

[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 Thr¹⁸³ and Tyr¹⁸⁵.

UniProt ID P45983 (Jnk1), P45984 (Jnk2)

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



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)