

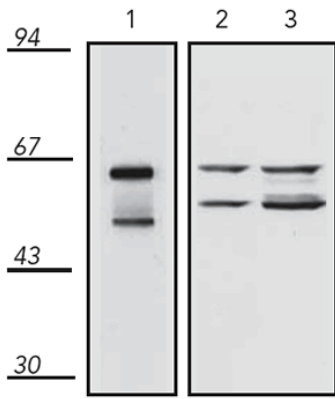
Cdc25A monoclonal antibody (DCS-120)

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](#).

Manuals, SDS & CofA

[View Online »](#)



← *cdc25A*

Western blot analysis of human HeLa cell lysate (1), mouse N1H/3T3 cell lysate (2), and rat Rat-1 cell lysate (3), probed with KAM-CC085.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cell division cycle 25A
Application	IP, WB
Clone	DCS-120
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
GenBank ID	AY137580
Gene/Protein Identifier	NM_001789 (RefSeq)
Host	Mouse
Immunogen	Recombinant human cdc25A.
Isotype	IgG2a
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (3µg/200-300µl)Western Blot (1-5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P30304
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)