

HEK293T host cell protein ELISA kit

Sensitive ELISA for the quantitative determination of host cell protein contamination in bulk products expressed in HEK293T expression systems.

The HEK293T host cell protein ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of HEK293T host cell protein contamination with results in 6 hours

Ordering Information

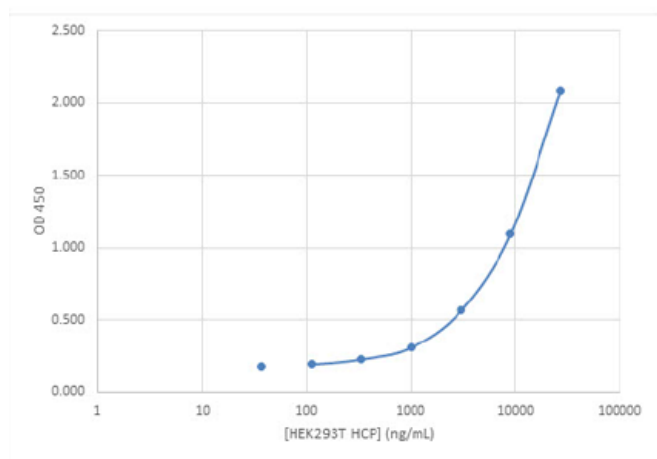
[Order Online »](#)

ENZ-KIT162-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of HEK293T host cell protein
- Low intra-assay variability allows for streamlined testing
- High throughput format with results in 6 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Standard curve of HEK293T host cell protein ELISA kit

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HEK293T HCP ELISA kit
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of contamination with HEK293T host cell protein.
Assay Time	6 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	1 x 96 well plate, protein standards, 1x dilution buffer, 10x PBS-T, reporting antibody, streptavidin-HRP conjugate, TMB substrate, stop solution, plate sealer
Sensitivity	37 ng/ml (37-27000 ng/ml)

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