

Anti-Phospho-MDM2 (Ser185)

Rabbit Polyclonal Antibody

Catalog # **M45-365R**

Lot # B3216-57

Cited Applications

ELISA, IP, WB

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

Specificity

Recognizes mouse MDM2 phosphorylated at Serine185

Cross Reactivity

Human, Mouse and Rat

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against synthesized peptide corresponding to amino acids 177-198 of mouse MDM2

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

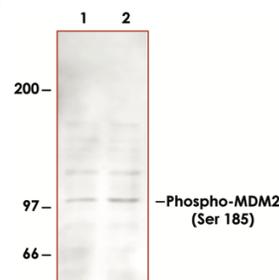
Scientific Background

MDM2 is a nuclear phosphoprotein that binds and inhibits transactivation by p53, as part of an autoregulatory negative feedback loop (1). Overexpression of the MDM2 gene product can lead to excessive inactivation of p53 thereby diminishing its tumor suppressor function. The inactivation of p53 is mediated by the E3 ubiquitin ligase activity of MDM2 which targets p53 for proteasomal degradation. MDM2 also affects the cell cycle, apoptosis, and tumorigenesis through interactions with other proteins, including retinoblastoma 1 and ribosomal protein L5 (2). Amplification of MDM2 is frequently observed in human sarcomas and this is consistent with the hypothesis that MDM2 binds to p53 which then leads to escape from p53-regulated growth control.

References:

1. Engel, T. et al: Elevated p53 and lower MDM2 expression in hippocampus from patients with intractable temporal lobe epilepsy. *Epilepsy Res.* 2007;77(2-3):151-6
2. Liu, G. et al: Genetic polymorphisms of MDM2, cumulative cigarette smoking and nonsmall cell lung cancer risk. *Int J Cancer.* 2008;122(4):915-8

Sample Data



Affinity Purified Anti-Phospho-MDM2 (Ser185) is shown to detect a 102 kDa band (arrow) corresponding to phosphorylated mouse MDM2 present in a 293T whole cell lysate. ~20 µg of lysate was loaded per lane for SDS-PAGE. Lane 1: Untreated cells, Lane 2: cells treated with IGF-1 (100ng/ml) for 20 min prior to harvest.

Anti-phospho-MDM2 (Ser185)

Rabbit Polyclonal Antibody

Catalog #	M45-365R
Lot #	B3216-57
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.