

# TEM2 (CT) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat TEM2 (CT)

### Instruction Manual

Catalog Number	PK-AB718-4361
Synonyms	TEM2 Antibody; Tumor endothelial marker 2, Ras homolog enriched in striatum, Rhes, RASD2
Description	Rhes, also known as tumor endothelial marker 2 (TEM2), is a small GTP-binding protein that is predominantly expressed in the striatal region of the brain. This protein belongs to the RASD subfamily of Ras-related GTP-binding protein superfamily and is thought to play a role in the normal development and function of the brain as mice lacking this gene showed increased anxiety levels and motor coordination deficits. Rhes was identified as TEM2 through analysis of genes whose expression was upregulated in tumor endothelium. Tumor endothelial markers 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.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TEM2 antibody was raised in rabbits against a 15 amino acid peptide near the carboxy terminus of the human TEM2.
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 <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TEM2 antibody can be used for detection of TEM2 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	Vargiu P, Morte B, Manzano J, et al. Thyroid hormone regulation of rhes, a novel RAS homolog gene expressed in the striatum. <i>Brain Res. Mol. Brain Res.</i> 2001; 94:1-8. Spano D, Branchi I, Rosica A, et al. Rhes is involved in striatal function. <i>Mol. Cell. Biol.</i> 2004; 24:5788-96. 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.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4361P; TEM2 Peptide Cat.No. PK-AB718-1320; Human Colon Tissue Lysate

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