

# Jnk2 monoclonal antibody (12C5)

The Stress-activated protein kinases alpha and gamma (SAPKalpha and gamma), also known as c-Jun NH2-terminal kinases 1 and 2 (Jnk1/2) are serine/ threonine protein kinases belonging to the MAP kinase family. The extracellular stimuli leading to cell stress include inflammatory cytokines (via Rho and Rac), growth factors, protein synthesis inhibitors and osmotic stress amongst others. Both Jnk1 and Jnk2 must be dually phosphorylated in the activation loop, at threonine 183 and tyrosine 185, to achieve full enzymatic activity.

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

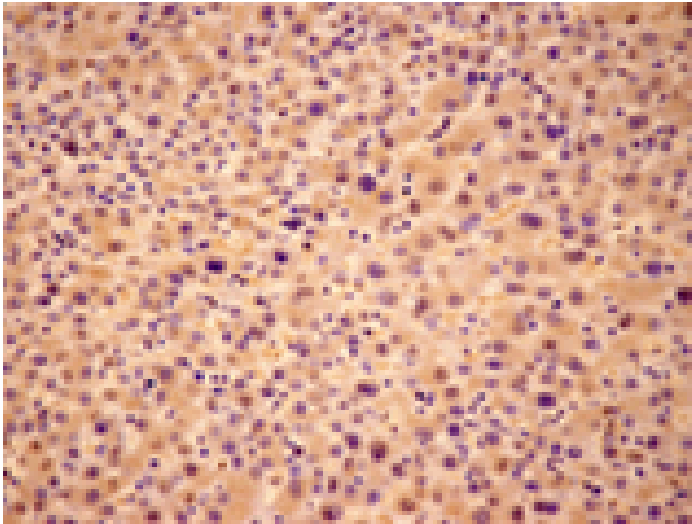
## Ordering Information

[Order Online »](#)

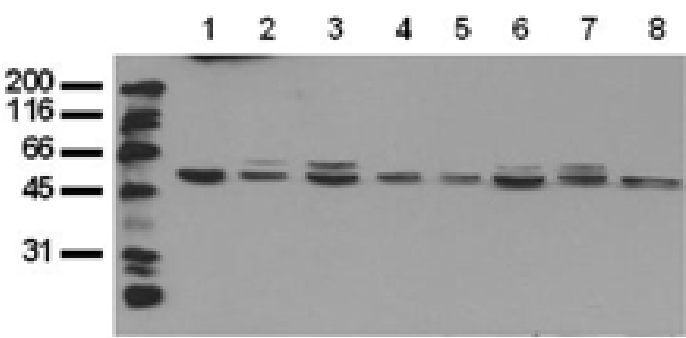
ADI-905-648-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human hepatoma tissue with Jnk2 mAb (12C5).



Western blot analysis of serum starved cancer cells A431 (1), A549 (2), SKOV3 (3), OVCAR5 (4), HaCaT (5), PC3 (6), HeLa (7), and HepG2 (8) probed with Jnk2 mAb (12C5).

# Handling & Storage

Use/Stability	Stable at -80°C up to 1 year, at 4°C up to 3 months.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	c-Jun N-terminal kinase 2, SAP kinase 1a, SAPK1a, MAP kinase 9, MAPK9
Application	ELISA, IHC, WB
Application Notes	Detects a band of ~54kDa by Western blot.
Clone	12C5
Formulation	Lyophilized from 1ml of 2x PBS containing 0.09% sodium azide, PEG, and sucrose.
Host	Mouse
Immunogen	Synthetic peptide corresponding to the sequence near the N-terminus of Jnk2.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	ELISA (0.05µg/ml)Immunohistochemistry (1:50)Western Blot (0.5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Reconstitution	Reconstitute with 1ml water (15 minutes at room temperature).
Species Reactivity	Dog, Human, Mouse

UniProt ID

P45984

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)