

Jnk3 monoclonal antibody (4G6)

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](#).

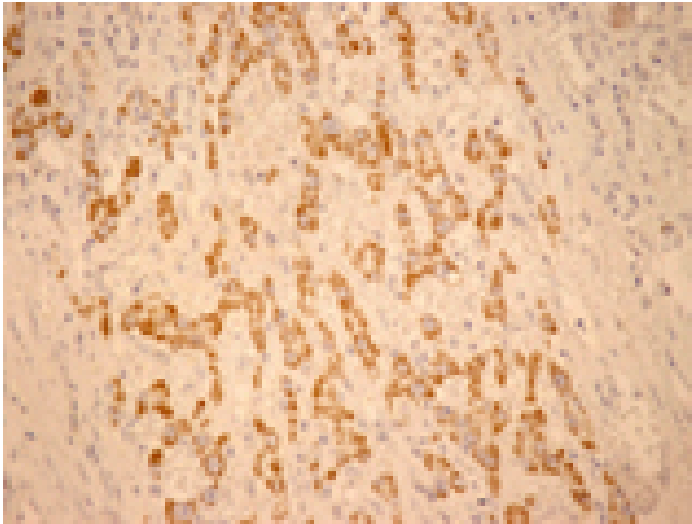
Ordering Information

[Order Online »](#)

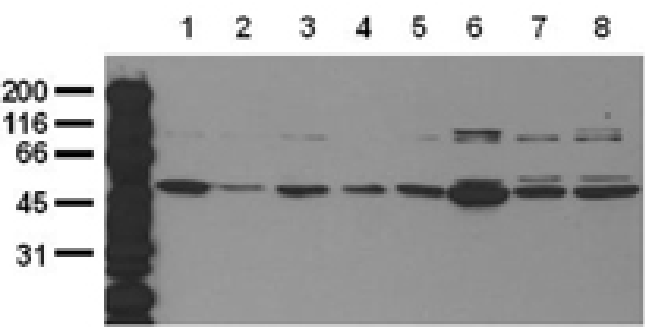
ADI-905-650-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human hepatoma tissue with Jnk3 mAb (4G6).



Western blot analysis of serum starved cancer cell A431 (1), A549 (2), SKOV3 (3), OVCAR5 (4), HaCaT (5), PC3 (6), HeLa (7), and HepG2 (8), probed with Jnk3 mAb (4G6).

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 3, SAP kinase 1b, SAPK1b, MAP kinase 10, MAPK10
Application	IHC, WB
Application Notes	Detects a band of ~54kDa by Western blot.
Clone	4G6
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 Jnk3.
Isotype	IgG1
Purity Detail	Thiophilic adsorption and size exclusion chromatography purified.
Recommendation Dilutions/Conditions	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	Human

UniProt ID

P53779

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com