GrpE (E. coli), (recombinant)

Recombinant GrpE (E. coli) is a 24 kDa cochaperone that assists DnaK/Hsp70 in protein folding, transport, disaggregation, and activation of partially folded proteins.

Greate an epitia fet & Darprotein which the encoded as a reprocistronic Escherichia coli gene (located at 56 min.) which is under the control of both substrates from the Dnak/Dna complex and s32 and s70 transcription factors. The GrpE protein is a so-called 63chaberone foecable mist on Swine Easters the Driak 7Hsp70) protein to effectively carry out DnaK-dependent chaperone activity (i.e. protein folding, protein transport, disaggregation of heat inactivated proteins, activation of mutant protein). This auxiliary role is critical to DnaK's chaperone biology since the grpE gene could not be deleted in a wild type background unless the host E. coli were previously adapted for deletions of the DnaK gene through the accumulation of extragenic suppressors. Also, in vitro these two proteins functionally interact; GrpE protein forms a salt resistant complex with DnaK protein, which is disrupted by ATP. It is suggested that DnaK and DnaJ assist protein folding by stabilizing an intermediate conformation; where as the GrpE protein, in an ATPdependent reaction release partially folded protein from the DnaK/J complex and transfer this immature protein to the GroEL/ES chaperone complex to finish the protein folding reaction. In some cases DnaK/GrpE could act independently of DnaJ chaperone protein: the "inactive" form of DnaA5 mutant protein is activated in the presence of the DnaK and GrpE proteins, and the presence of DnaJ inhibit this reaction. GrpE cochaperone protein assists DnaK chaperone during two crucial steps: binding to protein substrate and releasing DnaK chaperone from the complex with substrate.

Citations: 5

View Online »

Ordering Information

Order Online »

ADI-SPP-650-D	50µg
ADI-SPP-650-F	200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name HSP70 cofactor, Nucleotide exchange factor, HSP24, Heat

shock protein 24

Application Notes Western blot control.

Formulation Liquid. In 25mM HEPES, pH 7.6, containing 50mM

potassium chloride and 5.0% glycerol.

MW ~24kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Produced in E. coli. Source

UniProt ID P09372 (strain K12)

Last modified: October 9, 2025



eu@enzolifesciences.com