

# CSN8 (CT) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat CSN8 (COP9 signalosome complex subunit 8)

### Instruction Manual

Catalog Number	PK-AB718-4601
Synonyms	CSN8 Antibody; COP9 signalosome complex subunit 8, SGN8, JAB1-containing signalosome subunit 8
Description	The COP9 signalosome (CSN) is an evolutionarily conserved protein complex of the eight subunits that interacts with deubiquitinating enzymes and protein kinases and is highly homologous to the lid sub-complex of 26S proteasome. The CSN complex is an essential regulator of the ubiquitin conjugation pathway by mediating the deneddylation of the SCF-type E3 ligase complexes, which leads to a decrease in ubiquitin ligase activity of SCF-complexes such as SCF, CSA or DDB2. It is also involved in phosphorylation of p53, c-jun/JUN, ITPK1 and IRF8/ICSBP, possibly via its association with CK2 and PKD kinases. CSN8 encodes the smallest and the least conserved but first identified subunit of CSN. Recent studies show CSN8 is essential for Drosophila development and is essential for peripheral T cell homeostasis and antigen receptor-induced entry into the cell cycle from quiescence.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CSN8 antibody was raised in rabbits against a 15 amino acid peptide near the carboxy terminus of the human CSN8.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. CSN8 antibody can be used for detection of CSN8 by Western blot at 2 µg/mL.
Images	Available upon request.
References	Wei N and Deng XW. The COP9 signalosome. Annu. Rev. Cell Dev. Biol. 2003; 19:261-86. Bech-Otschir D, Seeger M, and Dubiel W. The COP9 signalosome: at the interface between signal transduction and ubiquitin-dependent proteolysis. J. Cell Sci. 2002; 115:467-73. Groisman R, Polanowska J, Kuraoka I, et al. The ubiquitin ligase activity in the DDB2 and regulated by the COP9 signalosome in response to DNA damage. Cell 2003; 113:357-67. Uhle S, Medalia O, Waldron R, et al. Protein kinase CK2 and protein kinase D are associated with the COP9 signalosome. EMBO J. 2003; 22:1302-12.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4601P; CSN8 Peptide Cat. No. PK-AB718-1304; Human Liver Tissue Lysate

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.