

# 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: 14

[View Online »](#)

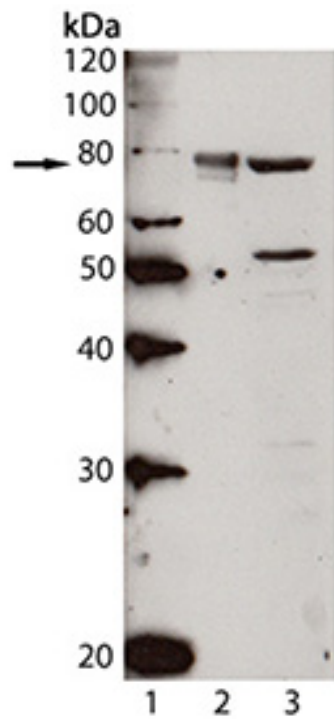
## Ordering Information

[Order Online »](#)

ADI-SPA-768-0050	50µg
ADI-SPA-768-0200	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Grp78, pAb (Prod. No. ADI-SPA-768): 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, Dog, Fungus, Guinea pig, Hamster, Human, Monkey, Mouse, Rabbit, Rat, Sheep, Xenopus
UniProt ID	P06761
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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)