

E. coli HCP ELISA Kit (host cell protein) ab240997

Overview

Product name	E. coli HCP ELISA Kit (host cell protein)
Detection method	Colorimetric
Assay type	Sandwich (quantitative)
Sensitivity	30 ng/ml
Range	3 ng/ml - 810 ng/ml
Assay time	3h 00m
Assay duration	Multiple steps standard assay
Product overview	<p><i>E. coli</i> HCP ELISA Kit (host cell protein) (ab240997) is designed to quantitatively measure host cell protein (HCP) contamination in bulk products expressed in <i>E. coli</i> expression systems.</p>

The provided 96-well plate has been pre-coated with Rabbit anti-*E. coli* HCP antibodies to capture *E. coli* proteins from biological samples. Following an incubation to allow capture of the *E. coli* protein by the antibodies on the plate, a second anti- *E. coli* HCP antibody, conjugated with biotin, is added and incubated to allow it to bind to the captured *E. coli* proteins. After incubation, the plate is washed and a Streptavidin-HRP conjugate is added and incubated for 30 min. The Streptavidin-HRP conjugate will be captured by any biotin labeled antibody bound to the plate. Following a wash step to remove unbound conjugate, TMB substrate is added and is converted by the captured HRP to a colored product in proportion to the amount of HCP bound to the plate. After a short incubation to allow color development, the reaction is stopped and the intensity of the generated color is detected in a spectrophotometer plate reader capable of measuring 450nm wavelength.

Platform	Pre-coated microplate (12 x 8 well strips)
----------	--

Properties

Storage instructions	Store at +4°C. Please refer to protocols.
----------------------	---

Components	1 x 96 tests
<i>E. coli</i> Protein Standard	1 x 600µl
10X PBST	1 x 30ml

Components	1 x 96 tests
5X Dilution Buffer	1 x 15ml
Coated Clear 96 Well Plates	1 unit
Plate Sealer	1 unit
Reporting Antibody	1 x 150µl
Stop Solution	1 x 15ml
Streptavidin-HRP Conjugate	1 x 400µl
TMB Substrate	1 x 15ml

**Relevance** Escherichia coli is a bacterium that is commonly found in the lower intestine of warm-blooded animals.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

#### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors