

CDKN2A polyclonal antibody

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

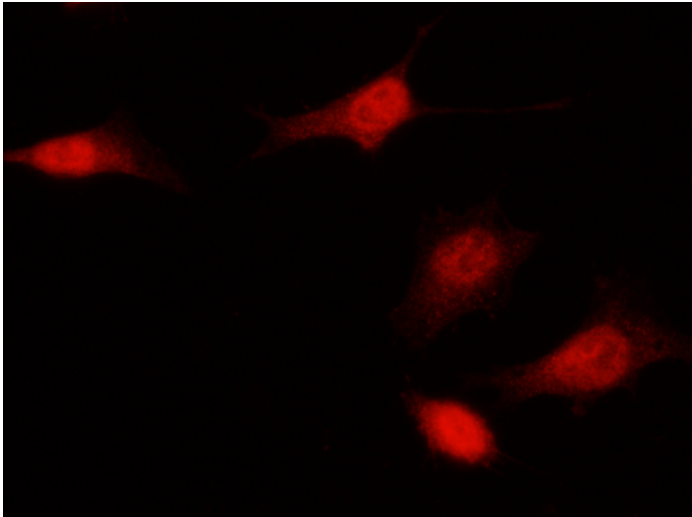
Ordering Information

[Order Online »](#)

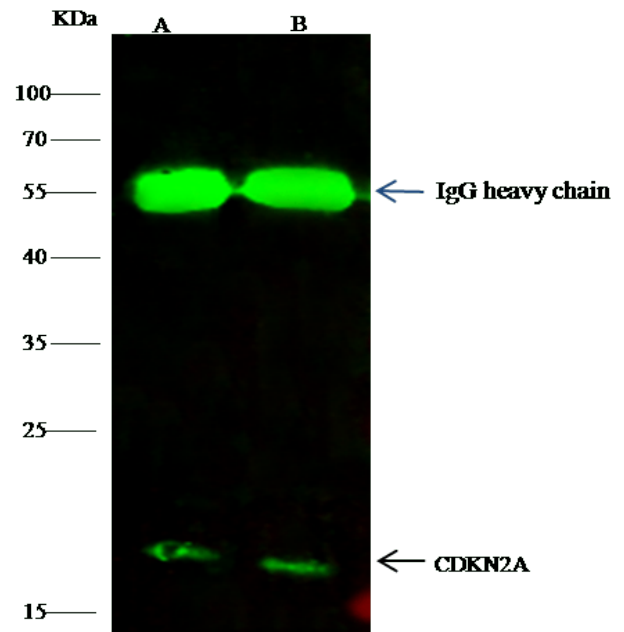
ENZ-ABS933-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of CDKN2A in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human CDKN2A polyclonal antibody (1:1000) at 4 °C overnight. Then cells were stained with the Alexa Fluor®594-conjugated Goat Anti-rabbit IgG secondary antibody (red). Positive staining was localized to cytoplasm and nucleus.



CDKN2A was immunoprecipitated using: Lane A: 0.5 mg HeLa Whole Cell Lysate. Lane B: 0.5 mg 293T Whole Cell Lysate. 2 µL anti-CDKN2A rabbit polyclonal antibody and 60 µg of Immomagnetic beads Protein G. Primary antibody: Anti-CDKN2A rabbit polyclonal antibody, at 1:200 dilution. Secondary antibody: DyLight 800-labeled antibody to rabbit IgG (H+L), at 1:5000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size: 17 kDa.



Anti-CDKN2A rabbit polyclonal antibody at 1:500 dilution. Lane A: Hela Whole Cell Lysate Lane B: 293T Whole Cell Lysate. Lysates/proteins at 30 µg per lane. Secondary Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:17 kDa.

Handling & Storage

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

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cyclin-dependent kinase inhibitor 2A, Tumor suppressor ARF
Application	ICC, IF, IP, WB
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the C-terminus of the human CDKN2A.
Isotype	IgG
Purity Detail	Protein A and antigen affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500-1:2'000)Immunocytochemistry (1:300-1:10'000)Immunofluorescence (1:300-1:10'000)Immunoprecipitation (1-4µl/mg of lysate)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
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)