Grp94 (canine), (recombinant)

Recombinant canine Grp94, an ER-resident Hsp90-family chaperone that assists in protein folding, assembly, and trafficking, and is upregulated under stress conditions

Green regulated protein 94) is an abundant resident endoplasmic reticulum (ER) lumenal stress protein, which together with cytosolic Hsp90 belongs to the Hsp90 family of molecular chaperones. Grp94 expression is upregulated by stress conditions such as glucose starvation and heat shock, which promote protein misfolding or unfolding. 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.

Citations: 1

View Online »

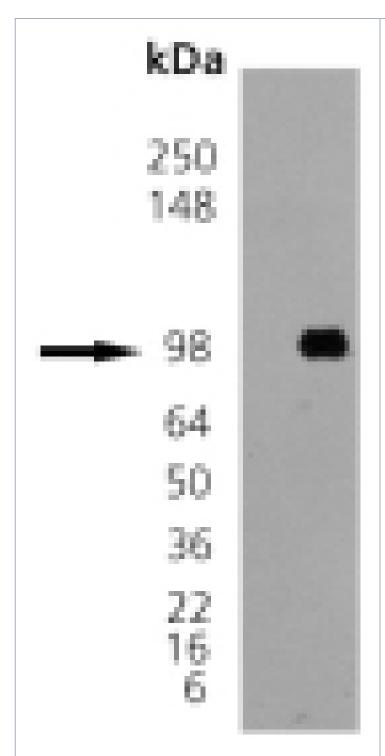
Ordering Information

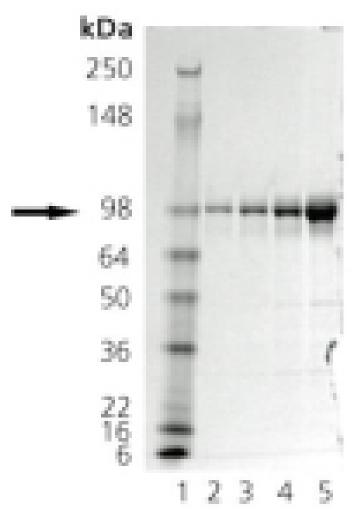
Order Online »

ADI-SPP-766-D	50µg
ADI-SPP-766-F	200µg

Manuals, SDS & CofA

View Online »





SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 0.5µg; Lane 3: 1µg; Lane 4: 2µg; Lane 5: 5µg of Grp94 (canine), (recombinant) detected by Coomassie Stain

Western Blot analysis: 100 ng of Grp94 (canine), (recombinant) probed with Prod. No. ADI-SPA-850.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tra1, Tumor rejection antigen 1, Hsp90B1, Gp96, Heat

shock protein 90 kDa beta member 1, Endoplasmin

Application Notes Western blot control.

Formulation Liquid. In 20mM MOPS, pH 7.5, containing 0.5mM EDTA.

MW ∼94kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in Sf21 insect cells. Produced in a baculovirus

expression system.

UniProt ID P41148

Last modified: October 9, 2025

