

# ERp72 (human), (recombinant) (His- tag)

Endoplasmic reticulum protein (ERp) 72 is a member of the protein disulfide isomerase (PDI) family that is localized to the endoplasmic reticulum. It contains three repeats of the thioredoxin-like regions that is postulated to represent three 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 glycosylated 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 »](#)

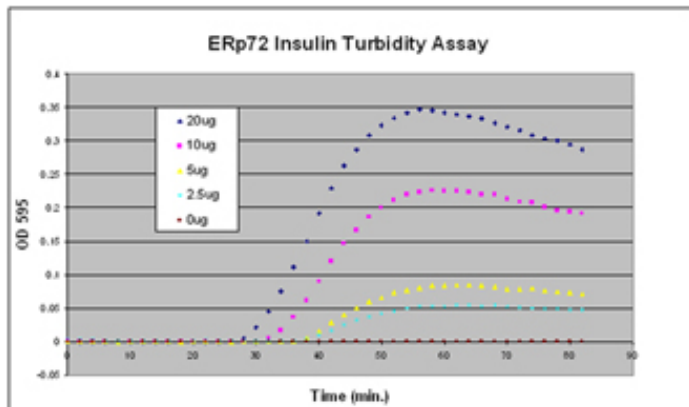
## Ordering Information

[Order Online »](#)

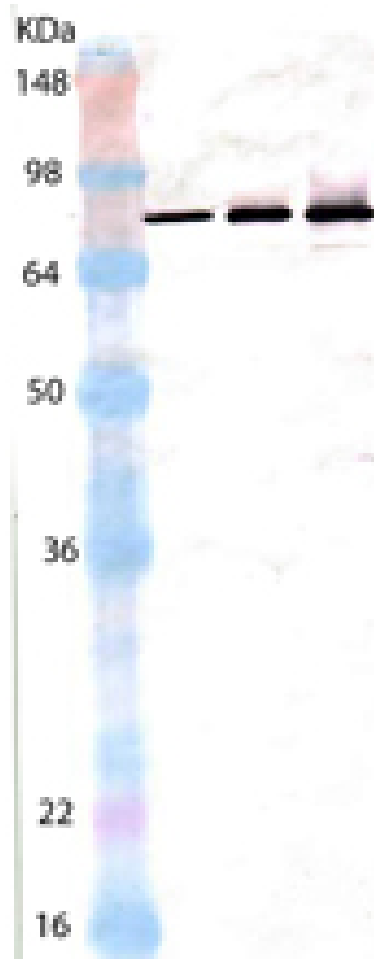
ADI-SPP-721-050	50µg
ADI-SPP-721-200	200µg

## Manuals, SDS & CofA

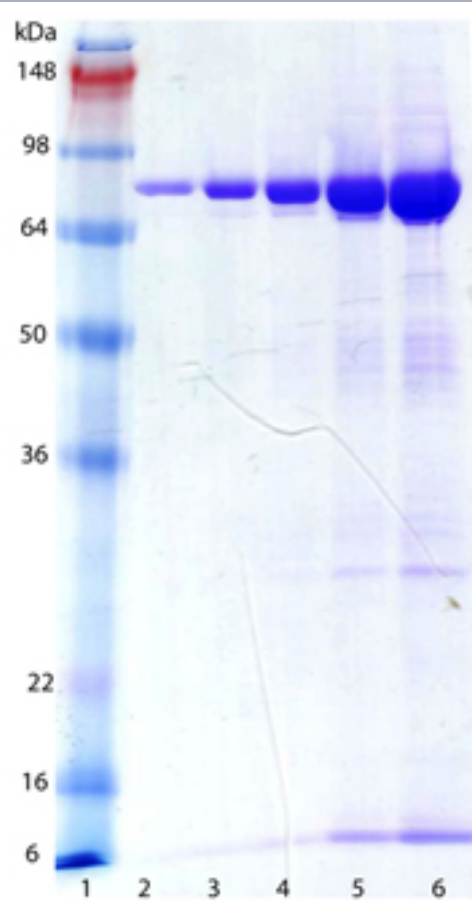
[View Online »](#)



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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)