

Proteasome 20S β 5 subunit (human) polyclonal antibody

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

Citations: 56

[View Online »](#)

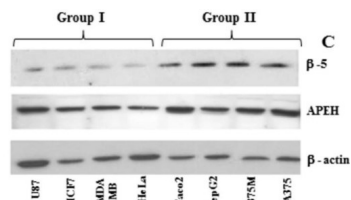
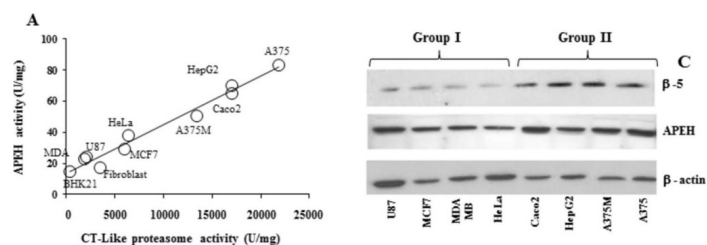
Ordering Information

[Order Online »](#)

BML-PW8895-0025	25 μ l
BML-PW8895-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Human cancer cell lines may be grouped according to the basal enzyme activities and expression levels of APEH and proteasome. Cells from eight human cancer lines (U87, MCF7, MDA-MB, HeLa, Caco2, HepG2, A375M, A375) and non-cancerous cells (BHK21, fibroblasts) were harvested at the pre-confluent stage and used for cytoplasmic or mRNA extract preparation. Basal APEH and proteasomal CT-like activities were measured in cytoplasmic extracts (A). The mRNA levels of APEH and β -5 subunit were evaluated by qRT-PCR and expressed as fold change in comparison to expressed levels in human fibroblast (B). Intracellular levels of β -5 and APEH were detected by immunoblotting (C upper panel). Typical Western blot was shown and data from three different analyses were normalized to the density of control protein (β -actin) and expressed as ratio over control (C lower panel). Results are presented as the mean values \pm SD of triplicate analyses from at least three different experiments.

*Significantly different ($P < 0.01$) from respective controls.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: RedOx status, proteasome and APEH: insights into anticancer mechanisms of t10,c12-conjugated linoleic acid isomer on A375 melanoma cells. *PLoS One* (2013)

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-5, Macropain ϵ chain, Proteasome chain 6, Proteasome subunit MB1, Proteasome subunit X
Application	IHC, WB
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Gene/Protein Identifier	PSMB5 (gene name)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 194-207 of human proteasome 20S (β 5 subunit).
Source	Purified from rabbit serum.
Species Reactivity	Human
Specificity	Recognizes human proteasome 20S β 5 subunit.
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	P28074
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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