

[pThr183/pTyr185]JNK1/2 polyclonal antibody

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA272-0020	20µl
BML-SA272-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	c-Jun N-terminal kinase 1/2, SAP kinase 1a/c, SAPK1a/c, MAP kinase 8/9, MAPK8/9
Application	IHC, WB
Crossreactivity	Does not cross-react with unphosphorylated JNK1&2.
Formulation	Liquid. In PBS containing BSA and sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human JNK1&2 phosphorylated at Thr183 and Tyr185. This region is conserved among many species, including mouse, rat, chicken, nematode and fly.
Species Reactivity	Human, Rat
Specificity	Recognizes JNK 1&2 dually phosphorylated at Thr ¹⁸³ and Tyr ¹⁸⁵ .
UniProt ID	P45983 (Jnk1), P45984 (Jnk2)
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