

DnaJ (E. coli), (recombinant)

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|--------------|------|
| ALX-201-1440 | 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 | Dry Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|---|
| Alternative Name | Heat shock protein 40, HSP40, Heat shock protein J |
| Formulation | Liquid. In 30mM HEPES/KOH, pH 8.2, containing 120mM KCl, 15mM NaCl and 4% glycerol. |
| Purity | ≥99% |
| Source | Produced in <i>E. coli</i> . |
| UniProt ID | P08622 |

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