Proteasome 19S Rpn12/S14 subunit (human) monoclonal antibody (P31-27)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-PW8835-0025	25µl
BML-PW8835-0100	100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 26S proteasome non-ATPase regulatory subunit 8

Application IHC, WB

Clone p31-27

Formulation Liquid. Containing 10mM sodium azide.

Gene/Protein Identifier PSMD8 (gene name)

Host Mouse

Immunogen Recombinant Rpn12 protein.

lsotype lgG1

Purity Detail Partially purified.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

Specificity Recognizes human proteasome 26S (19S regulator

complex) (Rpn12/S14 subunit).

UniProt ID P48556

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com