## Proteasome 19S Rpt1/S7 subunit (human) monoclonal antibody (MSS1-92)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

**Ordering Information** 

Order Online »

BML-PW9400-0025

25µl

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name 26S protease regulatory subunit 7, Proteasome 26S

subunit ATPase 2, MSS1

**Application** IP

Clone MSS1-92

Formulation Liquid. In TRIS buffer containing 10mM sodium azide.

**Host** Mouse

Immunogen Recombinant human Rpt1 protein.

lsotype lgG1

Purity Detail Ammonium sulphate precipitation and ion-exchanging

chromatography purified.

**Source** Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

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

complex.

UniProt ID P35998

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com