

Anti-RSL1D1

Rabbit Polyclonal Antibody

Catalog # **R76-363R**

Lot # B3216-71

Cited Applications

ELISA, IP, WB

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

Specificity

Recognizes the human PBK1 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to an internal sequence of human PBK1

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide.

Stability

1yr at -20°C from date of shipment

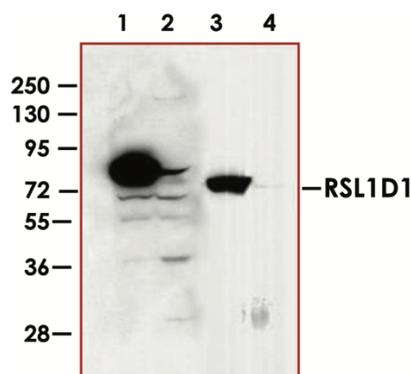
Scientific Background

This antibody is designed, produced, and validated as part of a collaboration between Rockland and the National Cancer Institute (NCI) and is suitable for Cancer, Immunology and Nuclear Signaling research. PBK1 protein (also known as Ribosomal L1 domain-containing protein 1, cellular senescence-inhibited gene protein, and CATX-11) was isolated from highly invasive first trimester trophoblast cells and has been proposed to regulate their naturally occurring invasive behavior (Huch et al., 1998). PBK1 was also found to be over-expressed in non-small-cell lung cancer (NSCLC) cells (Petroziello et al., 2004). A recent study suggests that PBK1 may up-regulate the urokinase-type plasminogen activator (uPA) gene, which plays an important role in cellular matrix degradation and activation of other protease systems involved in cell invasion (Tong et al., 2005).

References:

1. Huch, G., et al.: Identification of differentially expressed genes in human trophoblast cells by differential-display RT-PCR. Placenta. 1998; 19 (8): 557-567.
2. Petroziello, et al.: Suppression subtractive hybridization and expression profiling identifies a unique set of genes overexpressed in non-small-cell lung cancer. Oncogene. 2004; 23 (46): 7734-7745.
3. Tong, C., et al.: Identification of a novel nucleus protein involved in the regulation of urokinase in 95D cells. Acta Biochim. Biophys. Sin (Shanghai). 2005; 37 (5): 303-309.

Sample Data



Western blot using Anti-RSL1D1 antibody shows detection of over-expressed PBK1 in lysates from HeLa cells transfected with Flag-PBK1. Lanes 1 and 3 contain lysate from Flag-PBK1 transfected HeLa cells. Lanes 2 and 4 contain lysate from cells transfected with null vector. Lanes 1 and 2 were blotted with anti-Flag antibody. Lanes 3 and 4 were probed with a 1:500 dilution of anti-PBK1. The band at 75 kDa, indicated by the arrowhead, corresponds to PBK1.

Anti-RSL1D1

Rabbit Polyclonal Antibody

Catalog #	R76-363R
Lot #	B3216-71
Purification	Immunoaffinity 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.