

CHOP monoclonal antibody (9C8)

CHOP is a small nuclear protein that is capable of dimerizing with transcription factors C/EBP α and β . 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 »](#)

Ordering Information

[Order Online »](#)

ALX-804-551-C100	100 μ g
------------------	-------------

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 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](#).



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)