Proteasome 20S β2 subunit monoclonal antibody (MCP168)

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.

This antibody is covered by our Worry-Free Guarantee.

Citations: 15

View Online »

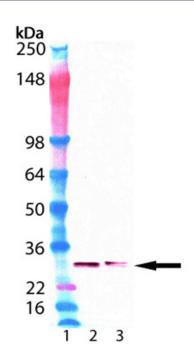
Ordering Information

Order Online »

BML-PW8145-0025	25µl
BML-PW8145-0100	100µl

Manuals, SDS & CofA

View Online »



Western Blot analysis of BML-PW8145: Lane 1: SeeBlueTMPlus2 MW Marker, Lane 2: HeLa cell lysate (Prod. No. ADI-LYC-HL100), Lane 3: Jurkat cell lysate (Prod. No. ADI-LYC-JK100).

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome subunit β type-7, Macropain chain Z

Application IHC, WB

Clone MCP168

Formulation Liquid. In PBS containing 10mM sodium azide.

Gene/Protein Identifier PSMB7 (gene name)

Host Mouse

Immunogen Dinitrophenylated proteasomes.

lsotype lgG1

Purity Detail Protein G-affinity purified ascites.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Yeast

Specificity Recognizes the β2 subunit of proteasome 20S.

UniProt ID Q99436

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com

Last modified: May 29, 2024