

E. coli host cell protein ELISA kit

Sensitive (30 ng/ml) ELISA for the quantitative determination of host cell protein contamination in bulk products expressed in *E. coli* expression systems.

The *E. coli* host cell protein ELISA kit is a complete, colorimetric, immunometric immunoassay kit for the quantitative determination of *E. coli* host cell protein contamination with results in 3 hours

Ordering Information

[Order Online »](#)

ENZ-KIT127-0001	96 wells
-----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of *E. coli* host cell protein, detecting as little as 30 ng/ml
- High throughput format with results in 3 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name E. coli HCP ELISA kit

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of contamination with *E.coli* 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 30 ng/ml (range 3-810 ng/ml)

Species Reactivity E. coli

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-
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)