[pSer^{32/36}]lκBα polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA412-0020

20µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Formulation Liquid. In PBS, pH 7.3, containing glycerol, BSA and

sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human

IκBα phosphorylated at Ser32 and Ser36.

Species Reactivity Human, Mouse, Porcine

Specificity Recognizes IκBα phosphorylated at Ser^{32/36}.

UniProt ID P25963

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

.

Last modified: May 29, 2024

