

COP9 signalosome

Csn6 subunit (human)

polyclonal antibody

The polyclonal antibody to Csn6 was generated by immunisation of rabbits with a synthetic peptide corresponding to residues [283-297] of the human Csn6 sequence. The antibody has been characterised by one-dimensional Western blotting. Vial contains a partially purified immunoglobulin preparation suspended phosphate-buffered saline containing 0.01M sodium azide. The COP9 signalosome, once defined as a repressor complex of light activated development in Arabidopsis, has recently been found in humans and is probably present in most multicellular organisms. The COP9 signalosome is closely related to the lid sub-complex of the 26S proteasome in structural composition and probably shares a common evolutionary ancestor. A multifaceted role of the COP9 signalosome in cell-signalling processes is hinted at by its associated novel kinase activity, as well as the involvement of its subunits in regulating multiple cell-signalling pathways and cell-cycle progression. The molecular genetic studies in Arabidopsis suggest that the complex functions as part of a highly conserved regulatory network, whose physiological rôle in animals has yet to be determined. The COP9 signalosome, also known as the the COP9 complex and JAB1-containing signalosome, is a conserved nuclear protein complex found in plants and animals. It is composed of eight distinct subunits, designated S1 to S8 (very recently re-designated Csn1-Csn8) according to the molecular weight of the mammalian complex subunits. Subunit composition and the subunit sequences are substantially conserved between the mammalian and the plant complexes, implying that the complex has a conserved cellular function in higher eukaryotic organisms. The role and function of Csn6 has yet to be determined although this subunit is known to interact with subunits Csn2 and Csn7 within the COP9 signalosome.

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

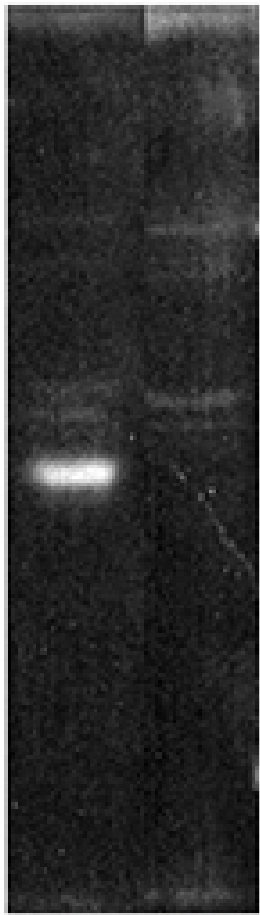
BML-PW8295-0025

25µl

BML-PW8295-0100	100µl
------------------------	-------

Manuals, SDS & CofA

[View Online »](#)



(a) **(b)**

Western blot analysis: Luminograph of a HeLa S3 cytosol preparation after SDS PAGE followed by blotting onto PVDF membrane and probing with (a) antibody BML-PW8295 and (b) antibody BML-PW8295 pre-absorbed with cognate peptide.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application Notes	Western Blot: The antibody has been characterised against both endogenous and purified signalosome preparations. Detects a band of ~33.5kDa.
--------------------------	--

Crossreactivity	Does not cross-react with other signalosome subunits.
------------------------	---

Formulation	Liquid. In PBS containing 10mM sodium azide.
--------------------	--

Host	Rabbit
------	--------

Immunogen	Synthetic peptide corresponding to aa 283-297 of human COP9 signalosome (Csn6 subunit).
------------------	---

Purity Detail	Partially purified immunoglobulin.
----------------------	------------------------------------

Recommendation Dilutions/Conditions	Western blot (an initial dilution of >1:1000 is recommended)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 Csn6 subunit of the COP9 signalosome.
--------------------	--

Technical Info / Product Notes

Immunoblotting – The antibody has been characterised against both endogenous and purified signalosome preparations. The antibody does not cross-react with other signalosome subunits. An initial dilution of >1:1000 is recommended.

Immunoprecipitation – No information currently available.

Immunohistochemistry – No information currently available.

UniProt ID

Q7L5N1

Worry-free Guarantee

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

.

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)