

# Human Cripto ELISA Kit ab245719

Recombinant SimpleStep ELISA<sup>®</sup>

[5 Images](#)

### Overview

**Product name** Human Cripto ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
protein	8			1.7%

**Sample type**

Cell culture supernatant, Serum, EDTA Plasma, Cit plasma

**Assay type**

Sandwich (quantitative)

**Sensitivity**

9.6 pg/ml

**Range**

26.56 pg/ml - 1700 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95	% - %
Serum	86	% - %
EDTA Plasma	86	% - %
Cit plasma	86	% - %

**Assay time**

1h 30m

**Assay duration**

One step assay

**Species reactivity**

**Reacts with:** Human

**Product overview**

Human Cripto ELISA Kit (ab245719) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Cripto protein in cell culture supernatant, cit plasma, edta plasma, and serum. It uses our proprietary SimpleStep ELISA<sup>®</sup> technology. Quantitate Human Cripto with 9.6 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.

**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 Cripto Capture Antibody	1 x 600µl
10X Human Cripto Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Human Cripto Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

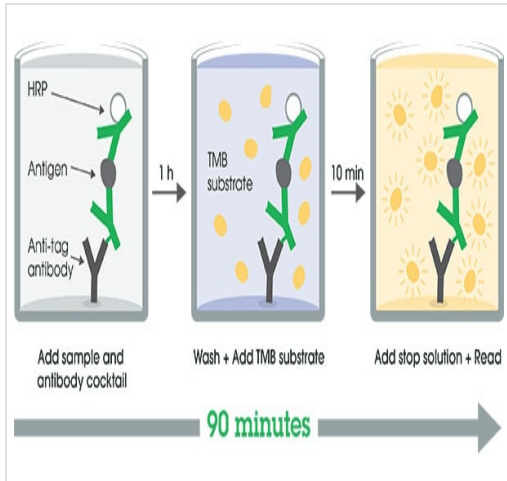
**Function** Could play a role in the determination of the epiblastic cells that subsequently give rise to the mesoderm.

**Tissue specificity** Preferentially expressed in gastric and colorectal carcinomas than in their normal counterparts.

**Sequence similarities** Contains 1 EGF-like domain.

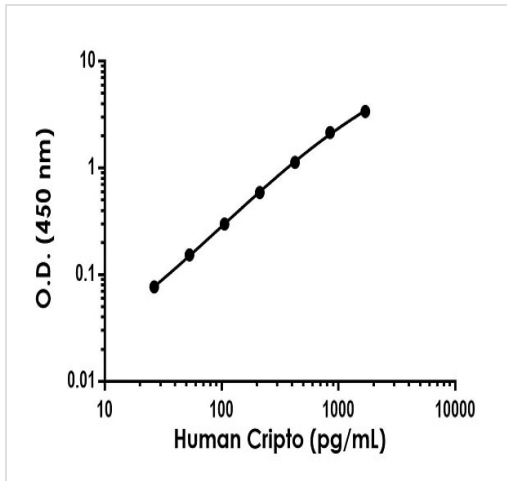
**Cellular localization** Cell membrane.

## Images



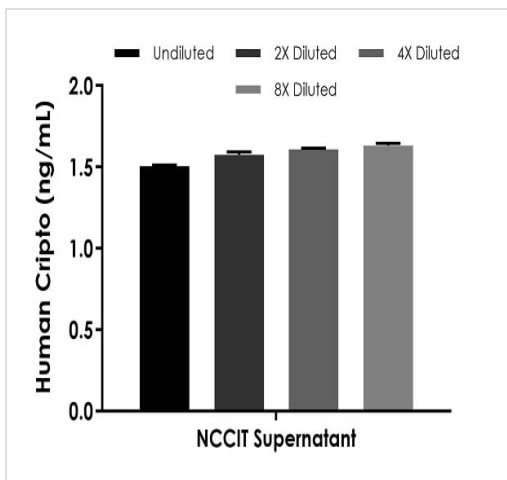
Other - Human Cripto ELISA Kit (ab245719)

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.



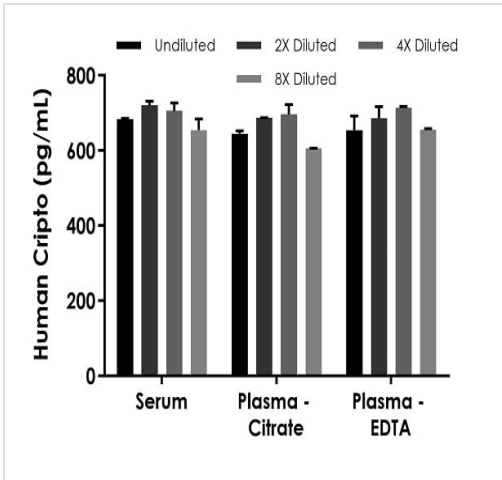
Example of human Cripto standard curve in Sample Diluent NS.

The Cripto 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 Cripto in human NCCIT supernatant samples.

The concentrations of Cripto were measured in duplicates, interpolated from the Cripto standard curves and corrected for sample dilution. Undiluted samples are as follows: NCCIT supernatant 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Cripto concentration was determined to be 1.58 ng/mL in NCCIT supernatant.



The concentrations of Cripto were measured in duplicates and interpolated from the Cripto standard curves. Undiluted samples are as follows: serum 12.5%, plasma (citrate) 12.5%, and plasma (EDTA) 12.5%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).

Interpolated concentrations of spiked Cripto in human serum and plasma samples.

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 Cripto ELISA Kit  
(ab245719)

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