

Chk1 monoclonal antibody (2G1D5)

The DNA damage checkpoint is a signal transduction pathway that has the ability to delay entry into mitosis. When DNA is damaged, the Check point kinases (CHKs) are known to act downstream of ATM to elicit appropriate responses such as cell cycle arrest. Chk1, for instance, is activated by ATM phosphorylation of serines 123, 178, 278, and 292 of the S phase-promoting cdc25A phosphatase and accelerates IR (ionizing radiation)induced degradation of cdc25A.

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

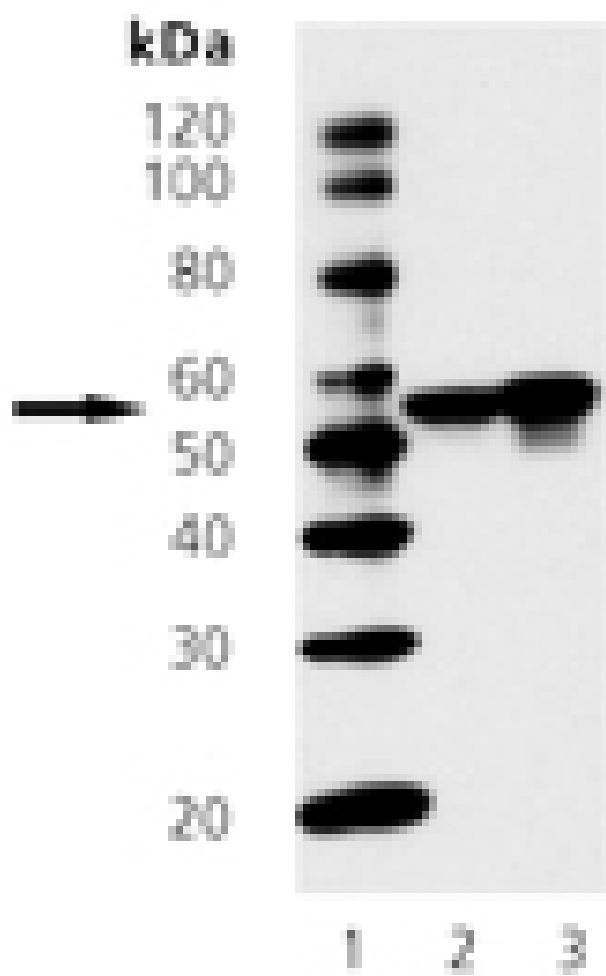
Ordering Information

[Order Online »](#)

ADI-905-770-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of CHK1 using HeLa and NIH/3T3 cell extracts. Lane 1: Molecular Weight Markers, Lane 2: HeLa, Lane 3: NIH/3T3.

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 Checkpoint kinase 1

Application WB

Application Notes Detects a band of ~55kDa by Western blot.

Clone 2G1D2

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

Host Mouse

Immunogen Recombinant human Chk1.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from mouse ascites.

Species Reactivity Human, Mouse, Rat

UniProt ID O14757

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)

This antibody is covered by our Worry-Free Guarantee.