

# PRODUCT INFORMATION

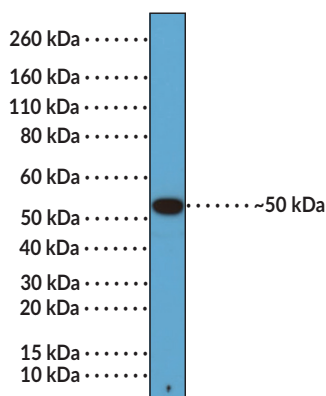
## $\alpha$ -Tubulin (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM113)

Item No. 32376

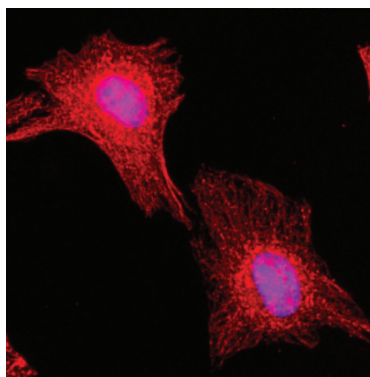
### Overview and Properties

<b>Contents:</b>	This vial contains 50 $\mu$ l of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	$\alpha$ -Tubulin Chain
<b>Immunogen:</b>	Peptide from the C-terminal region of $\alpha$ -tubulin
<b>Cross Reactivity:</b>	(+) $\alpha$ -Tubulin 1A, $\alpha$ -Tubulin 1B
<b>Species Reactivity:</b>	Species independent
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	$\geq$ 1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM113
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunocytochemistry (ICC) and Western blot (WB); the recommended starting dilution is 1:200 for ICC and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



WB of A431 cells using  $\alpha$ -Tubulin (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM113) at a dilution of 1:1,000.



Immunofluorescent labeling of HeLa cells with  $\alpha$ -Tubulin (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM113) at a 1:200 dilution (red). Nuclei have been labeled with DAPI (blue).

**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, 03/12/2024

**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

---

$\alpha$ -Tubulin is a cytoskeletal protein and constituent of microtubules, a cytoskeletal assembly that has roles in a variety of cellular processes, including cell motility, division, differentiation, and intracellular transport.<sup>1</sup>  $\alpha$ -Tubulin is highly conserved in eukaryotes and expressed in a cell- and isotype-specific manner.<sup>2</sup> There are eight human  $\alpha$ -tubulin isotypes that consist of an N-terminal domain, which binds GTP and is required for microtubule self-assembly, and a variable C-terminal tail, which contains interaction sites for microtubule-associated proteins (MAPs) and is subject to a variety of post-translational modifications that regulate microtubule function and stability.<sup>1,3-5</sup> Cayman's  $\alpha$ -Tubulin (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM113) can be used for immunocytochemistry (ICC) and Western blot (WB) applications. The antibody recognizes  $\alpha$ -tubulin at approximately 50 kDa.

## References

---

1. Gadadhar, S., Bodakuntla, S., Natarajan, K., *et al.* The tubulin code at a glance. *J. Cell Sci.* **130**(8), 1347-1353 (2017).
2. Binarová, P. and Tuszynski, J. Tubulin: Structure, functions and roles in disease. *Cells* **8**(10), 1294 (2019).
3. Vemu, A., Atherton, J., Spector, J.O., *et al.* Tubulin isoform composition tunes microtubule dynamics. *Mol. Biol. Cell* **28**(25), 3564-3572 (2017).
4. Valiron, O. New insights into microtubule elongation mechanisms. *Commun. Integr. Biol.* **4**(1), 10-13 (2011).
5. Zhang, F., Su, B., Wang, C., *et al.* Posttranslational modifications of  $\alpha$ -tubulin in Alzheimer disease. *Transl. Neurodegener.* **4**(9), (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