

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

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

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 Thr <sup>180</sup> and Tyr <sup>182</sup> .
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 <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)