



Glowing products for science

PSD95 Recombinant Monoclonal Mouse Antibody (rK28/43) – Biotium Choice

A recombinant mouse monoclonal antibody that recognizes PSD95. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.

Product Description

PSD95 Recombinant Monoclonal Mouse Antibody (rK28/43) is a recombinant mouse monoclonal antibody that recognizes PSD95. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for fluorescence microscopy. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 7 bright and photostable CF® Dyes, including [near-IR CF®740](#)
- Suitable for immunohistochemistry, immunofluorescence, and western

PSD-95, encoded by DLG4, is a MAGUK family scaffold protein (with PSD-93, SAP97, SAP102) that clusters membrane receptors and ion channels such as NMDA receptor subunits, Kv1 channels, and Caspr2. Containing PDZ, SH3, and GuK domains, PSD-95 localizes to postsynaptic densities, where it is essential for synapse formation and maintenance. PSD-95/DLG4 antibodies are commonly paired with presynaptic markers to visualize synapses by microscopy.

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable [CF® Dye](#) for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)



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

Product attributes

Antibody Number	P036
Antibody reactivity (target)	PSD95
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary
Host species	Mouse
Clonality	Recombinant Monoclonal
Clone	rK28/43
Isotype	IgG2a
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	1742
SwissProt	P78352
Immunogen	Human recombinant PSD95 fusion protein (amino acids 77-299 – PDZ domains 1 and 2)
Antibody target cellular localization	Cytoplasm
Verified antibody applications	IF (verified), IHC (frozen) (verified)
Species reactivity	Chicken, Human, Mouse, Non-human primate, Rat
Expected antibody applications	IHC (published for clone), WB (published for clone)
Antibody application notes	Immunofluorescence: 1-5 ug/mL; Optimal concentration to be determined by end-user.
Positive control	Brain
Shipping condition	Room temperature
Storage Conditions	Store at 2 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/2 mg/mL rBSA/0.05% azide
Antibody research areas	Neuroscience
Cell/tissue expression	Brain, Neurons
Filter by antibody application	IHC, FFPE (published), WB (published for clone), IF (verified), IHC, frozen (verified)
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant mouse IgG produced in Chinese Hamster Ovary (CHO) cell line

PSD95 Recombinant Monoclonal Mouse Antibody (rK28/43) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	500 uL	P036-488A-500UL	CF®488A Features
CF®568	562/584 nm	100 ug/mL	500 uL	P036-568-500UL	CF®568 Features
CF®594	593/615 nm	100 ug/mL	500 uL	P036-594-500UL	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	500 uL	P036-640R-500UL	CF®640R Features
CF®647	652/668 nm	100 ug/mL	500 uL	P036-647-500UL	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	500 uL	P036-680R-500UL	CF®680R Features
CF®740	742/767 nm	100 ug/mL	500 uL	P036-740-500UL	CF®740 Features

This datasheet was generated on December 24, 2025 at 05:13:11 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/psd95-recombinant-monoclonal-mouse-antibody-rk28-43-biotium-choice/>