

Product datasheet

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