

# DsbA (E. coli), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-201-153-C025 | 25µg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                    |                                                                                                                      |
|--------------------|----------------------------------------------------------------------------------------------------------------------|
| Handling           | For maximum product recovery after thawing, centrifuge the vial before opening the cap.<br>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 | Thiol:disulfide interchange protein DsbA |
| Formulation      | Liquid. In water.                        |
| Purity           | ≥99%                                     |
| Source           | Produced in <i>E. coli</i> .             |
| UniProt ID       | P0AEG4                                   |

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)