

PRODUCT NAME		
p53K372me1 polyclonal antibody		
Other names: TP53, P53, TRP53, LSF1		
Cat. No. C15410048 (pAb-048-050)	Type: Polyclonal ChIP-grade	Size: 50 µg/ 25 µl
Lot #: 001	Source: Rabbit	Concentration: 2.0 µg/µl

Description: Polyclonal antibody raised in rabbit against human p53 (tumor protein p53), methylated at K372, using a KLH-conjugated synthetic peptide.

Specificity: Human: positive
Other species: not tested

Applications	Suggested dilution	References
Western blotting	1:200	Fig 1
ChIP	5 µg/ChIP	Ref 1

Purity: Protein G purified polyclonal antibody in PBS containing 0.05% azide and 0.05% ProClin 300.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles.

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures

This antibody has been described in:

(1) Chuikov S, Kurash JK, Wilson JR, Xiao B, Justin N, Ivanov GS, McKinney K, Tempst P, Prives C, Gambelin SJ, Barlev NA and Reinberg D (2004) Regulation of p53 activity through lysine methylation. Nature. 432: 353-560.

Last data sheet update: April 22, 2010

Target description

The transcription factor p53 (UniProt/Swiss-Prot entry P04637) is a tumour suppressor that regulates the cellular response to diverse cellular stresses. Upon activation, p53 induces several target genes which leads to cell cycle arrest and DNA repair, or alternatively, to apoptosis. In unstressed cells, p53 is kept inactive by the ubiquitin ligase MDM2 which inhibits the activity and promotes the degradation. p53 is a short-lived protein which is activated and stabilized via different post-translational modifications. One of these modifications, methylation of lysine 372 has been shown to positively affect the stability of p53. Mutations in p53 are involved in a vast majority of human cancers.

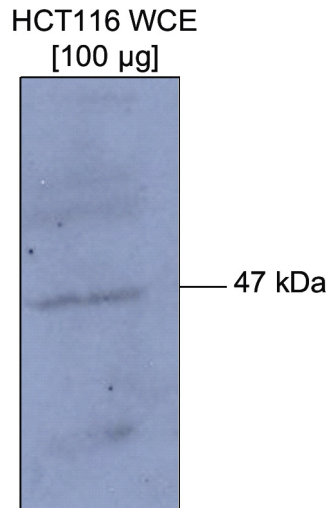


Figure 1

Western blot analysis using the Diagenode antibody directed against p53K372me1

Western blot was performed on 100 µg Colon carcinoma whole cell extract (HCT116 WCE) using the Diagenode antibody against p53K372me1 (Cat. No. pAb-048-050), diluted 1:200 in TBS-Tween containing 5% skimmed milk. The size and location of the protein of interest are indicated (Figure 1, kindly provided by Liu Xiangyu, Health Science Center, Peking University, China).