Proteasome 20S α5 subunit monoclonal antibody (MCP196)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

Ordering Information

Order Online »

BML-PW8125-0100

100µl

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 Proteasome subunit α type-5, Macropain ζ chain

Application IHC, WB

Clone MCP196

Formulation Liquid. In PBS containing 10mM sodium azide.

Gene/Protein Identifier PSMA5 (gene name)

Host Mouse

Immunogen Dinitrophenylated human placenta derived proteasomes.

Isotype lgG2a

Purity Detail Partially purified ascites.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rabbit, Rat

Specificity Recognizes the $\alpha 5$ subunit of the 20S proteasome.

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 P28066

Last modified: May 29, 2024



info-

eu@enzolifesciences.com