

CHO host cell protein ELISA kit

Sensitive (10 ng/ml) ELISA for the quantitative determination of host cell protein contamination in bulk products expressed in Chinese Hamster Ovary (CHO) expression systems.

The CHO host cell protein ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of Chinese Hamster Ovary (CHO) host cell protein contamination with results in 3 hours

Ordering Information

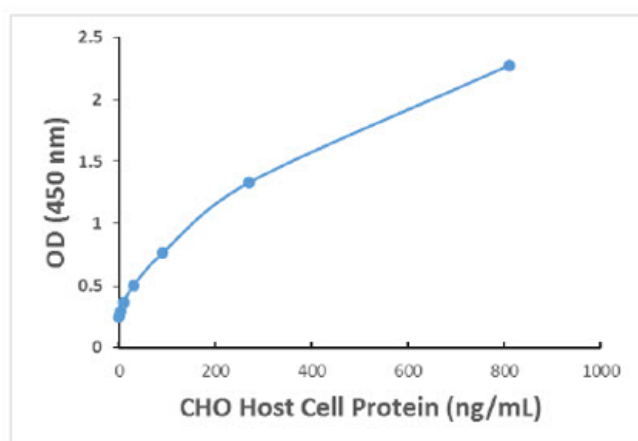
[Order Online »](#)

ENZ-KIT128-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of CHO host cell protein, detecting as little as 10 ng/ml
- Low intra-assay variability allows for streamlined testing
- High throughput format with results in 3 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis



Standard curve of CHO 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	CHO HCP ELISA kit
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of contamination with CHO host cell protein.
Assay Time	3 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	1 x 96 well plate, protein standards, 5x dilution buffer, 10x PBS-T, reporting antibody, streptavidin-HRP conjugate, TMB substrate, stop solution, plate sealer
Sensitivity	10 ng/ml (range 3-810 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-
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)