

# 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 »](#)

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

<b>Alternative Name</b>	26S proteasome non-ATPase regulatory subunit 4, Antisecretory factor 1, Multiubiquitin chain-binding protein
<b>Formulation</b>	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.
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Source</b>	GST-S5a (human Rpn10) is covalently coupled to agarose at a concentration of 0.5mg protein/ml of settled resin.
<b>UniProt ID</b>	P55036

**Last modified: May 29, 2024**



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)