

[pThr232]Aurora B polyclonal antibody

Aurora A, B, and C have emerged as key mitotic regulators of mammalian centrosome function, chromosome segregation and cytokinesis. Aurora B is one of several chromosomal passenger proteins involved in chromosome alignment at the metaphase plate and cytokinesis. Phosphorylation of a threonine residue in the Aurora activation loop (Thr232 in human Aurora B) is believed to be an essential regulatory mechanism in Aurora kinase activation.

This antibody is covered by our [Worry-Free Guarantee](#).

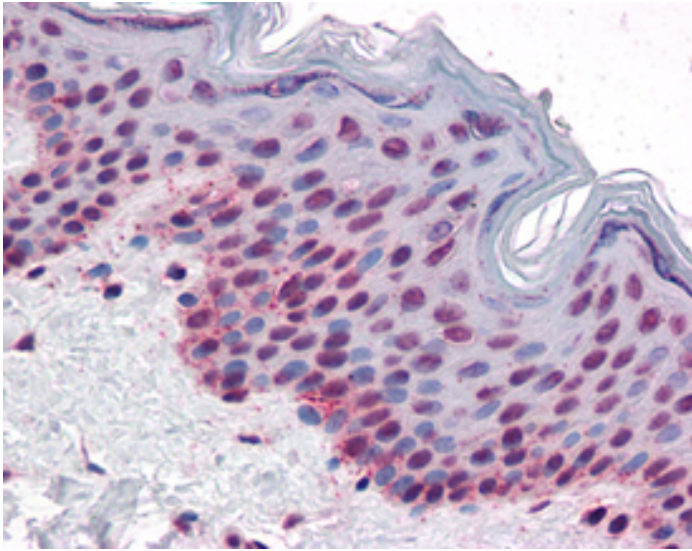
Ordering Information

[Order Online »](#)

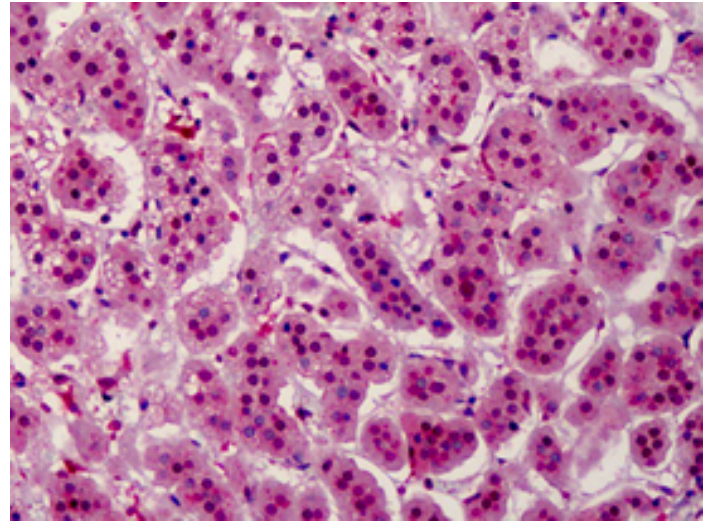
ADI-KAP-CC203-D	50µg
ADI-KAP-CC203-F	200µg

Manuals, SDS & CofA

[View Online »](#)



IHC Analysis: Human adrenal tissue stained with [pThr²³²]Aurora B, pAb at 5μg/ml.



IHC Analysis: Human skin tissue stained with [pThr²³²]Aurora B, pAb at 5μg/ml.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	IHC (PS), WB
Crossreactivity	Weakly cross-reacts with mouse, rat, bovine and pig Aurora B.
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	AAH80581
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human Aurora B phosphorylated at Thr232.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	Q96GD4
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)