

PRODUCT INFORMATION



Citrullinated GRP78 (human, recombinant)

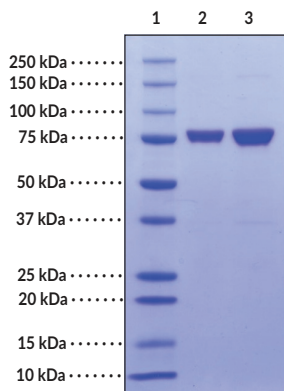
Item No. 25107

Overview and Properties

Synonyms:	BiP, Endoplasmic Reticulum Luminal Ca ²⁺ -Binding Protein GRP78, Glucose-Related Protein 78, Heat Shock Protein 5 (70 kDa), Hsp5 (70 kDa), Immunoglobulin Heavy Chain-Binding Protein
Source:	N-Terminal histidine-tagged human GRP78 purified from <i>E. coli</i> , citrullinated by PAD2
Amino Acids:	2-654
Uniprot No.:	P11021
Molecular Weight:	74.6 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	batch specific (≥90% estimated by SDS-PAGE)
Supplied in:	PBS, pH 7.4, with 10% glycerol
Protein	
Concentration:	batch specific mg/ml

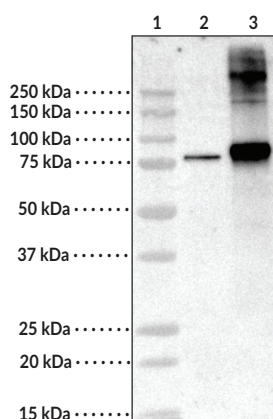
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: Citrullinated GRP78 (2 µg)
Lane 3: Citrullinated GRP78 (4 µg)

SDS-PAGE analysis of citrullinated GRP78. Citrullinated GRP78 at 2 µg (Lane 2) and 4 µg (Lane 3) stained with Coomassie on 4-20% SDS-PAGE.



Lane 1: MW Markers
Lane 2: GRP78
Lane 3: Citrullinated GRP78

Analysis of GRP78 citrullination. GRP78 and citrullinated GRP78 were reacted with Cayman's Citrulline-specific Probe-biotin (Item No. 17450) and detected using Streptavidin:HRP (Item No. 16747).

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, 02/18/2025

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) is a molecular chaperone that is ubiquitously expressed in the endoplasmic reticulum of mammalian cells.¹⁻³ GRP78 can be citrullinated at its 27 arginine residues by protein deiminases (PADs).⁴ The accumulation of citrullinated proteins *in vivo* leads to the production of anti-citrullinated protein antibodies (ACPAs) which perpetuate the inflammatory process.⁵ *In vitro*, ACPAs bind to citrullinated GRP78 expressed on the cell surface of peripheral blood mononuclear cells PMBCs and U937 cells leading to the production of TNF- α .^{6,7} In a mouse model of collagen-induced arthritis (CIA), anti-citrullinated GRP78 antibodies are found in the serum.⁴ Pre-immunization with citrullinated GRP78 prior to CIA induction shortens the time to joint inflammation and increases arthritis scores compared with non-citrullinated GRP78-immunized and non-immunized control mice. In autoimmune diseases such as rheumatoid arthritis, patient-derived serum contains higher levels of anti-citrullinated GRP78 antibodies than serum derived from patients with systemic lupus erythematosus and healthy controls.⁴

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. Shoda, H., Fujio, K., Shibuya, M., *et al.* Detection of autoantibodies to citrullinated BiP in rheumatoid arthritis patients and pro-inflammatory role of citrullinated BiP in collagen-induced arthritis. *Arthritis Res. Ther.* **13**(6), R191 (2011).
5. Kuhn, K.A., Kulik, L., Tomooka, B., *et al.* Antibodies against citrullinated proteins enhance tissue injury in experimental autoimmune arthritis. *J. Clin. Invest.* **116**(4), 961-973 (2006).
6. Lu, M.C., Lai, N.S., Yu, H.C., *et al.* Anti-citrullinated protein antibodies bind surface-expressed citrullinated Grp78 on monocyte/macrophages and stimulate tumor necrosis factor α production. *Arthritis Rheum.* **62**(5), 1213-1223 (2010).
7. Lu, M.C., Lai, N.S., Yin, W.Y., *et al.* Anti-citrullinated protein antibodies activated ERK1/2 and JNK mitogen-activated protein kinases via binding to surface-expressed citrullinated GRP78 on mononuclear cells. *J. Clin. Immunol.* **33**(3), 558-566 (2013).

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