

# CHOP monoclonal antibody (9C8)

CHOP is a small nuclear protein that is capable of dimerizing with transcription factors C/EBP  $\alpha$  and  $\beta$ . Once dimerized, this complex inhibits the normal binding and function of C/EBP to classical binding sites. Inversely, the C/EBP-CHOP dimer gains binding activity to other non-classical C/EBP stress related targets.

Under normal cellular conditions this protein is not expressed in detectable levels, but is highly unregulated during times of cellular/ER stress. Examples of CHOP inducing stress include: treatment with tunicamycin, nutrient starvation and reducing agents that interfere with the calcium flux across the ER membrane.

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

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Fig.1



## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** C/EBP homology protein, Growth arrest and DNA damage inducible protein 153, GADD153

**Application** ICC, IP, WB

**Application Notes** Detects a band of ~31kDa by Western blot.

**Clone** 9C8

**Formulation** Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium azide.

**Host** Mouse

**Immunogen** Recombinant mouse CHOP (C/EBP homology protein).

**Isotype** IgG2b

**Recommendation Dilutions/Conditions** Western Blot (1µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Mouse

**Specificity** Recognizes an N-terminal epitope.

**UniProt ID** P35639

**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)