



Anti-PERP

(TP53 apoptosis effector, PIGPC1, THW)

CATALOG NO.: 54078

BACKGROUND:

The p53 tumor-suppressor gene integrates numerous signals that control cell life and death. Several novel molecules involved in p53 network, including Chk2 (1), p53R2 (2), p53AIP1 (3), Noxa (4), PIDD (5), PID/MTA2 (6), MTBP (7) and PERP (8), were identified and their genes were cloned recently. PERP, also termed PIGPC1 and THW, is a plasma membrane protein (8-10). p53 binds to the promoter of PERP and transcriptionally activates PERP gene then the translated PERP protein mediates the p53 induced apoptosis (8). The expression of PERP causes cell death. PERP is a mediator of p53 induced apoptosis. PERP has sequence similarity to PMP-22/gas3 and is a new member of the PMP-22/gas3 family (8).

SOURCE AND REACTIVITY:

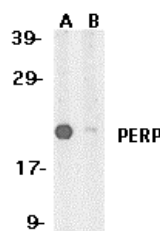
Rabbit anti-PERP polyclonal antibody was raised against a synthetic peptide corresponding to amino acids near the C-terminus of human PERP, which differ from the mouse sequence by three amino acids (7-9). Anti-PERP reacts with PERP at the molecular weight of 21 kDa on western blot. Species reactivity includes human, while others are not tested.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

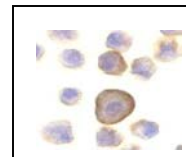
WB: 0.5 to 1 µg/ml.

Positive Control: A431 cell lysate



Western blot analysis of PERP expression in A431 whole cell lysates in the absence (A) and presence (B) of blocking peptide with anti-PERP at 1 µg/ml.

Immunocytochemistry of PERP in A431 cells with anti-PERP at 10 µg/ml.



This product is for in vitro research purposes only.

RELATED PRODUCTS:

A431 Cell Lysate, Catalog No. **29514**

Anti-CHK2 (NT), Catalog No. **54072**

Anti-P53R2 (NT), Catalog No. **54071**

STORAGE:

The antibody is supplied as purified IgG, in 1X PBS containing 0.02% sodium azide. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles..

REFERENCES:

1. Matsuoka, S. et al. *Science* **282**, 1893 (1998).
2. Tanaka, H. et al. *Nature* **404**, 42 (2000).
3. Oda, E. et al. *Science* **288**, 1053 (2000).
4. Oda, K. et al. *Cell* **102**, 849 (2000).
5. Lin, Y. et al. *Nat Genet.* **26**, 122 (2000).
6. Luo, J. et al. *Nature* **408**, 377 (2000).
7. Boyd, MT. et al. *J. Biol. Chem.* **275**, 31883 (2000).
8. Attardi, LD. Et al. *Genes Dev.* **14**, 704 (2000).
9. Goltsov, AA. et al. *GeneBank database*
10. Hildebrandt, T. et al. *GeneBank database*