# Grp94 monoclonal antibody (9G10) (PE conjugate)

Glucose-regulated protein 94 (Grp94, gp96), an abundant resident endoplasmic reticulum (ER) lumenal stress protein, belongs to the Hsp90 family of molecular chaperones along with cytosolic Hsp90. Grp94 and such other resident soluble proteins of the ER as the Ca 2+ binding protein subfamily (CaBP, CaBPI, CaBP2 and calreticulin) possess the C-terminal tetrapeptide Lys-Asp-Glu-Leu (KDEL), a sorting signal considered responsible for the retention of these proteins in the pre-Golgi compartments. Stress conditions such as glucose starvation and heat shock which promote protein misfolding or unfolding increase Grp94 expression. In addition to a homeostatic role in protein folding and assembly, Grp94 can function in the intracellular trafficking of peptides from the extracellular space to the MHC class I antigen processing pathway of antigen presentation cells. Grp94 and Hsp90 share high sequence identity and apparently identical adenosine nucleotide dependent modes of regulation, although previous data suggests that Hsp90 and Grp94 may differ in their nucleotide binding properties. The N-terminal domain of eukaryotic Hsp90 proteins contains a conserved adenosine nucleotide binding pocket which also serves as the binding site for the Hsp90 inhibitors geldanamycin and radicicol. However, the molecular basis for adenosine nucleotide-dependent regulation of Grp94 remains unclear. Data supports a ligand dependent regulation of Grp94 function, and suggests a model whereby Grp94 function is regulated through a ligand dependent conversion of Grp94 from an inactive to an active conformation.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

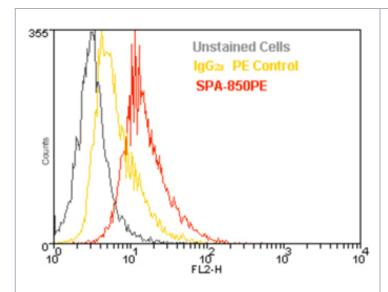
# **Ordering Information**

**Order Online** »

ADI-SPA-850PE-D	50µg
ADI-SPA-850PE-F	200µg

Manuals, SDS & CofA

View Online »



Flow cytometry analysis of  $10^6$  Jurkat cells stained using Grp94 mAb (9G10), PE Conjugate (ADI-SPA-850PE) at a concentration of  $10\mu g/ml$ .

## **Handling & Storage**

**Handling** Avoid vigorous centrifugation and vortexing. Do not freeze. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Endoplasmin, Tra1, Hsp90B1, Gp96

**Application** Flow Cytometry

**Application Notes**Detects a band of ~94kDa by Western blot.

**Clone** 9G10

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID M14772

**Host** Rat

**Immunogen** Native chicken Grp94.

**Isotype** IgG2a

Purity Detail Protein G affinity purified.

**Recommendation Dilutions/Conditions** Flow Cytometry (1:100)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from ascites.

**Species Reactivity**Bovine, Chicken, Dog, Guinea pig, Hamster, Human,

Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID P08110

Last modified: May 29, 2024



info-

eu@enzolifesciences.com