

Plk1 monoclonal antibody (3F8)

The polo-like kinases (PLKs) function in cell-cycle progression and in multiple stages of mitosis. PLK1 is activated by phosphorylation of Thr210 at the G2/M phase boundary. Active PLK1 participates in mitotic entry through activation of Cdc25C and nuclear import of cyclin B1, thereby activating the Cdc2/cyclin B kinase.

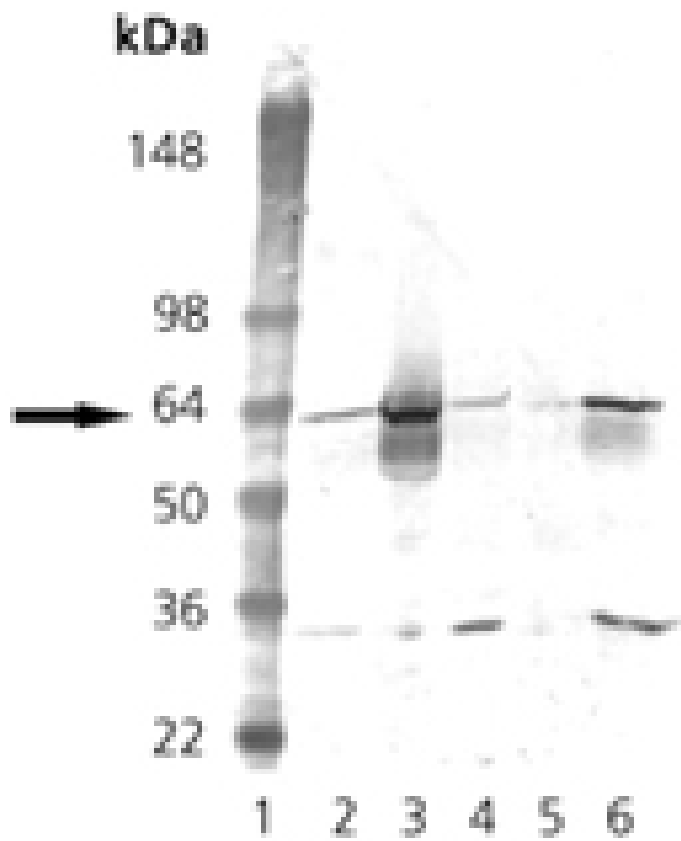
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of PLK1 mAb (3F8): Lane 1: MW marker, Lane 2: HeLa, Lane 3: Jurkat, Lane 4: PC-12, Lane 5: 3T3, Lane 6: HEK-293.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Polo-like kinase 1

Application WB

Application Notes Detects a band of ~62kDa by Western blot.

Clone 3F8

Crossreactivity Weakly cross-reacts with mouse Plk1.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Host Mouse

Immunogen Recombinant fragment from the middle of human PLK1.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Rat

UniProt ID P53350

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)