

Cdk1/cdc2 monoclonal antibody (5F6)

The cell division cycle (cdc) group of proteins consists of many protein classes, including cyclin dependent kinases (Cdks), phosphoprotein phosphatases, mitogen activated kinases, small GTPases, and cyclins. These proteins play a central role in regulating progression through the cell cycle through the formation and/or modification of key substrates. Cdc42 belongs to the Rho-family of GTPases, and promotes actin polymerization to form filopodia. Cdc25A is a dual-specificity phosphatase that acts on phosphorylated threonine and tyrosine residues of Cdks, leading to binding and activation of their partner cyclins.

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

Citations: 1

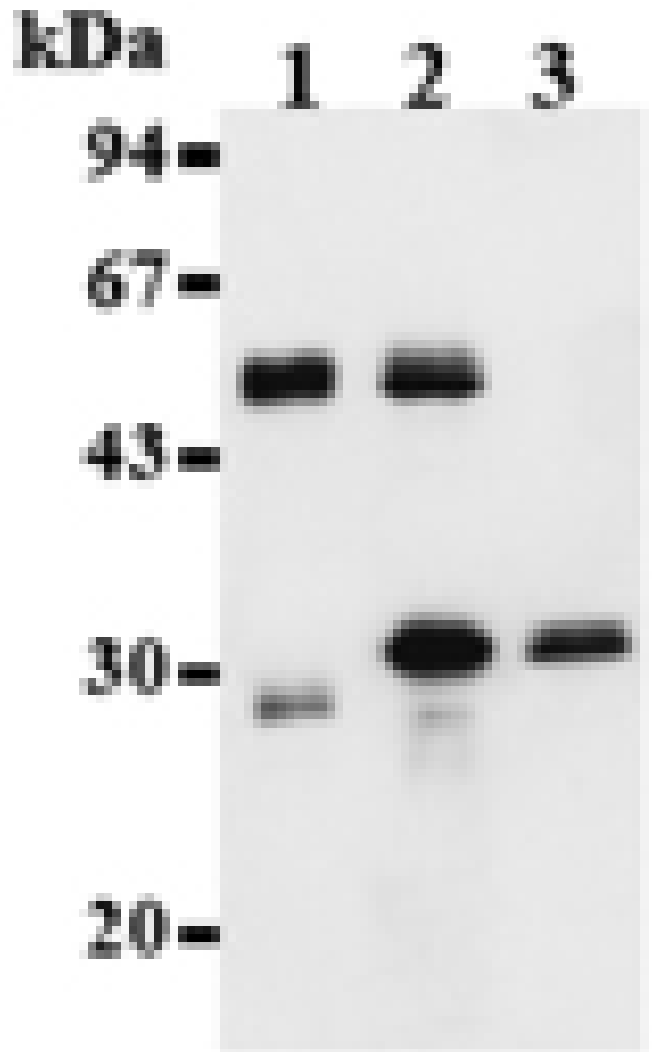
[View Online »](#)

Manuals, SDS & CofA

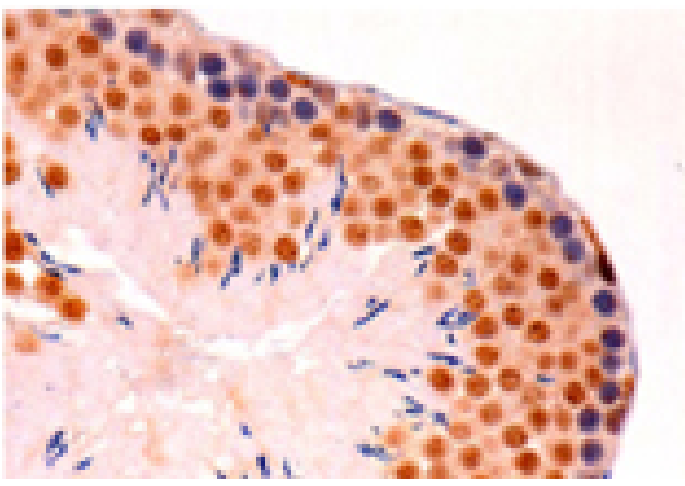
[View Online »](#)



Western blot analysis of human HeLa cell lysate probed with Cdk1/cdc2 mAb (5F6).



Immunoprecipitation analysis of Cdk1/cdc2 from A431 cells extract with normal mouse IgG (1) or Cdk1/cdc2 mAb (5F6) (2). After immunoprecipitation with the antibody, the immunocomplex was resolved by SDS-PAGE and immunoBlotted with Cdk1/cdc2 mAb (5F6). A431 crude lysate was resolved in lane 3.



Immunohistochemistry analysis of paraffin embedded section of rat testis using Cdk1/cdc2 mAb (5F6).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cyclin-dependent kinase 1, p34cdc2, Cell division cycle 2
Application	IHC, IP, WB
Application Notes	Detects a band of ~34kDa by Western blot.
Clone	5F6
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Mouse
Immunogen	Synthetic peptide corresponding to the sequence near the N-terminus of human Cdk1 (cdc2).
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (5µg/ml)Immunoprecipitation (2µg/200µl)Western Blot (1-5µg/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rabbit, Rat
UniProt ID	P06493
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



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)