

# SAPAP2 antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat SAPAP2 (SAP90/PSD-95-associated protein 2)

### Instruction Manual

Catalog Number	PK-AB718-4625
Synonyms	SAPAP2 Antibody: SAP90, PSD-95-associated protein 2, Disks large-associated protein 2, DLGAP2, DAP-2
Description	SAP90/PSD-95-associated protein 2 (SAPAP2, also known as DLGAP2) is a member of a protein family whose members specifically interact with PSD-95/SAP90, a membrane-associated guanylate kinase localized at postsynaptic density (PSD) in neuronal cells. Like the other SAPAP proteins, SAPAP2 is thought to be an adaptor protein that also interacts with different synaptic scaffolding proteins, cytoskeletal and signaling components, such as focal adhesion kinase (FAK) and proline-rich tyrosine kinase 2 (PYK2). SAPAP2 mRNA is targeted to cell bodies in a similar manner to SAPAP1 and -4, whereas SAPAP3 mRNA is detected mainly in cell bodies. At least three isoforms of SAPAP2 are known to exist. This SAPAP2 antibody will not cross-react with other SAPAP proteins.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	SAPAP2 antibody was raised in rabbits against a 16 amino acid peptide from near the center of human SAPAP2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	At least three isoforms of SAPAP2 are known to exist. This SAPAP2 antibody will not cross-react with other SAPAP 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.
	SAPAP2 antibody can be used for detection of SAPAP2 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	SAPAPs. A family of PSD-95/SAP90-associated proteins localized at postsynaptic density. <i>J. Biol. Chem.</i> 1997; 272:11943-51. Ranta S, Zhang Y, Ross B, et al. Positional cloning and characterization of the human DLGAP2 gene and its exclusion in progressive epilepsy with mental retardation. <i>Eur. J. Hum. Genet.</i> 2000; 8:381-4. Kindler S, Rehbein M, Classen B, et al. Distinct spatiotemporal expression of SAPAP transcripts in the developing rat brain: a novel dendritically localized mRNA. <i>Brain Res. Mol. Brain Res.</i> 2004; 126:14-21. Bongiorno-Borbone L, Kadare G, Benfenati F, et al. FAK and PYK2 interact with SAP/PSD-95-associated protein-3. <i>Biochem. Biophys. Res. Commun.</i> 2005; 337:641-6.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4625P; SAPAP2 Peptide Cat. No. PK-AB718-1284; L1210 Cell Lysate

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