

GrpE (E. coli), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-201-145-C025 | 25µg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---|
| Handling | For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. |
| Short Term Storage | +4°C |
| Long Term Storage | -80°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------|--|
| Alternative Name | HSP70 cofactor, Nucleotide exchange factor, HSP24, Heat shock protein 24 |
| Formulation | Liquid. In 30mM HEPES/KOH, pH 7.0, containing 10mM KCl and 5% glycerol. |
| Purity | ≥99% |
| Source | Produced in <i>E. coli</i> . |
| UniProt ID | P09372 (strain K12) |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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