

S5a (human) polyclonal antibody

S5a/Rpn10 is a major ubiquitin binding protein that binds preferentially to polyubiquitin chains. It is found as a non-ATPase subunit of the 19S regulatory complex of the 26S proteasome where it functions as a receptor for proteins destined for proteolytic degradation. In yeast, the homologue of S5a, Rpn10, is required for degradation of a subset of cellular proteins by the proteasome. Unlike other proteasome subunits, S5a exists predominantly as a free protein in the cytosol (i.e. not incorporated into the proteasome), though its precise role in the cytosol has yet to be fully elucidated. S5a contains two approximately 15 amino acid ubiquitin interacting motifs (UIM) which are responsible for its binding to polyubiquitin chains.

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

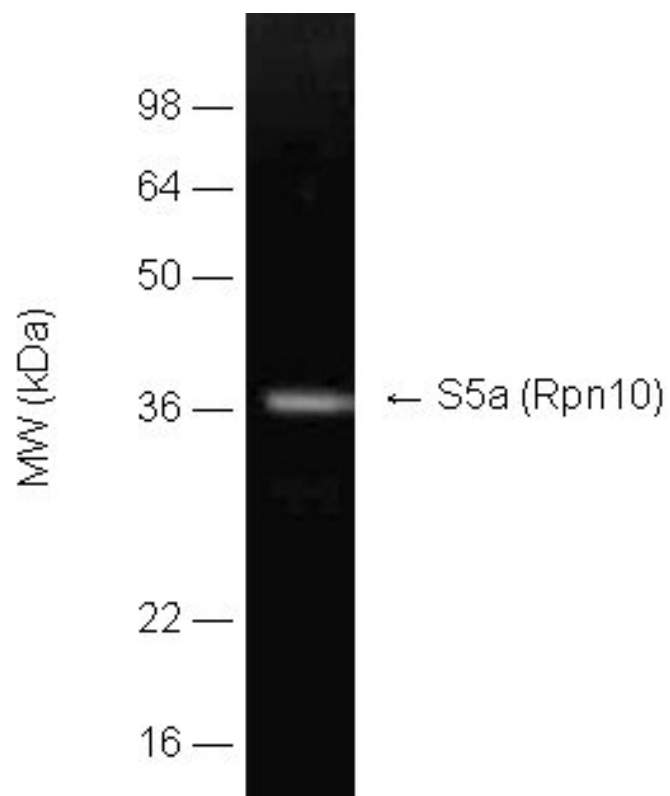
Ordering Information

[Order Online »](#)

BML-PW1045-0025	25µl
BML-PW1045-0100	100µl

Manuals, SDS & CofA

[View Online »](#)



Detection of S5a (Rpn10) 19S proteasome regulator subunit in purified 26S proteasome sample (50ng, BML-PW9310) by Western blotting using BML-PW1045 at 1:1000 dilution.

Handling & Storage

Handling	Store unopened vial at -20°C until required for use. Store diluted antibody at 2-4°C and use within 1 month. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Rpn10, MCB1, PSMD4, 26S proteasome regulatory subunit S5A, Mbp1, Sun1, p54, Antisecretory factor 1 (ASF1)
Application	WB
Formulation	Liquid. In PBS containing 0.01% sodium azide.
Host	Rabbit
Immunogen	Full-length recombinant GST-tagged human S5a (Accession number: P55036).
Purity Detail	Purified by immunodepletion of GST-reactive antibodies followed by Protein A purification.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Human
Specificity	Recognizes the Rpn10/S5a subunit of the 19S regulator complex.
UniProt ID	P55036



Worry-Free Guarantee

ENZO LIFE SCIENCES

Phone: 800.942.0430
info-
usa@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

European Sales Office
ENZO LIFE SCIENCES

(ELS) AG
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg

Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France

Phone: +33 1 72 440 1335
info-
fr@enzolifesciences.com

Germany

Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland

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