

DnaJ (E. coli), (recombinant)

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-144-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	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