## 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

eu@enzolifesciences.com