

## Product datasheet

# Human Colipase ELISA Kit ab230937

SimpleStep ELISA<sup>®</sup>

[6 Images](#)

### Overview

**Product name** Human Colipase ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
EDTA plasma	5			4.9%

Inter-assay

Sample	n	Mean	SD	CV%
EDTA plasma	3			11.1%

**Sample type**

Cell culture supernatant, Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma, Tissue Homogenate

**Assay type**

Sandwich (quantitative)

**Sensitivity**

3 pg/ml

**Range**

5 pg/ml - 300 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	103	102% - 104%
Urine	90	86% - 94%
Serum	98	97% - 99%
Hep Plasma	108	101% - 121%
EDTA Plasma	111	104% - 124%

Sample type	Average %	Range
Cit plasma	111	104% - 120%
Tissue Homogenate	102	100% - 103%

**Assay time**

1h 30m

**Assay duration**

One step assay

**Species reactivity**

**Reacts with:** Human

**Does not react with:** Cow

**Product overview**

Colipase *in vitro* SimpleStep ELISA<sup>®</sup> (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Colipase protein in humanserum, plasma, urine, cell culture supernatant, and cell culture extract samples.

The SimpleStep ELISA<sup>®</sup> employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

Colipase is a small protein used to anchor pancreatic lipase, thereby increasing its hydrophobicity and preventing lipase from elution by bile salts.

**Platform**

Pre-coated microplate (12 x 8 well strips)

**Properties**

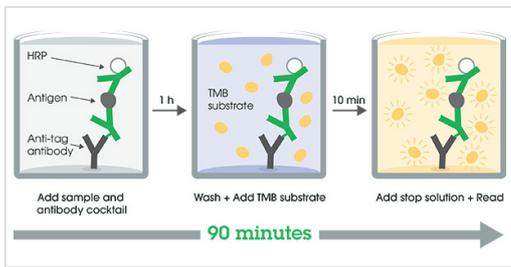
**Storage instructions**

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Human Colipase Capture Antibody	1 x 600µl
10X Human Colipase Detector Antibody	1 x 600µl
10X Wash Buffer PT ( <a href="#">ab206977</a> )	1 x 20ml
50X Cell Extraction Enhancer Solution ( <a href="#">ab193971</a> )	1 x 1ml
5X Cell Extraction Buffer PTR ( <a href="#">ab193970</a> )	1 x 10ml
Antibody Diluent 4BI	1 x 6ml
Human Colipase Lyophilized Recombinant Protein	2 vials

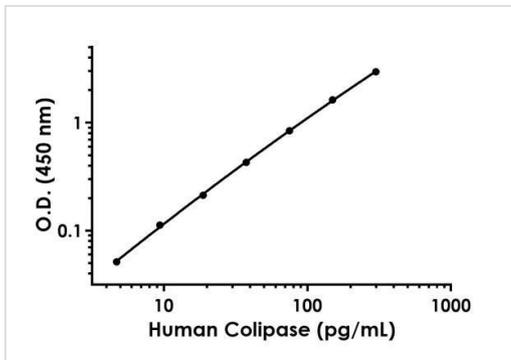
Components	1 x 96 tests
Plate Seals	1 unit
Sample Diluent NS	1 x 100ml
SimpleStep Pre-Coated 96-Well Microplate ( <a href="#">ab206978</a> )	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

**Images**



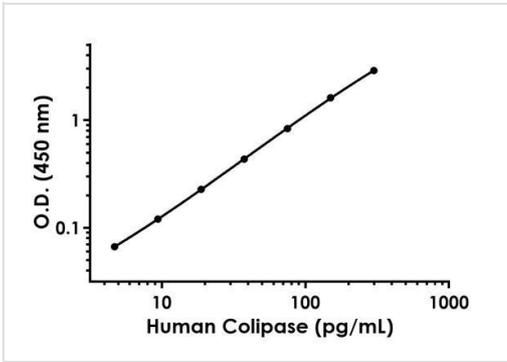
Other - Human Colipase ELISA Kit ([ab230937](#))

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.



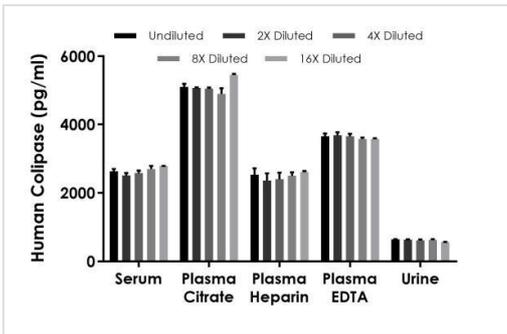
Example of human Colipase standard curve in Sample Diluent NS

Background-subtracted data values (mean +/- SD) are graphed.



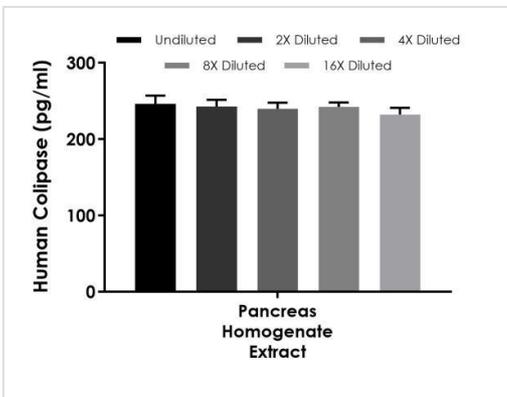
Example of human Colipase standard curve in 1X Cell Extraction Buffer PTR

Background-subtracted data values (mean +/- SD) are graphed.



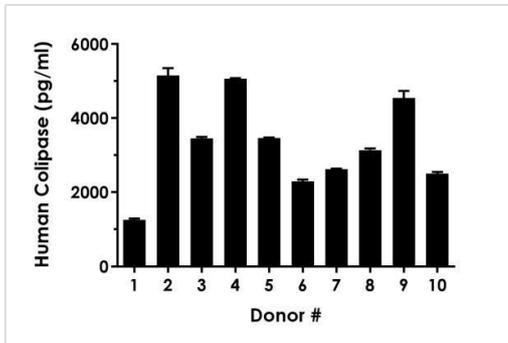
Interpolated concentrations of native Colipase in human serum, plasma, and urine samples

The concentrations of Colipase were measured in duplicates, interpolated from the Colipase standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 10%, plasma (citrate) 3.1%, plasma (heparin) 10%, plasma (EDTA) 6.3%, and urine 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Colipase concentration was determined to be 2684 pg/mL in serum, 5049 pg/mL in plasma (citrate), 2526 pg/ml in plasma (heparin), 3666 pg/ml in plasma (EDTA), and 613 pg/mL in urine.



Interpolated concentrations of native Colipase in human pancreas homogenate extract based on a 0.04 µg/mL extract load

The concentrations of Colipase were measured in duplicate and interpolated from the Colipase standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Colipase concentration was determined to be 243 pg/mL in human pancreas homogenate extract at this loading.



Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Colipase concentration was determined to be 3348 pg/mL with a range of 1229 – 5295 pg/mL.

Serum from ten individual healthy human male donors was measured in duplicate

**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