

### Human LECT2 ELISA Kit ab269554

リコンビナント SimpleStep ELISA

画像数 7

#### 製品の概要

製品名 Human LECT2 ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Serum	8			4.7%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Serum	3			7.4%

サンプルの種類 Serum, Tissue Extracts, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 0.706 ng/ml

検出範囲 1.56 ng/ml - 100 ng/ml

添加回収試験 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Serum	94	91% - 96%
Tissue Extracts	120	120% - 120%
Hep Plasma	101	99% - 103%
EDTA Plasma	115	113% - 118%
Cit plasma	114	111% - 117%

全工程の試験時間 1h 30m

ステップ One step assay

## 種交差性

交差種: Human

非交差種: Mouse, Rat, Cow

## 製品の概要

Human LECT2 ELISA Kit (ab269554) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of LECT2 protein in hep plasma, serum, cit plasma, tissue extracts, and edta plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human LECT2 with 0.706 ng/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.

## 特記事項

Leukocyte cell derived chemotaxin 2 (LECT2) is a highly conserved protein that functions with major development in liver regeneration, immune modulation, bone growth, neuronal development, cancer, and several more in pathophysiology. Studies have shown LECT2 to be a positive regulator of chondrocyte proliferation. Moreover, the liver is considered to be the source of LECT2 due to exceptionally high levels of mRNA. However, expression in these cells is extremely low or undetectable suggesting that hepatocytes secrete LECT2 instantly after production.

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.

## 試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

## 製品の特性

### 保存方法

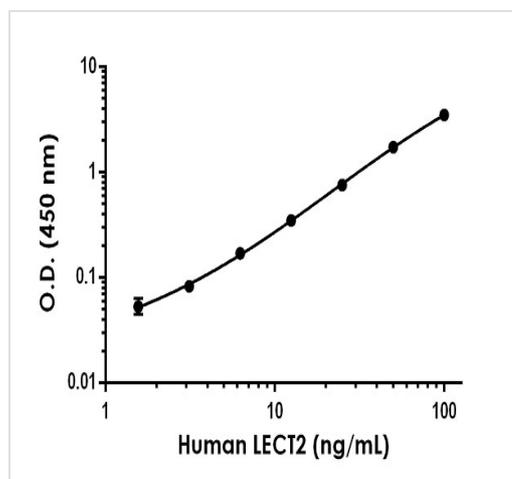
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Human LECT2 Capture Antibody	1 x 600µl
10X Human LECT2 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml

内容	1 x 96 tests
Antibody Diluent 4BI	1 x 6ml
Human LECT2 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

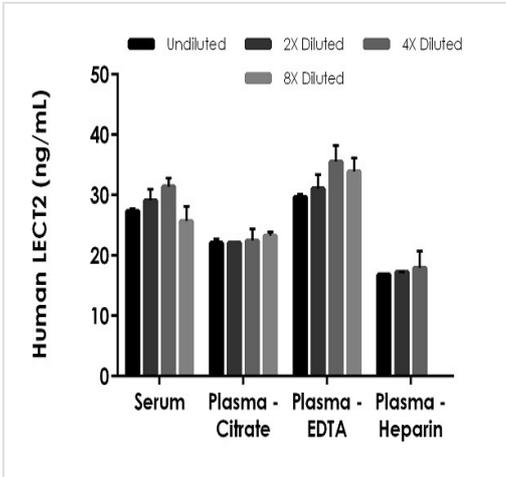
<b>機能</b>	Has a neutrophil chemotactic activity. Also a positive regulator of chondrocyte proliferation.
<b>組織特異性</b>	Highly expressed in adult and fetal liver and weakly in testis. Not expressed in bone marrow.
<b>配列類似性</b>	Belongs to the LECT2/MIM-1 family.
<b>細胞内局在</b>	Cytoplasm. Secreted.

## 画像



The LECT2 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.

