kappa, rVHH (N2507.MFAB) – MiniMab™

Conjugation	Ex/Em	Conc.	STORM¹ compatibility	Catalog No.	Dye Features
<u>CF®488A</u>	490/516 nm	100 ug/mL	Yes	N005-488A-500UL	CF®488A Features
CF®498	498/519 nm	100 ug/mL	Yes	N005-498-500UL	
CF®568	562/584 nm	100 ug/mL	Yes	N005-568-500UL	CF®568 Features
CF®583R	585/609 nm	100 ug/mL	Yes	N005-583R-500UL	CF®583R Features
CF®647	652/668 nm	100 ug/mL	Yes	N005-647-500UL	CF®647 Features
CF®660C	667/685 nm	100 ug/mL	Yes	N005-660C-500UL	CF®660C Features
CF®680	681/698 nm	100 ug/mL	Yes	N005-680-500UL	CF®680 Features

¹ 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 6, 2025 at 06:36:02 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/alpaca-anti-mouse-fab-kappa-rvhh-n2507-mfab-minimab/