Grp78/BiP (hamster), (recombinant)

Recombinant hamster Grp78/BiP, a 78 kDa ER chaperone that ensures proper protein folding, assembly, and glycosylation, while maintaining cellular homeostasis and

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 »

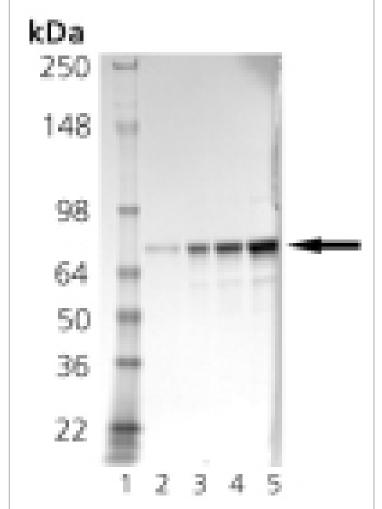
Ordering Information

Order Online »

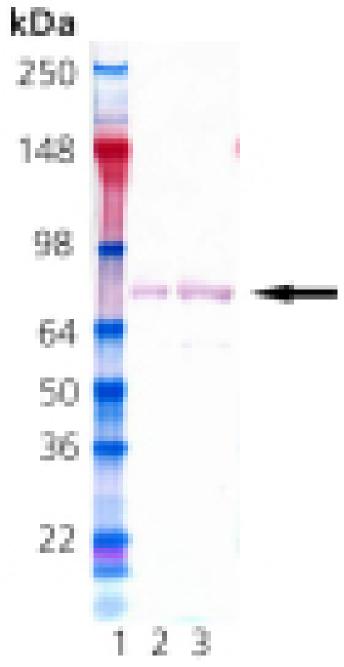
ADI-SPP-765-D	50µg
ADI-SPP-765-F	200μg

Manuals, SDS & CofA

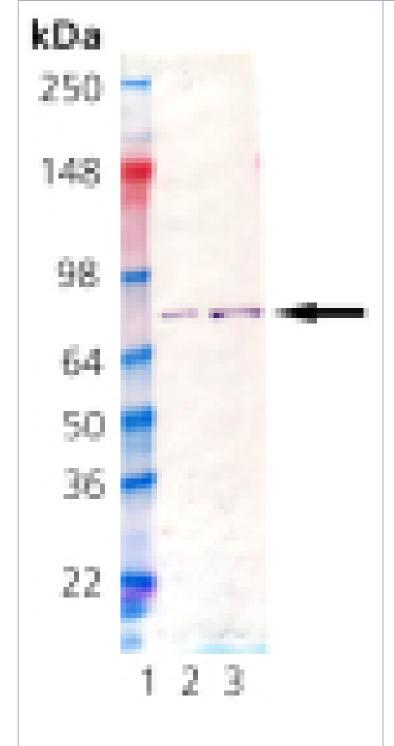
View Online »



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 0.5µg, Lane 3: 1µg, Lane 4: 2µg, Lane 5: 5µg of Recombinant Grp78 Hamster Protein.



Western Blot analysis: Lane 1: MW marker, Lane 2 and 3: Recombinant Grp78 Hamster Protein probed with anti-Recombinant Grp78 (Bip) pAb.



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

Last modified: October 9, 2025



eu@enzolifesciences.com