



Glowing products for science

PSD95 Recombinant Alpaca VHH (SdAb2601.PSD95) - MiniMab™

Recombinant Alpaca VHH (SdAb2601.PSD95) targeting PSD95. Part of our MiniMab™ SdAb series that have been engineered for optimal conjugate performance.



Product Description

PSD95 Recombinant Alpaca VHH (SdAb2601.PSD95) recognizes PSD95 (also known as PSD-95, SAP-90, or DLG4), a MAGUK family protein critical for building a dynamic scaffold structure at postsynaptic sites. 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 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

PSD95 functions as an organizing centrum of receptors, ion channels, kinases, and other relevant players in achieving synaptic plasticity and molecular changes during long-term potentiation. This recombinant alpaca VHH is cross-reactive with human, mouse, and rat PSD-95

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.

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

Product attributes

Antibody number	N009
Antibody reactivity (target)	PSD95
Antibody type	MiniMab™ SdAb (VHH), Primary
Host species	Alpaca
Clonality	Recombinant single-domain antibody
Clone	SdAb2601.PSD95
Isotype	VHH
Molecular weight	95-110 kDa
Synonyms	discs large homolog 4; Disks large homolog 4; FLJ97752; FLJ98574; post synaptic density; post-synaptic density protein 95; Postsynaptic density protein 95; PSD-95; PSD-95 alpha 2b; PSD-95 beta; SAP-90; Synapse-associated protein 90; synapse-associated protein SAP90; Tax interaction protein 15
Human gene symbol	DLG4
Entrez gene ID	13385
Immunogen	Recombinant Mouse PSD95 protein (PDZ domains 1 and 2)
Antibody target cellular localization	Cytoplasm
Verified antibody applications	IF (verified)
Species reactivity	Human, Mouse, Rat
Antibody application notes	Immunofluorescence: 0.1 ug/mL; Optimal concentration to be determined by end-user.
Positive control	Brain, Retina
Shipping condition	Room temperature
Storage Conditions	Store conjugates at 2 °C to 8 °C, Protect fluorescent conjugates 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
Product origin	Recombinant alpaca VHH produced in E.coli, Recombinant BSA produced in Chinese hamster ovary cells
Antibody research areas	Neuroscience
Cell/tissue expression	Brain, Neurons

PSD95 Recombinant Alpaca VHH (SdAb2601.PSD95) – MiniMab™

Conjugation	Ex/Em	Conc.	STORM ¹ compatibility	Catalog No.	Dye Features
CF@488A	490/516 nm	100 ug/mL	Yes	N009-488A-200UL	CF@488A Features
CF@498	498/519 nm	100 ug/mL	Yes	N009-498-200UL	
CF@568	562/584 nm	100 ug/mL	Yes	N009-568-200UL	CF@568 Features
CF@583R	585/609 nm	100 ug/mL	Yes	N009-583R-200UL	CF@583R Features
CF@647	652/668 nm	100 ug/mL	Yes	N009-647-200UL	CF@647 Features
CF@660C	667/685 nm	100 ug/mL	Yes	N009-660C-200UL	CF@660C Features
CF@680	681/698 nm	100 ug/mL	Yes	N009-680-200UL	CF@680 Features
CF@740	742/767 nm	100 ug/mL	No	N009-740-200UL	CF@740 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 September 18, 2026 at 05:46:23 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/psd95-recombinant-alpaca-vhh-sdab2601-psd95-minimab/>