## Proteasome 20S α6 subunit monoclonal antibody (MCP106)

Manuals, SDS & CofA

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

Ordering Information

Order Online »

BML-PW9390-0001

1ml

**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** PROS30, Proteasome subunit α type-1, Macropain subunit

C2, Proteasome v chain

**Application** WB

Clone MCP106

**Formulation** Liquid. Tissue culture supernatant containing 10mM

sodium azide.

Gene/Protein Identifier PSMA1 (gene name)

Host Mouse

Human placenta-derived proteasomes. **Immunogen** 

Isotype lgG1

Source Purified from hybridoma tissue culture supernatant.

Human, Mouse, Rabbit **Species Reactivity** 

**Technical Info / Product Notes** Various systems for the nomenclature of the proteasome

> subunits have been established. This may be a source of confusion as the system on UniProt differs from "standard" nomenclature as described in the literature. The UniProt ID and Gene Name will help to clearly identify the proteins.

**UniProt ID** P25786

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com