

# Jnk1 monoclonal antibody (5D10)

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](#).

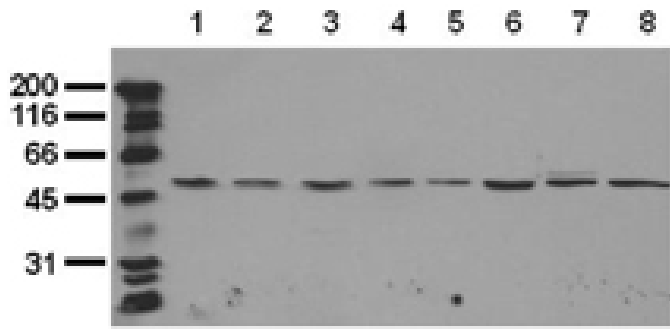
## Ordering Information

[Order Online »](#)

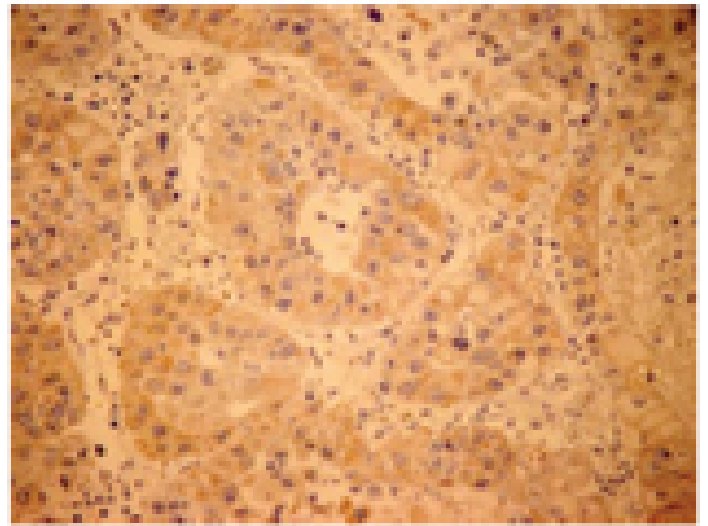
ADI-905-649-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



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 Jnk1 mAb (5D10).



Immunohistochemistry analysis of human hepatoma tissue with Jnk1 mAb (5D10).

## 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 1, SAP kinase 1c, SAPK1c, MAP kinase 8, MAPK8

**Application** ELISA, IHC, WB

**Application Notes** Detects a band of ~45kDa by Western blot.

**Clone** 5D10

**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 Jnk1.

**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, Rat

**UniProt ID** P45983



ENZO LIFE SCIENCES,

Worry Free Guarantee  
Phone: 800.942.0430

info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office

(ELS) AG  
Phone: +41 61 926 8989

info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands

Phone: +32 3 466 0420

info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France

Phone: +33 1 42 44 65 65

info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany

Phone: +49 7621 5500 526

info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland

Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900

info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)

**This antibody is covered by our Worry-Free Guarantee**