[pThr²⁴⁶]PRAS40 (human) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA360-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proline-rich Akt substrate 40

Application WB

Formulation Liquid. In PBS, pH 7.3, containing BSA and sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human

PRAS40 (proline-rich Akt substrate 40) phosphorylated at

Thr246.

Species Reactivity Human

Specificity Recognizes PRAS40 phosphorylated at pThr²⁴⁶.

UniProt ID Q96B36

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

•

Last modified: May 29, 2024

