

E. coli host cell protein standards

Quantitative protein standards for Western blotting internal controls or ELISA standards.

E. coli host cell protein standards are native proteins that can be used as an internal control for the Western blot analysis of CHO Host Cell Proteins or as standards in an ELISA.

Ordering Information

[Order Online »](#)

ENZ-PRT121-0025	25µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance

Clear liquid.

Application Notes

Protein standards for the *E. coli* host cell protein ELISA kit or for Western blot analysis.

Formulation

Phosphate-buffered solution.

Source

Purified from *E. coli* host cells.

Last modified: December 1, 2025



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-
fr@enzolifesciences.com](mailto:info-fr@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)