

# PTK7 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat PTK7 (NT)

## Instruction Manual

Catalog Number	PK-AB718-4299
Synonyms	PTK7 Antibody: Protein tyrosine Kinase 7, colon carcinoma kinase-4, CCK-4
Description	Protein-tyrosine kinases (PTKs) play important roles in regulating cell proliferation and differentiation during development. One member of the PTK family, PTK7, has been suggested to regulate the planar cell polarity (PCP) pathway in vertebrates and may play a role in neural convergent extension and neural tube closure. PTK7 has also been implicated in the development of cancer. Loss of PTK7 expression was seen in several melanoma cell lines and biopsies. Conversely, high-throughput analysis of acute myeloid leukemia samples showed an increased level of PTK7 expression compared to normal bone marrow and purified CD34+ cells. Multiple isoforms of PTK7 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	PTK7 antibody was raised in rabbits against a 15 amino acid peptide from near the amino terminus of human PTK7.
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. PTK7 antibody can be used for detection of PTK7 by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL.
Images	Available upon request.
References	Park SK, Lee HS and Lee ST. Characterization of the human full-length PTK7 cDNA encoding a receptor protein kinase-like molecule closely related to chick KLG. <i>J. Biochem.</i> 1996; 119:235-9. Lu X, Borchers AG, Jolicoeur C, et al. PTK7/CCK-4 is a novel regulator of planar cell polarity in vertebrates. <i>Nature</i> 2004; 430:93-8. Easty DJ, Mitchell PJ, Patel K, et al. Loss of expression of receptor tyrosine kinase family genes PTK7 and SEK in metastatic melanoma. <i>Int. J. Cancer</i> 1997; 1061-5. Muller-Tidow C, Schwable J, Steffen B, et al. High-throughput analysis of genome-wide receptor tyrosine kinase expression in human cancers identifies potential novel drug targets. <i>Clin. Cancer Res.</i> 2004; 10:1241-9.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4299P; PTK7 Peptide Cat.No. PK-AB718-1320; Human Colon Tissue Lysate

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