

# HSP40 (human), (recombinant) (His- tag)

## Ordering Information

[Order Online »](#)

|                  |       |
|------------------|-------|
| ALX-201-274-C100 | 100µg |
|------------------|-------|

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** SDS-PAGE of HSP40 (human) (recombinant) (His) (Prod. No. ALX-201-274).

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Heat shock protein 40, Hdj1, DnaJ protein homolog 1

**Concentration** 1mg/ml

**Endotoxin Content** <1 EU/µg protein (determined by LAL method).

**Formulation** 0.2µm filtered solution. In PBS, pH 7.4, containing 10% glycerol

**MW** ~40.2kDa

**Purity** ≥90% (SDS-PAGE)

**Source** Produced in *E.coli* and fused to a His-tag.

**UniProt ID** P25685

**Last modified: May 29, 2024**



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)