

Product datasheet

Human IDO ELISA Kit ab245710

Recombinant SimpleStep ELISA®

★★★★★ [1 Abreviews](#) [2 References](#) [4 Images](#)

Overview

Product name Human IDO ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
cell extract	21			3%

Inter-assay

Sample	n	Mean	SD	CV%
cell extract	3			3.6%

Sample type Cell culture extracts

Assay type Sandwich (quantitative)

Sensitivity 13 pg/ml

Range 78.125 pg/ml - 5000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture extracts	93	81% - 100%

Assay time 1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Human

Product overview

Human IDO ELISA Kit (ab245710) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IDO protein in cell culture extracts. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human IDO with 13 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This

approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

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	10 x 96 tests
10X Human IDO Capture Antibody	1 x 600µl	10 x 600µl
10X Human IDO Detector Antibody	1 x 600µl	10 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml	10 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml	2 x 50ml
Antibody Diluent 5BI	1 x 6ml	10 x 6ml
Human IDO Lyophilized Recombinant Protein	2 vials	10 x 2 vials
Plate Seals	1 unit	10 units
Sample Diluent NS (ab193972)	1 x 12ml	0 x 0ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	10 units
Stop Solution	1 x 12ml	1 x 120ml
TMB Development Solution	1 x 12ml	1 x 120ml

Function Catalyzes the cleavage of the pyrrol ring of tryptophan and incorporates both atoms of a molecule of oxygen.

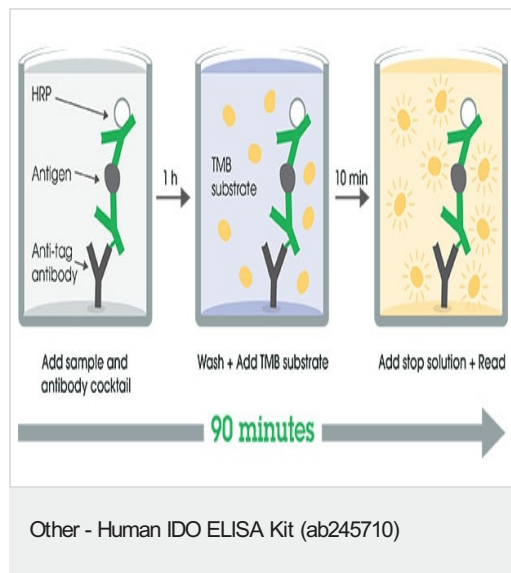
Pathway

Amino-acid degradation; L-tryptophan degradation via kynurenine pathway; L-kynurenine from L-tryptophan: step 1/2.

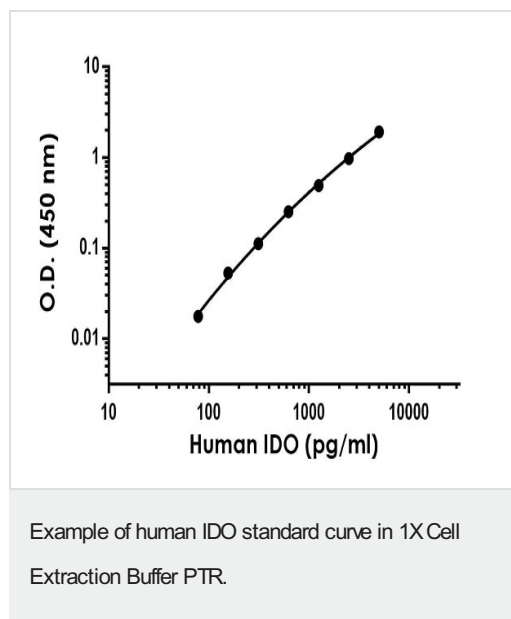
Sequence similarities

Belongs to the indoleamine 2,3-dioxygenase family.

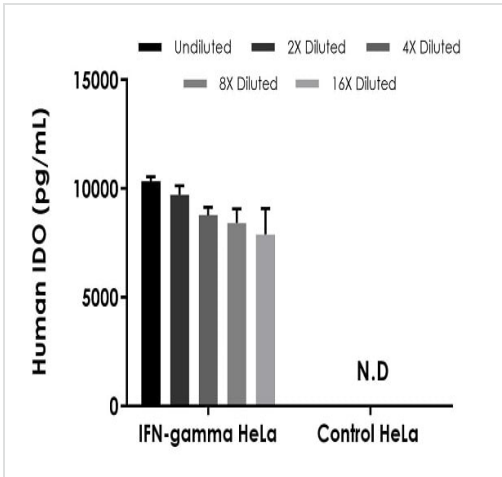
Images



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



The IDO standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native IDO in IFN stimulated HeLa cell extracts based on a 20 µg/mL extract load.

The concentrations of IDO were measured in duplicate and interpolated from the IDO standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IDO concentration was determined to be 9033 pg/mL. IDO was not detected in naive HeLa cells with a 20 µg/mL extract load.

Powered by recombinant antibodies

Research with confidence
Consistent and reproducible results

Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Human IDO ELISA Kit
(ab245710)

To learn more about the advantages of recombinant antibodies see [here](#).

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