Proteasome 19S Rpn10/S5a subunit (human), (recombinant) (GST-

the 26S proteasome, S5a, recognises and binds multi-ubiquitinylated proteins containing chains of at least four ubiquitin moieties. Purified recombinant S5a retains the ability to bind multi-ubiquitinylated 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-ubiquitinylated proteins containing chains of at least four ubiquitin moieties.

Citations: 1

View Online »

Ordering Information

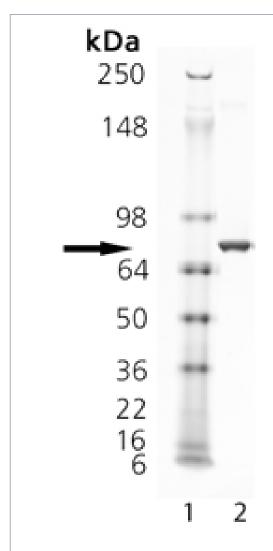
Order Online »

BML-UW8465-0100

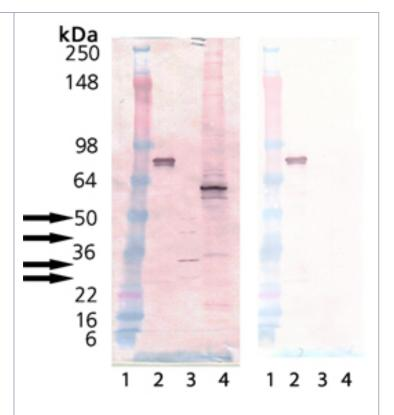
100µg

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 μ g/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 NotesUseful 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

Last modified: May 29, 2024

