

Rim3 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse Rim3 (NT)

Instruction Manual

Catalog Number	PK-AB718-4471
Synonyms	Rim3 Antibody: Rab3-interacting molecule 3, regulating synaptic membrane exocytosis 3, RIM3 gamma, RIMS3, NIM3
Description	Rab3-interacting molecules (RIMs) are synaptic proteins necessary for neural transmission and plasticity. While both Rim1 and Rim 2 are thought to be effector proteins for Rab3, binding to Rab3 on synaptic vesicles in a GTP-dependent manner, less is known of Rim3. Expression of Rim3 in PC12 cells induced a significant increase in calcium-triggered exocytosis, with no appreciable change in the baseline release, suggesting that it plays a role in the regulation of exocytosis. Rim3 protein localizes primarily to neuronal dendrites and the postsynaptic densities, as opposed to Rim1 which is found in presynapse locations, indicating that Rim3 may contribute to synapse transmission and plasticity. This antibody is predicted to have no cross-reactivity to other Rim proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rim3 antibody was raised in rabbits against a 17 amino acid peptide from near the amino terminus of human Rim3.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
Specificity	This antibody is predicted to have no cross-reactivity to other Rim proteins.
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Rim3 antibody can be used for detection of Rim3 by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 5 µg/mL.
Images	Available upon request.
References	Wang Y, Sugita S, and Sudhof TC. The RIM/NIM family of neuronal C2 domain proteins: interactions with Rab3 and a new class of Src homology 3 domain proteins. <i>J. Biol. Chem.</i> 2000; 275:20033-44. Liang F, Zhang B, Tang J, et al. RIM3gamma is a postsynaptic protein in the rat central nervous system. <i>J. Comp. Neurol.</i> 2007; 503:501-10.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4471P; Rim3 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.