

# TEM4 (IN) antibody (pAb)

## Rabbit Anti-Human TEM4 (IN)

### Instruction Manual

Catalog Number	PK-AB718-4367
Synonyms	TEM4 Antibody; Tumor endothelial marker 4, ARHGEF17, Rho guanine nucleotide exchange factor 17, p164RHOGEF
Description	Rho GTPases, which are activated by specific guanine-nucleotide exchange factors (GEFs), play pivotal roles in several cellular functions. TEM4, encoding a protein containing 1510 amino acids, contains a RhoGEF-specific Dbl homology (DH) domain but lacks their typical pleckstrin homology domain. TEM4 is a Rho-specific GEF with novel structural and regulatory properties and predominant expression in the heart. It couples tyrosine kinase signals with the activation of the rho/rac GTPases, thus leading to cell differentiation and/or proliferation. Elevated levels of TEM4, TEM5, TEM6, TEM7 and TEM7R were also raised in breast cancer tissues. TEM4 could also prove to be useful targets therapeutically.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TEM4 antibody was raised in rabbits against an 18 amino acid peptide near the center of human TEM4.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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. TEM4 antibody can be used for detection of TEM4 by Western blot at 1 - 2 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	Ruemenapp U, Freichel-Blomquist A, Wittinghofer B, et al. A mammalian Rho-specific guanine-nucleotide exchange factor (p164-RhoGEF) without a pleckstrin homology domain. <i>Biochem. J.</i> 2002; 366:721-8. 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. 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 &amp; Experimental Metastasis</i> 2004; 21:31-7. Nanda A and St Croix B. Tumor endothelial markers: new targets for cancer therapy. <i>Curr. Opin. Oncol.</i> 2004; 16:44-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4367P; TEM4 Peptide Cat.No. PK-AB718-1205; Jurkat Cell Lysate

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