

# Proteasome 19S

## Rpn10/S5a subunit

### (human),

### (recombinant) (GST-tag)

A subunit of the 26S proteasome, S5a, recognises and binds multi-ubiquitinated proteins containing chains of at least four ubiquitin moieties. Purified recombinant S5a retains the ability to bind multi-ubiquitinated proteins in isolation, when immobilised on glutathione-Sepharose, and after being eluted from glutathione-Sepharose, run on SDS-PAGE, and blotted onto nitrocellulose. Human S5a sequence, with an N-terminal fusion of glutathione S-transferase (*Schistosoma japonicum*), was expressed in *E. coli* with GST tag at amino terminus.

S5a (RPN10), a subunit of the 19S regulator of the 26S proteasome, binds multi-ubiquitinated proteins containing chains of at least four ubiquitin moieties.

Citations: 1

[View Online »](#)

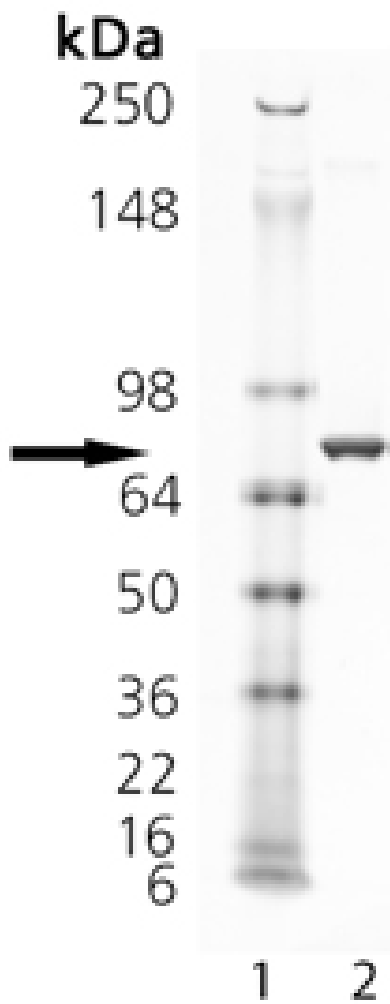
### Ordering Information

[Order Online »](#)

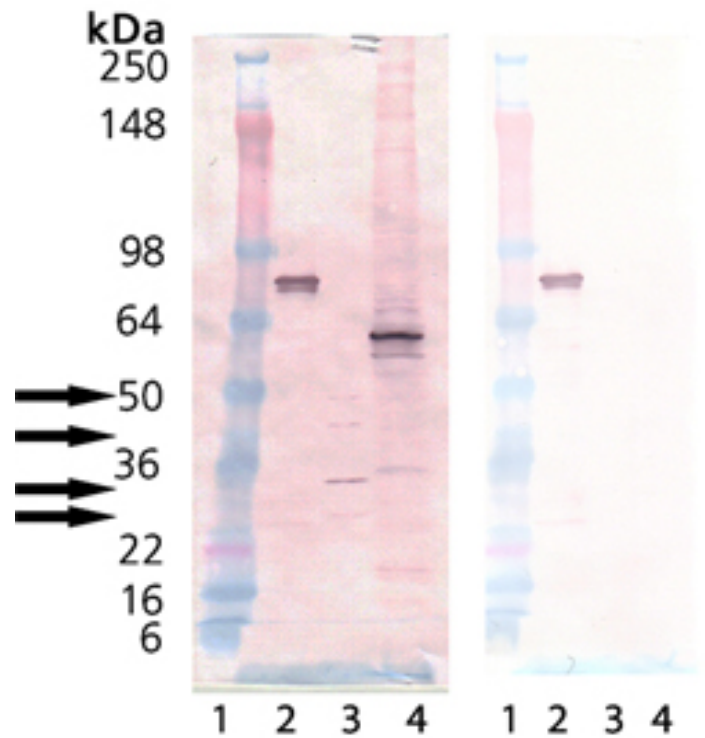
BML-UW8465-0100	100µg
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis of Rpn10/S5a. Lane 1: MW Marker, Lane 2: 1 µg.



Western Blot Analysis of Rpn10/S5a (Prod. No. BML-UW8465). Lane 1: MW Marker, Lane 2: Rpn10/S5a, Lane 3: Multiubiquitin chains (Prod. No. BML-UW8860), Lane 4: HeLa lysate (Prod. No. ADI-LYC-HL100). The blot on the left was probed with 10 µg/ml Rpn10, right with blocking buffer only. Rpn10 was detected by probing with GST polyclonal antibodies.

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Activity** Bound to mutliubiquitin chains of length >4 when conjugated to agarose (Prod. No. BML-UW8635).

**Alternative Name** 26S proteasome non-ATPase regulatory subunit 4, Antisecretory factor 1, Multiubiquitin chain-binding protein

**Appearance** Colorless liquid.

**Application Notes** Useful for the isolation of multi-ubiquitinylated proteins.

**Formulation** Liquid. In 20mM TRIS, pH 7.5, containing 500mM sodium chloride.

**MW** ~78kDa (observed)

**Purity** ≥95% (SDS-PAGE)

**Purity Detail** Purified by multi-step chromatography.

**Source** Produced in *E. coli*.

**UniProt ID** P55036



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)