

# HEK293T Host Cell Protein Standards

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

HEK293T host cell protein standards are native proteins that can be used as an internal control for the Western blot analysis of HEK293T Host Cell Proteins or as standards in an ELISA.

## Ordering Information

[Order Online »](#)

ENZ-PRT311-0025	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      Used as internal control for the Western blot analysis of HEK293T host cell proteins or development of HCP ELISA

Source      Purified from HEK293T 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)