

Blm10

(*Saccharomyces cerevisiae*) polyclonal antibody

Blm10 (mammalian homolog PA200) is a large 246kDa regulatory protein that binds to the 20S proteasome core particle (CP). Blm10 recognizes pre-activated open gate CP which is assumed to exist in equilibrium with inactive closed gate CP. Blm10 distinguishes gate conformations, regulates CP activation, and has been shown to be involved in CP maturation.

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

Citations: 5

[View Online »](#)

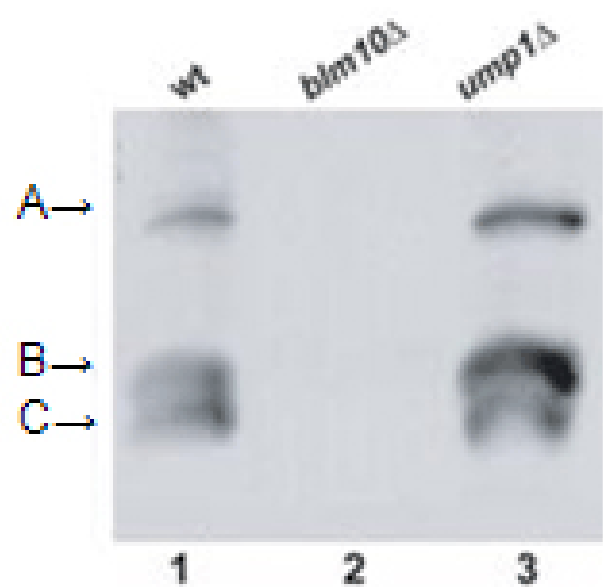
Ordering Information

[Order Online »](#)

BML-PW0570-0025	25µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C. Dilute with PBS, pH 7.2-7.4, containing 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store diluted antibody at +4°C, do not freeze and use within 1 month.
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	Blm3, PA200 homolog
Application	WB
Formulation	Liquid. Neat serum diluted in PBS containing 0.01% sodium azide.
Host	Rabbit
Immunogen	Recombinant N-terminal protein fragment (aa 1-157) of <i>Saccharomyces cerevisiae</i> Blm10.
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.
Species Reactivity	<i>Saccharomyces cerevisiae</i>
UniProt ID	P43583

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-fr@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

Last modified: May 29, 2024