

PRODUCT INFORMATION

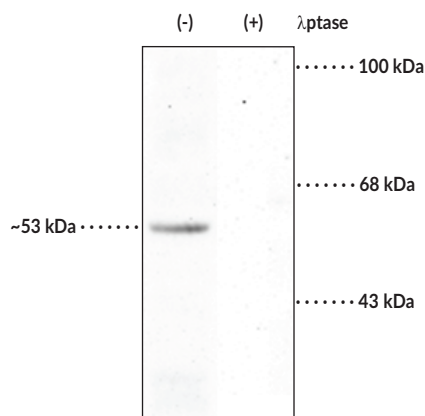
p53 (Phospho-Ser³⁹²) Polyclonal Antibody

Item No. 10004807

Overview and Properties

Contents:	This vial contains 10 µg of affinity-purified IgG polyclonal antibody.
Synonyms:	Antigen NY-CO-13, Cellular Tumor Antigen p53, Phosphoprotein p53, Transformation-related Protein 53, Tumor Protein p53, Tumor Suppressor p53
Immunogen:	Synthetic peptide corresponding to amino acid residues surrounding phospho-Ser ³⁹² of human p53 conjugated to keyhole limpet hemocyanin (KLH)
Molecular Weight:	~53 kDa
Cross Reactivity:	(+) Human, rat p53 (phospho-Ser ³⁹²)
Species Reactivity:	(+) Human, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	100 µl of 10 mM HEPES, pH 7.5, with 50% glycerol, 0.1% BSA, and 150 mM sodium chloride
Host:	Rabbit
Isotype:	IgG
Applications:	Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



(-): WB of rat brain nuclear fraction lysate showing specific immunolabeling of the ~53 kDa p53 phosphorylated at Ser³⁹².
 (+): Phosphospecificity is shown where the immunolabeling is completely eliminated by blot treatment with λ. phosphatase (λ.ptase, 1,200 units for 30 minutes).

WARNING
 THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
 This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
 Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/21/2021

CAYMAN CHEMICAL
 1180 EAST ELLSWORTH RD
 ANN ARBOR, MI 48108 · USA
 PHONE: [800] 364-9897
 [734] 971-3335
 FAX: [734] 971-3640
 CUSTSERV@CAYMANCHEM.COM
 WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

p53 is a transcription factor and tumor suppressor encoded by *TP53* in humans with roles in cell cycle arrest, apoptosis, senescence, differentiation, and DNA repair.^{1,2} It is composed of two N-terminal intrinsically disordered transactivation domains (TADs), a proline-rich domain (PRD), a DNA-binding domain (DBD), a tetramerization domain (TD), and an intrinsically disordered C-terminal domain.¹ p53 is ubiquitously expressed, with protein levels increasing under conditions of extra- or intracellular stress, such as DNA damage, oncogene activation, oxidative stress, or hypoxia.³ It is activated and translocated to the nucleus in response to cellular stress *via* post-translational modifications, such as phosphorylation, methylation, and acetylation, where it binds p53 consensus DNA-binding elements and regulates transcription of its target genes in a cell-, tissue-, and stress signal type-specific manner.^{1,3} p53 is phosphorylated at Ser³⁹² by IκB kinase β (IKKβ) in glutamine-deficient cancer cells.⁴ Loss-of-function and/or gain-of-function mutations in *TP53* occur in approximately 50% of human cancers.² Cayman's p53 (Phospho-Ser³⁹²) Polyclonal Antibody can be used for Western blot (WB) applications. This antibody recognizes p53 at ~53 kDa from human and rat samples.

References

1. Liu, Y., Tavana, O., and Gu, W. p53 Modifications: Exquisite decorations of the powerful guardian. *J. Mol. Cell Biol.* **11**(7), 564-577 (2019).
2. Stein, Y., Rotter, V., and Aloni-Grinstein, R. Gain-of-function mutant p53: All the roads lead to tumorigenesis. *Int. J. Mol. Sci.* **20**(24), 6197 (2019).
3. Yue, X., Zhao, Y., Xu, Y., *et al.* Mutant p53 in cancer: Accumulation, gain-of-function and therapy. *J. Mol. Biol.* **429**(11), 1595-1606 (2017).
4. Gabra, M.B.I., Yang, Y., Lowman, X.H., *et al.* IKKβ activates p53 to promote cancer cell adaptation to glutamine deprivation. *Oncogenesis* **7**(11), 93 (2018).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM