

Proteasome 20S β 5t subunit (human) monoclonal antibody (b8-17)

When cells are stimulated with pro-inflammatory cytokines, most of their constitutively expressed proteasomes are replaced with immunoproteasomes, which increase the production of peptides for presentation on MHC class I molecules. In addition, cortical thymic epithelial cells selectively express a type of proteasome known as the thymoproteasome, in which the β 5i immunoproteasome subunit has been replaced by the thymus specific proteasome subunit β 5t. The thymoproteasome is required for the positive selection of thymocytes and shaping the T cell receptor repertoire of cytotoxic T lymphocytes. In addition mice deficient in β 5t display defective development of CD8(+) T cells in the thymus suggesting a key role for β 5t and the thymoproteasome in generating MHC class I-restricted CD8(+) T cells.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BML-PW1020-0025	25 μ l
BML-PW1020-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)

MW (kDa)

40 —

36 —

21 —



← Precursor

← β5t

Western blot showing detection of 20S proteasome subunit β5t in C57/BL6 mouse thymus lysate by BML-PW1020 at 1:1000 dilution.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Proteasome subunit β type-11
Application	WB
Clone	b8-17
Formulation	Liquid. In PBS containing 0.01% sodium azide.
Gene/Protein Identifier	PSMB11 (gene name)
Host	Mouse
Immunogen	Recombinant human 20S proteasome β 5t subunit.
Isotype	IgG1
Purity Detail	Ion exchange purified.
Recommendation Dilutions/Conditions	Western Blot (1:500-1:1,000)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 β 5t subunit of the 20S proteasome.

Technical Info / Product Notes	The hybridoma secreting the antibody clone b8-17 was generated by fusion of splenocytes from Balb/cmice which had received repeated immunisation with recombinant human β 5t subunit protein (accession number: A5LHX3). Immunoglobulin was purified from tissue culture supernatant by ion exchange chromatography.
UniProt ID	A5LHX3
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com