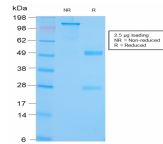


# GRP94 / HSP90B1(Glucose-Regulated Protein 94) (Endoplasmic Reticulum Marker) Monoclonal Rabbit Antibody (HSP90B1/3168R)



## Product Description

Recognizes a protein of 94 kDa, which is identified as the glucose-regulated protein 94 (grp94) and also tumor rejection antigen (gp96). Grp94 shows a high degree of sequence homology with the heat shock protein 90 (hsp90). This MAb is highly specific to grp94 and shows minimal cross-reaction with other members of the HSP90 family. Grp's are a class of proteins unresponsive to heat shock and are induced by glucose deprivation. Grp94 has been studied as a prognostic factor in breast cancer. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

## Product attributes

Antibody number	#3168
Antibody reactivity (target)	GRP94, HSP90B1
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	HSP90B1/3168R
Isotype	IgG
Molecular weight	94 kDa
Synonyms	Glucose Regulated Protein 94; 94kDa glucose-regulated protein; Endoplasmic; Endothelial cell (HBMEC) glycoprotein (ECGP); Heat shock protein 90kDa beta member 1 (HSP90B1); Stress inducible tumor rejection antigen GP96; TRA1; Tumor rejection antigen-1 (gp96)
Human gene symbol	HSP90B1
Entrez gene ID	7184
SwissProt	P14625
Unigene	192374
Immunogen	Recombinant full-length human HSP90B1 protein
Verified antibody applications	Flow (intracellular) (verified), IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Chicken, Cow, Dog, Guinea pig, Hamster, Horse, Human, Monkey, Mouse, Pig, Rabbit, Rat, Sheep, Xenopus laevis
Positive control	HepG2, U20S cells. About 50% of breast tumors express grp94 protein.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Tumor expression	Breast cancer

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.