

Alpaca Anti-Mouse IgG2a/b, rVHH (N2603.M2AB)- MiniMab™



VHH single-domain antibody that binds to mouse-IgG2a and IgG2b, part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.

Product Description

Alpaca Anti-Mouse IgG2a/b, rVHH (N2603.M2AB) is a secondary antibody that binds to mouse-IgG2a and IgG2b subtypes with high affinity. 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.

Product attributes

| | |
|-------------------------------------|--|
| Antibody number | N010 |
| Antibody type | MiniMab™ SdAb (VHH), Secondary |
| Clonality | Recombinant single-domain antibody |
| Host species | Alpaca |
| Clone | N2603.M2AB |
| Isotype | VHH |
| Antibody reactivity (target) | Mouse IgG2a, Mouse IgG2b |
| Synonyms | Mouse IgG2; Mouse IgG, Fc; Fc region specific; alpaca monoclonal Nanobody |
| Species reactivity | Mouse |
| Verified antibody applications | IF (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 IgG2a/b, rVHH (N2603.M2AB)- MiniMab™

| Conjugation | Ex/Em | Conc. | STORM ¹ compatibility | Catalog No. | Dye Features |
|-------------------------|------------|-----------|----------------------------------|---------------------------------|----------------------------------|
| CF@405S | 404/431 nm | 100 ug/mL | No | N010-405S-500UL | CF@405S Features |
| CF@488A | 490/516 nm | 100 ug/mL | Yes | N010-488A-500UL | CF@488A Features |
| CF@568 | 562/584 nm | 100 ug/mL | Yes | N010-568-500UL | CF@568 Features |
| CF@583R | 585/609 nm | 100 ug/mL | Yes | N010-583R-500UL | CF@583R Features |
| CF@594 | 593/615 nm | 100 ug/mL | No | N010-594-500UL | CF@594 Features |
| CF@640R | 642/663 nm | 100 ug/mL | No | N010-640R-500UL | CF@640R Features |
| CF@647 | 652/668 nm | 100 ug/mL | Yes | N010-647-500UL | CF@647 Features |
| CF@660R | 662/682 nm | 100 ug/mL | No | N010-660R-500UL | CF@660R Features |
| CF@680 | 681/698 nm | 100 ug/mL | Yes | N010-680-500UL | CF@680 Features |
| CF@680R | 680/701 nm | 100 ug/mL | Yes | N010-680R-500UL | CF@680R Features |
| CF@740 | 742/767 nm | 100 ug/mL | No | N010-740-500UL | CF@740 Features |
| CF@770 | 770/797 nm | 100 ug/mL | No | N010-770-500UL | CF@770 Features |

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