

TEM7 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat TEM7 (IN)

Instruction Manual

Catalog Number	PK-AB718-4341
Synonyms	TEM7 Antibody; Tumor endothelial marker 7, TEM3, plexin domain-containing protein 1, PLXDC1
Description	Tumor endothelial marker 7 (TEM7) was identified through serial analysis of gene expression on endothelial cells isolated from human normal and malignant colorectal tissues. Further experiments verified TEM7 was highly expressed in the endothelium of numerous other cancer types including breast, lung and brain tumors. At least four isoforms of TEM7 are known to exist; these include intracellular, secreted, and membrane-bound forms. A homologous protein, TEM7R (also known as PLXDC2), acts as a binding partner to TEM7 and is also abundantly expressed in the endothelium of malignant colorectal cancer but is absent or rare in normal colon mucosa. High expression of TEM7 is associated with metastasis and poor survival of patients with osteogenic sarcoma.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TEM7 antibody was raised in rabbits against a 16 amino acid peptide from near the center of human TEM7.
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. TEM7 antibody can be used for detection of TEM7 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	St. Croix B, Rago C, Velculescu V, et al. Genes expressed in human tumor endothelium. Science 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. Cancer Res. 2004; 64:8507-11. Carson-Weber EB, Watkins DN, Nanda A, et al. Cell surface tumor epithelial markers are conserved in mice and humans. Cancer Res. 2001; 61:6649-55. Fuchs B, Mahlum E, Halder C, et al. High expression of tumor endothelial marker 7 is associated with metastasis and poor survival of patients with osteogenic sarcoma. Gene 2007; 399:137-43.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4341P; TEM7 Peptide Cat.No. PK-AB718-1464; Rat Liver Tissue Lysate

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