

Anti-PTK7

Rabbit Polyclonal Antibody

Catalog # P95-63R

Lot # O2361-45

Cited Applications

E, WB, IHC, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the PTK7 protein

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

PTK7 antibody was raised against a 18 amino acid synthetic peptide from near the carboxy terminus of human PTK7

Formulation

PBS + 0.02% sodium azide.

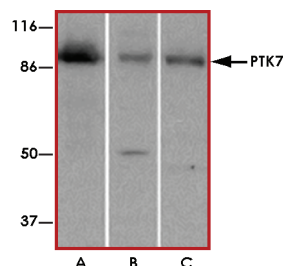
Stability

1yr at -20°C from date of shipment.

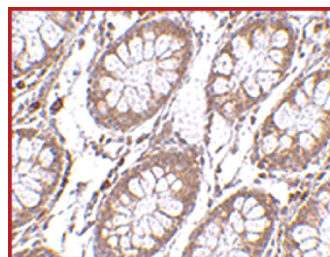
References

1. Park SK et al: Characterization of the human full-length PTK7 cDNA encoding a receptor protein kinase-like molecule closely related to chick KLG. J. Biochem.1996; 119:235-9.
2. Lu X et al: PTK7/CCK-4 is a novel regulator of planar cell polarity in vertebrates. Nature2004; 430:93-8.
3. Easty DJ et al: Loss of expression of receptor tyrosine kinase family genes PTK7 and SEK in metastatic melanoma. Int. J. Cancer1997; 1061-5.
4. Muller-Tidow C et al: High-throughput analysis of genome-wide receptor tyrosine kinase expression in human cancers identifies potential novel drug targets. Clin. Cancer Res.2004; 10:1241-9.

Sample Data



Western blot analysis of PTK7 in (A) human colon, (B) mouse kidney and (C) rat liver tissue lysate with PTK7 antibody at 1 µg/mL.



Immunohistochemistry of PTK7 in human colon tissue with PTK7 antibody at 2.5 µg/mL.

Anti-PTK7

Rabbit Polyclonal Antibody

Catalog Number

P95-63R

Specific Lot Number

O2361-45

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.