

# Proteasome activator 11S $\beta$ subunit polyclonal antibody

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW8240-0025	25 $\mu$ l
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Proteasome activator complex subunit 2, PA28 $\beta$
Application	IHC, WB
Formulation	Liquid. Antiserum containing 10mM sodium azide.
GenBank ID	2351198 (mouse)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 15-31 of mouse PA28 $\beta$ .
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes the $\beta$ subunit of proteasome activator 11S.
UniProt ID	P97372 (mouse), Q9UL46 (human)
Worry-free Guarantee	<b>This antibody is covered by our <a href="#">Worry-Free Guarantee</a></b>

Last modified: May 29, 2024



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)