

Jnk1 monoclonal antibody (1C2)

JNK1 is a member of the MAP kinase family. MAP kinases act as an integration point for multiple biochemical signals, and are involved in a wide variety of cellular processes such as proliferation, differentiation, transcription regulation and development. JNK1 is activated by various cell stimuli, and targets specific transcription factors, and thus mediates immediate-early gene expression in response to cell stimuli. The activation of JNK1 by tumor-necrosis factor alpha (TNF-alpha) is found to be required for TNF-alpha induced apoptosis. JNK1 is also involved in UV radiation induced apoptosis, which is thought to be related to cytochrom c-mediated cell death pathway. Studies of the mouse counterpart of JNK1 gene suggested that it play a key role in T cell proliferation, apoptosis and differentiation. Four alternatively spliced transcript variants encoding distinct isoforms have been reported.

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

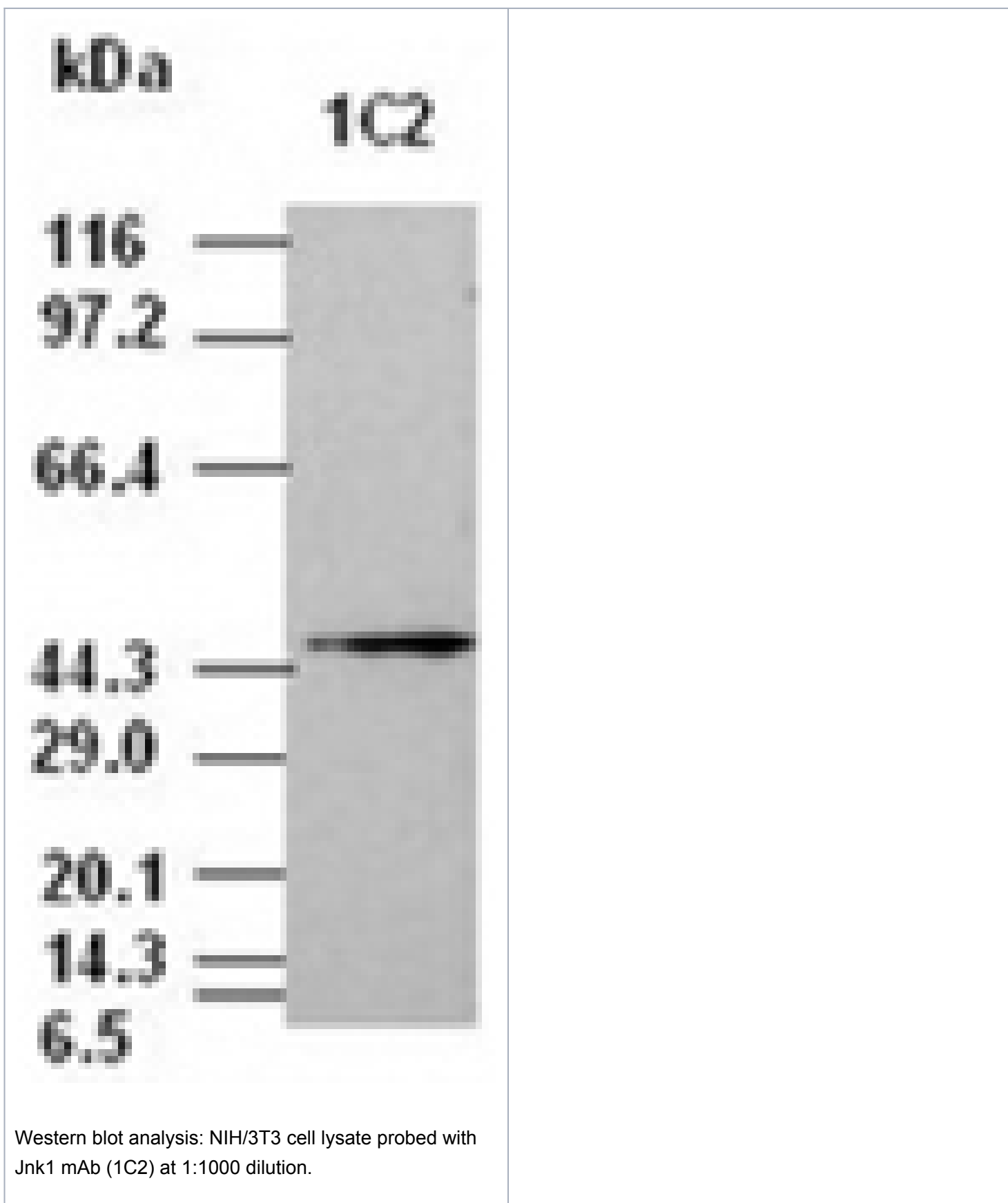
Ordering Information

[Order Online »](#)

ADI-905-862-100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name c-Jun N-terminal kinase 1, SAP kinase 1c, SAPK1c, MAP kinase 8, MAPK8

Application WB

Application Notes Detects a band of ~45kDa (full length) by Western blot.

Clone 1C2

Formulation Liquid. In PBS, pH 7.3, containing 1% BSA, 50% glycerol and 0.02% sodium azide.

Host Mouse

Immunogen Recombinant human Jnk1 (aa 1-384).

Isotype IgG2a

Purity Detail Protein G affinity purified.

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

Species Reactivity Human, Mouse

UniProt ID P45983

Worry-free Guarantee This antibody is covered by our [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-
fr@enzolifesciences.com](mailto:info-fr@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)