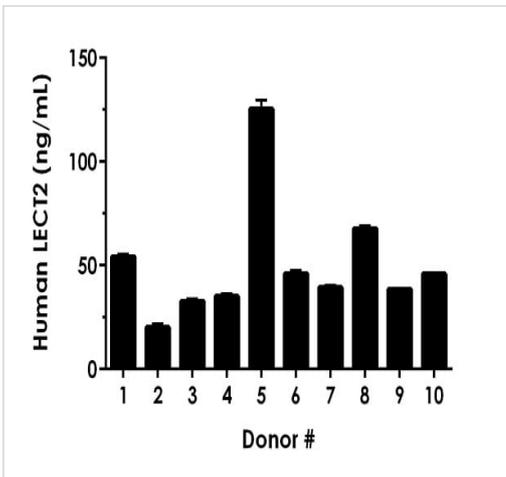
Example of human LECT2 standard curve in Sample Diluent NS.



Interpolated concentrations of native LECT2 in human serum and plasma samples.

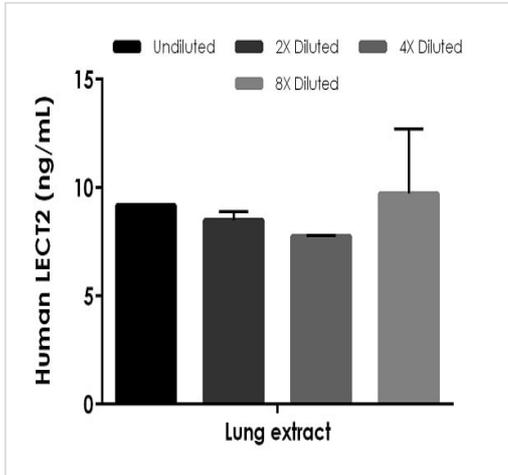
The concentrations of LECT2 were measured in duplicates, interpolated from the LECT2 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 25%, plasma-citrate 25%, plasma-EDTA 25%, and plasma-heparin 25%.

The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean LECT2 concentration was determined to be 28.4 ng/mL in serum, 22.5 ng/mL in plasma-citrate, 32.5 ng/mL in plasma-EDTA, and 13.0 ng/mL in plasma-heparin.



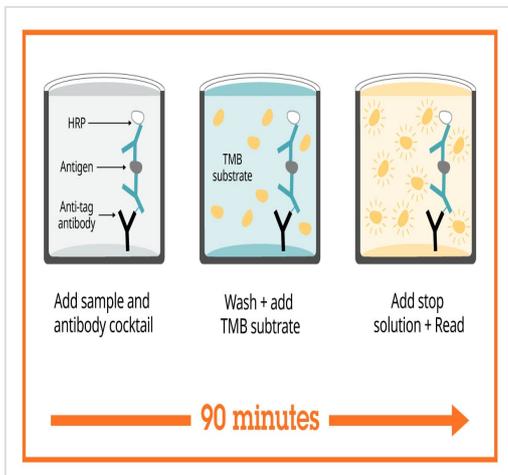
Serum from ten individual healthy human female donors were measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean LECT2 concentration was determined to be 51 ng/mL with a range of 20 - 125 ng/mL.



Interpolated concentrations of native LECT2 in human lung extract based on a 1,000 µg/mL extract load.

The concentrations of LECT2 were measured in duplicate and interpolated from the LECT2 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean LECT2 concentration was determined to be 8.8 ng/mL in lung extract.



Sandwich ELISA - Human LECT2 ELISA Kit (ab269554)

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.

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 LECT2 ELISA Kit  
(ab269554)

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

Get more done with  
SimpleStep ELISA



**Easy to use**  
Single-wash 90-minute protocol



**Flexible**  
Matched antibody pairs available



**Precision antibodies**  
High sensitivity, specificity and reproducibility



**Scalable**  
Now in 10-pack and 384-well formats

Sandwich ELISA - Human LECT2 ELISA Kit  
(ab269554)

To learn more about the advantages of SimpleStep ELISA<sup>®</sup> kits 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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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