

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: 11

[View Online »](#)

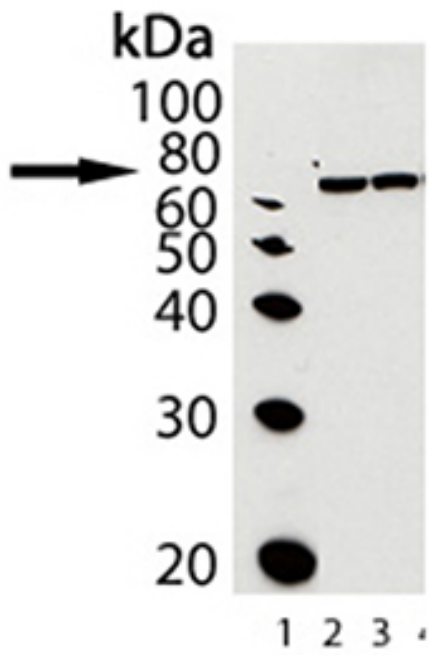
Ordering Information

[Order Online »](#)

ADI-SPA-769-0200	200µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Grp78, pAb (Prod. No. ADI-SPA-769): Lane 1: MW marker, Lane2: Grp78/BiP (hamster), (recombinant) (Prod. No. ADI-SPP-765), Lane3: Liver (rat), (microsome extract) (Prod. No. ADI-LYT-RM100)

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, Glucose Regulated Protein 78, Hsp70-5, Heat shock protein 70 family

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.

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)