

# Grp78/BiP (hamster), (recombinant)

Grp78, a 78 kDa glucose-regulated protein also known as BiP or immunoglobulin heavy chain binding protein, is a stress-response protein 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. Grp78 is critical for maintenance of cell homeostasis and the prevention of apoptosis. Grp78 protein levels provide a reliable biomarker of hypoglycemia as well as serve a neuroprotective function in neurons exposed to glutamate and oxidative stress. Reduced Grp78 levels occur in the brains of Alzheimer's Disease patients, and decreased Grp78 expression appears to be associated with missense mutations in the human presenilin-1 (PS1) gene. The induction of the Grp78 protein is associated with the development of drug-resistance to anti-tumor drugs.

Citations: 7

[View Online »](#)

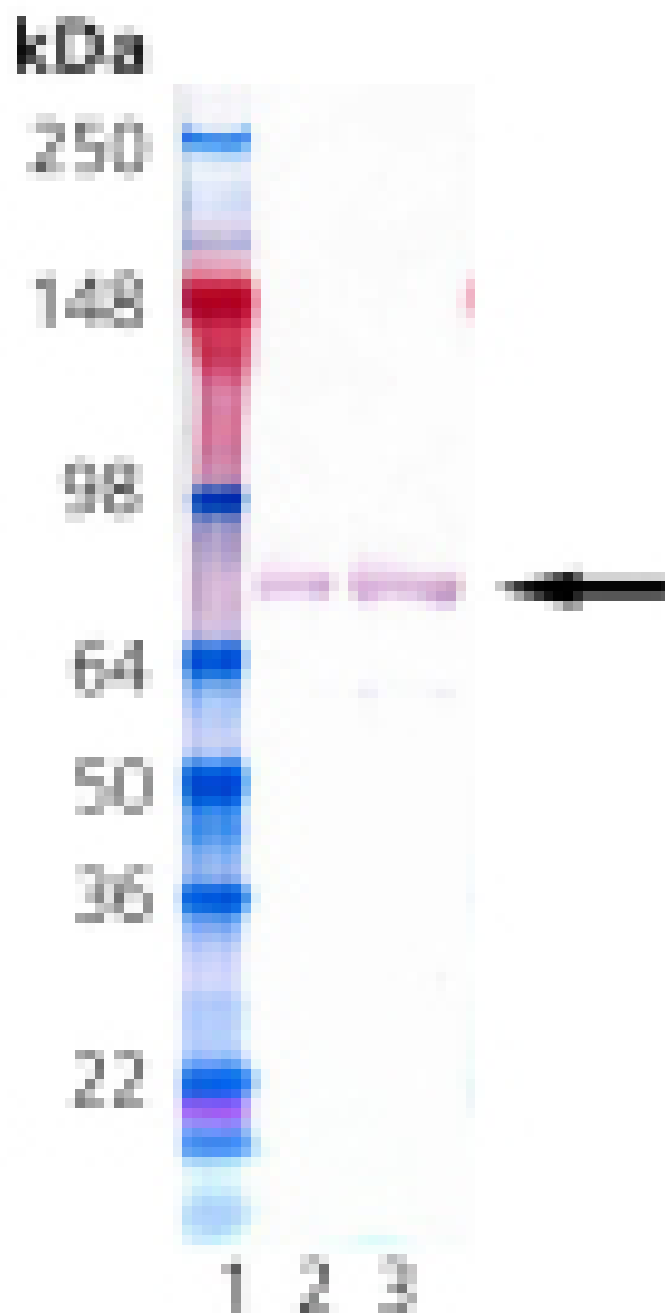
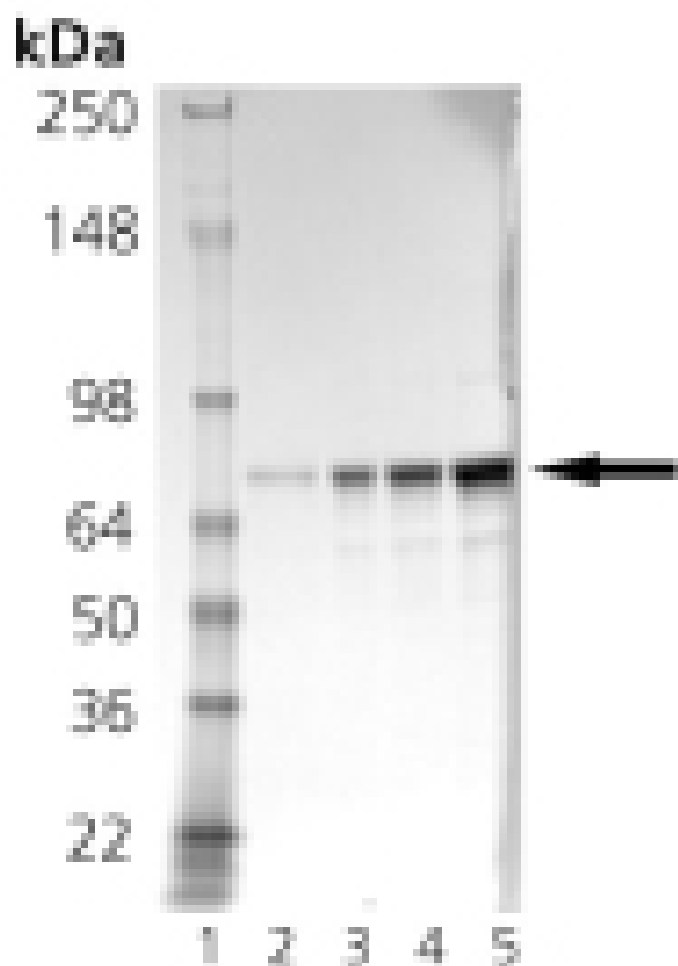
## Ordering Information

