PI31 polyclonal antibody

Recombinant PI31 has been shown to compete with the proteasome regulators PA28 and 19S PA700 for binding to 20S, indicating that PI31 binds the 20S α -rings and may function by hindering substrate access to the 20S catalytic channel.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

Order Online »

BML-PW9710-0025	25µl
BML-PW9710-0100	100μΙ

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome inhibitor 31

Application WB

Formulation Liquid. In PBS containing 10mM sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 241-271 of C-

terminal of human PI31.

Species Reactivity Human, Mouse

UniProt ID Q92530

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

•

Last modified: May 29, 2024

