

TEM5 (IN) antibody (pAb)

Rabbit Anti-Human/Rat TEM5 (IN)

Instruction Manual

Catalog Number	PK-AB718-4371
Synonyms	TEM5 Antibody; Tumor endothelial marker 5, G protein-coupled receptor 124, GPR124
Description	Tumor endothelial markers (TEMs) are significantly up-regulated during angiogenesis and neoangiogenesis that are crucial for the growth of solid tumors. TEMs localized on the cell surface and conserved across species are of particular interest for future development of anti-angiogenic therapies. These include TEMs such as TEM1, TEM5, TEM7 and TEM8. TEM5 is a member of the adhesion family of G protein coupled receptors and is localized on the surface of endothelial cells. TEM5 is a seven-pass transmembrane receptor, unlike TEM1, TEM7 and TEM8 which span the membrane once. TEM5 is abundantly expressed in tumor vessels, heart, placenta, ovary, small intestine, and colon. Proteolytically processed soluble TEM5 mediates endothelial cell survival during angiogenesis by linking integrin to glycosaminoglycans.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TEM5 antibody was raised in rabbits against a 14 amino acid peptide near the center of human TEM5.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, 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. TEM5 antibody can be used for detection of TEM5 by Western blot at 2 and 4 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	Carson-Walter EB, Watkins DN, Nanda A, et al. Cell surface tumor endothelial markers are conserved in mice and humans. <i>Cancer Res.</i> 2001; 61:6649-55. Yamamoto Y, Irie K, Nanda A, et al. Direct binding of the human homologue of the Drosophila disc large tumor suppressor gene to seven-pass transmembrane proteins, tumor endothelial marker 5 (TEM5), and a novel TEM5-like protein. <i>Oncogene</i> 2004; 23:3889-97. Davies G, Cunnick GH, Mansel RE, et al. Levels of expression of endothelial markers specific to tumour-associated endothelial cells and their correlation with prognosis in patients with breast cancer. <i>Clinical & Experimental Metastasis</i> 2004; 21:31-7. Vallon M and Essler M. Proteolytically processed soluble tumor endothelial marker (TEM) 5 mediates endothelial cell survival during angiogenesis by linking integrin alpha(v)beta3 to glycosaminoglycans. <i>J. Biol. Chem.</i> 2006; 281:34179-88.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4371P; TEM5 Peptide Cat.No. PK-AB718-1465; Rat Kidney Tissue Lysate

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