

PRODUCT INFORMATION

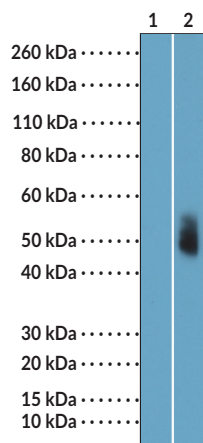
Aurora A (Phospho-Thr²⁸⁸)/Aurora B (Phospho-Thr²³²)/Aurora C (Phospho-Thr¹⁹⁸) Rabbit Monoclonal Antibody (Clone RM454)

Item No. 35904

Overview and Properties

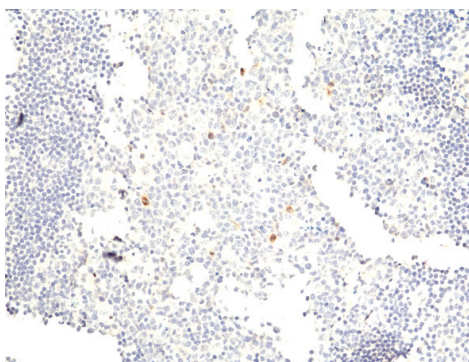
Contents:	This vial contains 100 µl protein A-affinity purified monoclonal antibody.
Immunogen:	A phospho-peptide corresponding to Phospho-Aurora A (Thr ²⁸⁸)/Aurora B (Thr ²³²)/Aurora C (Thr ¹⁹⁸)
Cross Reactivity:	(+) Aurora A (Thr ²⁸⁸)/Aurora B (Thr ²³²)/Aurora C (Thr ¹⁹⁸); (-) Aurora A/Aurora B/Aurora C without phosphorylation at Aurora A (Thr ²⁸⁸)/Aurora B (Thr ²³²)/Aurora C (Thr ¹⁹⁸)
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:	RM454
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:50-1:200 for IHC and 1:1,000-1:3,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: HeLa cell lysates untreated
Lane 2: HeLa cell lysates treated

WB of HeLa cell lysates untreated or treated with nocodazole and calyculin A using Aurora A (Phospho-Thr²⁸⁸)/Aurora B (Phospho-Thr²³²)/Aurora C (Phospho-Thr¹⁹⁸) Rabbit Monoclonal Antibody (Clone RM454) at a dilution of 1:2,500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using Aurora A (Phospho-Thr²⁸⁸)/Aurora B (Phospho-Thr²³²)/Aurora C (Phospho-Thr¹⁹⁸) Rabbit Monoclonal Antibody (Clone RM454) at a dilution of 1:50.

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, 01/30/2024

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

The Aurora kinases, Aurora A, -B, and -C, are a family of serine/threonine kinases encoded by distinct genes, *AURKA*, *AURKB*, and *AURKC*, respectively, that have roles in mitosis and meiosis.^{1,2} They are composed of an ATP-binding N-terminal domain, a highly conserved catalytic domain, and a short C-terminal domain that contains an activation loop and mediates cofactor binding.¹ Each Aurora kinase contains a key threonine residue in its C-terminal activation loop, threonine 288 (Thr²⁸⁸), Thr²³², and Thr¹⁹⁸ in Aurora A, -B, and -C, respectively, that is autophosphorylated upon cofactor binding, inducing a conformational change that mediates kinase activation.^{1,2} Aurora A and Aurora B have roles in mitosis, whereas Aurora C regulates meiosis.² Aurora A and Aurora B are constitutively expressed in mitotic cells, where Aurora A localizes to the spindle poles and Aurora B localizes to the central spindle.² Aurora A has roles in mitotic entry, centrosome maturation and separation, and spindle assembly, whereas Aurora B regulates chromosome condensation and alignment, spindle assembly, and cytokinesis. Aurora C is expressed in germ cells during meiosis, localizes to the chromosome midbody, and is required for spermatogenesis.^{2,3} Aurora A, -B, and -C are overexpressed in a variety of cancers, including breast, colorectal, ovarian, and cervical cancers, and the pan-Aurora kinase inhibitor AMG 900 (Item No. 19176) inhibits tumor growth in an HCT116 mouse xenograft model.^{3,4} Cayman's Aurora A (Phospho-Thr²⁸⁸)/Aurora B (Phospho-Thr²³²)/Aurora C (Phospho-Thr¹⁹⁸) Rabbit Monoclonal Antibody (Clone RM454) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Willems, E., Dedobbeleer, M., Digregorio, M., *et al.* The functional diversity of Aurora kinases: A comprehensive review. *Cell Div.* **13**, 7 (2018).
2. Goldenson, B. and Crispino, J.D. The aurora kinases in cell cycle and leukemia. *Oncogene* **34(5)**, 537-545 (2015).
3. Tang, A., Gao, K., Chu, L., *et al.* Aurora kinases: Novel therapy targets in cancers. *Oncotarget* **8(14)**, 23937-23954 (2017).
4. Payton, M., Bush, T.L., Chung, G., *et al.* Preclinical evaluation of AMG 900, a novel potent and highly selective pan-aurora kinase inhibitor with activity in taxane-resistant tumor cell lines. *Cancer Res.* **70(23)**, 9846-9854 (2010).

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