

# 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
BML-PW8240-0100	100 $\mu$ 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, PA28β

**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 PA28β.

**Species Reactivity**      Human, Mouse, Rat

**Specificity**      Recognizes the β subunit of proteasome activator 11S.

**UniProt ID**      P97372 (mouse), Q9UL46 (human)

**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)