

---

## PARD3 Polyclonal Antibody

(Catalog # A51053)

---

### Background

Adapter protein involved in asymmetrical cell division and cell polarization processes. Seems to play a central role in the formation of epithelial tight junctions. Targets the phosphatase PTEN to cell junctions. Involved in Schwann cell peripheral myelination. Association with PARD6B may prevent the interaction of PARD3 with F11R/JAM1, thereby preventing tight junction assembly. The PARD6-PARD3 complex links GTP-bound Rho small GTPases to atypical protein kinase C proteins. Required for establishment of neuronal polarity and normal axon formation in cultured hippocampal neurons.

### Description

PARD3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q8TEW0

### Purification

Protein G purified

### Immunogen

Recombinant Human Partitioning defective 3 homolog protein (1068-1356AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

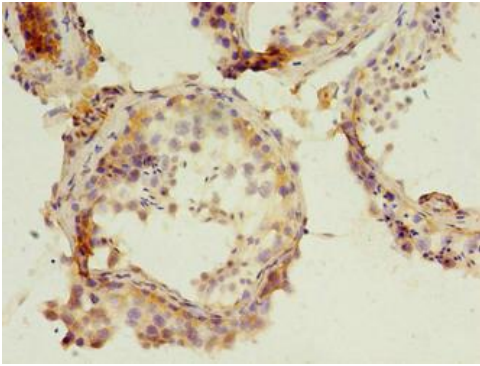
### Alternative Names

Partitioning defective 3 homolog, PAR-3, PARD-3, Atypical PKC isotype-specific-interacting protein, ASIP, CTCL tumor antigen se2-5, PAR3-alpha, PARD3, PAR3, PAR3A

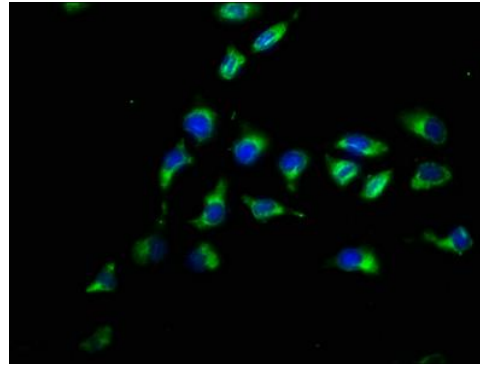
### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:100-1:1000, IF:1:50-1:500

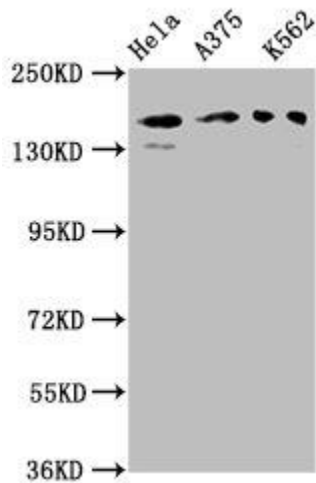
---



Immunohistochemistry of paraffin-embedded human testis tissue using PARD3 Antibody at dilution of 1:100



Immunofluorescence staining of A549 cells with PARD3 Antibody at 1:200, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



**Western Blot**

Positive WB detected in: HeLa whole cell lysate, A375 whole cell lysate, K562 whole cell lysate

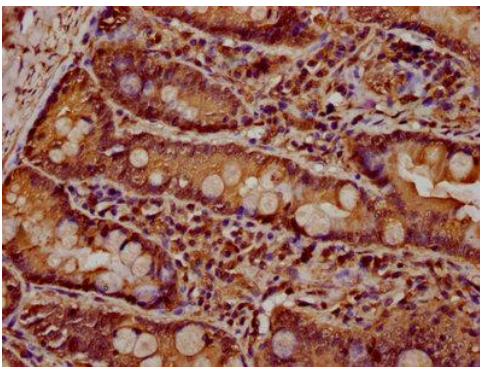
All lanes: PARD3 antibody at 3ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 152, 142, 147, 139, 114, 109, 110, 150 kDa

Observed band size: 152 kDa



IHC image of PARD3 Antibody diluted at 1:600 and staining in paraffin-embedded human small intestine tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.