

# [pThr183/pTyr185]JNK1/2 polyclonal antibody

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	c-Jun N-terminal kinase 1/2, SAP kinase 1a/c, SAPK1a/c, MAP kinase 8/9, MAPK8/9
<b>Application</b>	IHC, WB
<b>Crossreactivity</b>	Does not cross-react with unphosphorylated JNK1&2.
<b>Formulation</b>	Liquid. In PBS containing BSA and sodium azide.
<b>Host</b>	Rabbit
<b>Immunogen</b>	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.
<b>Species Reactivity</b>	Human, Rat
<b>Specificity</b>	Recognizes JNK 1&2 dually phosphorylated at Thr <sup>183</sup> and Tyr <sup>185</sup> .
<b>UniProt ID</b>	P45983 (Jnk1), P45984 (Jnk2)
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)