

Plxdc2 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Plxdc2 (NT)

Instruction Manual

Catalog Number	PK-AB718-4417
Synonyms	Plxdc2 Antibody: Plexin domain-containing protein 2, Tumor endothelial marker 7-related protein, TEM7R
Description	Plxdc2, also known as Tumor endothelial marker 7-related (TEM7R) encodes a protein with 57% amino acid identity to TEM7, the most abundant tumor endothelial marker. Plxdc2 is strongly expressed in the endothelial cells of the tumor stroma, but not in the endothelial cells of normal colonic tissue. Plxdc2 is also expressed at high levels in vessels of some normal tissues, with highest expression in muscle and lung. Plxdc2 and TEM7 may be important for tumor angiogenesis in humans. Cortactin was identified as a protein capable of binding to the extracellular region of both TEM7 and Plxdc2, and may provide new opportunities for the delivery of therapeutic and imaging agents to the vessels of solid tumors.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Plxdc2 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human Plxdc2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
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. Plxdc2 antibody can be used for detection of Plxdc2 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Carson-Weber EB, Watkins DN, Nanda A, et al. Cell surface tumor epithelial markers are conserved in mice and humans. <i>Cancer Res.</i> 2001; 61:6649-55. Nabda A and St. Croix Bl. Tumor endothelial markers: new targets for cancer therapy. <i>Curr Opin Oncol.</i> 2004; 16:44-9. St. Croix B, Rago C, Velculescu V, et al. Genes expressed in human tumor endothelium. <i>Science</i> 2000; 289:1197-202. Nanda A, Buckhaults P, Seaman S, et al. Identification of a binding partner for the endothelial cell surface proteins TEM7 and TEM7R. <i>Cancer Res.</i> 2004; 64:8507-11.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4417P; Plxdc2 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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