

Proteasome 20S α 2 subunit monoclonal antibody (MCP21) (agarose immobilized)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW8335-0500	0.5ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Store unopened vial at 4°C until required for use. DO NOT FREEZE.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome subunit α type-2, Macropain subunit C3

Application WB

Clone MCP21

Formulation Liquid. In PBS containing 0.01M sodium azide.

Gene/Protein Identifier PSMA2 (gene name)

Host Mouse

Immunogen Human placenta derived proteasomes.

Isotype IgG1

Source Purified from mouse ascites.

Species Reactivity Bovine, Human, Rabbit

Specificity Recognizes the $\alpha 2$ subunit of the 20S proteasome.

UniProt ID P25787

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

.