Proteasome 19S Rpn8/S12 subunit (human) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-PW8180-0025

25µl

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 26S proteasome non-ATPase regulatory subunit 7

Application IHC, WB

Formulation Liquid. In PBS containing 10mM sodium azide.

Host Sheep

Immunogen Recombinant human proteasome 26S (19S regulator

complex) (Rpn8/S12 subunit) (aa 1-205)

Species Reactivity Human

Specificity Recognizes Rpn8/S12 subunit of the 19S regulator

complex.

UniProt ID P51665

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

•

Last modified: May 29, 2024

