

PRODUCT INFORMATION



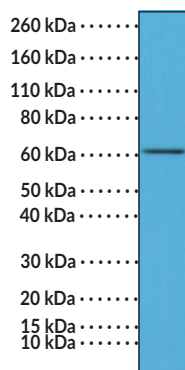
NF- κ B (p65) Rabbit Monoclonal Antibody (Clone RM273)

Item No. 32223

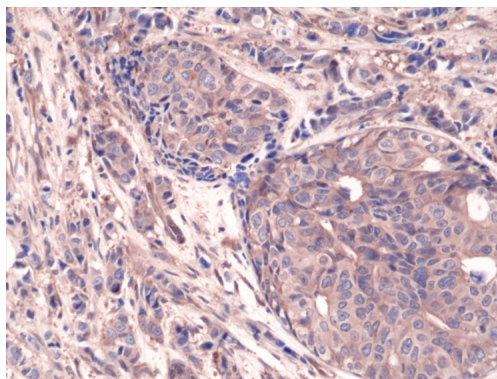
Overview and Properties

Contents:	This vial contains 100 μ l of protein A-affinity purified monoclonal antibody.
Synonyms:	Nuclear Factor NF- κ B p65 Subunit, Transcription Factor p65
Immunogen:	Peptide from the C-terminal region of NF- κ B (p65)
Cross Reactivity:	(+) NF- κ B
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:	RM273
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:500-1:1,500 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 MCF-7 cells. MCF-7 cells were subjected to WB using a 1:500 dilution of NF- κ B (p65) Rabbit Monoclonal Antibody (Clone RM273).



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast cancer tissue using a 1:1,250 dilution of NF- κ B (p65) Rabbit Monoclonal Antibody (Clone RM273).

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, 10/16/2020

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

PRODUCT INFORMATION



Description

NF- κ B p65 is a ubiquitously expressed transcription factor that is a subunit of the NF- κ B complex and is encoded by the *RELA* gene in humans.¹ It is composed of an N-terminal Rel homology domain, which contains the nuclear localization signal (NLS), and mediates dimerization, nuclear localization, and DNA and protein interactions, and two C-terminal transactivation domains that are subject to a variety of post-translational modifications and regulate the transcriptional activity of p65.^{1,2} NF- κ B p65 regulates the expression of a large number of genes in response to inflammatory and environmental cues that play critical roles in innate and adaptive immunity and cellular differentiation.² Silencing of *Rela* induces tumor cell apoptosis in a murine Lewis lung carcinoma model, and *RELA* silencing in THP-1 monocytes decreases secreted levels of IL-1 β and TNF- α induced by LPS.^{3,4} Genome-wide deletion of *Rela* in mice is embryonic lethal.⁵ NF- κ B p65 is overexpressed in the inflamed joints of patients with rheumatoid arthritis, and naïve CD4 T cells isolated from the whole blood of patients with multiple sclerosis have increased phosphorylation of NF- κ B p65.^{6,7} Cayman's NF- κ B (p65) Rabbit Monoclonal Antibody (Clone RM273) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Peterson, J.M., Bakkar, N., and Guttridge, D.C. NF- κ B signaling in skeletal muscle health and disease. *Curr. Top. Dev. Biol.* Pavlath, G.K., editor, Elsevier (2011).
2. Giridharan, S. and Srinivasan, M. Mechanisms of NF- κ B p65 and strategies for therapeutic manipulation. *J. Inflamm. Res.* **11**, 407-419 (2018).
3. Qu, Y., Zhang, X. and Wu, R. Knockdown of NF- κ B p65 subunit expression suppresses growth of nude mouse lung tumor cell xenografts by activation of Bax apoptotic pathway. *Neoplasma* **62(1)**, 34-40 (2015).
4. Wu, C., Zhao, J., Zhu, G., et al. siRNA directed against NF- κ B inhibits mononuclear macrophage cells releasing proinflammatory cytokines *in vitro*. *Mol. Med. Rep.* **16(6)**, 9060-9066 (2017).
5. Beg, A.A. and Baltimore, D. An essential role for NF- κ B in preventing TNF- α -induced cell death. *Science* **274(5288)**, 782-784 (1996).
6. Makarov, S.S. NF- κ B in rheumatoid arthritis: A pivotal regulator of inflammation, hyperplasia, and tissue destruction. *Arthritis Res.* **3(4)**, 200-206 (2001).
7. Housley, W.J., Fernandez, S.D., Vera, K., et al. Genetic variants associated with autoimmunity drive NF κ B signaling and responses to inflammatory stimuli. *Sci. Transl. Med.* **7(291)**, 291ra93 (2015).

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