

DsbC (E. coli), (recombinant)

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-268-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

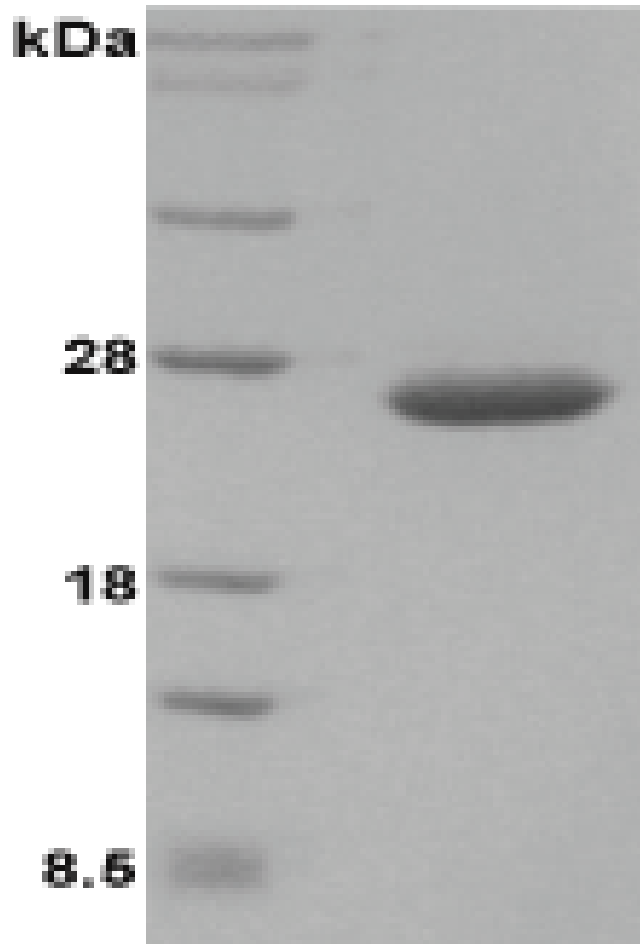


Figure: SDS-PAGE of DsbC (*E. coli*) (recombinant)
(Prod. No. [ALX-201-268](#)).",DsbC (E. coli)

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 Thiol:disulfide interchange protein DsbC

Concentration 1mg/ml

Formulation Liquid. In 20mM TRIS, pH 7.5, containing 2mM EDTA.

MW ~23.5kDa

Purity ≥95% (SDS-PAGE)

Source Produced in *E. coli*.

UniProt ID P0AEG6

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)