

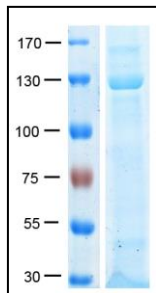
Rpn2 Protein (26S Proteasome Regulatory Subunit 2)

Cat. # PS120

Background: The 26S proteasome is the central protease in ubiquitin-mediated degradation of cellular proteins. It consists of a 20S catalytic core capped by one or two 19S regulatory particles. Rpn1, Rpn2, and Rpn10 are located at the base of the 19S. Rpn1 and Rpn2 are the largest proteasomal subunits (110 and 104 kDa, respectively). Deletion of either subunit is lethal, and mutations in yeast result in impaired proteasome function, accumulation of polyubiquitinated proteins, and improper nuclear proteasome localization. Recent findings based on the silencing of Rpn2 in cancer cells indicate that this protein is a potential key therapeutic target for drug-resistant tumors.

Product Information

Quantity:	50 µg
Molecular weight:	104.2kDa
Tag:	His ₆
Buffer:	20 mM sodium phosphate pH 7.4, 500 mM NaCl, 200 mM imidazole, 1mM β – mercaptoethanol.
Source:	Yeast recombinant, expressed in <i>E. coli</i> . Accession number P32565.
Storage:	-80°C. Avoid repeated freeze/thaw cycles



Recombinant Rpn2 protein (His₆ tag) was expressed in *E. coli* and purified on the Ni-NTA resin using a standard protocol. 5µg of purified protein was resolved on 10% PAGE and stained with Coomassie blue.

References

1. Saeki, Y. et al (2002) Identification of ubiquitin-like protein-binding subunits of the 26S proteasome. *Biochem. Biophys. Res. Commun.*
2. Wendler P. et al. (2004) The bipartite nuclear localization sequence of Rpn2 is required for nuclear import of proteasomal base complexes via karyopherin alpha and proteasome functions. *J Biol Chem* 279(36):37751-62
3. Kimura Y. et al. (2003) N-Terminal modifications of the 19S regulatory particle subunits of the yeast proteasome. *Arch Biochem Biophys* 409(2):341-8.
4. Gomez TA. et al (2011) Identification of a functional docking site in the Rpn1 LRR domain for the UBA-UBL domain protein Ddi1. *BMC Biol.* May 31;9:33.
5. He J, et al. (2012) The structure of the 26S proteasome subunit Rpn2 reveals its PC repeat domain as a closed toroid of two concentric α -helical rings. *Structure* 20(3):513-21.
6. Rosenzweig R. et al (2012) Rpn1 and Rpn2 coordinate ubiquitin processing factors at proteasome. *J Biol Chem.* Apr 27;287(18):14659-71.
7. Huber EM and Groll M. (2012) The 19S cap puzzle: a new jigsaw piece. *Structure.* 20(3):387-8
8. Lander GC. et al. (2012) Complete subunit architecture of the proteasome regulatory particle. *Nature.* Jan 11;482(7384):186-91.