

# PRODUCT INFORMATION



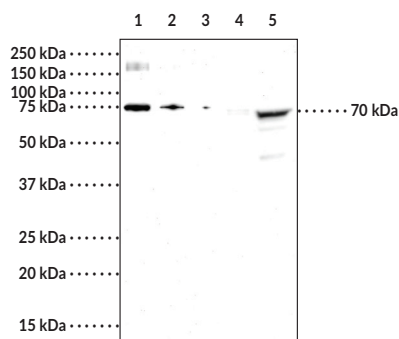
## GRP78 Polyclonal Antibody

Item No. 24533

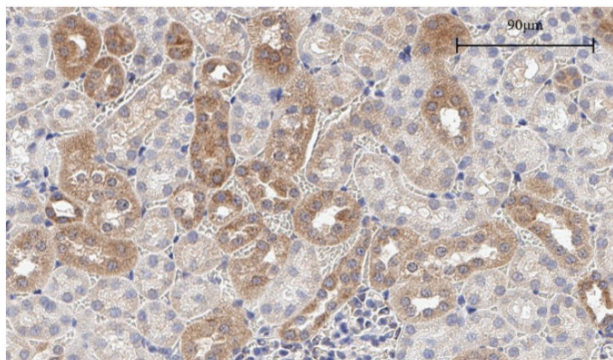
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µg of protein A-purified GRP78 polyclonal antibody
<b>Synonyms:</b>	BiP, Endoplasmic Reticulum Luminal Ca <sup>2+</sup> -Binding Protein GRP78, 78 kDa Glucose-Regulated Protein, Glucose-Regulated Protein 78, Heat Shock 70 kDa Protein 5, HspA5, Immunoglobulin Heavy Chain-Binding Protein
<b>Immunogen:</b>	Human recombinant GRP78 protein (full length)
<b>Species Reactivity:</b>	(+) Human
<b>Uniprot No.:</b>	P11021
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	ELISA, Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1 µg/ml for ELISA and 5 µg/ml for IHC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: GRP78 Recombinant Protein (0.02 µg)  
Lane 2: GRP78 Recombinant Protein (0.005 µg)  
Lane 3: GRP78 Recombinant Protein (0.001 µg)  
Lane 4: Hsp70 (HspA1A) Recombinant Protein (0.1 µg) [negative control]  
Lane 5: HeLa Heat Shock Cell Lysate (50 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human kidney tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with GRP78 Polyclonal Antibody (Item No. 24533) at a 1:200 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/22/2023

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

Glucose-regulated protein 78 kDa (GRP78), also known as heat shock 70 kDa protein 5 (HspA5) and immunoglobulin heavy chain-binding protein (BiP), is a glucose-regulated protein that is constitutively expressed in the lumen of the endoplasmic reticulum (ER).<sup>1-3</sup> It is composed of two functional domains, an N-terminal nucleotide-binding domain that contains an ATP catalytic site and a C-terminal substrate binding domain that binds hydrophobic polypeptides.<sup>4</sup> GRP78 functions as a molecular chaperone, assisting in the translocation of polypeptides from the cytosol into the ER, folding of nascent polypeptides, as well as refolding and preventing aggregation of misfolded proteins. It also plays a role in the ER-assisted degradation (ERAD) and unfolded protein response (UPR) pathways.<sup>5,6</sup> GRP78 chaperone activity is driven by an ATPase cycle that is regulated by ER-localized DnaJ-like protein co-factors and nuclear exchange factors.<sup>7,8</sup> Expression of GRP78 is upregulated in response to ER stress caused by viral and bacterial infections as well as various cancers.<sup>9</sup> ER stress can also promote extracellular secretion of GRP78 leading to its anti-inflammatory functions in immunity.<sup>10</sup> Cayman's GRP78 Polyclonal Antibody can be used for Western blot and ELISA applications. The antibody recognizes GRP78 at ~72 kDa from human samples.

## References

1. Vogel, J.P., Misra, L.M., and Rose, M.D. Loss of BiP/GRP78 function blocks translocation of secretory proteins in yeast. *J. Cell. Biol.* **110(6)**, 1885-1895 (1990).
2. Simons, J.F., Ferro-Novick, S., Rose, M.D., et al. BiP/Kar2p serves as a molecular chaperone during carboxypeptidase Y folding in yeast. *J. Cell. Biol.* **130(1)**, 41-49 (1995).
3. Mayer, M.P. and Bukau, B. Hsp70 chaperones: Cellular functions and molecular mechanism. *Cell Mol. Life Sci.* **62(6)**, 670-684 (2005).
4. Yang, J., Nune, M., Zong, Y., et al. Close and allosteric opening of the polypeptide-binding site in a human Hsp70 chaperone BiP. *Structure* **23(12)**, 2191-2203 (2015).
5. Plemper, R.K., Böhmler, S., Bordallo, J., et al. Mutant analysis links the translocon and BiP to retrograde protein transport for ER degradation. *Nature* **388(6645)**, 891-895 (1997).
6. Okamura, K., Kimata, Y., Higashio, H., et al. Dissociation of Kar2p/BiP from an ER sensory molecule, Ire1p, triggers the unfolded protein response in yeast. *Biochem. Biophys. Res. Commun.* **279(2)**, 445-450 (2000).
7. Szabo, A., Langer, T., Schröder, H., et al. The ATP hydrolysis-dependent reaction cycle of the *Escherichia coli* Hsp70 system DnaK, DnaJ, and GrpE. *Proc. Natl. Acad. Sci. USA* **91(22)**, 10345-10349 (1994).
8. Behnke, J., Feige, M.J., and Hendershot, L.M. BiP and its nucleotide exchange factors Grp170 and Sil1: Mechanisms of action and biological functions. *J. Mol. Biol.* **42(7)**, 1589-1608 (2015).
9. Booth, L., Roberts, J.L., Cash, D.R., et al. GRP78/BiP/HSPA5/Dna K is a universal therapeutic target for human disease. *J. Cell. Physiol.* **230(7)**, 1661-1676 (2015).
10. Shields, A.M., Panayi, G.S., and Corrigan, V.M. Resolution-associated molecular patterns (RAMP): RAMParts defending immunological homeostasis. *Clin. Exp. Immunol.* **165(3)**, 292-300 (2011).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM