



EAGLE
BIOSCIENCES

**Anti-Glucose Regulated Protein
(Grp) 75 (N-Terminus)**

Product Number: GRP00-A200

Aliquot: 200 μ L

Storage: -70° C

PRODUCT SPECIFICATION SHEET

Contents:	Polyclonal rabbit antiserum to human Glucose Regulated Protein (Grp) 75
Immunogen:	N-terminal fragment of human Grp 75 (amino acids 47-69)
Specificity:	Human and mouse Grp 75
Application:	Western blot analysis, chemiluminescent detection at 1-2 minute exposure.
Quantity:	200 μ L rabbit antiserum
Working Dilution:	A dilution of 1:500 for Western blots will yield a positive result with Grp 75. Secondary antibody used is anti-rabbit IgG peroxidase diluted at 1:100,000 (Cat# 605220; Boehringer-Mannheim; Indianapolis, IN)
Storage:	Stable for several weeks at 4°C, several months at -20°C, and indefinitely at -70°C. Avoid repeated freeze/thaw cycles.

References:

1. Wadhwa, R., Kaul, S.C., Mitsui, Y., *Cell Struct. Function*, **19** : 1-10 (1994)
2. Kim, H.T., Nelson, Clayberger, Sanjanwala, M., Sklar, J., Krensky, A.M., *J. Immunol.*, **154** : 1614-1623 (1995)
3. Mizzen, L.A., Chang, C., Garrels, J.I., Welch, W.J., *Biol Chem.*, **264** : 20664-20675 (1989)
4. Webster, T.J., Naylor, D.J., Hartman, D.J., Hoj, P.B., Hoogenraad, N.J., *DNA and Cell Biol.*, **13**: 1213-1220 (1994)

Product Developed and Manufactured in the USA

EAGLE BIOSCIENCES, INC.
20A NW Blvd., Suite 112
Nashua, NH 03063

Phone: 617-419-2019
Fax: 617-419-1110
Email: info@EagleBio.com
www.EagleBio.com