

Grp78/BiP polyclonal antibody

Grp78 (Glucose Regulated 78kDa Protein), is also known as BiP or Immunoglobulin heavy chain binding protein. It is a stressresponse protein which is induced by agents or conditions that adversely affect endoplasmic reticulum (ER) function. This protein is essential for the proper glycosylation, folding, and assembly of many membrane bound and secreted proteins as well as for maintenance of cell homeostasis and the prevention of apoptosis.

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

Citations: 51

[View Online »](#)

Ordering Information

[Order Online »](#)

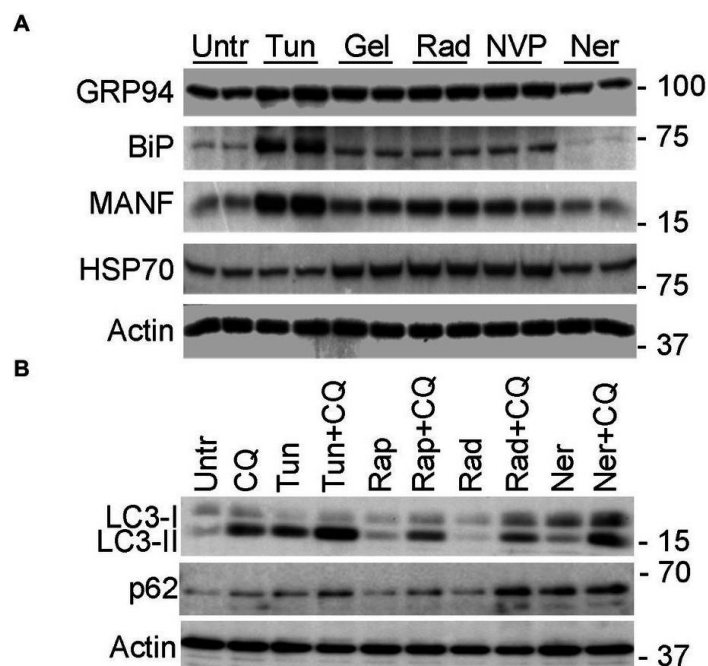
ADI-SPA-826-D	50µg
ADI-SPA-826-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis: Lane 1: MW marker, Lane 2: 150 ng of Grp78 (BiP) recombinant protein (Prod. No. ADI-SPP-765), Lane 3: 20 μg of CHO-K1 cell lysate, Lane 4: 20 μg of mouse liver microsomal tissue extract (Prod. No. ADI-LYT-MM100).



Induction of UPR (A,C) and autophagy genes (B,D) by L1000-ER/UPR/autophagy drugs. Cultured GECs were incubated for 24 h with drugs (10 μM concentrations). In panel B, drugs were added together with CQ, as indicated. Lysates were immunoblotted, as indicated. (A,B) Representative immunoblots. (C–E) Densitometric quantification (the expression of each protein is normalized to actin). * $p < 0.05$, ** $p < 0.01$, + $p < 0.001$, ++ $p < 0.0001$ vs. untreated. Mean \pm SE, 3 experiments performed in duplicate. (E) Autophagosomal flux, as determined by the level of LC3-II in the presence of chloroquine (densitometric quantification of specific lanes in panel B). * $p < 0.05$, ** $p < 0.01$, vs. chloroquine. Tunicamycin and rapamycin are stimulators of the UPR or autophagy (positive controls). Untr, untreated; Tunic, tunicamycin; Gel, geldanamycin; Rad, radicicol; NVP, NVP-AUY922; Ner, neratinib; Rap, rapamycin; CQ, chloroquine.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Analysis of gene expression and use of connectivity mapping to identify drugs for treatment of human glomerulopathies. *Front Med (Lausanne)* (2023)

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	HspA5, HSP70-5, Glucose-regulated protein 78, Heat shock 70kDa protein 5
Application	WB
Application Notes	Detects a band of ~78kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M14050
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of rat Grp78 (BiP). The sequence is completely conserved in mouse and hamster.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, ECL)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	Bovine, Fungus, Hamster, Monkey, Mouse, Rabbit, Rat, Xenopus
UniProt ID	P06761

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)