

COP9 signalosome (human), (purified)

Originally identified in *Arabidopsis* as a suppressor of light-dependent development and is essential for development in plants, and is itself both a target for kinase activity and associates with and coordinates activity of kinases. Although the CSN complex and the 'lid' of the 19S regulator of the 26S proteasome share significant structural similarities, their functions seem to be different and neither protein-protein interactions nor associated activities described for the CSN have been identified for the 19S 'lid'. The CSN appears to be a platform connecting signalling with degradation by the 26S proteasome.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW9425-0020	20µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CSN

Formulation Liquid. In 20mM TRIS, pH 7.2, containing 10% glycerol, 50mM KCl and 1mM mercaptoethanol.

Source Isolated from human erythrocytes.



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)