

# CENP-A monoclonal antibody (3-19)

Centromere protein A (CENP-A), a 17-19 kDa centromere-specific histone variant, is 62% identical to the carboxy-terminal domain of histone H3. In the presence of DNA, CENP-A forms an octameric complex with histones H4, H2A, and H2B. CENP-A defines active centromere regions by forming centromere-specific nucleosomes on which kinetochores are assembled. CENP-A resides specifically in the inner plate of the kinetochore, and is essential for kinetochore targeting of CENP-C and other kinetochore components. CENP-A is required for nucleosomal packaging of centromeric DNA at interphase.

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

Citations: 35

[View Online »](#)

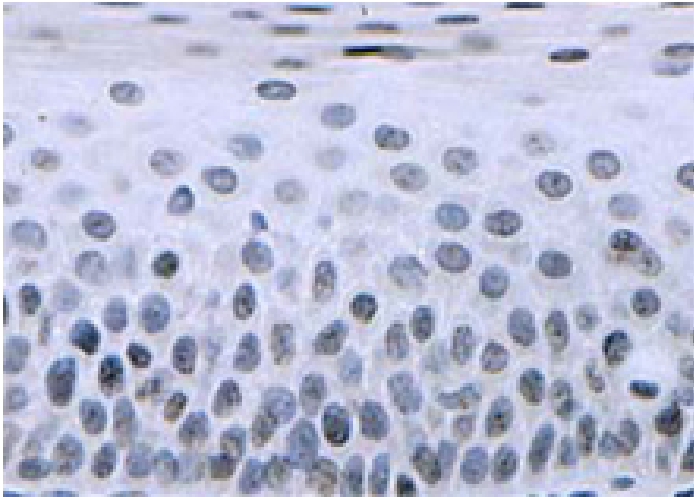
## Ordering Information

[Order Online »](#)

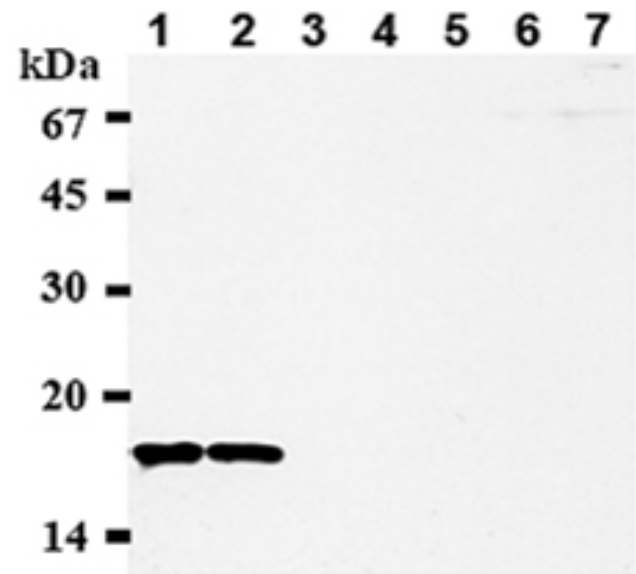
ADI-KAM-CC006-E	100µg
-----------------	-------

## Manuals, SDS & CofA

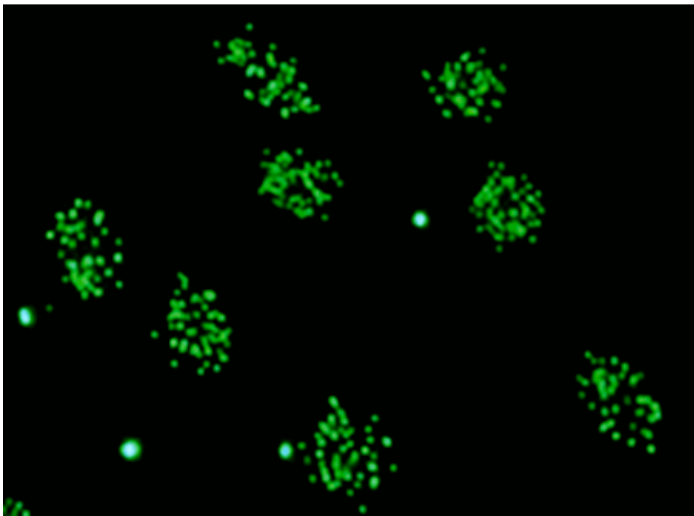
[View Online »](#)



Immunohistochemical detection of CENP-A on paraffin embedded section of a squamous epithelium of human tonsil.



Western blot analysis of Jurkat (1), Raji (2), PC12 (3), Rat-1 (4), WR19L (5), C2C12 (6), and NIH3T3 (7) cell lysates probed with CENP-A mAb (3-19).



Immunocytochemical detection of CENP-A on acetone fixed HEp-II with CENP-A mAb (3-19).

# Handling & Storage

Long Term Storage                    -20°C

Shipping                                Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Centromere protein A
Application	ICC, IHC, WB
Application Notes	Detects bands of ~18kDa by Western blot.
Clone	3-19
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Mouse
Immunogen	Synthetic peptide corresponding to a portion of human CENP-A .
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (10µg/ml)Immunohistochemistry (5µg/ml)Western Blot (1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P49450
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)