Proteasome 20S β7 subunit (human) monoclonal antibody (MCP444)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-PW9150-0025	25µl
BML-PW9150-0100	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 β type-4, PROS-26, Macropain β

chain

Application IP, WB

Clone MCP444

Formulation Liquid. Ascites in PBS, containing 0.01 M sodium azide.

Gene/Protein Identifier PSMB4 (gene name)

Host Mouse

Immunogen Human placenta derived proteasomes.

lgG1

Purity Detail Purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

Specificity Recognizes the β7 subunit of the 20S proteasome.

Technical Info / Product NotesVarious 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.

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com