

# 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](#).

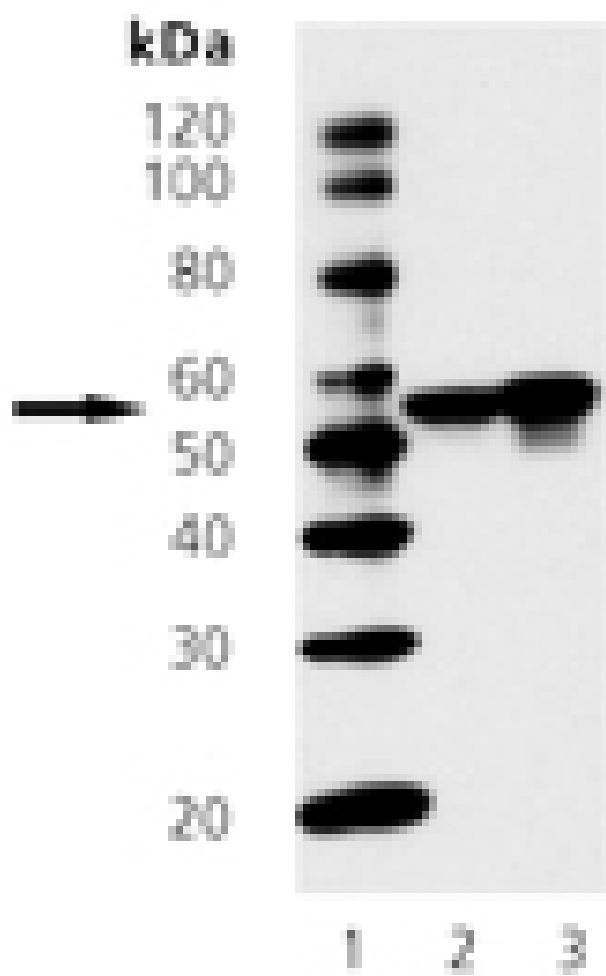
## Ordering Information

[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 Dilutions/Conditions	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	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)