

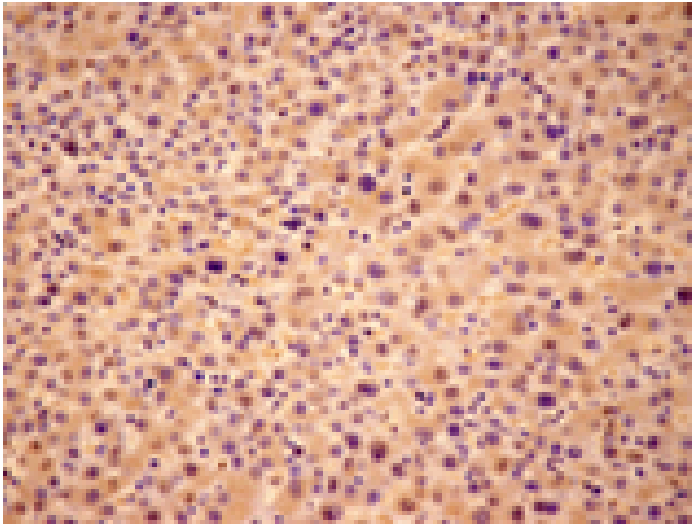
Jnk2 monoclonal antibody (12C5)

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.

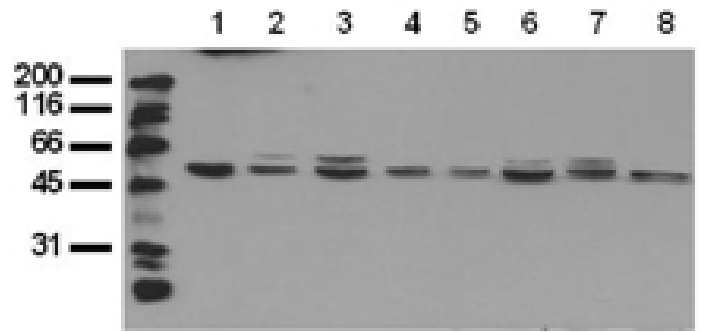
This antibody is covered by our [Worry-Free Guarantee](#).

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)