Proteasome 20S β2 subunit monoclonal antibody (MCP165)

The hybridoma secreting the antibody to subunit $\beta 2$ was generated by fusion of splenocytes from Balb/c mice which had received repeated immunization with dinitrophenylated proteasomes. The antibody (clone MCP165) has been characterised by one- and two dimensional Western blotting. The antibody has not been demonstrated to immunoprecipitate the native 20S proteasome. The immunoglobulin subclass is lgG_1 .

This antibody is covered by our Worry-Free Guarantee.

Citations: 11

View Online »

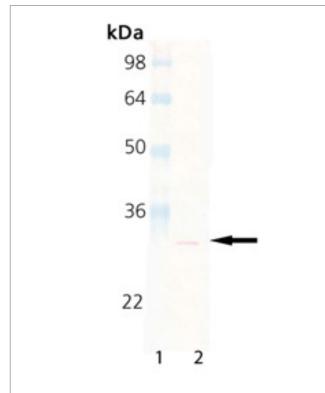
Ordering Information

Order Online »

BML-PW9300-0025	25µI
BML-PW9300-0100	100μΙ

Manuals, SDS & CofA

View Online »



Western Blot Analysis of Proteasome 20S β2 subunit, mAb (MCP165), (Prod. No. BML-PW9300): Lane 1: MW Marker, Lane 2: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100)

Handling & Storage

Use/Stability Store unopened vial at -20°C until required for use. Aliquot undiluted antibody into

smaller volumes (not less than 10µl) prior to freezing if appropriate. The use of high quality 'antiserum-grade' plastic or glass vials is recommended. Store diluted antibody at

2-4°C (do not freeze) and use within 1 month.

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-7, Macropain chain Z

Application IHC, WB

Application Notes Immunoprecipitation: Does not immunoprecipitate intact

proteasomes from HeLa cell preparations.

Immunohistochemistry: Has been shown to have

applicability in immunohistochemical applications using

human tissue.

Western blot: Single dimension SDS-PAGE of a HeLa cell lysate followed by Western blotting gives a single band

with a relative molecular weight of ~28kDa.

Clone MCP165

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

Gene/Protein Identifier PSMB7 (gene name)

Host Mouse

Immunogen Dinitrophenylated proteasomes.

lsotype lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions

Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source

Purified from hybridoma tissue culture supernatant.

Species Reactivity

Human, Mouse

Specificity

Recognizes the β2 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

Q99436 (human), P70195 (mouse)

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

