

# CHO host cell protein standards

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

CHO 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-PRT122-0025	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 CHO host cell protein ELISA kit or for Western blot analysis.

**Formulation** Liquid. In PBS only.

**Source** Purified from Chinese Hamster Ovary (CHO) host cells.



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)