

PRODUCT INFORMATION



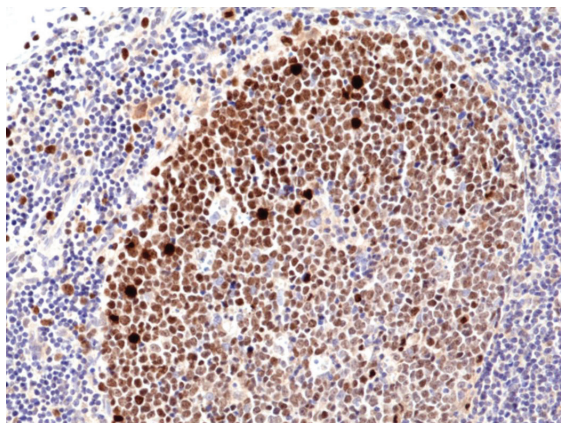
Ki-67 Rabbit Monoclonal Antibody (Clone RM360)

Item No. 32309

Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Antigen Ki-67, Marker of Proliferation Ki-67
Immunogen:	Peptide corresponding to the internal region of human Ki-67
Cross Reactivity:	(+) Ki-67
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:	RM360
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemical staining of formalin-fixed and paraffin-embedded tonsil tissue using Ki-67 Rabbit Monoclonal Antibody (Clone RM360) at a dilution of 1:1,000.

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

Ki-67 is a non-histone nuclear and nucleolar protein and marker of cell proliferation encoded by *MKI67* in humans.^{1,2} It is composed of an N-terminal forkhead-associated domain, a protein phosphatase 1-binding domain, a central conserved domain, 16 concatenated direct repeats, and six C-terminal leucine-arginine rich motif repeats.² Ki-67 expression and localization is cell cycle-specific. It is localized to the perinuclear region during G₁, the fibrillary component of the nucleolus during interphase, and the perichromosomal layer during chromatin condensation, with Ki-67 levels increasing from late G₁ phase to S phase and peaking at mitosis.¹ Ki-67 is associated with proliferative activity of intrinsic cell populations in tumors and is commonly used as a marker of tumor aggressiveness in patients with breast, lung, prostate, cervical, or CNS cancers.³ Ki-67 is also commonly used as a marker of adult neurogenesis.⁴ Cayman's Ki-67 Rabbit Monoclonal Antibody (Clone RM360) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Menon, S.S., Guruvayoorappan, C., Sakthivel, K.M., *et al.* Ki-67 protein as a tumour proliferation marker. *Clin. Chim. Acta* **491**, 39-45 (2019).
2. Sun, X. and Kaufman, P.D. Ki-67: More than a proliferation marker. *Chromosoma* **127(2)**, 175-186 (2018).
3. Li, L.T., Jiang, G., Chen, Q., *et al.* Ki67 is a promising molecular target in the diagnosis of cancer (review). *Mol. Med. Rep.* **11(3)**, 1566-1572 (2015).
4. Kee, N., Sivalingam, S., and Wojtowicz, J.M. The utility of Ki-67 and BrdU as proliferative markers of adult neurogenesis. *J. Neurosci. Methods* **115(1)**, 97-105 (2002).

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