

PRODUCT INFORMATION



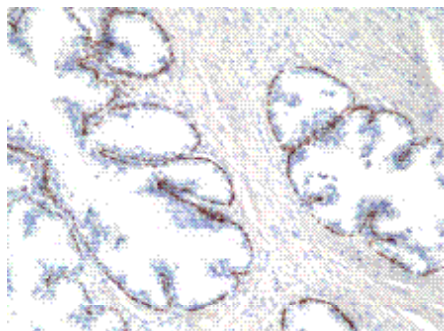
Δ Np63 (N-Term) Rabbit Monoclonal Antibody (RM392)

Item No. 32316

Overview and Properties

Contents:	This vial contains 100 μ l of protein A-affinity purified monoclonal antibody.
Synonym:	p40
Immunogen:	Peptide from the N-terminal region of human Δ Np63
Cross Reactivity:	(+) Δ Np63
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM392
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC); the recommended starting dilution is 1:50-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemical staining of formalin-fixed and paraffin-embedded human prostate tissue using Cayman's Δ Np63 (N-Term) Rabbit Monoclonal Antibody (RM392) at a dilution of 1:200.

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, 02/21/2024

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

p63 is a transcription factor and member of the p53 gene family encoded by *TP63* in humans.^{1,2} Alternative promoter use and splicing of *TP63* mRNA produces several variants, including α , β , γ , δ , and ϵ isoforms in full-length or N-terminally truncated forms known as TAp63 and Δ Np63, respectively. Δ Np63 is expressed in kidney, spleen, thymus, and epithelia but not the heart, liver, testes, or brain and is localized to the cytosol as a tetramer.² It is activated in response to various signaling pathways, including EGFR, Wnt, and NF-kB, and has roles in cell survival, senescence, stemness, and metabolism, as well as oncogenesis. Δ Np63 levels are positively correlated with poor response to chemotherapeutics in various cancers. N- and C-terminal mutations in Δ Np63 are associated with epidermal defects and skin fragility in patients with ankyloblepharon ectodermal dysplasia and clefting (AEC).^{2,3} Cayman's Δ Np63 (N-Term) Rabbit Monoclonal Antibody (RM392) can be used for immunohistochemistry (IHC).

References

1. Soares, E. and Zhou, H. Master regulatory role of p63 in epidermal development and disease. *Cell. Mol. Life Sci.* **75(7)**, 1179-1190 (2017).
2. Fisher, M.L., Balinth, S., and Mills, A.A. p63-related signaling at a glance. *J. Cell Sci.* **133(17)**, jcs228015 (2020).
3. Koster, M.I., Marinari, B., Payne, A.S., et al. Δ Np63 knockdown mice: A mouse model for AEC syndrome. *Am. J. Med. Genet. A.* **149A(9)**, 1942-1947 (2009).

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