ERp72 (human), (recombinant) (Histag)

Recombinant human ERp72 (His-tag), an ERresident protein disulfide isomerase that

Enclosive min reticulum protein (FRP) 72 is a member of the protein disulfide isomerase (PDI) family that is localized to the endoplasmic synthesized glycoproteins and acts as a reticulum. It contains three repeats of the thioredoxin-like regions that is posturated to represent in earlier independently acting catalytic domains (CGHC) of PDI activity. ERp72 contains the sequence at its carboxyl terminus which serves as its ER retention signal. Together with other ER resident proteins such as BiP, GRP94, and PDI, they serve as the molecular chaperones for proper folding of newly translocated and glycolsylated proteins such as thyroglobulin (Tg) and human chorionic gonadotropin (hCG)-b. ERp72 expression is regulated by the level of misfolding proteins in the ER, as the amount of ERp72 increased in response to epithelial ischemia, a condition that perturbs the maturation of secretory proteins.

Citations: 1

View Online »

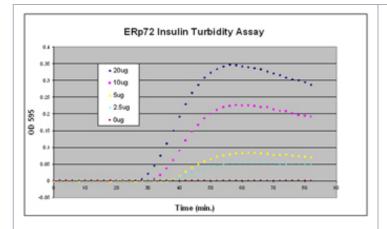
Ordering Information

Order Online »

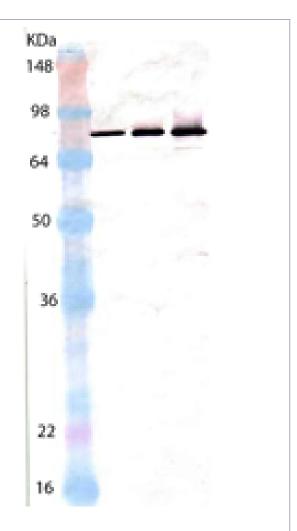
ADI-SPP-721-050	50µg
ADI-SPP-721-200	200µg

Manuals, SDS & CofA

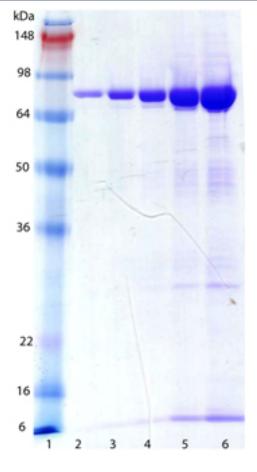
View Online »



Activity Assay: Varying amounts of ERp72 were incubated with insulin in the presence of DTT. Activity is measured as an increase in absorbance over time.



Western blot analysis of ERp72: Lane 1: MW marker, Lane 2: 25ng, Lane 3: 50ng, Lane 4: 100ng.



SDS PAGE analysis of ERp72: Lane 1: MW Marker, Lane 2: 0.25ug, Lane 3: 0.5ug, Lane 4: 1.0ug, Lane 5:

2.5ug, Lane 6: 5.0ug.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PDIA4, Protein disulfide isomerase A4

Application Notes Western blot control.

Formulation Liquid. In 50mM TRIS, pH 7.5, containing 150mM sodium

chloride and 1mM EDTA.

MW ~68kDa

Purity ≥85% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in E. coli. Human ERp72 is fused at the N-

terminus to a His-tag.

UniProt ID P13667

Last modified: October 9, 2025

