

# [pThr<sup>180</sup>/pTyr<sup>182</sup>]p38 polyclonal antibody

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

## Ordering Information

[Order Online »](#)

BML-SA266-0020	20µl
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application      WB

Crossreactivity      Does not cross-react with unphosphorylated forms of p38.

Formulation      Liquid. In PBS containing BSA and sodium azide.

Host      Rabbit

Immunogen      Synthetic peptide corresponding to a portion of human p38 phosphorylated at Thr180 and Tyr182.

Species Reactivity      Carp, Dog, Human, Mouse, Rat

Specificity      Recognizes p38 isoforms p38 $\alpha$ , p38 $\beta$ , and p38 $\gamma$  phosphorylated dually at Thr<sup>180</sup> and Tyr<sup>182</sup>.

UniProt ID      Q16539

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)