

Histone Recombinant Alpaca VHH (SdAb2511.HISTONE), MiniMab™



VHH nanobody that binds to histone H2A-H2B heterodimers, part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.

Product Description

Histone Recombinant Alpaca VHH (SdAb2511.HISTONE) binds to histone H2A-H2B heterodimers and has broad species reactivity. 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 nanobody has been validated for immunofluorescence microscopy 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

Histone Recombinant Alpaca VHH (SdAb2511.HISTONE) labels chromatin throughout the cell cycle by specifically binding histone H2A-H2B heterodimers, enabling visualization of interphase chromatin, chromatin condensation, and mitotic chromosomes.

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.

Histone Recombinant Alpaca VHH (SdAb2511.HISTONE), MiniMab™

Conjugation	Ex/Em	Conc.	STORM ¹ compatibility	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	Yes	N008-488A-1ML	CF®488A Features
CF®568	562/584 nm	100 ug/mL	Yes	N008-568-1ML	CF®568 Features
CF®583R	585/609 nm	100 ug/mL	Yes	N008-583R-1ML	CF®583R Features
CF®594	593/615 nm	100 ug/mL	Yes	N008-594-1ML	CF®594 Features
CF®647	652/668 nm	100 ug/mL	Yes	N008-647-1ML	CF®647 Features
CF®660C	667/685 nm	100 ug/mL	Yes	N008-660C-1ML	CF®660C Features
CF®680	681/698 nm	100 ug/mL	Yes	N008-680-1ML	CF®680 Features
CF®740	742/767 nm	100 ug/mL	No	N008-740-1ML	CF®740 Features

NANOBODY is a registered trademark of ABLYNX.

This datasheet was generated on January 29, 2026 at 11:24:59 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/histone-recombinant-alpaca-vhh-sdab2511-histone-minimab/>

Product attributes

Antibody number	N008
Antibody type	Biotium Choice Primary, MiniMab™ SdAb (VHH), Primary
Clonality	Recombinant single-domain antibody
Host species	Alpaca
Clone	SdAb2511.HISTONE
Isotype	VHH
Antibody reactivity (target)	Histone H2A/H2B
Synonyms	H2A histone family member, Histone H2A-H2B heterodimers, Histone-Binding-Protein, Histone-Nanobody, and Histone-single domain antibody (sdAb)
Species reactivity	Broad reactivity
Verified antibody applications	IF (verified)
Positive control	HeLa cells.
Antibody application notes	Immunofluorescence: 0.1 ug/mL; Optimal concentration to be determined by end-user.
Antibody research areas	Cancer
Antibody/conjugate formulation	Conjugates: PBS/0.1% rBSA/0.05% azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Shipping condition	Room temperature
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