Proteasome 19S Rpn7/S10a subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

Order Online »

BML-PW8225-0025	25μΙ
BML-PW8225-0100	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 26S proteasome non-ATPase regulatory subunit 6, Breast

cancer-associated protein SGA-113M, Phosphonoformate

immuno-associated protein 4, Proteasome regulatory

particle subunit p44S10

Application IP, WB

Formulation Liquid. In PBS containing 10mM sodium azide.

GenBank ID D14663

Host Rabbit

Immunogen Recombinant full length human proteasome 26S (19S

regulator complex) (Rpn7/S10a subunit).

Species Reactivity Human, Yeast

Specificity Recognizes the Rpt7/S10a subunit of the 19S regulator

complex.

UniProt ID Q15008

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

.

Last modified: May 29, 2024

