

---

## PSMB2 Polyclonal Antibody

(Catalog # A52680)

---

### Background

The proteasome is a multicatalytic proteinase complex with a highly ordered ring shaped 20S core structure. The core structure is composed of 4 rings of 28 non identical subunits; 2 rings are composed of 7 alpha subunits and 2 rings are composed of 7 beta subunits. Proteasomes are distributed throughout eukaryotic cells at a high concentration and cleave peptides in an ATP/ubiquitin dependent process in a non-lysosomal pathway. An essential function of a modified proteasome, the immunoproteasome, is the processing of class I MHC peptides. PSMB2 is a member of the proteasome B-type family, also known as the T1B family, that is a 20S core beta subunit. The proteasome is a multicatalytic proteinase complex with a highly ordered ring shaped 20S core structure. The core structure is composed of 4 rings of 28 non identical subunits; 2 rings are composed of 7 alpha subunits and 2 rings are composed of 7 beta subunits. Proteasomes are distributed throughout eukaryotic cells at a high concentration and cleave peptides in an ATP/ubiquitin dependent process in a non-lysosomal pathway. An essential function of a modified proteasome, the immunoproteasome, is the processing of class I MHC peptides. This gene encodes a member of the proteasome B type family, also known as the T1B family, that is a 20S core beta subunit. Publication Note: This RefSeq record includes a subset of the publications that are available for this gene. Please see the Entrez Gene record to access additional publications.

### Description

PSMB2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300. 50% Glycerol, 0.01M PBS, pH 7.4.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

P49721

### Purification

Protein G purified

### Immunogen

Recombinant Human Proteasome subunit beta type-2 protein (1-201AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

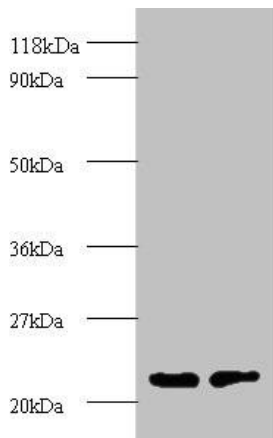
### Alternative Names

Macropain subunit C7-I, Multicatalytic endopeptidase complex subunit C7-I, Proteasome component C7-I

### Application

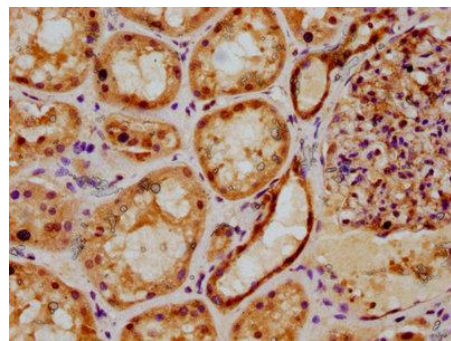
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

---

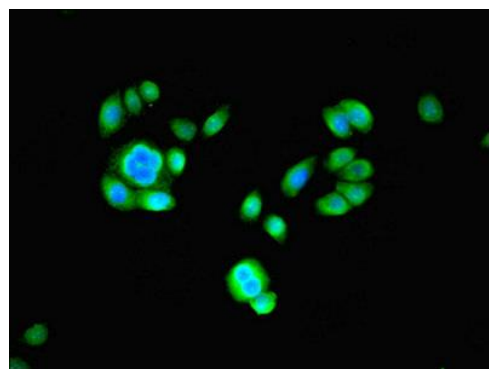


#### Western blot

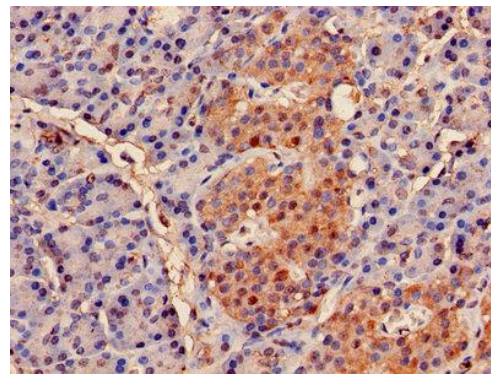
All lanes: PSMB2 Polyclonal Antibody at 2ug/ml  
 Lane 1: EC109 whole cell lysate  
 Lane 2: 293T whole cell lysate  
 Secondary  
 Goat polyclonal to Rabbit IgG at 1/15000 dilution  
 Predicted band size: 22.9kDa  
 Observed band size: 22.9kDa



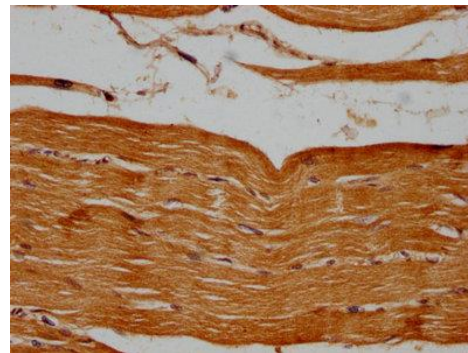
IHC image of PSMB2 Polyclonal Antibody diluted at 1:500 and staining in paraffin-embedded human kidney tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescent analysis of PC3 cells using PSMB2 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunohistochemistry of paraffin-embedded human pancreatic tissue using PSMB2 Polyclonal Antibody at dilution of 1:100



IHC image of PSMB2 Polyclonal Antibody diluted at 1:500 and staining in paraffin-embedded human skeletal muscle tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.