## Proteasome 20S β2i subunit polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

BML-PW8210-0100

100µl

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 subunit beta type-10, Macropain chain

subunit MECI-1

**Application** IHC, WB

**Formulation** Liquid. Antiserum containing 10mM sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to aa 44-73 of human

proteasome 20S (β2i subunit).

Species Reactivity Human

**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 P40306

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

\_

Last modified: May 29, 2024

