[pThr¹⁸⁰/pTyr¹⁸²]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 p38α, p38β, and p38γ

phosphorylated dually at Thr¹⁸⁰ and Tyr¹⁸².

UniProt ID Q16539

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



uk@enzolifesciences.com