Proteasome 20S α2 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.

lsotype lgG1

Source Purified from mouse ascites.

Species Reactivity Bovine, Human, Rabbit

Specificity Recognizes the α2 subunit of the 20S proteasome.

UniProt ID P25787

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

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com