## Proteasome 19S Rpt1/S7 subunit (yeast) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

BML-PW8255-0100

100µl

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** Dilute to working strength with phosphate buffered saline pH 7.2-7.4 and 1% normal

goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store diluted antibody

at 2-4°C and use within 1 month.

Handling Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name CIM5, 26S protease regulatory subunit 7 homolog

**Application** WB

**Application Notes**Detects a band of ~52kDa by Western blot.

**Formulation** Liquid. Neat serum.

GenBank ID 464862

**Host** Rabbit

Immunogen Recombinant protein corresponding to the N-terminal 100

amino acids of YTA3.

**Recommendation Dilutions/Conditions** Western Blot (1:10,000-1:50,000)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Yeast

Specificity Recognizes the Rpt1/S7 subunit of the 19S regulator

complex.

UniProt ID P33299

Last modified: May 29, 2024



info-

eu@enzolifesciences.com