

Proteasome 19S

Rpn10/S5a subunit

(human),

(recombinant) (GST-tag) (agarose immobilized)

Rpn10/S5a, a subunit of the 19S regulator of the 26S proteasome, binds multi-ubiquitylated proteins containing chains of at least four ubiquitin moieties. Rpn10/S5a-agarose retains this ability and may be used to identify and purify multi-ubiquitylated proteins.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW8635-0500	0.5ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Formulation Liquid. The resin is supplied as a slurry in 50mM TRIS, pH 7.5, containing 150mM sodium chloride, 50% glycerol, 1mM DTT and 0.01% sodium azide.

Purity ≥95% (SDS-PAGE)

Source GST-S5a (human Rpn10) is covalently coupled to agarose at a concentration of 0.5mg protein/ml of settled resin.

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-
be@enzolifesciences.com](mailto:info-be@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)