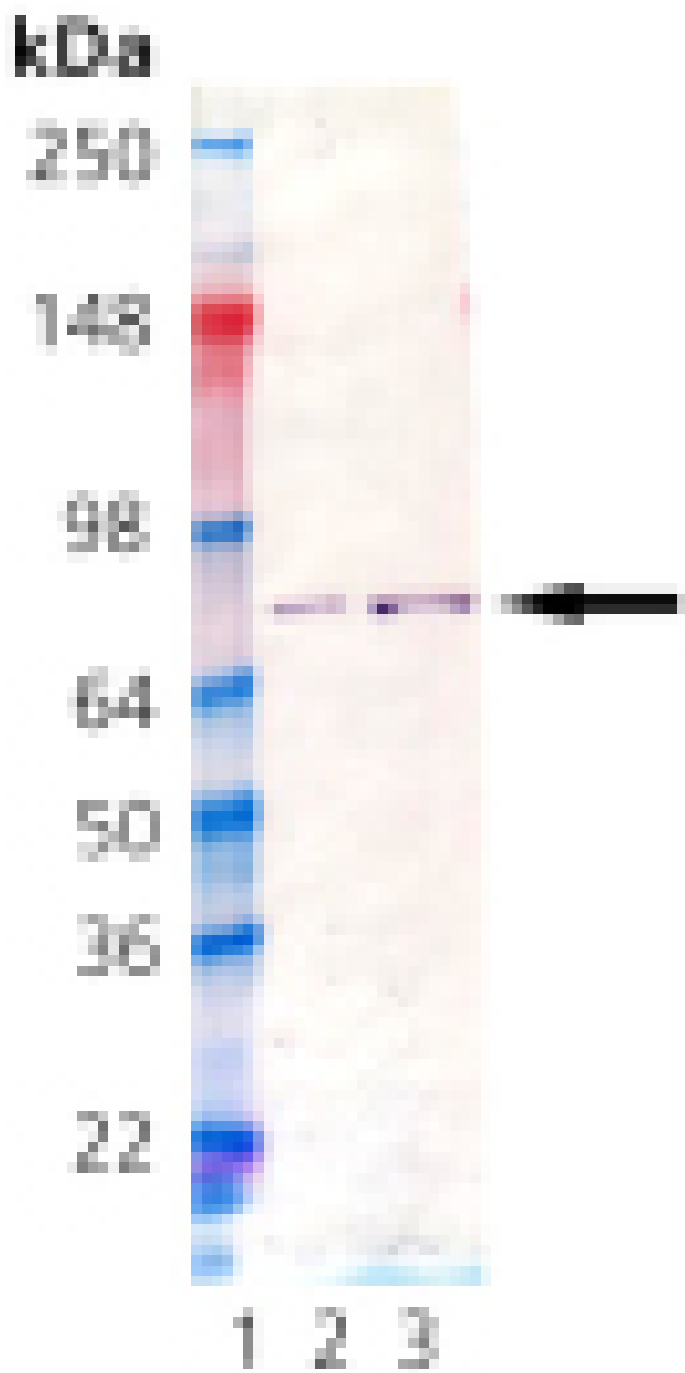
[Order Online »](#)

ADI-SPP-765-D	50µg
ADI-SPP-765-F	200µg

## Manuals, SDS & CofA

[View Online »](#)





Western Blot analysis: Lane 1: MWM, Lane 2 and 3: Recombinant Grp78 Hamster Protein probed with Anti-KDEL (Grp78, Grp94) mAb.

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Alternative Name**      HspA5, Hsp70-5, Heat shock 70 kDa protein 5

**Application Notes**      Western blot control.

**Formulation**      Liquid. In 50mM Tris, pH 7.5, containing 150mM NaCl, 10% glycerol and 1mM DTT

**MW**      ~78kDa

**Purity**      ≥80% (SDS-PAGE; Western blot)

**Purity Detail**      Purified by multi-step chromatography.

**Source**      Produced in *E. coli*.

**UniProt ID**      P07823



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)