

PRODUCT INFORMATION



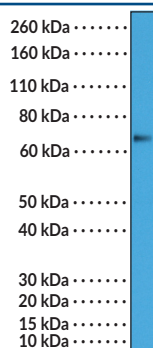
p63 Rabbit Monoclonal Antibody (Clone RM383)

Item No. 32314

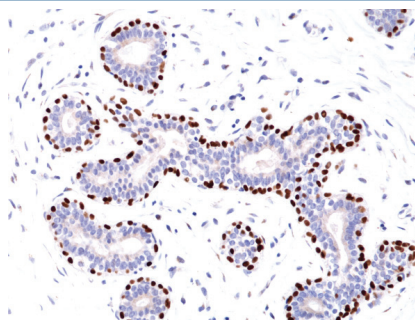
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	TP73L, Tumor Protein p63, Tumor Protein p73-like
Immunogen:	Peptide corresponding to residues near the C-terminus of human p63
Cross Reactivity:	(+) p63
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM383
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:500-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

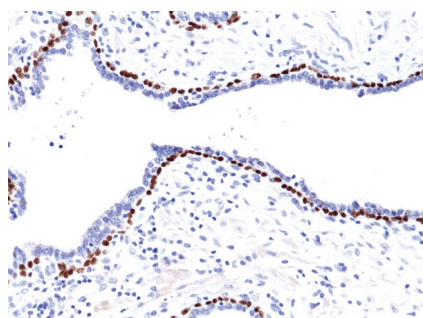
Images



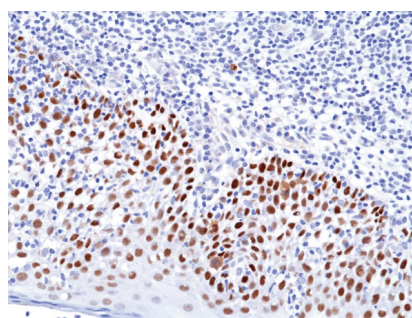
WB of A431 cell lysates using p63 Rabbit Monoclonal Antibody (Clone RM383) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast tissue using p63 Rabbit Monoclonal Antibody (Clone RM383) at a 1:100 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human prostate tissue using p63 Rabbit Monoclonal Antibody (Clone RM383) at a 1:100 dilution.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using p63 Rabbit Monoclonal Antibody (Clone RM383) at a 1:100 dilution.

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, 02/22/2024

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 with roles in epidermal development and lineage specification, as well as cell proliferation, death, senescence, and differentiation.^{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, which exhibit distinct functions, tissue expression, and cellular localization. Mutations in *TP63* are associated with various developmental disorders, including ectrodactyly ectodermal dysplasia-clefting syndrome, split-hand/foot malformations, Rapp-Hodgkin syndrome, and primary ovarian insufficiency. Cayman's p63 Rabbit Monoclonal Antibody (Clone RM383) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

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).

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